

Agenda

17:00 Welcome

Lydia Gan

Marketing Manager,
Southeast Asia & Taiwan,
Life Sciences Solutions,
Thermo Fisher Scientific

17:05 Speeding up your cell therapy
workflow: Next-generation
immunoassay

Ryan Lim

Field Application Scientist
Immunoassay & Cell Therapy, Thermo
Fisher Scientific

17:35 When setting up qPCR reactions
become a matter of minutes

Angeline Kan

Product Manager,
Liquid Handling Solution, Thermo Fisher
Scientific

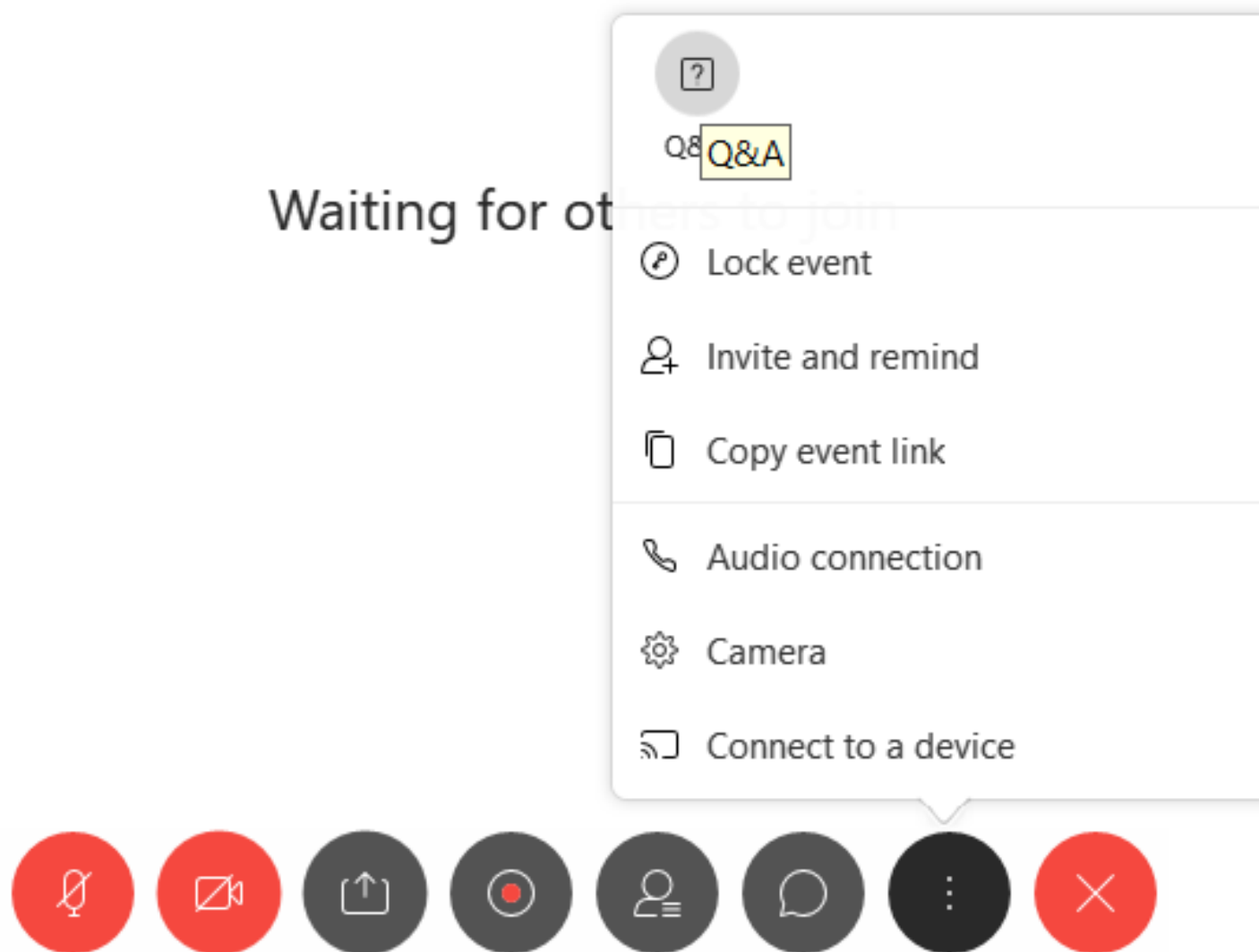
17:45 Q&A

17:55 Closing



Q&A

If you have any questions that you would like to pose, please access the Q&A section as shown here





ThermoFisher
SCIENTIFIC

Speeding up your cell therapy workflow: Next generation Immunoassay

Ryan Lim; PhD

Field Application Scientist (Immunoassay & Cell Therapy)
(Asia Pacific and Japan)

A Unique Customer Value Proposition for Pharma and Biotech

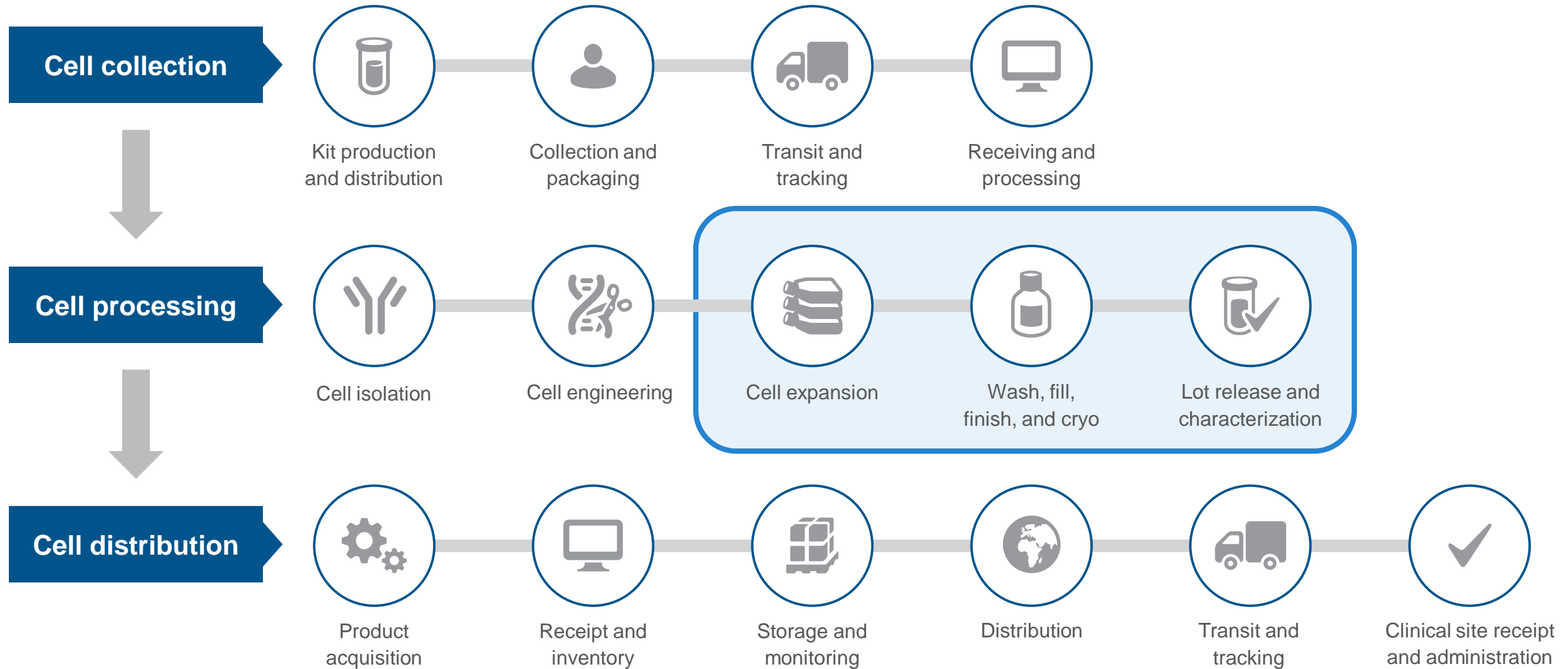
Leading life sciences offering

Leading contract development and manufacturing organization



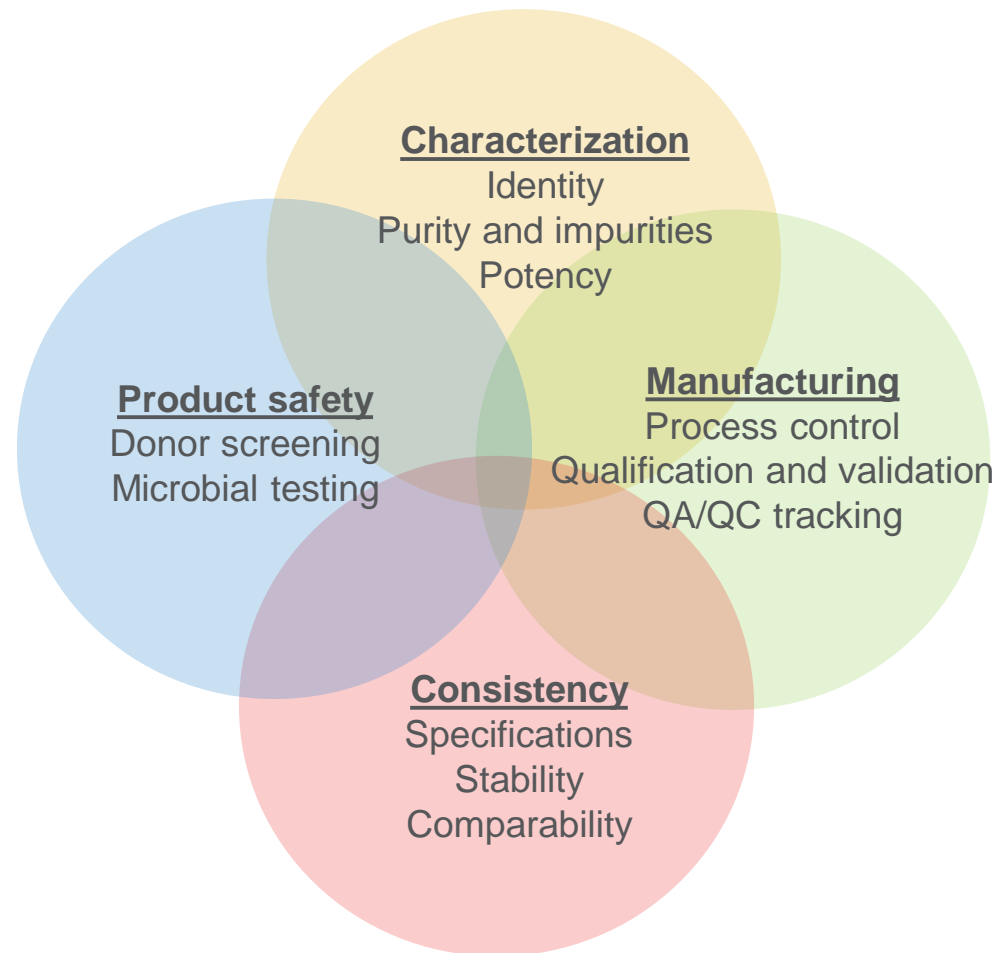
Partnering to support our customers—from discovery to molecule to medicine

Cell and Gene Therapy Workflows



Safety, Characterization, Manufacturing and Consistency are one

Critical factors that cannot be considered separately



- Product Safety
- Characterization
- Manufacturing
- Consistency

It is not possible to consider one aspect on its own.

What does the US FDA Guidelines recommend regarding ancillary reagents

PRODUCT MANUFACTURING AND CHARACTERIZATION INFORMATION

Ancillary Reagents/ Raw materials

- Reagents can affect the safety, potency, and purity of the final product, especially by introducing adventitious agents.
- Things required to declare for reagent:
 - Concentration of the reagent at the manufacturing step at which it is used; vendor/supplier; source; reagent quality; COA or cross-reference letters
- Qualification Program
 - If not FDA-approved or cleared, additional testing may be needed
 - includes safety testing (sterility, endotoxin, mycoplasma, and adventitious agents), functional analysis, purity testing, and assays (e.g., residual solvent testing) to demonstrate absence of potentially harmful substances.
 - Extent of testing depend on how the specific reagent is used in the manufacturing process.
- Determination of Removal of Reagents from Final Product
 - test the final product for residual manufacturing reagents with known or potential toxicities and describe the test procedures you use to detect residual levels of these reagents in the final product

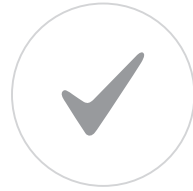
Our solutions



Cell Therapy Systems (CTS) Products—Designed for Cell Therapy Manufacturing

Gibco™ CTS™ products provide you with a proven choice so you can transition your cell therapy to the clinic with confidence.

- 20 years of cGMP manufacturing cell therapy products
- Extensive safety testing panel and FDA drug master files
- Proven use in FDA-approved cell therapies and over 100 clinical trials



cGMP-compliant manufacturing

- Manufactured in conformity with **cGMP** for medical devices, 21 CFR Part 820
- **FDA-registered** manufacturing site with an ISO 13485–certified quality management system



Testing and documentation

- Traceability documentation—including **Drug Master Files (DMFs)** and certificates of origin
- **Extensive safety testing**—including sterility, endotoxin, adventitious agent, and mycoplasma on applicable products



Proven use

- Used in **FDA-approved CAR-T therapies**^[1,2] and the first FDA-approved therapeutic cancer vaccine ^[3]
- Used in **over 100 clinical trials**



1. Tangying Lily Lu et al. (2016) A rapid cell expansion process for production of engineered the autologous CAR-T cell therapies. *Hum Gene Ther Methods*.
2. <http://thermofisher.mediaroom.com/2017-08-30-First-FDA-Approved-Cell-Therapy-for-Leukemia-Utilizes-Thermo-Fisher-Scientifics-CTS-Dynabeads-Technology>
3. Madan RA et al. (2011) Sipuleucel-T: harbinger of a new age of therapeutics for prostate cancer. *Expert Rev Vaccines*.

Cell Therapy Media and reagents --- Solutions for your cell type

Workflow Solutions Product Selection Guide

Cell type	T cell	MSC	PSC	NSC	HSC
Media/ supplements	<ul style="list-style-type: none"> Gibco™ CTS™ OpTmizer™ T-Cell Expansion SFM Gibco™ CTS™ AIM-V™ Medium Gibco™ CTS™ Immune Cell SR 	<ul style="list-style-type: none"> Gibco™ CTS™ StemPro™ MSC SFM Gibco™ StemPro™ MSC SFM XenoFree Gibco™ MesenPRO R3™ Medium 	<ul style="list-style-type: none"> Gibco™ CTS™ Essential 8™ Medium Gibco™ CTS™ KnockOut™ SR XenoFree Medium Gibco™ CTS™ KnockOut™ DMEM 	<ul style="list-style-type: none"> Gibco™ CTS™ B-27™ Supplement, XenoFree Gibco™ CTS™ KnockOut™ DMEM/F-12 Gibco™ CTS™ N-2 Supplement Gibco™ CultureOne™ Supplement (100X) 	<ul style="list-style-type: none"> Gibco™ StemPro™-34 SFM
Growth factors/cytokines	<ul style="list-style-type: none"> Gibco™ CTS™ IL2 (Interleukin 2) Recombinant Human Protein Gibco™ CTS™ IL4 (Interleukin 4) Recombinant Human Protein Gibco™ CTS™ IL7 (Interleukin 7) Recombinant Human Protein Gibco™ IL15 Recombinant Human Protein Gibco™ IL21 Recombinant Human Protein Gibco™ CTS™ GM-CSF Recombinant Human Protein Gibco™ CTS™ TNF Recombinant Human Protein 	<ul style="list-style-type: none"> Gibco™ CTS™ FGF-Basic Full Length Recombinant Human Protein Gibco™ CTS™ TGFB1 Recombinant Human Protein 	<ul style="list-style-type: none"> Gibco™ CTS™ FGF-Basic Full Length Recombinant Human Protein Gibco™ CTS™ TGFB1 Recombinant Human Protein Gibco™ CTS™ SCF (C-Kit Ligand) Recombinant Human Protein Gibco™ CTS™ FLT3 Ligand Recombinant Human Protein Gibco™ BMP4 Recombinant Human Protein Gibco™ EGF Recombinant Human Protein Gibco™ Activin A Recombinant Human Protein 	<ul style="list-style-type: none"> Gibco™ CTS™ FGF-Basic Full Length Recombinant Human Protein Gibco™ EGF Recombinant Human Protein 	<ul style="list-style-type: none"> Gibco™ CTS™ IL2 (Interleukin 2) Recombinant Human Protein Gibco™ IL3 Recombinant Human Protein Gibco™ CTS™ IL4 (Interleukin 4) Recombinant Human Protein Gibco™ IL5 Recombinant Human Protein Gibco™ CTS™ IL6 (Interleukin 6) Recombinant Human Protein Gibco™ CTS™ IL7 (Interleukin 7) Recombinant Human Protein Gibco™ CTS™ SCF (C-Kit Ligand) Recombinant Human Protein Gibco™ CTS™ GM-CSF Recombinant Human Protein Gibco™ CTS™ FLT3 Ligand Recombinant Human Protein Gibco™ TPO (Thrombopoietin), Recombinant Human Protein Gibco™ M-CSF Recombinant Human Protein
Extracellular matrix	<ul style="list-style-type: none"> NA 	<ul style="list-style-type: none"> Gibco™ CTS™ CELLstart™ Substrate 	<ul style="list-style-type: none"> Gibco™ CTS™ Vitronectin (VTN-N) Recombinant Human Protein Gibco™ CTS™ CELLstart™ Substrate Gibco™ rhLaminin-521 	<ul style="list-style-type: none"> Gibco™ CTS™ CELLstart™ Substrate 	<ul style="list-style-type: none"> NA
Reagents	<ul style="list-style-type: none"> Gibco™ CTS™ Dynabeads™ CD3/CD28 Gibco™ CTS™ DPBS Gibco™ CTS™ GlutaMAX™-I Supplement Gibco™ L-Glutamine Gibco™ LV-MAX™ Lentiviral Production System 	<ul style="list-style-type: none"> Gibco™ L-Glutamine Gibco™ StemPro™ Osteogenesis Differentiation Kit Gibco™ StemPro™ Chondrogenesis Differentiation Kit Gibco™ StemPro™ Adipogenesis Differentiation Kit Gibco™ CTS™ DPBS Gibco™ CTS™ GlutaMAX™-I Supplement Gibco™ CTS™ TrypLE™ Select Enzyme Invitrogen™ Lipofectamine™ Stem Transfection Reagent 	<ul style="list-style-type: none"> Invitrogen™ CTS™ CytoTune™-iPS 2.1 Sendai Reprogramming Kit Gibco™ PSC Cardiomyocyte Differentiation Kit Gibco™ PSC Dopaminergic Neuron Differentiation Kit Gibco™ PSC Definitive Endoderm Induction Kit Gibco™ CTS™ DPBS Gibco™ CTS™ GlutaMAX™-I Supplement Gibco™ CTS™ TrypLE™ Select Enzyme Invitrogen™ Lipofectamine™ Stem Transfection Reagent 	<ul style="list-style-type: none"> Gibco™ CTS™ DPBS Gibco™ CTS™ GlutaMAX™-I Supplement Gibco™ CTS™ TrypLE™ Select Enzyme Invitrogen™ Lipofectamine™ Stem Transfection Reagent 	<ul style="list-style-type: none"> Invitrogen™ Dynabeads™ CD34 Positive Isolation Kit Gibco™ CTS™ DPBS Gibco™ CTS™ GlutaMAX™-I Supplement
Preservation	<ul style="list-style-type: none"> Gibco™ CTS™ Synth-a-Freeze™ Medium 	<ul style="list-style-type: none"> Gibco™ CTS™ Synth-a-Freeze™ Medium 	<ul style="list-style-type: none"> Gibco™ CTS™ Synth-a-Freeze™ Medium Gibco™ PSC Cryopreservation Kit 	<ul style="list-style-type: none"> Gibco™ CTS™ Synth-a-Freeze™ Medium 	<ul style="list-style-type: none"> Gibco™ CTS™ Synth-a-Freeze™ Medium



Regarding efficacy: Potency assay

What is potency

Purpose for Potency

- Potency measurements are a necessary part of product characterization testing, comparability studies, and stability protocols, which are used to establish that a consistently manufactured product is administered during all phases of clinical investigation.
- No single test can predict clinical efficacy, hence demonstrate by “substantial evidence”
- Need to develop a release assay to measure your product’s potency.

What should be measured for Potency

Product Characterization

- Recommend that you measure a wide range of product attributes in addition to the tests used for routine lot release

Mechanism of Action

- Ideally have MOA, but due to complexity, there is some leeway, but all attempt must first be made

1. FDA Potency Tests for Cellular and Gene Therapy Products (Final Guidance for Industry): JANUARY 2011
2. International Society for Cellular Therapy perspective on immune functional assays for mesenchymal stromal cells as potency release criterion for advanced phase clinical trials. **Cytotherapy**. 2016 Feb;18(2):151-9

Potency Assay Development

- Recommended to start early as it:
 - Demonstrate product activity, quality and consistency throughout product development
 - Generate a collection of data to support specifications for lot release
 - Provide a basis for assessing manufacturing changes
 - Evaluate product stability
 - Recognize technical problems or reasons a different assay might be preferable
- If you do **not provide sufficient assurance** of potency of product lots to be used in your pivotal trial(s), your trial may be considered “deficient in design to meet its stated objectives” and may be placed on clinical hold (21 CFR 312.42(b)(2)(ii)).

Potency Assay Validation

- Product meets prescribed requirements of potency (21 CFR 601.2), which requires that you validate your potency assay (21 CFR 211.165(e)).

Accuracy

Specificity

Linearity and Range

System Suitability

Robustness

Precision (Repeatability, intermediate precision and reproducibility)

Cell Therapy cell type and the cytokine connection

MSC

MSC Immunomodulatory Potential

- By affecting B-cells, Dendritic cells, Monocytes/macrophages, NK cells and T-cells
 - Effect have been attributed to its secretion of cytokine

Table 1 Cytokines secreted by mesenchymal stem cells and the corresponding target cells²

Cytokines secreted by MSCs	Target cells
IL-10	Mph, Neu, DCs, Th1, Tregs, Tr1, tumor cells
IL-6	Neu, Mo, DCs, B, Th2, Tregs, Th17, CD8+FoxP3+
TGFβ	Mph, NK, DCs, B, T, Tregs
Chemokines	Neu, Mo, NK, Eo, Baso, DCs, Ly
CCL-2/MCP-1	Mph, EC, PL, Th2, Th17
CCL-5/RANTES	Neu, Mo, DCs, Th1, Tregs, CD8+FoxP3+
IDO	Mo, DCs, B, T, Tregs
VEGF	DCs, EC, Th1, Th17, Tregs
ICAM	T, MSCs
PGE2	Mph, Mo, NK, DCs, T, Tr1

NK

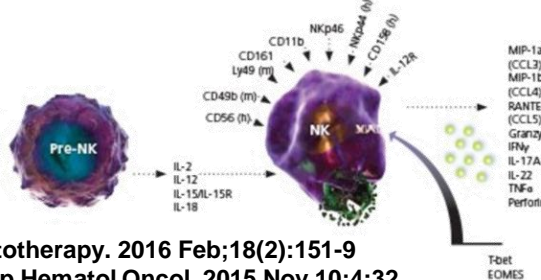
NK Cell

Cytokine-Induced
NK (CIK)-NKT

Cytokines T
Lymphocyte (CTL)

Critical Factors for efficacy to study

- Release of cytotoxic granules to mediate cell apoptosis
 - Perforin; granzyme
- Secreted cytokines and chemokines
 - IFN-gamma, TNF-alpha, IL-4, IL-5, IL-10, IL-17A, and IL-22
- Death-Ligand/death receptor-mediated cell apoptosis
 - TNF/TNF-R1; FasL/Fas; TRAIL/TRAIL-R1/TRAIL-R2



1. Chapter 20 - Cytotoxic T Cells , **Immunology for Pharmacy 2012 Pages 162-168**
2. Secretion of immunoregulatory cytokines by mesenchymal stem cells **World J Stem Cells. 2014 Nov 26; 6(5): 552–570.**
3. Immunomodulation by Mesenchymal Stem Cells (MSCs): Mechanisms of Action of Living, Apoptotic, and Dead MSCs **Front Immunol. 2019 Jun 4;10:1191**
4. Cell-based immunotherapy with cytokine-induced killer (CIK) cells: From preparation and testing to clinical application. **Hum Vaccin Immunother. 2017 Jun; 13(6): 1379–1387.**
5. International Society for Cellular Therapy perspective on immune functional assays for mesenchymal stromal cells as potency release criterion for advanced phase clinical trials. **Cytotherapy. 2016 Feb;18(2):151-9**
6. The combination of dendritic cells-cytotoxic T lymphocytes/cytokine-induced killer (DC-CTL/CIK) therapy exerts immune and clinical responses in patients with malignant tumors. **Exp Hematol Oncol. 2015 Nov 10;4:32**

Examples of reviewing cytokine signatures for CART-T immunotherapy

[Sci Transl Med](#). 2011 Aug 10;3(95):95ra73. doi: 10.1126/scitranslmed.3002842.

T cells with chimeric antigen receptors have potent antitumor effects and can establish memory in patients with advanced leukemia.

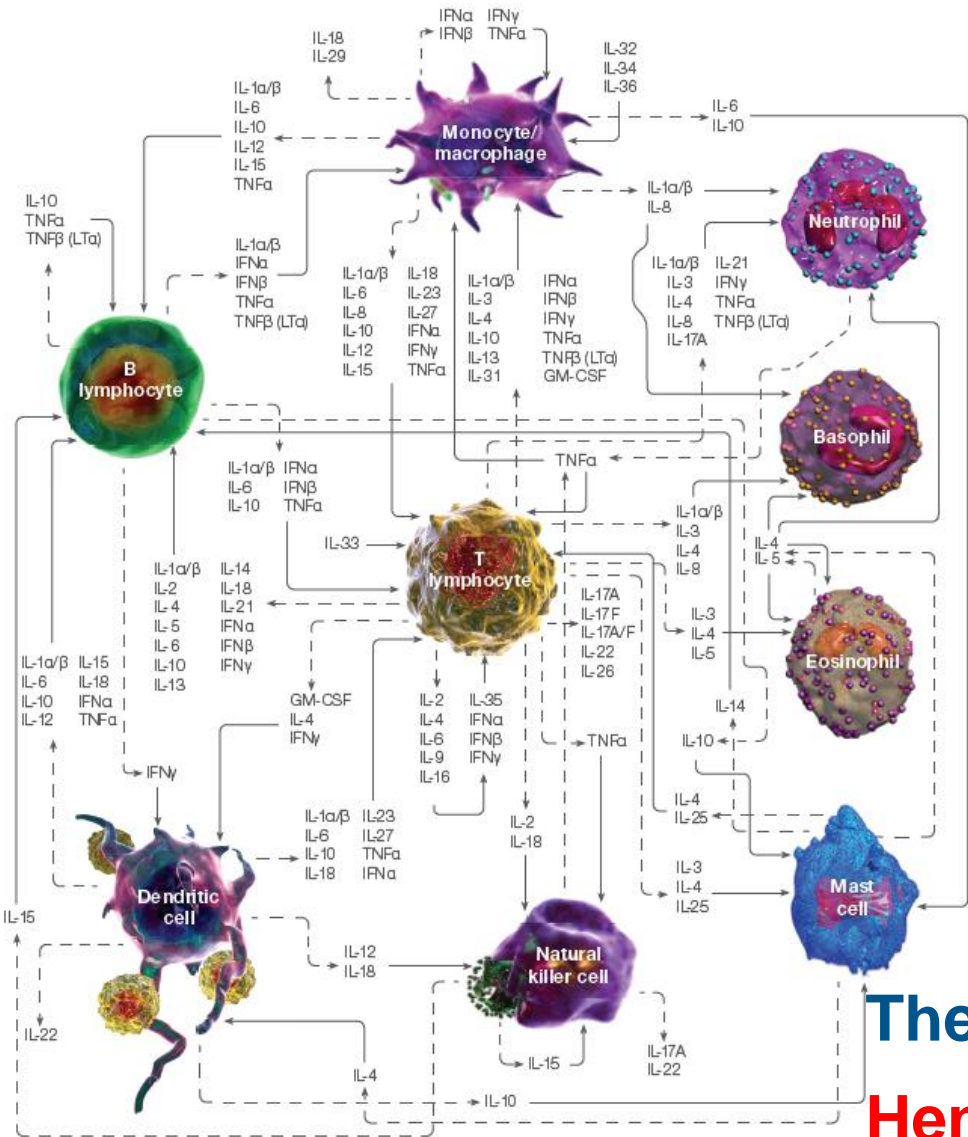
[Kalos M¹](#), [Levine BL](#), [Porter DL](#), [Katz S](#), [Grupp SA](#), [Bagg A](#), [June CH](#).

This study investigated 30 different cytokines using a 30-Plex panel over multiple timepoints. 11 cytokines showed significant changes post-transfusion.

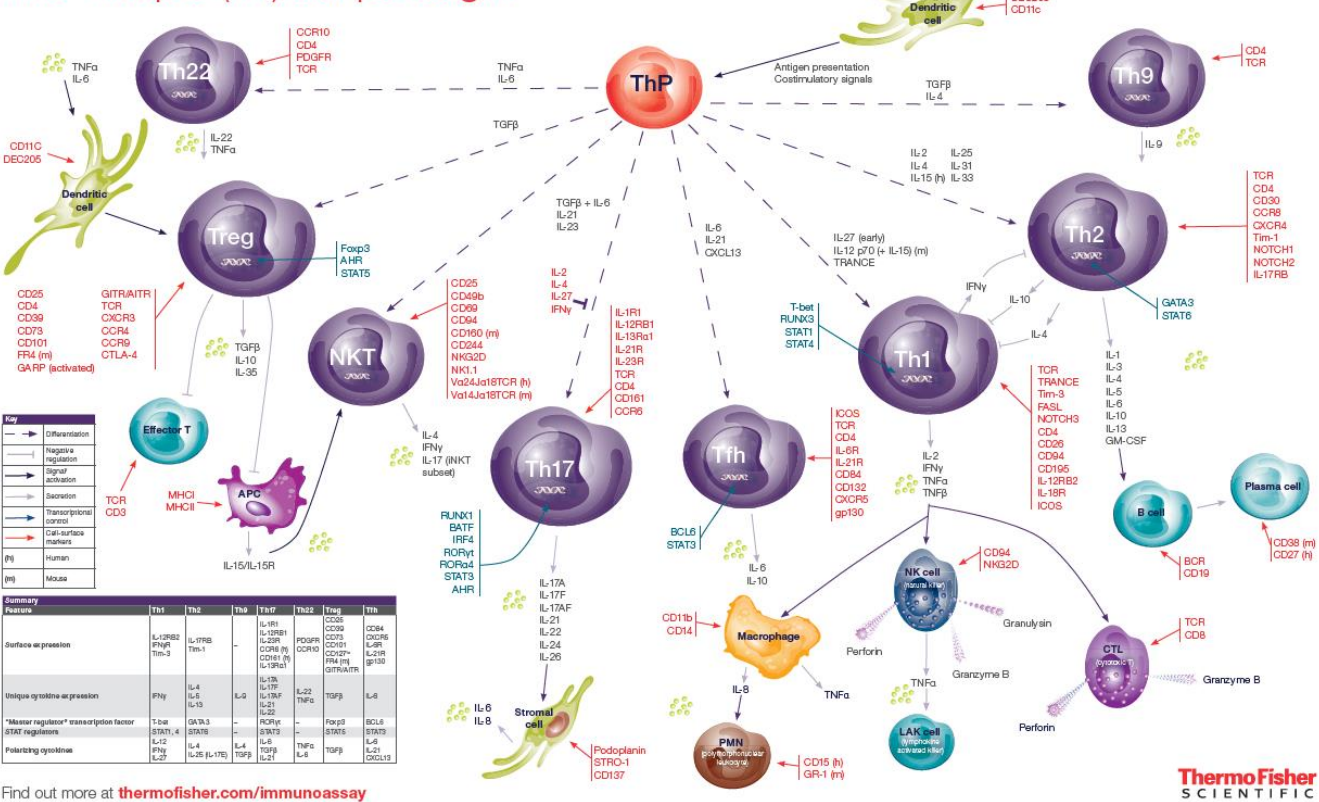
Excerpt from YESCARTA™ (axicabtagene ciloleucel) Package Insert

“After YESCARTA infusion, pharmacodynamic responses were evaluated over a 4-week interval by measuring transient elevation of cytokines, chemokines and other molecules in blood. Levels of cytokines and chemokines such as IL-6, IL-8, IL-10, IL-15, TNF- α , IFN- γ , and sIL2R α were analyzed. Peak elevation was observed within the first 14 days after infusion, and levels generally returned to baseline within 28 days. “

Cytokine Network



The T helper (Th) cell paradigm



There are more than just one cytokine in play
Hence, conventional methods would not be effective

Our Solutions



Direct and consistent: (Improving Gold-standard technology)

Challenges of ...



Consistent Data

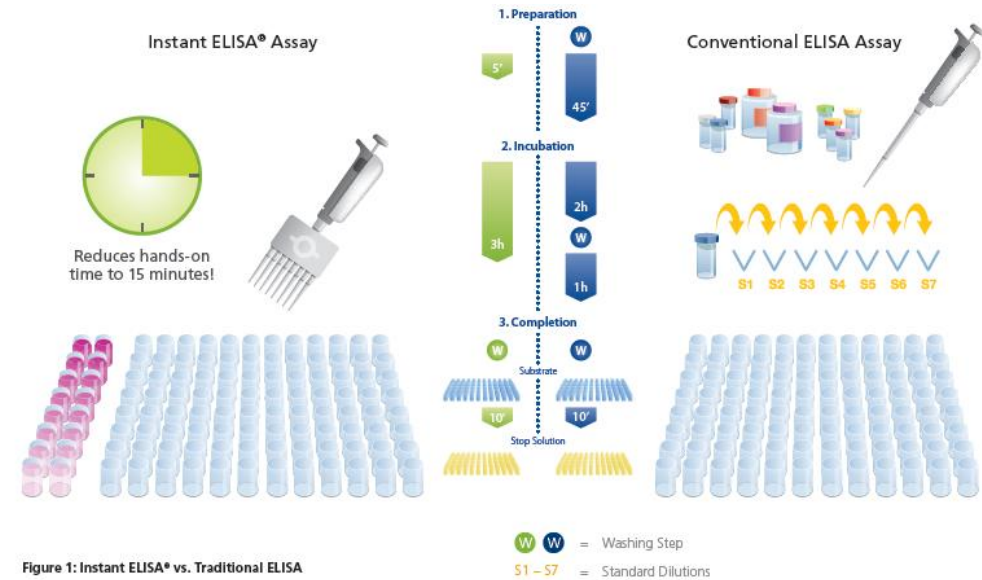
- Variability due to multiple steps
- Need direct method



Time-to-results

- Long testing cycles delays lot release

Innovation: Instant™ ELISA Kit



- Increase consistency of data by reducing human error
- Time saving due to single wash

Instant™ ELISA Kit

17 Steps for Conventional ELISA

- | | |
|----|--|
| 01 | Washing of coated plate |
| 02 | Reconstitution of standard protein |
| 03 | Addition of sample diluent to standard wells |
| 04 | Titration of standard curve |
| 05 | Addition of sample diluent |
| 06 | Sample addition |
| 07 | Dilution of biotin conjugate |
| 08 | Addition of biotin conjugate |
| 09 | Incubation |
| 10 | Preparation of streptavidin-HRP |
| 11 | Washing step |
| 12 | Addition of streptavidin-HRP |
| 13 | Incubation |
| 14 | Washing step |
| 15 | Addition of TMB substrate |
| 16 | Addition of stop solution |
| 17 | Calculation of results |

7 Steps for Instant ELISA

- | | |
|----|---------------------------|
| 01 | Rehydration of plate |
| 02 | Sample addition |
| 03 | Incubation |
| 04 | Washing step |
| 05 | Addition of TMB substrate |
| 06 | Addition of stop solution |
| 07 | Calculation of results |

Greater consistency in shorter time

- Standard curve is pre-made in lyophilized form.

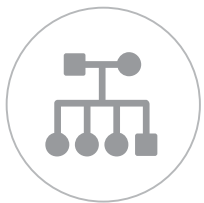
- Single wash step only

- Antibodies are pre-mixed

Lesser hands-on time = Lesser chance of error

Solving Challenges — (The solution we want)

Challenges of ...



Product complexity

- Requires multiple targets
- Validation required



Time-to-results

- Long testing cycles delays lot release



Trust in data

- Accuracy, precision, sensitivity, range and robustness
- Simple and straight forward

Innovation: ProcartaPlex (Multiplex Immunoassay)



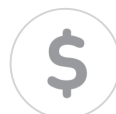
Specific like a sandwich ELISA (antibody pair)



Analysis of up to 80 analytes (50 with the Luminex® MAGPIX® system) in a single sample



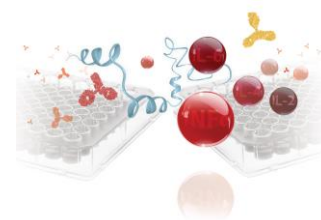
Low sample volume needed—25 µL for plasma or serum, or 50 µL for cell culture supernatant



Save cost— price per target/sample decrease with more protein targets



Save time— run equivalent of multiple ELISAs in 1 day (4-5hrs)



Secreted proteins simultaneously measured using ProcartaPlex Cytokine Assay.

Bead-Based Assay Technology

Invitrogen™ ProcartaPlex™ immunoassays apply Luminex® xMAP® technology

80 sets of superparamagnetic 6.5 µm microspheres

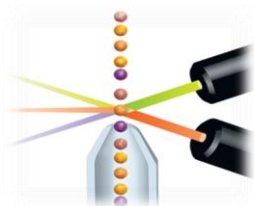
Each bead set is associated with a specific analyte—in essence, giving rise to multiple different immunoassays in a single well.

Beads are dyed using a particular bead-to-dye ratio that gives each a particular signature.

Each bead color is assigned to an analyte.

A different target can be bound to each unique bead type.

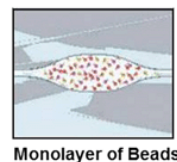
Microsphere suspension



Detects intensity of phycoerythrin (PE) with higher **quantitation**

Detects bead with higher **specificity** (analyte)

LED-based analysis



0.5 sec dwell time

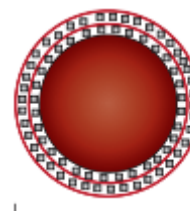
Magnetic Capture

Interrogate label with Green LED (525 nm)

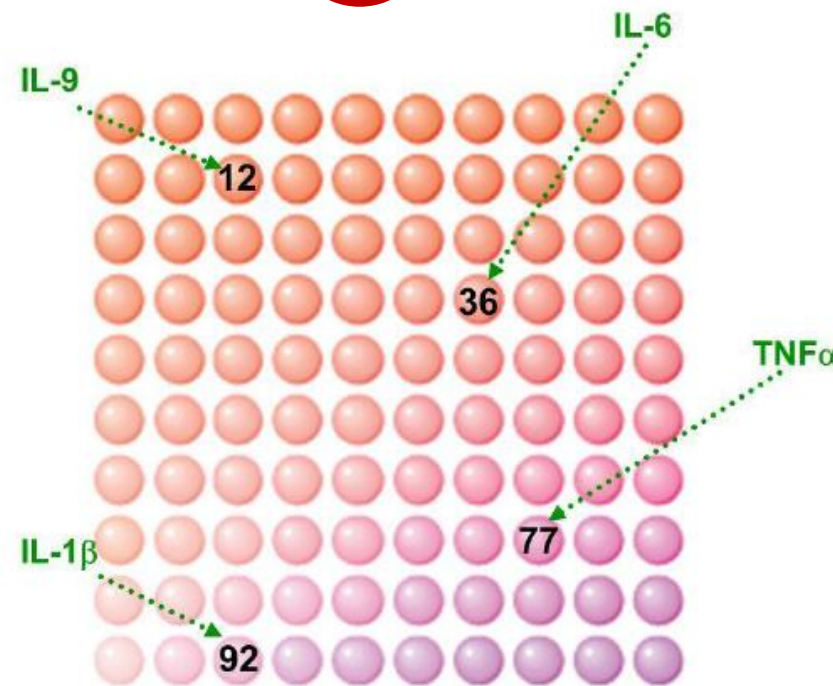
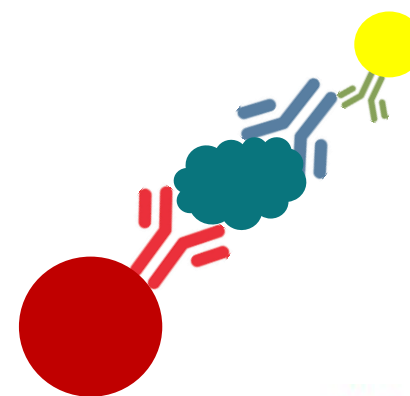
Identify and quantify with CCD imager

Interrogate label with Red LED (635 nm)

Magnetic Microsphere



6.5 microns



Procartaplex VS ELISA



IL-8



IFN-γ



TNF-α



GM-CSF

Traditional ELISA

Procartaplex



Save Sample



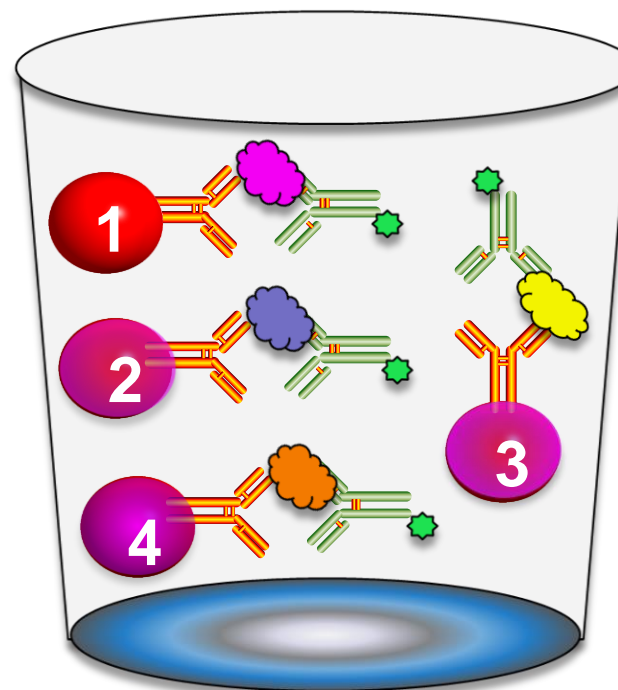
Save Money



Save Time



Increase accuracy



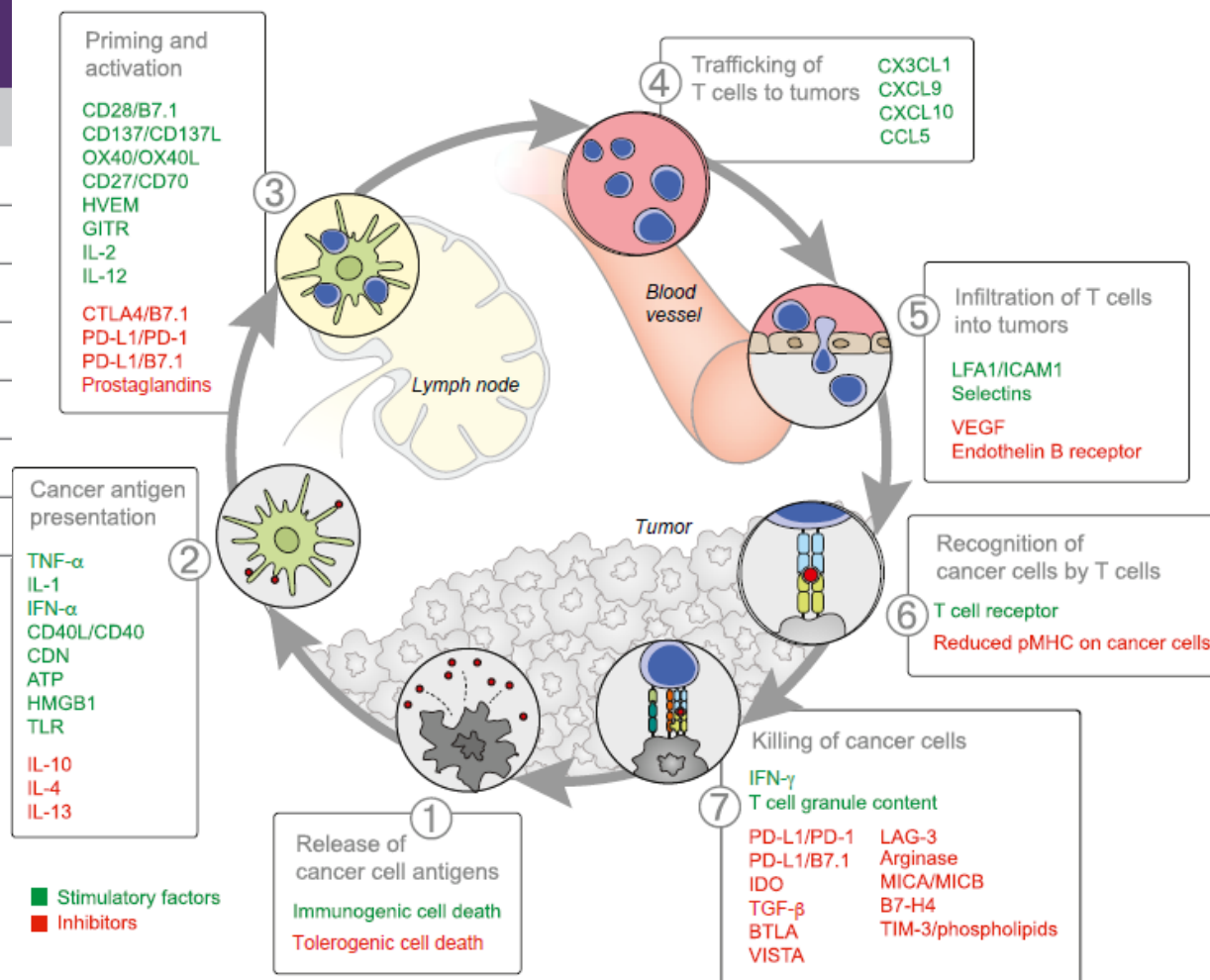
25-50uL sample/well

Immuno-panels

Immuno-Oncology Checkpoint 14-Plex Human ProcartaPlex Panel 1 Cat. No. EPX14A-15803-901

Preconfigured targets

CD28	PD-1
CD80 (B7-1)	PD-L1
CD137	PD-L2
CD27	IDO
HVEM	BTLA
GITR	LAG-3
CTLA4	TIM-3



Immuno-Oncology Checkpoint 14-Plex Human ProcartaPlex Panel 2 Cat. No. EPX140-15815-901

Preconfigured targets

Arginase-1	Siglec 7
CD112 (nectin-2)	Siglec 9
CD155 (PVR)	sMICA
CD73 (NT5E)	sMICB
CD96 (tactile)	sULBP 1
E-cadherin	sULBP 3
Perforin	sULBP 4

ProcartaPlex Custom Panel Configurator and Selection Tool

www.thermofisher.com/order/luminex/

The panel is created by single click buttons and will turn blue when selected.

Species →

1. What species of protein do you need to detect?

Human Mouse Rat Porcine Canine

Non Human Primate

Assay →

2. What panel type do you want?

ProcartaPlex High Sensitivity ProcartaPlex AB Isotyping

Coagulation Apoptosis

Sample →

3. What sample type do you want?

Serum/Plasma Other

Instrument →

4. What instrument type do you want?

MAGPIX Luminex100/200 Flexmap3D

5. Select your analytes:
Please select one or more protein targets.

4-1BBL Elafin IL-2R MCP-2 (CCL8) RANTES (CCL5)

Adiponectin Emmprin IL-3 MCP-3 (CCL7) RBP4

Analyte selection
Blue: analyte selected
Red: analyte not compatible

Your selection list:

1. 4-1BBL (43)
2. IL-2R (22)
3. IL-3 (73)
4. IL-4 (20)

Next Step

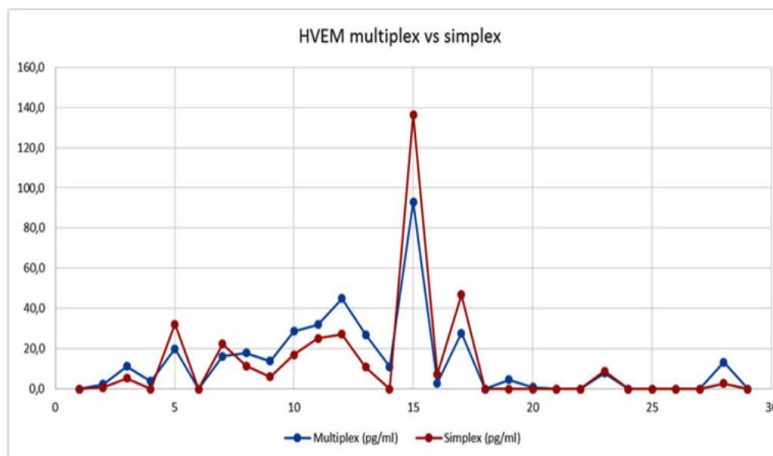
ProcartaPlex: Analyte specificity and Scalability with Consistency

Cross-reactivity testing of analytes in preconfigured panels

		Cross-reactivity in %													
		BTLA	GITR	HVEM	IDO	LAG-3	PD-1	PD-L1	PD-L2	TIM-3	CD28	CD80	CD137	CD27	CD152
Each individual recombinant protein spiked in S1 concentration.	BTLA	--	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%
	GITR	0,0%	--	0,0%	0,0%	0,0%	0,0%	2,9%	0,0%	4,4%	0,0%	0,0%	0,0%	0,0%	0,0%
	HVEM	0,0%	0,0%	--	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%
	IDO	0,0%	0,0%	0,0%	--	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%
	LAG-3	0,2%	1,3%	0,3%	0,9%	--	0,8%	1,8%	2,7%	9,8%	0,8%	0,2%	0,8%	0,9%	1,3%
	PD-1	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%
	PD-L1	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%
	PD-L2	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%
	TIM-3	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%	0,0%	0,0%	0,0%	0,0%
	CD28	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%	0,0%	0,0%	0,0%
	CD80	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%	0,0%	0,0%
	CD137	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%	0,0%
	CD27	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--	0,0%
	CD152	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	0,0%	--

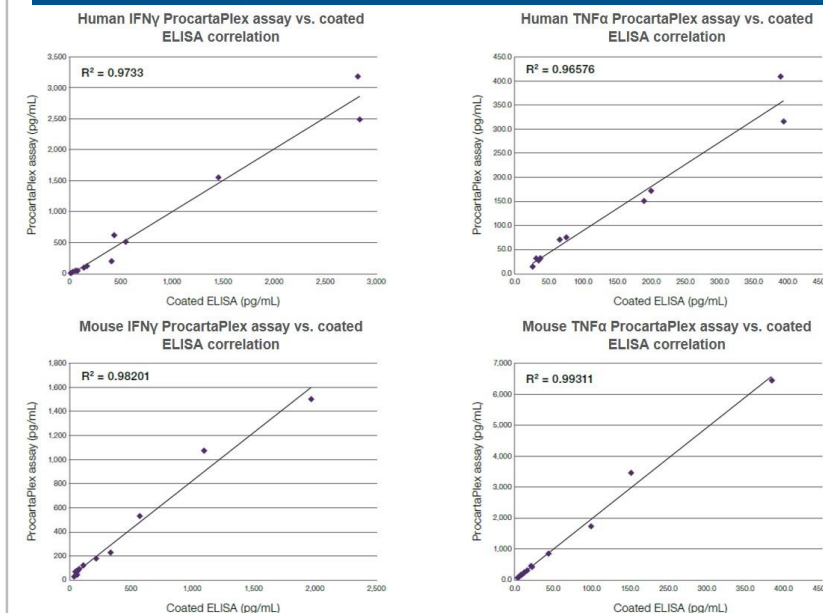
- Measuring a matrix spiked with known concentrations of a variety of different but related analytes
- Measuring a single recombinant protein within a mixture of different bead conjugates

Scalability with consistency



- We use the same antibodies between bigger panels and smaller panel
- Allow for translation of data

Data mappable to ELISA



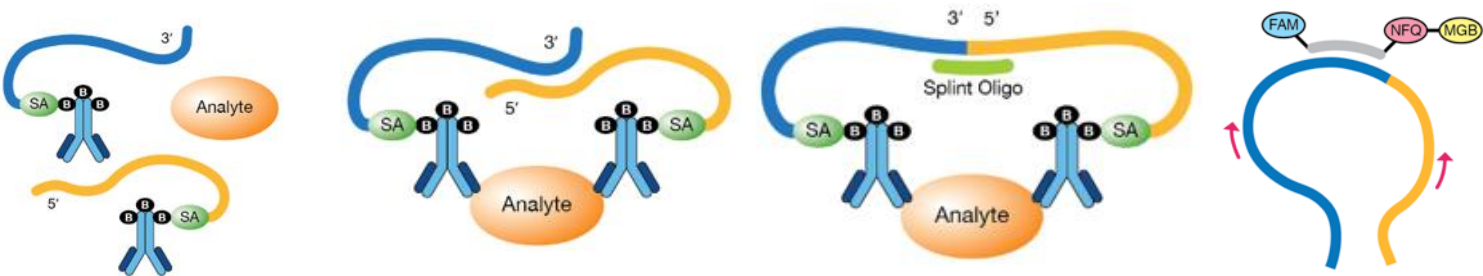
- Switch easily from ProcataPlex assays to ELISAs with reliable results
- Most of our ProcataPlex assays use the same antibody pairs as our ELISAs`

Introducing ProQuantum Immunoassays – a Leap in Next Gen Innovation

A simple, yet powerful platform for target-specific protein quantitation

High-sensitivity	Detect low levels of proteins (fg/ml) with greater sensitivity than traditional methods like ELISA
Small sample consumption	Use 2-5 µL of sample (e.g., 2 µL vs 75 µL for triplicate wells with other methods)
Fast, easy workflow	Homogeneous (no wash) assay can provide sample to answer in 2 hours
Broad dynamic range	Can reach 5 logs, minimizing sample dilutions to ensure falling within the range
No proprietary instrument to purchase	Runs on any real-time PCR instrument
Includes intuitive cloud-based software	For robust data analysis and statistical group wise comparison

Specificity of matched antibody pairs
+
Sensitivity and broad dynamic range of qPCR



Standard curve wizard

Standard curve wizard interface showing input fields for assay parameters and a visual representation of the standard curve.

Plate setup wizard

Plate setup wizard interface showing a 96-well plate layout and options for sample distribution.

Customized lab instructions

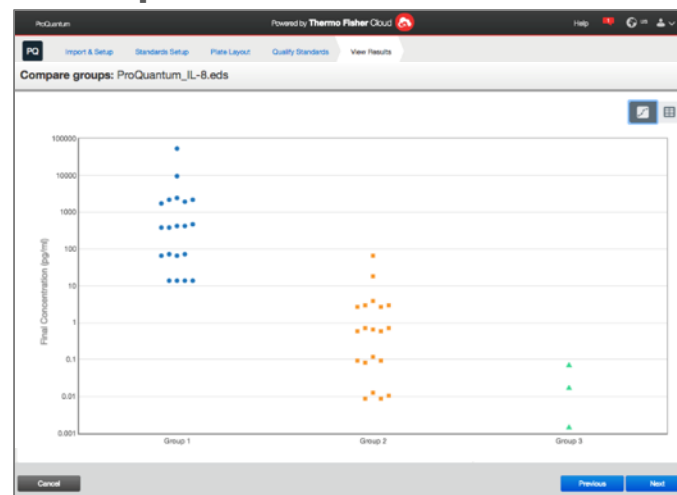
Customized lab instructions interface showing step-by-step protocols for PCR setup and reaction mixture preparation.

Sample analyte concentrations

Sample analyte concentrations interface showing a table of results and a graph of the standard curve.

Well	ID	ACT	Measured concentration	%CV	%Rec
A2,B2,C2	S2	14.233	1000.000	7.430	91.300
A3,B3,C3	S3	13.300	200.000	6.780	100.000
A4,B4,C4	S4	11.533	40.000	5.820	100.300
A5,B5,C5	S5	9.800	8.000	14.017	78.300
A6,B6,C6	S6	8.400	1.600	5.326	128.000
A7,B7,C7	S7	6.487	0.320	9.123	111.667
A8,B8,C8	S8	4.600	0.064	5.812	95.300
A9,B9,C9	S9	2.900	0.013	7.098	78.300
A10,B10,C10	S10	2.233	0.003	45.476	128.300
D1	U1	15.467	5328.225	-	-
D2	U2	14.267	991.990	-	-
D3	U3	13.167	243.243	-	-
D4	U4	11.367	42.269	-	-
D5	U5	9.867	8.552	-	-

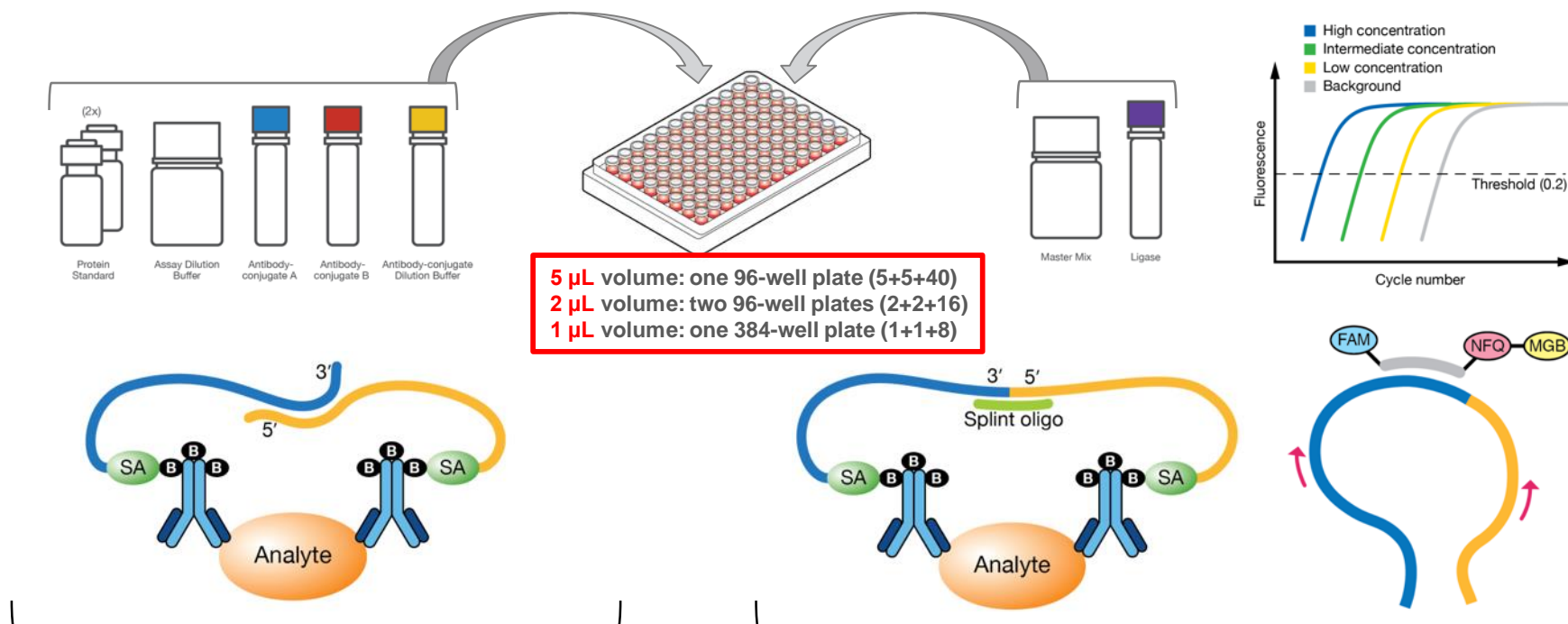
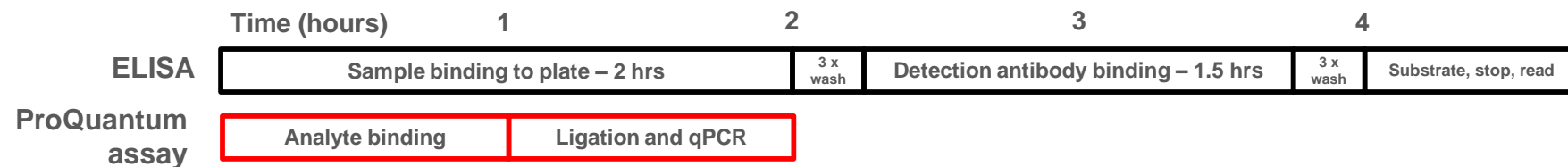
Group wise statistics



Powered by **Thermo Fisher Cloud**

- Guided assay setup is lab notebook / e-notebook friendly
- Export data GraphPad Prism software
- Instrument run monitoring app
- Compatible with **ALL** real-time PCR instruments
- Linked to online training resources

ProQuantum High-sensitivity Immunoassays are Rapid and Robust



Step 1: analyte binding

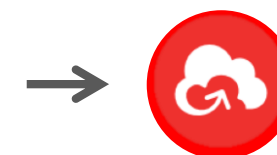
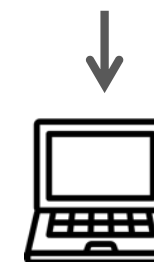
- 5 µL sample + 5 µL antibodies
- 1 hr incubation at room temperature

Step 2: ligation and qPCR readout

- Add 40 µL ligase and qPCR master mix
- In-instrument ligation and 40 cycles qPCR

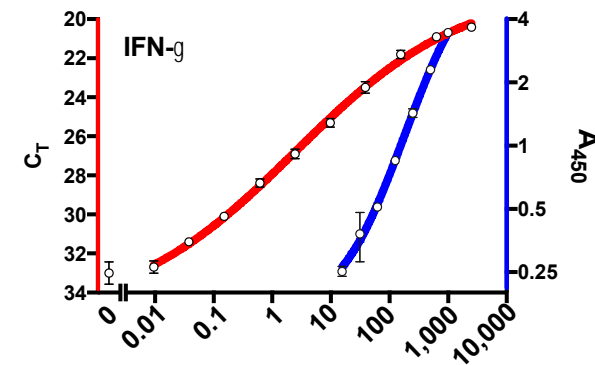
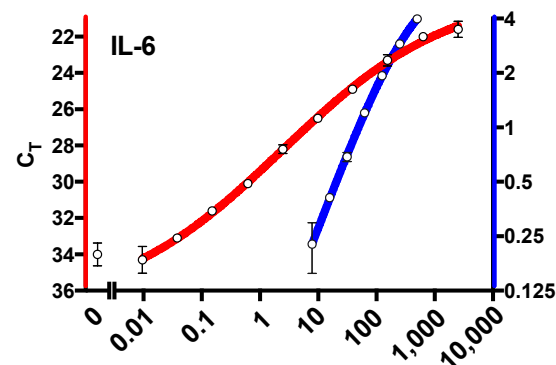
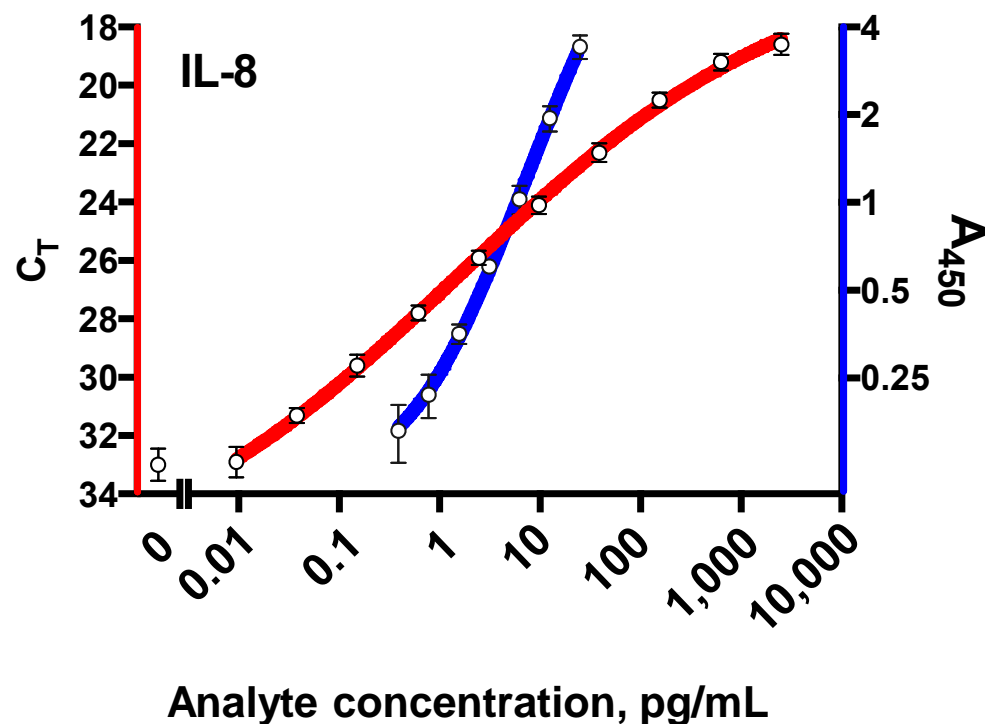


Transfer .EDS file or .CSV file with Ct values in plate map orientation



Superior Sensitivity and Dynamic Range Using Only a Fraction of Sample Volume

● ProQuantum immunoassay ■ ELISA



Assay characteristics and performance include:

- ✓ Serum spike-and-recovery: 80% - 120%
- ✓ Plasma spike-and-recovery: 70% - 130%
- ✓ Parallelism and linearity of dilution: >0.9
- ✓ Protein standard calibration to NIBSC or leading ELISA
- ✓ Inter- and intra-assay reproducibility: <15% CV of calculated concentrations (in linear range of assay)
- ✓ Analyte cross-reactivity: <10%
- ✓ Benchmarking to ELISA

Our Targets list

HUMAN					MOUSE	
Adiponectin	IFN- α	IL-18	IL-13	MIP-1 α	GM-CSF	LAP/TGF β
CEA	IFN γ	IL-21	IL-17A	MIP-1 β	Granzyme B	MCP-1
cMET	IgA	IL-4	IL-17F	SAA	IFN α	TNF α
EGF	IgE	IL-5	IP-10	Tau pT181	IFN γ	VEGF
Eotaxin	IgG4	IL-6	LAP/TGF β	Tau Total	IL-1 α	
EPO	IgM	IL-7	Leptin	TNF α	IL-2	
Flt-3 Ligand	IL-1 α	IL-8	MCP-1	TNF- β	IL-4	
Fractalkine	IL-1 β	IL-10	MCP-2	Troponin I	IL-5	
GM-CSF	IL-2	IL-12p40	MCP-3	VEGF	IL-6	
GRO- α	IL-3	IL-12p70	MDC		IL-17A	

Coming soon	
HUMAN	MOUSE
IL-2R	IFN β
IL-22	IL-22

www.thermofisher.com/ProQuantum

Discovering elevated cytokines in severe cases

Lancet. 2020 Feb 15;395(10223):497-506. doi: 10.1016/S0140-6736(20)30183-5. Epub 2020 Jan 24.

Clinical features of patients infected with 2019 novel coronavirus in Wuhan, China.

Huang C¹, Wang Y², Li X³, Ren L⁴, Zhao J⁵, Hu Y⁶, Zhang L¹, Fan G⁷, Xu J⁸, Gu X⁷, Cheng Z⁹, Yu T¹, Xia J¹, Wei Y¹, Wu W¹, Xie X¹, Yin W⁶, Li H², Liu M¹⁰, Xiao Y⁴, Gao H¹¹, Guo L⁴, Xie J⁵, Wang G¹², Jiang B³, Gao Z¹³, Jin Q⁴, Wang J¹⁴, Cao B¹⁵.

✚ Author information

Erratum in

Department of Error. [Lancet. 2020]

- Researchers observed that in patient with a severe response **IL2, IL7, IL10, GSCF, IP10, MCP1, MIP1A and TNF-α**

Lung injury from cytokine storm

Nat Rev Immunol. 2020 Apr 6. doi: 10.1038/s41577-020-0305-6. [Epub ahead of print]

In the eye of the COVID-19 cytokine storm.

Vaninov N¹.

✚ Author information

1 Sinai Immunology Review Project, Icahn School of Medicine at Mount Sinai, New York, NY, USA. sinai.immunology@gmail.com.

- Cytokine storm shown to be a cause for the lung damage

Clinical trials using hMSCs to reduce cytokine levels

- As of 20th April 2020 there is current 45 cell therapy trials for Covid-19¹
- Of which a vast majority uses hMSC¹
- Recent references

Aging and Disease

[www.aginganddisease.org](http://dx.doi.org/10.14336/AD.2020.0228)

Volume 11, Number 2, 216-228, April 2020

<http://dx.doi.org/10.14336/AD.2020.0228>

Original Article

Transplantation of ACE2⁺ Mesenchymal Stem Cells Improves the Outcome of Patients with COVID-19 Pneumonia

Zikuan Leng^{1,5,6}, Rongjia Zhu^{1,6}, Wei Hou^{1,6}, Yingmei Feng^{1,6}, Yanlei Yang¹, Qin Han², Guangliang Shan², Fanyan Meng¹, Dongshu Du¹, Shibua Wang², Junfen Fan¹, Wenjing Wang², Luchan Deng², Hongbo Shi², Hongjun Li¹, Zhongjie Hu¹, Fengchun Zhang¹, Jinning Gao¹, Hongjian Liu^{1,7}, Xiaoxia Li¹, Yangyang Zhao¹, Kan Yin¹, Xijing He², Zhengchao Gao¹, Yibin Wang¹, Bo Yang¹, Ronghua Jin^{1,8}, Ilya Stambler^{9,10,11}, Lee Wei Lim^{9,10,12}, Huanxing Su^{9,10,13}, Alexey Moskalev^{9,10,14}, Antonio Cano^{9,10,15}, Sasanka Chakrabarti¹⁶, Kyung-Jin Min^{9,10,17}, Georgina Ellison-Hughes^{9,10,18}, Calogero Caruso^{9,10,19}, Kunlin Jin^{9,10,20}, Robert Chunhua Zhao^{1,2,9,10}

¹School of Life Sciences, Shanghai University, Shanghai, China. ²Institute of Basic Medical Sciences Chinese Academy of Medical Sciences, School of Basic Medicine Peking Union Medical College, Beijing, China. ³Beijing YouAn Hospital, Capital Medical University, Beijing, China. ⁴Department of Rheumatology and Clinical Immunology, Peking Union Medical College Hospital, Chinese Academy of Medical Sciences and Peking Union Medical College, Beijing, China. ⁵Department of Orthopaedics, the First Affiliated Hospital of Zhengzhou University, Zhengzhou, China. ⁶Institute of Stem Cell and Regeneration Medicine, School of Basic Medicine, Qingdao University, Shandong, China. ⁷Department of Orthopaedics, the Second Affiliated Hospital of Xi'an Jiaotong University, Xi'an, China. ⁸Department of Neurosurgery, the First Affiliated Hospital of Zhengzhou University, Zhengzhou, China. ⁹The Executive Committee on Anti-aging and Disease Prevention in the framework of Science and Technology, Pharmacology and Medicine Themes under an Interactive Atlas along the Silk Roads, UNESCO, Paris, France. ¹⁰International Society on Aging and Disease, Fort Worth, Texas, USA. ¹¹The Geriatric Medical Center "Shmuel Harofe", Beer Yaakov, affiliated to Sackler School of Medicine, Tel-Aviv University, Tel-Aviv, Israel. ¹²School of Biomedical Sciences, Li Ka Shing Faculty of Medicine, University of Hong Kong, Hong Kong, China. ¹³Institute of Chinese Medical Science, University of Macau, Macau, China. ¹⁴Institute of Biology, Komi Science Center of Russian Academy of Sciences, Syktyvkar, Russia. ¹⁵Department of Pediatrics, Obstetrics and Gynecology, University of Valencia, Valencia, Spain. ¹⁶Maharishi Markandeshwar Deemed University, Mullana-Ambala, India. ¹⁷Department of Biological Sciences, Inha University, Incheon, South Korea. ¹⁸Faculty of Life Sciences & Medicine, King's College London, London, UK. ¹⁹Department of Biomedicine, Neuroscience and Advanced Diagnostics, University of Palermo, Palermo, Italy. ²⁰University of North Texas Health Science Center, Fort Worth, TX76107, USA.

[Received February 25, 2020; Revised February 28, 2020; Accepted February 28, 2020]

Clinical remission of a critically ill COVID-19 patient treated by human umbilical cord mesenchymal stem cells

Bing Liang^{1,*}, Junhui Chen^{2,*}, Tao Li^{3,*}, Haiying Wu^{4,*}, Wenjie Yang^{1,*}, Yanjiao Li⁵, Jianchun Li¹, Congtao Yu², Fangang Nie¹, Zhaoxia Ma², Mingxi Yang¹, Panrong Nie⁶, Yanfeng Gao^{2,7,*}, Chuanyun Qian^{4,*}, Min Hu^{2,5,8,*}

¹ Department of Critical Care Medicine, Baoshan People's Hospital, Baoshan 678000, China

² Intervention and Cell Therapy Center, Peking University Shenzhen Hospital, Shenzhen 518035, China

³ Yunnan Yasheng Medical Technology Co., Ltd., Kunming 650021, China

⁴ Emergency Department of the First Affiliated Hospital of Kunming Medical University, EICU/MICU, Kunming 650032, China

⁵ Yunnan Key Laboratory for Basic Research on Bone and Joint Diseases & Yunnan Stem Cell Translational Research Center, Kunming University, Kunming 650214, China

⁶ Department of Neonatology, Baoshan People's Hospital, Baoshan, 678000, China

⁷ School of Pharmaceutical Sciences (Shenzhen), Sun Yat-sen University, Shenzhen 517108, China

⁸ Yunnan Jici Institute for Regenerative Medicine co., Ltd., Kunming 650106, China

* These authors contribute equally.

* Correspondence to: E-mail: huminyun@163.com (Hu M.); qianchuanyun@126.com (Qian C.); yfgao@hotmail.com (Gao Y.)

1. <https://celltrials.org/>

How Thermo Fisher can partner you with MSC workflow for COVID-19

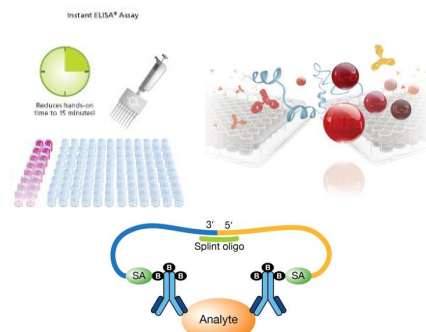
MSC Expansion



Wash and fill



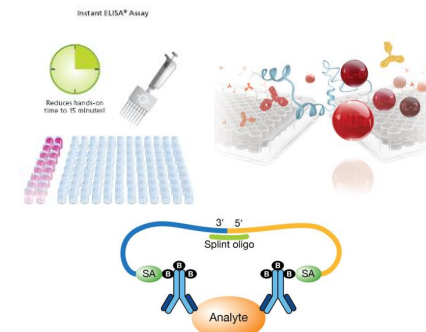
Potency testing



Purity and Safety testing



Patient severity classification



Cell culture plastics

- ❖ Nunc™ TripleFlask™ Treated Cell Culture Flasks
- ❖ Cell Factory or bioreactor

Reagents

- ❖ StemPro™ MSC SFM XenoFree
- ❖ CTS™ GlutaMAX™-I
- ❖ CTS™ CELLstart™
- ❖ CTS™ TrypLE™
- ❖ CTS™ TGFB1/FGF2 Recombinant Human Protein

- Wash, media exchange, formulate, fill
 - ❖ Rotea™ counterflow centrifugation

Cytokine testing with

- ❖ Instant ELISA
- ❖ ProcartaPlex
- ❖ ProQuantum

Safety and purity testing

- ❖ MycoSEQ™ Mycoplasma Detection Assay
- ❖ ProQuantum
- ❖ TaqMan Gene Expression Arrays
- ❖ Attune NxT for cellular analysis
- ❖ PIERCE LAL endotoxin kit

Cytokine testing with

- ❖ Instant ELISA
- ❖ ProcartaPlex
- ❖ ProQuantum

Complementary RUO and CTS Products for MSC expansion

Seamless transition from basic research to clinical applications

- Our Research Use Only (RUO) portfolio supports cost-effective basic research
- Transition to Gibco™ Cell Therapy Systems (CTS™) products as you translate into clinical research

RUO	CTS
KnockOut DMEM/F-12	CTS KnockOut DMEM/F-12
KnockOut DMEM	CTS KnockOut DMEM
GlutaMAX Supplement	CTS GlutaMAX-I Supplement
TGF-β 1 Recombinant Human Protein	CTS TGF-β 1 Recombinant Human Protein
Synth-a-Freeze Medium	CTS Synth-a-Freeze Medium
TrypLE Select Enzyme	CTS TrypLE Select Enzyme
DPBS	CTS DPBS

Cell Processing System Specifically Developed for Cell Therapy Manufacturing



Gibco™ Rotea™ Counterflow centrifuge

Key benefits

- ✓ **Process flexibility**—user-programmable software enables you to create and optimize a broad range of protocols for cell separation, washing, and concentration
- ✓ **High cell recovery and viability**—gentle fluidized bed supports low shear processing, enabling >95% cell recovery with no loss in cell viability
- ✓ **Low output volumes**—proprietary technology that can deliver as little as 5 mL of concentrate
- ✓ **Appropriate for research through commercial manufacturing**—closed single-use kit; processing speed and compact size enable seamless commercial scale-up

The content provided herein may relate to products that have not been officially released and is subject to change without notice.

Process Flexibility- A Multi-purpose System for Cell Therapy Manufacturing

Currently suitable for wash/concentration between cell therapy processing workflow steps

Cell therapy processing



Thaw cells



Cell isolation and activation



Cell engineering



Cell expansion



Fill, finish, and cryopreservation



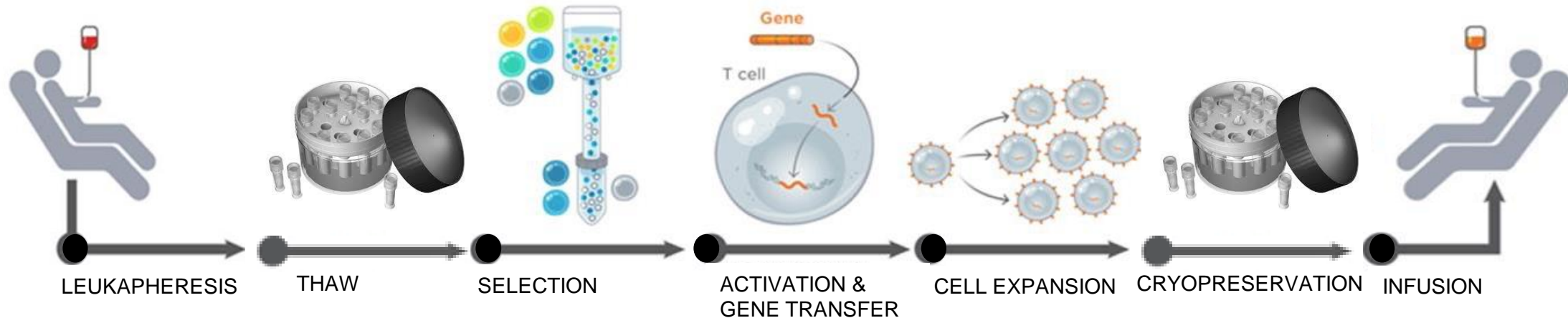
Lot release and characterization

Future applications can include isolation

Isolate PBMC from apheresis product

Isolate individual immune cell types (eg: T or NK cells)

Isolate engineered cells (eg: CAR-T, CAR-NK)



The Complete Gibco™ Rotea™ System

A complete, closed cell processing system



Instrument

- Compact
- Multipurpose
- Easily integrates into your process



Single-use kit

- Sterile, closed kit
- Flexible input and output ports



Software

- Intuitive, user-programmable interface
- Preset protocols available
- Supports GMP manufacturing

The content provided herein may relate to products that have not been officially released and is subject to change without notice.

Highly flexible instrument and software with sterile, single-use kit that scales with you from research through commercial manufacturing

Takeaway message

START IT RIGHT AND EARLY

- Using the right starting material and reagents to prevent future issues
- Consider potency early, know what you want to test

CYTOKINE EXPRESSION IS CRITICAL FOR MULTIPLE ASPECTS

- The need to measure and establish cytokine signature for potency and safety

MAINTAIN CONSISTENCY WITH A CHEAPER, SHORTER AND FASTER WORKFLOW

- Shorter work flow, higher throughput





Thermo Fisher Scientific and its affiliates are not endorsing, recommending, or promoting any use or application of products presented by third parties during this event. Information and materials presented or provided by third parties are provided as-is and without warranty of any kind, including regarding intellectual property rights and reported results. Parties presenting images, text, and materials represent they have the right to do so.

Intended use of the products mentioned in this presentation vary. For specific intended use statements please refer to the product label. © 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

The world leader in serving science

Q&A

