

Validation & Assay Performance Summary



CellSensor[®] :Myc-*bla* HCT116 Cell Line

Cat. no. K1467

CellSensor[®] Cell-Based Assay Validation Packet

This cell-based assay has been thoroughly tested and validated by Invitrogen and is suitable for immediate use in a screening application. The following information illustrates the high level of assay testing completed and the validation of assay performance under optimized conditions.

Pathway Description

Increased wild-type MYC expression occurs frequently in human cancers. Myc up-regulation occurs as a consequence of activation of one or more signaling pathways that induce MYC expression and function as a regulator of gene transcription. These include MAPK, PI3K and Wnt- β -catenin pathways. The target genes regulated by Myc are involved in the many biological activities attributed to Myc, including growth, transformation, proliferation and angiogenesis.

HCT116 is a colon cancer cell line which expresses a mutated form of β -catenin. This form of β -catenin leads to the accumulation of β -catenin and constitutive activation of downstream genes such as MYC.

Cell Line Description

CellSensor[®] Myc-*bla* HCT116 contains a beta-lactamase reporter gene under the control of Myc binding sequences. The construct was transduced into HCT116 cells by lentivirus. This cell line is a clonal population isolated by flow cytometry. It has been validated for cell plating density and DMSO tolerance. The signaling pathway has been validated using RNAi against c-MYC and ICG-001, an inhibitor of the wnt- β -catenin pathway. The expression of the mutated version of β -catenin in HCT116 cells results in the constitutive activation of beta-lactamase in this CellSensor[®] line, which can be knocked down by Myc RNAi and ICG-001.

Validation Summary

Testing and validation of this assay was evaluated in a 384-well format using LiveBLazer™-FRET B/G Substrate.

1. Primary agonist dose response under optimized conditions (n=3)

Z'-Factor	= 0.66
Response Ratio	= 3.4
Recommended cell no. cells/well	= 8000
Recommended [DMSO]	= 0.5-1%
Recommended compound incubation time	= 24 hours

2. Alternate Stimuli

n.a.

3. Stealth™ RNAi Testing

See below

4. Small molecule inhibitor Testing

See below

5. Cell culture and maintenance

See Cell Culture and Maintenance Section and Table 1

Assay Testing Summary

6. Assay performance with variable cell number

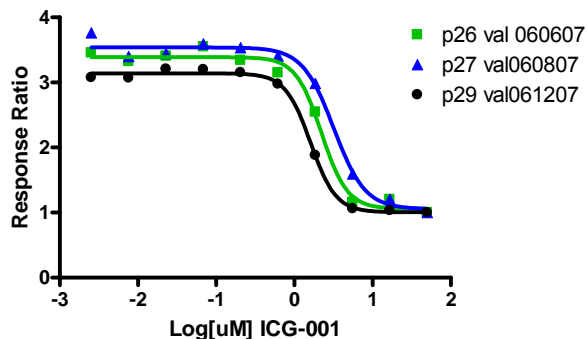
7. Assay performance with variable substrate loading time

8. Assay performance with variable DMSO concentration

9. Assay performance with variable compound incubation time

Determination of assay window

Figure 1 — Myc-*bla* HCT-116 response to ICG-001 under optimized conditions



	p26 val 060607	p27 val060807	p29 val061207
BOTTOM	1.061	1.054	1.006
TOP	3.388	3.538	3.139
LOGEC50	0.3461	0.5036	0.2072
HILLSLOPE	-2.638	-2.177	-2.684
EC50	2.219	3.189	1.611

Myc-*bla* HCT116 cells (8000 cells/well) were assayed on three separate days represented by the three curves shown on the graph. Cells were plated the day prior to the assay in a 384-well format and treated with indicated amount of ICG-001 in the presence of 0.1% DMSO in Assay Medium for 24 hours. Cells were then loaded with LiveBLazer™-FRET B/G Substrate for 120 minutes. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the response ratio was calculated as the 460/530 Emission Ratio of untreated cells divided by the 460/530 Emission Ratio of cells treated with 50 μ M ICG-001. The response ratio was plotted for the indicated treatment (n=16 for each data point).

Target validation with RNAi

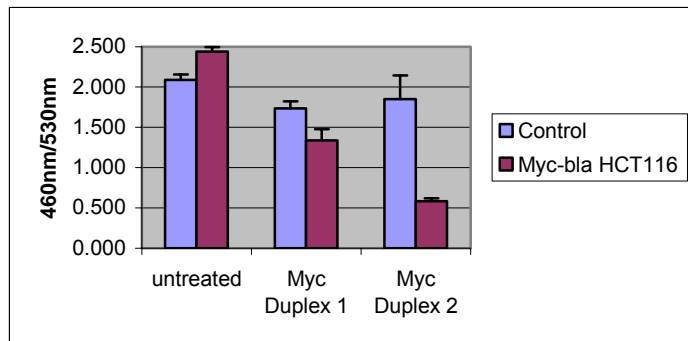


Figure 2 — Myc-*bla* HCT116 response to treatment with Myc RNAi

Myc-*bla* HCT116 or control cells (10000 cells/well) were plated the day prior to the transfection in a 96-well format in growth medium. The cells were transfected with the indicated RNAi duplexes (20 nM final concentration, Invitrogen, #12936-50) using Lipofectamine™RNAiMax (Invitrogen, #13778-075) according to manufacturer instructions, and incubated for 72 hours with the RNAi. Cells were then loaded with LiveBLazer™-FRET B/G Substrate for 120 minutes. Emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the 460/530 Ratios plotted for each treatment (n=4 for each data point).

Cell Culture and Maintenance

Thaw cells in Growth Medium without Blastocidin and culture them in Growth Medium with Blastocidin. Pass or feed cells at least twice a week and maintain them in a 37°C/5% CO₂ incubator. Maintain cells between 10% and 85% confluency. Do not allow cells to reach confluence.

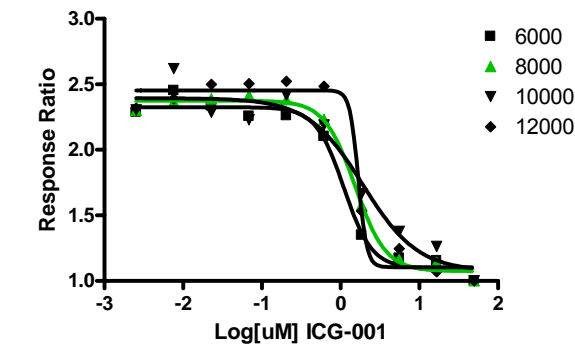
Note: We recommend passing cells for three passages after thawing before using them in the beta-lactamase assay. For more detailed cell growth and maintenance directions, please refer to protocol.

Table 1 – Cell Culture and Maintenance

Component	Growth Medium	Assay Medium	Freezing Medium
McCoy's 5A Medium	90%	—	—
OPTI-MEM	—	90%	—
Dialyzed FBS Do Not Substitute!	10%	0.5%	—
NEAA	—	0.1 mM	—
HEPES (pH 7.3)	—	10 mM	—
Sodium Pyruvate	—	1 mM	—
Penicillin (antibiotic)	100 U/ml	100 U/ml	—
Streptomycin (antibiotic)	100 µg/ml	100 µg/ml	—
Blasticidin (antibiotic)	5 µg/ml	—	—
Recovery™ Cell Culture Freezing Medium	—	—	100%

Assay Performance with Variable Cell Number

Figure 3 — ICG-001 inhibition with different plating cell numbers/well

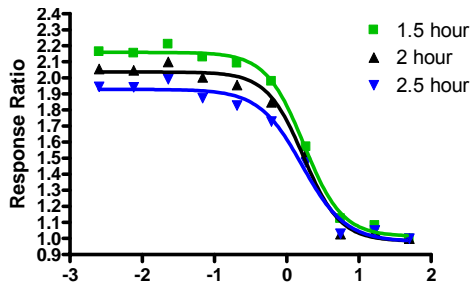


	6000	8000	10000	12000
BOTTOM	1.097	1.074	1.072	1.105
TOP	2.325	2.375	2.394	2.453
LOGEC50	0.04235	0.1848	0.2848	0.2263
HILLSLOPE	-2.432	-2.227	-1.240	-8.016
EC50	1.102	1.530	1.927	1.684

Myc-*bla* HCT116 cells were plated the day prior to the assay at the indicated number of cells/well in a 384-well format in growth medium and treated with indicated amount of ICG-001 in Assay Medium for 24 hours. Cells were then loaded with LiveBLazer™-FRET B/G Substrate for 120 minutes. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the response ratio was calculated as the 460/530 Emission Ratio of untreated cells divided by the 460/530 Emission Ratio of cells treated with 50 μ M ICG-001. The response ratio was plotted for the indicated treatment (n=8 for each data point).

Assay Performance with Variable Substrate Loading Time

Figure 4 — ICG-001 inhibition with various substrate loading times

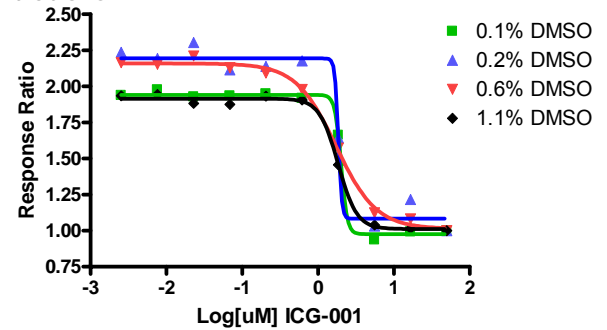


	1.5 hour	2 hour	2.5 hour
BOTTOM	1.013	0.9840	0.9808
TOP	2.159	2.037	1.928
LOGEC50	0.2380	0.2328	0.2150
HILLSLOPE	-1.641	-1.703	-1.456
EC50	1.730	1.709	1.641

Myc-*bla* HCT116 cells were plated the day prior to the assay at 8000 cells/well in a 384-well format in growth medium and treated with indicated amount of ICG-001 in Assay Medium for 24 hours. Cells were then loaded with LiveBLazer™-FRET B/G Substrate for 1.5, 2 and 2.5 hours. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the response ratio was calculated as the 460/530 Emission Ratio of untreated cells divided by the 460/530 Emission Ratio of cells treated with 50 μ M ICG-001. The response ratio was plotted for the indicated treatment (n=8 for each data point).

Assay Performance with Variable DMSO Concentration

Figure 5 — ICG-001 inhibition with various DMSO concentrations

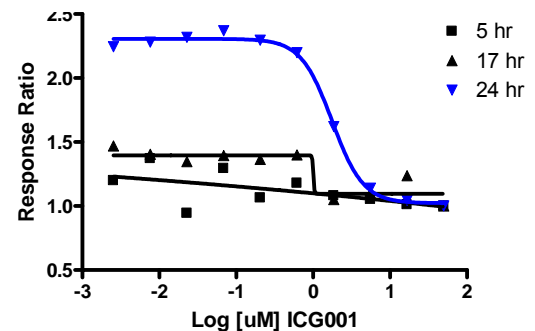


	0.1% DMSO	0.2% DMSO	0.6% DMSO	1.1% DMSO
BOTTOM	0.9756	1.083	1.013	1.011
TOP	1.941	2.194	2.159	1.915
LOGEC50	0.3042	0.2680	0.2380	0.2646
HILLSLOPE	-10.74	-19.22	-1.641	-3.513
EC50	2.015	1.853	1.730	1.839

Myc-*bla* HCT116 cells (8000 cells/well) were plated the day prior to the assay in a 384-well format and treated with indicated amount of ICG-001 in the presence of indicated concentrations of DMSO in Assay Medium for 24 hours. Cells were then loaded with LiveBLazer™-FRET B/G Substrate for 120 minutes. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the response ratio was calculated as the 460/530 Emission Ratio of untreated cells divided by the 460/530 Emission Ratio of cells treated with 50 μ M ICG-001. The response ratio was plotted for the indicated treatment (n=8 for each data point).

Compound Incubation Time

Figure 6 — Compound incubation time



	5 hr	17 hr	24 hr
BOTTOM	-1663	1.097	1.022
TOP	1.692	1.396	2.306
LOGEC50	82.95	0.003941	0.2471
HILLSLOPE	-0.04157	-64.78	-2.150
EC50	Value too large	1.009	1.766

Myc-*bla* HCT116 cells were plated at 8000 cells/well in a 384-well format in growth medium and treated with indicated amount of ICG-001 in Assay Medium for 5, 17 and 24 hours. Cells were then loaded with LiveBLazer™-FRET B/G Substrate for 2 hours. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and the response ratio was calculated as the 460/530 Emission Ratio of untreated cells divided by the 460/530 Emission Ratio of cells treated with 50 μ M ICG-001. The response ratio was plotted for the indicated treatment (n=8 for each data point).

|