

Human Aggregated Beta Amyloid ELISA Kit

Catalog Number KHB3491 (96 tests)

Pub. No. MAN0014907 Rev. 3.0 (31)

CAUTION! This kit contains materials with small quantities of sodium azide. Sodium azide reacts with lead and copper plumbing to form explosive metal azides. Upon disposal, flush drains with a large volume of water to prevent azide accumulation. Avoid ingestion and contact with eyes, skin and mucous membranes. In case of contact, rinse affected area with plenty of water. Observe all federal, state, and local regulations for disposal.

Note: For safety and biohazard guidelines, see the "Safety" appendix in the *ELISA Technical Guide* (Pub. no. MAN0006706). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Product description

The Invitrogen™ Human Aggregated Beta Amyloid ELISA Kit is a solid-phase sandwich Enzyme-Linked Immunosorbent Assay (ELISA). This assay is designed to detect and quantify the level of aggregated beta amyloid (A β) in human tissue homogenates, ventricular fluid, CSF, tissue culture supernatant, and buffered solution. The assay recognizes both natural and recombinant human aggregated A β , and does not cross-react with mouse or rat.

Extracellular amyloid plaques are composed primarily of beta amyloid. Associated with the extracellular plaques are activated microglia and astrocytes. An important neurotoxic form of A β is an oligomer, composed of approximately 12 subunits, with a composite molecular weight of 54 kDa. This oligomeric form of A β (also known as Amyloid Derived Diffusible Ligand or ADDL) can be separated from fibrillar and protofibrillar forms of aggregated beta amyloid by high speed centrifugation or by size exclusion methods.

Contents and storage

Upon receipt, store the kit at 2°C to 8°C.

Contents	Cat. No. KHB3491 (96 tests)
Hu Aggregated A β Standard, lyophilized; contains 0.1% sodium azide.	2 vials
Standard Diluent Buffer; contains 0.1% sodium azide; red dye ^[1]	60 mL
Antibody Coated Plate, 96-well plate	1 plate
Hu Aggregated A β Biotin Conjugate; contains 0.1% sodium azide; blue dye ^[1]	11 mL
Streptavidin-HRP (100X); contains 3.3 mM thymol	0.125 mL
HRP Diluent; contains 3.3 mM thymol; yellow dye ^[1]	25 mL
Wash Buffer Concentrate (25X)	100 mL
Stabilized Chromogen, Tetramethylbenzidine (TMB)	25 mL
Stop Solution	25 mL
Plate Covers, adhesive strips	3

^[1] In order to help our customers avoid any mistakes in pipetting the ELISAs, we provide colored Standard Diluent Buffer, Detection Antibody, and HRP Diluent to help monitor the addition of solutions to the reaction wells. This does not in any way interfere with the test results.

Materials required but not supplied

- Distilled or deionized water
- Calibrated adjustable precision pipettes and glass or plastic tubes for diluting solutions; beakers, flask and cylinders for preparation of reagents
- Microtiter plate reader with software capable of measurement at or near 450 nm
- Plate washer—automated or manual (squirt bottle, manifold dispenser, or equivalent)

Before you begin

IMPORTANT! Reagents are lot-specific. Do not mix or interchange different reagent lots from various kit lots.

- Review the **Procedural guidelines** and **Plate washing directions** in the *ELISA Technical Guide* available at thermofisher.com.
- Allow reagents to reach room temperature before use. Mix to redissolve any precipitated salts.

Prepare 1X Wash Buffer

1. Dilute 16 mL of Wash Buffer Concentrate (25X) with 384 mL of deionized or distilled water. Label as 1X Wash Buffer.
2. Store the concentrate and 1X Wash Buffer in the refrigerator. Use the diluted buffer within 14 days.

Prepare Tissue Extraction Buffer

Note: See the *ELISA Technical Guide* for detailed information on preparing Tissue Extraction Buffer.

1. Prepare 5 mL of Tissue Extraction Buffer.
Tissue Extraction Buffer consists of 25 mM Tris (pH 7.4), and 150 mM NaCl.
2. Immediately before use, add 1.46 μ M pepstatin A, 0.154 μ M aprotinin, 2.03 μ M leupeptin, 0.5 mM AEBSF, and 0.29 mM PMSF.

If desired, supplement the buffer with phosphatase inhibitors: 0.05 mM fenvalerate, 0.05 mM cantharidin, 1 mM Na₃VO₄, 1 mM Na₄P₂O₇, and 50 mM NaF.

Prepare aggregated A β sample (liquid)

Due to shared epitopes, the following analytes may cause interference with this assay: APP, monomeric A β 40, monomeric A β 42, protofibrillar A β , and fibrillar A β . Sample preparation should therefore be carefully considered when using this assay.

1. Collect samples in pyrogen/endotoxin-free polypropylene tubes.

Note: Aggregated A β is sensitive to freeze-thaw cycles. Avoid multiple freeze-thaw cycles of frozen samples. Thaw completely and mix well prior to analysis.

2. Clarify samples by centrifugation prior to analysis.
 - a. Centrifuge at $14,000 \times g$ for 10 minutes to minimize fibrils in aggregated A β -containing samples.
 - b. Centrifuge at $100,000 \times g$ for 1 hour at 4°C to minimize fibrils and protofibrils.

Size exclusion methods, such as gel permeation chromatography or ultrafiltration, can also improve assay performance.

Prepare tissue homogenate

1. Determine the mass of the tissue sample in a microcentrifuge tube.
2. Add $10 \times$ the tissue mass of Tissue Extraction Buffer, then homogenize thoroughly.
3. Sonicate the homogenized tissue sample (2 blasts, 10 seconds each).
4. Centrifuge the homogenate at $100,000 \times g$ for 1 hour at 4°C .
5. Carefully transfer the clear supernatant fraction into clean microcentrifuge tubes on ice.

The supernatant fraction contains TBS-soluble aggregates of A β .

Pre-dilute samples

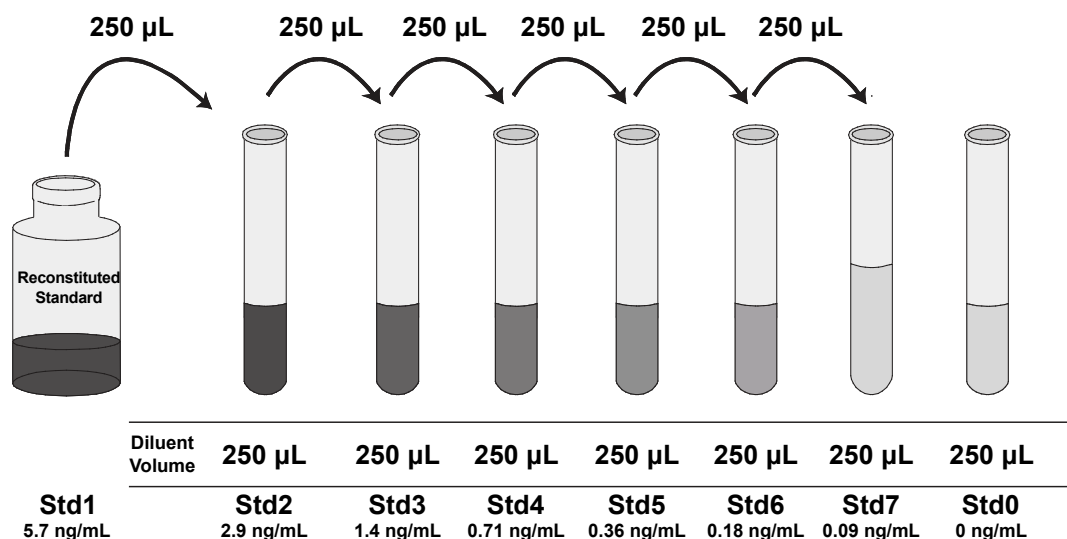
Sample concentrations should be within the range of the standard curve. Because conditions may vary, each investigator should determine the optimal dilution for each application.

- Perform sample dilutions with Standard Diluent Buffer.
- Dilute samples prepared in Tissue Extraction Buffer 1:2 or greater in Standard Diluent Buffer (e.g., 25 μL sample into 25 μL buffer).

Dilute standards

Note: Use glass or plastic tubes for diluting standards.

1. Reconstitute Hu Aggregated A β Standard to 5.7 ng/mL with Standard Diluent Buffer. Refer to the standard vial label for instructions. Swirl or mix gently and allow the contents to sit for 10 minutes to ensure complete reconstitution. Label as 5.7 ng/mL human aggregated A β . **Use the standard within 1 hour of reconstitution.**
2. Add 250 μL Standard Diluent Buffer to each of 7 tubes labeled as follows: 2.9, 1.4, 0.71, 0.36, 0.18, 0.09, and 0 ng/mL human aggregated A β .
3. Make serial dilutions of the standard as shown in the following dilution diagram. Mix thoroughly between steps.
4. Remaining reconstituted standard should be discarded or frozen in aliquots at -80°C for further use. Standard can be frozen and thawed one time only without loss of immunoreactivity.



Prepare 1X Streptavidin-HRP solution

Note: Prepare 1X Streptavidin-HRP within 15 minutes of usage.

The Streptavidin-HRP (100X) is in 50% glycerol, which is viscous. To ensure accurate dilution:

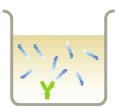



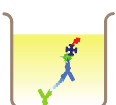
1. For each 8-well strip used in the assay, pipet 10 μL Streptavidin-HRP (100X) solution, wipe the pipette tip with clean absorbent paper to remove any excess solution, and dispense the solution into a tube containing 1 mL of Streptavidin-HRP Diluent. Mix thoroughly.
2. Return the unused Streptavidin-HRP (100X) solution to the refrigerator.

Perform Assay (Total assay time: 4 hours)

IMPORTANT! Perform a standard curve with each assay.

- Allow all components to reach room temperature before use. Mix all liquid reagents prior to use.
- Determine the number of 8-well strips required for the assay. Insert the strips in the frames for use. Re-bag any unused strips and frames, and store at 2°C to 8°C for future use.



1 Bind antigen 	1.1. Add 100 µL of standards, controls, or samples (see “Pre-dilute samples” on page 2) to the appropriate wells. Leave the wells for chromogen blanks empty. 1.2. Cover the plate with a plate cover and incubate for 2 hours at room temperature. 1.3. Thoroughly aspirate the solution and wash wells 4 times with 1X Wash Buffer.
2 Add Biotin Conjugate 	2.1. Add 100 µL Hu Aggregated Aβ Biotin Conjugate solution into each well except the chromogen blanks. 2.2. Cover the plate with plate cover and incubate for 1 hour at room temperature. 2.3. Thoroughly aspirate the solution and wash wells 4 times with 1X Wash Buffer.
3 Add Streptavidin-HRP 	3.1. Add 100 µL 1X Streptavidin-HRP solution (see page 2) into each well except the chromogen blanks. 3.2. Cover the plate with a plate cover and incubate for 30 minutes at room temperature. 3.3. Thoroughly aspirate the solution from the wells and wash wells 4 times with 1X Wash Buffer.
4 Add Stabilized Chromogen 	4.1. Add 100 µL Stabilized Chromogen to each well. The substrate solution begins to turn blue. 4.2. Incubate for 30 minutes at room temperature in the dark. Note: TMB should not touch aluminum foil or other metals.
5 Add Stop Solution 	Add 100 µL Stop Solution to each well. Tap the side of the plate to mix. The solution in the wells changes from blue to yellow.

Read the plate and generate the standard curve

- Read the absorbance at 450 nm. Read the plate within 2 hours after adding the Stop Solution.
- Use curve-fitting software to generate the standard curve. A 4 parameter algorithm provides the best standard curve fit. Optimally, the background absorbance may be subtracted from all data points, including standards, unknowns and controls, prior to plotting.
- Read the concentrations for unknown samples and controls from the standard curve. Multiply value(s) obtained for sample(s) by the appropriate factor to correct for the sample dilution.

Note: Dilute samples producing signals greater than the upper limit of the standard curve in Standard Diluent Buffer and reanalyze. Multiply the concentration by the appropriate dilution factor.

Performance characteristics

Standard curve example

The following data were obtained for the various standards over the range of 0 to 5.7 ng/mL human aggregated Aβ.

Standard Human Aggregated Aβ (ng/mL)	Optical Density (450 nm)
5.7	3.09
2.9	2.22
1.4	1.46
0.71	0.57
0.36	0.37
0.18	0.26
0.09	0.17
0	0.09

Inter-assay precision

Samples were assayed 48 times in multiple assays to determine precision between assays.

Parameters	Sample 1	Sample 2	Sample 3
Mean (ng/mL)	2.82	1.36	0.33
Standard Deviation	0.16	0.12	0.03
% Coefficient of Variation	5.74	8.99	7.61

Intra-assay precision

Samples of known human aggregated Aβ concentrations were assayed in replicates of 16 to determine precision within an assay.

Parameters	Sample 1	Sample 2	Sample 3
Mean (ng/mL)	2.84	1.48	0.32
Standard Deviation	0.19	0.09	0.03
% Coefficient of Variation	6.68	6.29	9.26

Linearity of dilution

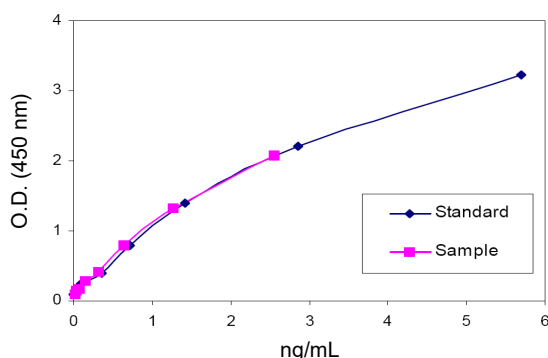
Aggregated A β was diluted in Standard Diluent Buffer over the range of the assay and measured. Linear regression analysis of sample values versus the expected concentrations yielded a correlation coefficient of 0.99.

Dilution	Measured (ng/mL)	Expected	
		Conc. (ng/mL)	%
Neat	2.55	2.55	100
1/2	1.35	1.28	106
1/4	0.737	0.638	115
1/8	0.344	0.319	108
1/16	0.214	0.159	134
1/32	0.088	0.079	110

Parallelism

Aggregated A β peptide was serially diluted in Standard Diluent Buffer. The optical density of each dilution was plotted against the human aggregated A β standard curve. The standard accurately reflects aggregated human aggregated A β content in a sample.

Parallelism of Human Aggregated Beta Amyloid ELISA



Sensitivity

The analytical sensitivity of the assay is <0.01 ng/mL human aggregated A β . This was determined by adding two standard deviations to the mean O.D. obtained when the zero standard was assayed 30 times.

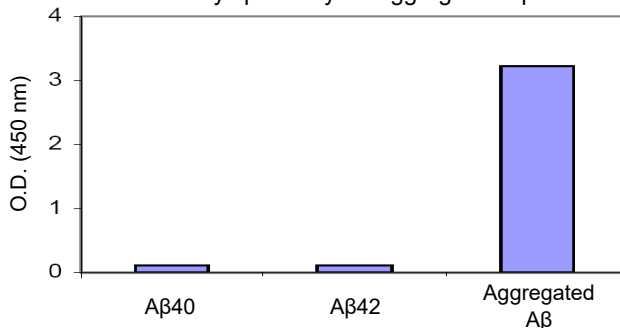
Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

Specificity

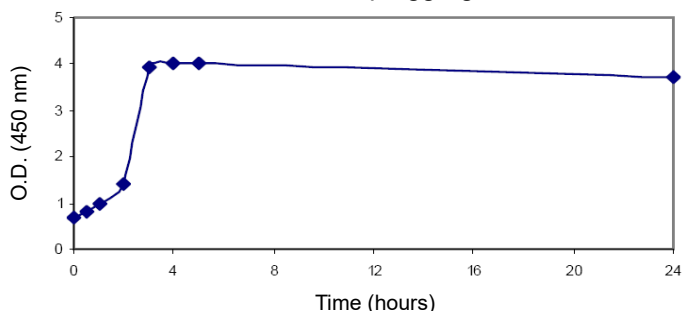
The specificity of the Human Aggregated Beta Amyloid ELISA Kit for human aggregated A β was determined by measuring samples containing A β 40 peptide, A β 42 peptide, and aggregated A β . The data confirms that the assay recognizes the aggregated form of A β , but not the nonaggregated A β 40 or A β 42 peptides.

Assay specificity for aggregated A β



Human A β 40 peptide was treated with 100% HFIP at 1 mg/mL for 1 hour then dried with nitrogen gas. The dried peptide was resuspended with water and incubated at 37°C. Samples were collected over a 24 hour period and analyzed with the Human Aggregated Beta Amyloid ELISA Kit. A β aggregation appeared to reach maximal level at 4 hours post-reconstitution.

Induction of A β aggregation



Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN0014907

Revision	Date	Description
3.0 (31)	1 July 2022	<ul style="list-style-type: none">In "Standard curve example" and "Linearity of dilution", the concentration units of human aggregated Aβ were corrected to ng/mL.The symbols definition table was replaced with a reference to thermofisher.com/symbols-definition.
2.0 (30)	17 August 2017	Baseline for this revision history. Updated to the latest Thermo Fisher template, with associated updates to logos, trademarks, legal, and style.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.