



LEGEND MAX™ β -Amyloid x-42 ELISA Kit Components and Protocol

Kit Components

• Capture Antibody Coated Plate	1 Stripwell plate
• 1-42 Standard	(2) 20 μ g vial
• 5X Wash Buffer	125mL
• Standard Diluent	2 x 3mL
• 2X Incubation Buffer	12mL
• HRP Detection Antibody	40 μ L
• TMB Substrate	22mL
• Plate Sealer	2
• 1.5mL Microcentrifuge Tubes	20

LEGEND MAX™ β -Amyloid x-42 Protocol

- Equilibrate all kit components to room temperature before use.
- The 96 well plate for this kit is in stripwell format and can be used in more than one experiment. The calculations below are based on running a full plate, please adjust accordingly if running a partial plate.

I. Preparation of Standard Intermediates (refer to Table 1)

- Label (2) 1.5mL microcentrifuge tubes as intermediate #1 & 2. (use enclosed tubes)
- Add 990 μ L of standard diluent to intermediate tubes #1 & 2.
- Reconstitute one 20 μ g vial of 1-42 standard with 80 μ L of Standard Diluent. **Mix well by inversion. Do not vortex.** Concentration will be 250 μ g/mL.
- Incubate 30 minutes at room temperature.

Note: During 30 minute incubation; proceed thru appropriate sections below (e.g. sections II, III and V).

- After 30 minutes incubation, mix well by inversion, do not vortex.




Once reconstituted, standard **must** be used within the same day.

- Add 10 μ L from the vial of reconstituted 1-42 standard to 990 μ L of standard diluent in intermediate tube #1. Mix well by inversion, do not vortex.
- Remove 10 μ L from intermediate #1 tube and add to 990 μ L of standard diluent in intermediate #2 tube. Mix well by inversion, do not vortex.
- The final concentration of standard in intermediate tube #2 will be 25ng/mL.

TABLE 1.

Tube Number	Volume of Standard	Volume of Standard Diluent (μ L)	Concentration (ng/mL)
Reconstituted Standard	0	80	250,000
Intermediate #1	10 μ L of reconstituted standard	990	2,500
Intermediate #2	10 μ L of Intermediate #1	990	25

LEGEND MAX™ β -Amyloid x-42
ELISA Kit with Pre-coated Plate 
Catalog No.: 842401 (Previously
Covance Cat. No. SIG-38956)

II. Preparation of 1X Incubation Buffer

- I. Label a 50mL centrifuge tube as “1X Incubation Buffer”
- J. Dilute 2X Incubation buffer to 1X by adding 10mL of 2X incubation buffer to 10mL of lab grade water* in the 50mL tube labeled “1X Incubation Buffer”.
- K. Mix well by vortexing. This will be diluent for the standard curve and samples.

*Note: Lab grade filtered water such as injection grade, cell culture grade, Reverse Osmosis De-Ionization (RODI).

III. Preparation of 1X Wash Buffer

- A. Label a 1L container “1X Wash buffer”.
- B. Dilute 5X wash buffer to 1X for use. Mix 125mL of 5X Wash buffer with 500mL of lab grade water for a total volume of 625mL.

IV. Preparation of Standard Curve (Refer to Table 2)

- A. Label (8) 1.5mL microcentrifuge tubes as #1-8 (use enclosed tubes).
- B. Aliquot 240 μ L of the 1X incubation buffer (made previously in step II) to each of the standard curve tubes (#2-8) and 490 μ L to tube #1 (standard curve top point).
- C. Remove 10 μ L from intermediate #2 and add to 490 μ L 1X incubation buffer in tube #1 (this will be the top point of the standard curve, final concentration will be 250pg/mL).
- D. Mix well by inversion, do not vortex.
- E. Continue making 1.8 fold serial dilutions by adding 300 μ L of the previous dilution to 240 μ L of 1X incubation buffer in tubes #2-7. Mix well by inversion between each dilution.

TABLE 2

Tube Number	Volume of Standard (μ L)	Volume 1X Incubation Buffer	Intermediate Concentration (pg/mL)
1	10 μ L intermediate #2	490	500.0
2	300 μ L of #1	240	277.8
3	300 μ L of #2	240	154.3
4	300 μ L of #3	240	85.7
5	300 μ L of #4	240	47.6
6	300 μ L of #5	240	26.5
7	300 μ L of #6	240	14.7
8	0	240	0

V. Sample Preparation

- A. Dilute samples in 1X incubation buffer to 2X concentration. **Mix well by inversion.**

For example, if the final sample dilution should be 1:10, dilute the sample 1:5 in 1X incubation buffer. The sample will then be diluted 1:2 in step VII for a final dilution factor of 1:10.

LEGEND MAX™ β -Amyloid x-42
 ELISA Kit with Pre-coated Plate
 Catalog No.: 842401 (Previously
 Covance Cat. No. SIG-38956)



- B. Run samples in duplicate or triplicate.

Note: All sample matrices will perform differently in the kit. It is important that you determine the ideal dilution for your particular sample type. It is good practice to run 2-3 dilutions per sample to ensure at least 1 dilution falls within the range of the standard curve. For most sample types, a good starting dilution would be 1:5-1:10. Due to the format of the assay, samples are not able to run neat.

VI. Preparation of diluted HRP Detection Antibody

- A. Label a 15mL tube as “Diluted HRP Detection Antibody”
- B. Add 6mL of 1X Incubation buffer to the tube labeled “Diluted HRP Detection Antibody”
- C. Add 6 μ L of HRP Detection Antibody and mix well by vortexing

VII. Running the Assay

Day 1

- A. Remove plate from foil pouch. Remove extra stripwells if necessary and store in resealable foil pouch at 2-8°C until use.
- B. Add 300 μ L per well of 1X Wash Buffer.
- C. Dump out wash buffer and pat dry on paper towels.
- D. Add 50 μ L of each standard to the plate in duplicate or triplicate. Follow the plate layout outlined in Table 3 below.

Note: Wells E1-E3 contain the zero or blank sample (Std #8)

- E. Add 50 μ L of each sample to the plate in duplicate or triplicate.
- F. Add 50 μ L per well of diluted HRP detection antibody to all wells.
- G. Cover plate with plate sealer.
- H. Mix the plate gently on a plate shaker for one minute.
- I. Incubate overnight at 2-8°C.

NOTE: Once diluted with 50 μ L of diluted HRP detection antibody, the final standard curve concentrations will be outlined in Table 4. Please use these concentrations to generate your standard curve.

TABLE 3.

	1	2	3	4	5	6	7	8	9	10	11	12
A	Std 1	Std 1	Std 1	Sample 1	Sample 1							
B	Std 2	Std 2	Std 2	Sample 2	Sample 2							
C	Std 3	Std 3	Std 3	Sample 3	Sample 3							
D	Std 4	Std 4	Std 4	Sample 4	Sample 4							
E	Std 8	Std 8	Std 8	Sample 5	Sample 5							
F	Std 5	Std 5	Std 5	Sample 6	Sample 6							
G	Std 6	Std 6	Std 6	Sample 7	Sample 7							
H	Std 7	Std 7	Std 7	Sample 8	Sample 8							





Day 2

- A. Remove plate from refrigerator and dump contents.
- B. Wash plate by adding 300 μ L of 1X Wash Buffer per well.
- C. Dump out 1X Wash Buffer and pat dry.
- D. Repeat steps B-C 4 more times for a total of 5 washes.
- E. Add 200 μ L TMB substrate to each well.

Note: If running multiple plates, stagger the addition of substrate so that all plates aren't ready for reading simultaneously.

- F. Incubate 40-50 minutes at room temperature in the dark.
- G. Read Plate at 620nm.

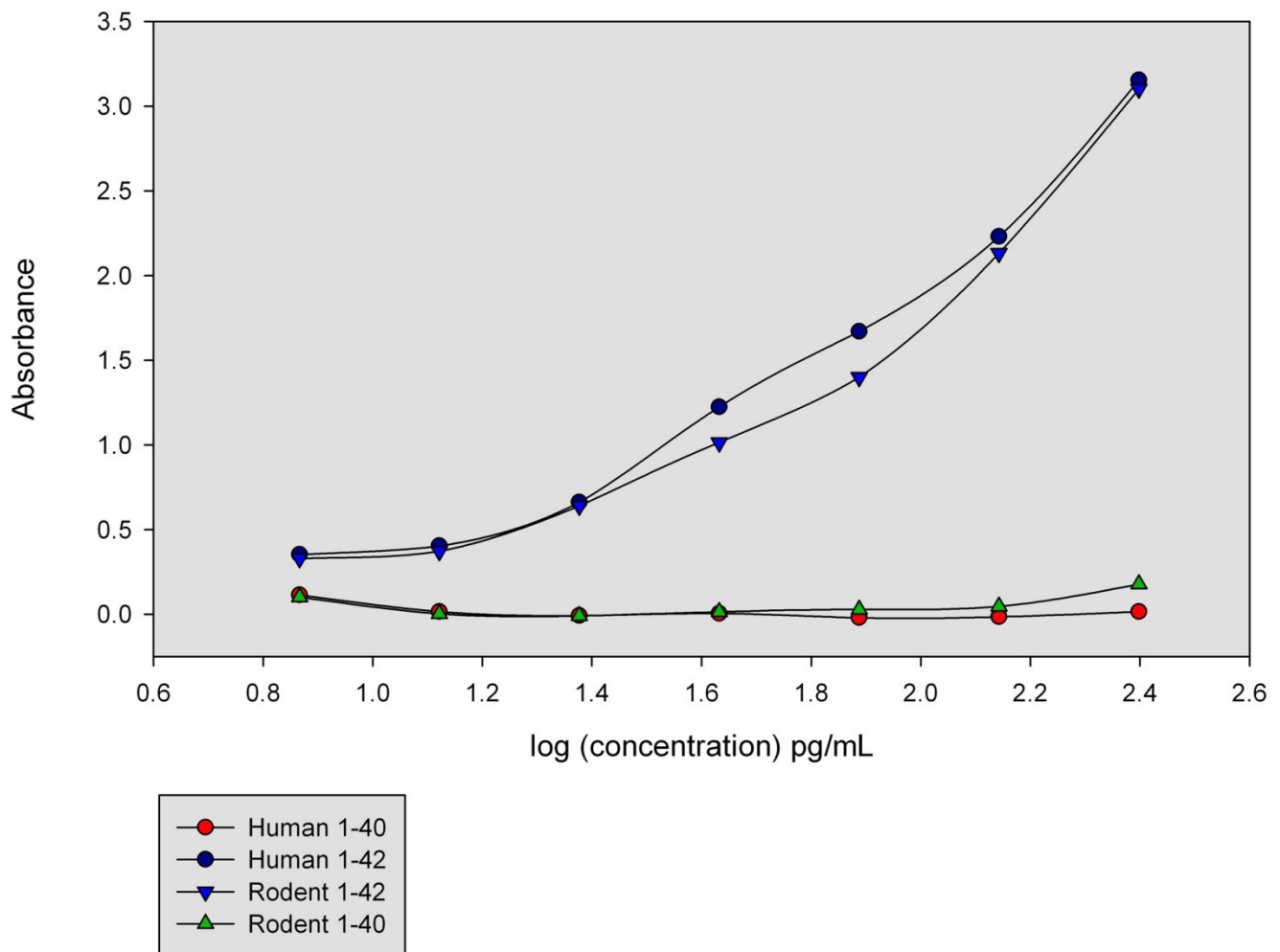
TABLE 4.

Standard#	Final Concentration (pg/mL)
1	250.0
2	138.9
3	77.2
4	42.9
5	23.8
6	13.2
7	7.4
8	0





4-Parameter Standard Curve





Preparation of Brain Samples for LEGEND MAX™ β -Amyloid x-42 ELISA

Procedure

I. Soluble and insoluble tissue fractionation.

This protocol generates soluble and insoluble subcellular fractions for analysis of A β partitioning in tissues.

- A. Homogenize in TBS with protease inhibitors (Pierce sells excellent solid tablets for use with homogenization buffers) at 5 mLs per 1 g tissue. Teflon/glass homogenizer with 5-6 passes on ice. EDTA¹ (2 mM) can be a useful addition to homogenization buffers (see notes below).
- B. Spin for 20 minutes x 350,000 g (or equivalent)
- C. Remove supernatant (S1 or soluble fraction) - usually contains < 5 % of the total brain A β pool in young mice and decreases with age.
- D. Resuspend pellet (agitate to break up) with: (use only one method below)
 - i. Detergent² - 1% triton OR 0.6% SDS (see notes) in TBS/inhibitor buffer using same volume as homogenization step (15 minutes incubation).
 - ii. 70% Formic acid², 40-50% of homogenization volume (incubate 30 minutes).
- E. Spin a second time (20 minutes x 350,000 g or equivalent) and remove supernatant - detergent or formic acid extract fraction (E1).
- F. OPTIONAL - if detergent extract was prepared above, then remaining pellet can be further extracted with formic acid - 70% formic acid, half homogenization volume, spin down 10 minutes in microfuge full speed (10 minutes x 14,000 g). Remove formic acid extract (E2).
- G. S1 and E1 fractions are best assayed immediately but can be stored at -20°C (including non-neutralized formic acid samples)

II. Whole tissue extraction

This protocol generates a single fraction containing the total A β tissue load.

- A. Homogenize tissue as above in TBS with protease inhibitors and EDTA.
- B. To homogenate add concentrated Triton X-100, SDS (to final 1 or 0.6% respectively) or formic acid (to final 70%). Agitate by pipetting up and down to mix. Detergents and formic acid are added after tissue homogenization to reduce foaming and other potential problems.
- C. Spin for 20 minutes x 350,000 g (or equivalent).
- D. Remove supernatant for assay (whole brain extract).




NB High salt and biomolecule concentrations in formic acid whole tissue extracts can sometimes result in visible precipitants forming in later steps. These samples retain strong A β signal but we have not established to what degree signal attenuated under these conditions.






Notes on Preparation and Assaying of Samples

¹EDTA in homogenization buffer helps solubilize A β and increases recovery of A β in the soluble fraction by as much as 50% for mild β -amyloid loaded brain (Cherny, 1999, JBC). HOWEVER, it is unclear if EDTA-solubilized material is free in solution or a precipitate in vivo. Mice at different ages/stages give different soluble/chelator yields (presumably because of shifts in quantity/quality of chelator-soluble A β forms with age). Hence, caution should be exercised if experiments are comparing A β levels in S1/EDTA fractions ("soluble" A β) from different aged mice.

²Choice of type of detergent or formic acid will depend on experimental circumstance.

-  Non-ionic Triton X-100 extracts a portion of insoluble A β . As with chelators, the portion of insoluble A β extracted by non-ionic detergents is dependent on age/stage of mice. For early stage of β -amyloid deposition (i.e. mice >12 months) up to 80% of total mouse brain A β is extracted by Triton X-100. Proportion drops rapidly 20% or less in old animals (i.e. mice >18 months) with heavy mature β -amyloid deposition. BioLegend A β ELISA's (and most other ELISA's) are compatible with non-ionic detergents.
-  The ionic detergent SDS gives more complete A β extraction from insoluble pellets (>90% with young and \approx 50% for older mice). HOWEVER, SDS is not compatible with many ELISA systems. BioLegend A β kits are resistant to SDS and samples with \leq 0.2% SDS can be used with minimal loss in sensitivity (< 10%). Hence, E1 fractions with SDS should have a final in assay dilution of at least 3-fold.
-  Early protocols used formic acid to extract A β from AD brain. For mature β -amyloid deposits in human and old mice brain the highest extraction yields are still observed with this buffer. HOWEVER, a difficult multi-step process is required to prepare formic acid fractions for assay (both E1 and whole tissue extracts). Normalization between samples is also problematic.

Steps to prepare formic acid extracts for A β assay;



-  The bulk of the formic acid buffer needs to be evaporated off (down to 10-20% of original vol) with a nitrogen stream. If dried completely formic acid brain samples form a plastic-like pellet that is highly resistant to resolubilization.
-  Aliquots of the remaining material must be neutralize (to pH 7) with 5 M NaOH in 1M Tris. Indicator should be added to the neutralization buffer to signal when pH 7 is reached.
-  Assay immediately - DO NOT FREEZE neutralized samples (precipitates of highly enriched insoluble hydrophobic species will form that interfere with assay signal).

Care should be taken to have equivalent pH for all samples. Neutralized samples should have a final in assay dilution of at least 10- fold. Normalizing between samples is problematic. Residual formic acid interferes with most protein assays and protein levels may not be proportional to starting materials in any case. Careful attention to maintaining equivalent sample volumes is often the most practical approach.








TBS = Tris, pH 7.4 (50 mM) + NaCl (150 mM)



Helpful Hints

-  Washing is critical in this assay. To ensure proper washing, a plate washer is highly recommended for all washes. Pipette washing is not vigorous enough and can result in high background and poor sensitivity.
-  Beta Amyloid is a highly hydrophobic protein and is very sticky. Mixing of the standards and samples is also a key step. Mix tubes by quickly inverting or shaking them for 10-15 seconds per tube before removing material for the next serial dilution. This will help ensure good linearity with your standard curve and samples. Also, minimize the number of times a pipette goes in and out of the standard tubes, this will reduce any loss of peptide.

Frequently Asked Questions

-  I am getting very high background and my standard curve is not linear at the low end. What is causing this?
This is typically a washing issue. Inadequate washing will lead to high background and poor sensitivity at the low end of the curve. To fix the issue, either use a plate washer or squirt bottle for all washing steps. Pipette washing will not be vigorous enough. This should help reduce the background and in turn give better sensitivity at the low end.
-  I have a plateau at the top end of my standard curve, what could be causing this?
Typically this is caused by not preparing your standard curve from intermediate #2 and using the reconstituted standard instead. The concentration is too high to be detected by the reader and will cause a plateau of the first few points on the standard curve.
-  My samples are undetectable in the assay, what is causing this?
Generally there are 2 causes. First, ensure that you added the diluted HRP detection antibody to the plate prior to the overnight incubation. Omission of this step will result in no signal. Another reason could be due to very low levels of A Beta in your samples. Due to the limitations of the kit, the samples must be diluted a minimum of 1:2 in the assay and cannot be run undiluted.
-  What are the recommended dilutions for my samples?
This is going to vary based on your specific samples. We recommend running 3 different dilutions for each sample, which helps ensure one of the dilutions will fall within the range of the standard curve. If you are testing a large number of samples, consider doing a trial run with a subset of samples to optimize the dilution factors.
-  What types of samples will work with this assay?
Currently, this assay works with TCS, plasma, CSF and tissue homogenate samples. It is not recommended for use with serum samples.
-  What type of curve fit should I use ?
We recommend using a 4 parameter curve fit to generate the standard curve for this assay.
-  What if I have further questions?
Please call our technical support at 1.877.273.3103 or send an e-mail to tech@biolegend.com. We will be more than happy to answer any questions you may have.

Customer Service

Tel: 1.877.246.5343 (US & Canada)