

AlphaLISA Amyloid beta 1-40 HS Detection Kit

Product number: AL373C/F

Caution: For Laboratory Use. A research product for research purposes only.

Contents

	Page
Product Information.....	2
Quality Control.....	2
Analyte of Interest.....	3
Description of the AlphaLISA Assay	3
Precautions.....	3
Kit content: Reagents and Materials.....	4
Recommendations.....	5
Assay Procedure.....	5
Data Analysis.....	8
Assay Performance Characteristics.....	9
CSF Experiments.....	11
Troubleshooting Guide.....	11

Product Information

Application: This kit is designed for the quantitative determination of Amyloid beta 1-40 (A β 40) in cerebrospinal fluid and cell culture supernatants using a homogeneous AlphaLISA assay (no wash steps). The assay shows negligible cross-reactivity with other forms of Amyloid beta (i.e. A β 1-38 and A β 1-42).

Sensitivity: Lower Detection Limit (LDL): 42.2 pg/mL
Lower Limit of Quantification (LLOQ): 89.0 pg/mL
EC₅₀: 9.1 ng/mL

Dynamic range: Kit designed to detect [A β 40] between: 42.2 – 10,000 pg/mL (Figure 1).

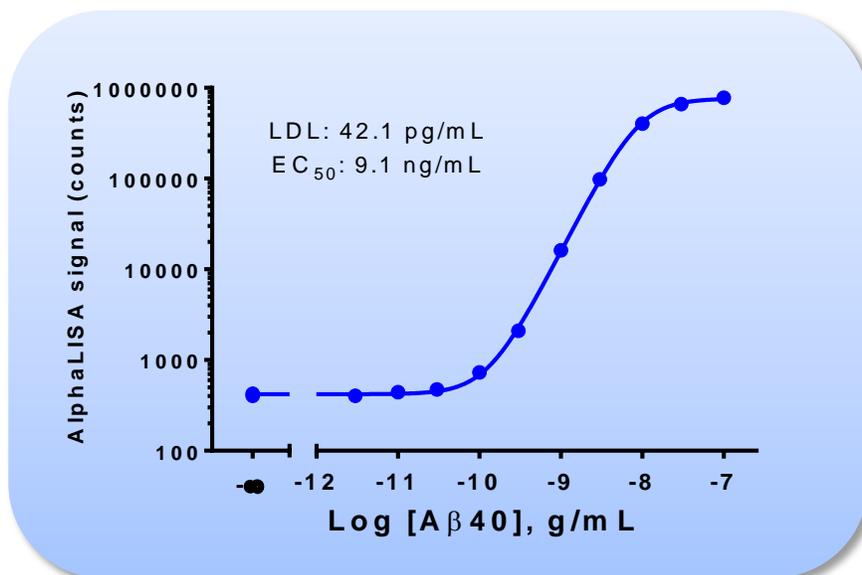


Figure 1. Typical sensitivity curve in AlphaLISA Immunoassay Buffer. The data was generated using a white Optiplate™-384 microplate and the EnVision® Multilabel Plate Reader 2103 with Alpha option.

Storage: Store kit in the dark at +4°C. **Store analyte at -20°C.** Once the **A β 1-40 analyte** has been reconstituted, the analyte is stable for up to 6 weeks at 4°C. For long term storage, aliquot and store at -20°C. Limit the number of freeze-thaw cycles.

Stability: This kit is stable for at least 6 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

Quality Control

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals, EC₅₀ and LDL were measured on the EnVision Multilabel Plate Reader with Alpha option using the protocol described in this technical data sheet. We certify that these results meet our quality release criteria. Maximum counts may vary between bead lots and the instrument used, with no impact on LDL measurement.

Analyte of Interest

Amyloid beta (A β) is a short peptide derived from a transmembrane protein, the amyloid precursor protein (APP), by proteolysis. The β - and γ -secretases cleave the respective N- and C-terminal ends of the A β sequence, liberating the A β peptide from APP. A β 40 is the major species of A β produced by neurons and other cells, and accounts for over 70% of total A β produced, while the remainder is typically composed A β 42 and other size variants ranging from 36 to 43 AA in length. Interestingly, it has been observed by some that the A β 40/A β 42 ratio increased in cerebrospinal fluid of AD patients. A β 40 could be an important biomarker for detecting Alzheimer's disease progression. This AlphaLISA kit has been designed to detect A β 40 in cell culture supernatants and cerebrospinal fluid.

Description of the AlphaLISA Assay

AlphaLISA technology allows for the detection of molecules of interest in buffer, cell culture media, serum and plasma in a highly sensitive, quantitative, reproducible and user-friendly mode. In an AlphaLISA assay, a Biotinylated Anti-Analyte Antibody binds to the Streptavidin-coated Alpha Donor beads, while another Anti-Analyte Antibody is conjugated to AlphaLISA Acceptor beads. In the presence of the analyte, the beads come into close proximity. The excitation of the Donor beads provokes the release of singlet oxygen molecules that triggers a cascade of energy transfers in the Acceptor beads, resulting in a sharp peak of light emission at 615 nm (Figure 2). Combining this assay with an AlphaPLEX 645- or AlphaPLEX 545 - based kit will allow the quantification of 2 (or more) analytes in the same well.

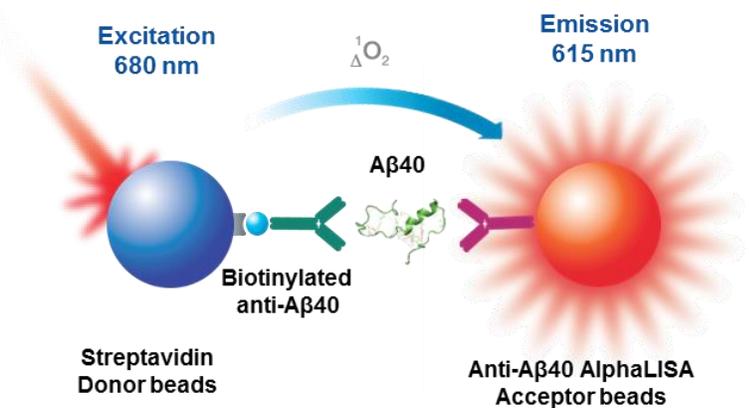


Figure 2. AlphaLISA assay principle.

Precautions

- The AlphaScreen[®] Donor beads are light-sensitive. All the other assay reagents can be used under normal light conditions. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters (preferred) or Roscolux filters #389 from Rosco) can be applied to light fixtures.
- All blood components and biological materials should be handled as potentially hazardous. The analyte included in this kit is from a source.
- Some analytes are present in saliva. Take precautionary measures to avoid contamination of the reagent solutions.
- The Biotinylated Anti-Analyte Antibody contains sodium azide. Contact with skin or inhalation should be avoided.

Kit Content: Reagents and Materials

Kit components	AL373C (500 assay points ^{***})	AL373F (5000 assay points ^{***})
AlphaLISA Anti-A β 40 Acceptor beads stored in PBS, 0.05% Proclin-300, pH 7.2	50 μ L @ 5 mg/mL (1 brown tube, <u>white</u> cap)	500 μ L @ 5 mg/mL (1 brown tube, <u>white</u> cap)
Streptavidin (SA)-coated Donor beads stored in 25 mM HEPES, 100 mM NaCl, 0.05% Proclin-300, pH 7.4	200 μ L @ 5 mg/mL (1 brown tube, <u>black</u> cap)	2 x 1 mL @ 5 mg/mL (2 brown tubes, <u>black</u> caps)
Biotinylated Antibody Anti-A β 40 stored in PBS, 0.1% Tween-20, 0.05% NaN ₃ , pH 7.4	50 μ L @ 500 nM (1 tube, <u>black</u> cap)	500 μ L @ 500 nM (1 tube, <u>black</u> cap)
Amyloid- β 1-40 peptide*	20 μ g, HFIP Film * (1 tube, <u>clear</u> cap)	20 μ g, HFIP Film * (1 tube, <u>clear</u> cap)
AlphaLISA Immunoassay Buffer (10X) **	10 mL, 1 small bottle	100 mL, 1 large bottle

* Please note that the Amyloid- β 1-40 analyte film may not be visible to the naked eye. One vial contains an amount of A β 40 sufficient for performing 100 standard curves. Additional vials can be ordered separately (cat # AL373S).

** Extra buffer can be ordered separately (cat # AL000C: 10 mL, cat # AL000F: 100 mL).

*** The number of assay points is based on an assay volume of 50 μ L in 96- or 384-well assay plates using the kit components at the recommended concentrations.

Sodium azide should **not** be added to the stock reagents. High concentrations of sodium azide (> 0.001 % final in the assay) might decrease the AlphaLISA signal. Note that sodium azide from the Biotinylated Antibody stock solution will not interfere with the AlphaLISA signal (0.0001% final in the assay).

Specific additional required reagents and materials:

The following materials are recommended:

Item	Suggested source	Catalog #
TopSeal™-A Adhesive Sealing Film	PerkinElmer Inc.	6050195
EnVision®-Alpha Reader	PerkinElmer Inc.	-
Dimethyl Sulfoxide	Sigma	D2650

Recommendations

- The volume indicated on each tube is guaranteed for single pipetting. Multiple pipetting of the reagents may reduce the theoretical amount left in the tube. To minimize loss when pipetting beads, it is preferable not to pre-wet the tip.
- Centrifuge all tubes (including lyophilized analyte) before use to improve recovery of content (2000g, 10-15 sec).
- Re-suspend all reagents by vortexing before use.
- Use Milli-Q[®] grade H₂O (18 MΩ•cm) to dilute 10X AlphaLISA Immunoassay Buffer.
- When diluting the standard or samples, change tips between each standard or sample dilution. When loading reagents in the assay microplate, change tips between each standard or sample addition and after each set of reagents.
- When reagents are added to the microplate, make sure the liquids are at the bottom of the well.
- Small volumes may be prone to evaporation. It is recommended to cover microplates with TopSeal-A Adhesive Sealing Films to reduce evaporation during incubation. Microplates can be read with the TopSeal-A Film.
- AlphaLISA signal is detected using an EnVision Multilabel Reader 2103 equipped with the Alpha option using the following settings: Total Measurement Time: 550 ms, Laser 680 nm Excitation Time: 180 ms, Mirror: D640as (Barcode# 444), Emission Filter: Wavelength 570nm, bandwidth: 100nm, Transmittance 75%, (Barcode# 244).
- AlphaLISA signal will vary with temperature and incubation time. For consistent results, identical incubation times and temperature should be used for each plate.
- The standard curves shown in this technical data sheet are provided for information only. A standard curve must be generated for each experiment. The standard curve should be performed in AlphaLISA Immunoassay buffer.

Assay Procedure

IMPORTANT: PLEASE READ THE RECOMMENDATIONS BELOW BEFORE USE

- The protocol described below is an **example** for generating one standard curve in a 50 µL final assay volume (48 wells, triplicate determinations) and 452 samples. The protocols also include testing samples in 384 well plates. If different amounts of samples are tested, the volumes of all reagents must be adjusted accordingly, as shown in the table below. ****These calculations do not include excess reagents to account for losses during transfer of solutions or dead volumes.*
- The standard dilution protocol is provided for information only. As needed, the number of replicates or the range of concentrations covered can be modified.
- Use of four background points in triplicate (12 wells) is recommended when LDL/LLOQ is calculated. One background point in triplicate (3 wells) can be used when LDL/LLOQ is not calculated.

		Volume				
Format	# of data points	Final	Sample	AlphaLISA beads / Biotin Antibody MIX	SA-Donor beads	Plate recommendation
AL373C	250	100 µL	10 µL	10 µL	80 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
	500	50 µL	5 µL	5 µL	40 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate™-384 (cat # 6005350)
	1 250	20 µL	2 µL	2 µL	16 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate™-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	2 500	10 µL	1 µL	1 µL	8 µL	Light gray AlphaPlate-1536 (cat # 6004350)
AL373F	5 000	50 µL	5 µL	5 µL	40 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290) Light gray AlphaPlate-384 (cat # 6005350)
	12 500	20 µL	2 µL	2 µL	16 µL	Light gray AlphaPlate-384 (cat # 6005350) ProxiPlate-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	25 000	10 µL	1 µL	1 µL	8 µL	Light gray AlphaPlate-1536 (cat # 6004350)

Protocol for A β 40 Total AlphaLISA Assay

2 Step Protocol – Dilution of standards in 1X AlphaLISA Immunoassay Buffer or cell culture medium. The protocol described below is for one standard curve (48 wells) and 452 sample wells. *If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly.*

Steps for Preparing Reagents

- 1) Preparation of 1X AlphaLISA Immunoassay Buffer:
Add 5 mL of 10X AlphaLISA Immunoassay Buffer to 45 mL H₂O.
- 2) Preparation of A β 40 analyte standard dilutions (Do NOT vortex analyte!):
 1. Reconstitute 20 μ g **A β 40-HFIP film** gently with 10 μ L of Dimethyl sulfoxide (DMSO; Sigma-Aldrich D2650) by swirling the tip at the bottom of the tube and repeatedly pipetting up and down for 10 – 15 seconds. **Do not vortex the analyte!** Dilute dissolved A β 40 film to 20 μ g/mL with 990 μ L of 1X AlphaLISA Immunoassay Buffer
 - a. Store on ice immediately after reconstitution. The reconstituted **A β 40** analyte film is stable up to 6 weeks at 4°C
 - b. Prepare standard dilutions as follows in 1X AlphaLISA Immunoassay Buffer or cell culture medium (change tip between each standard dilution):

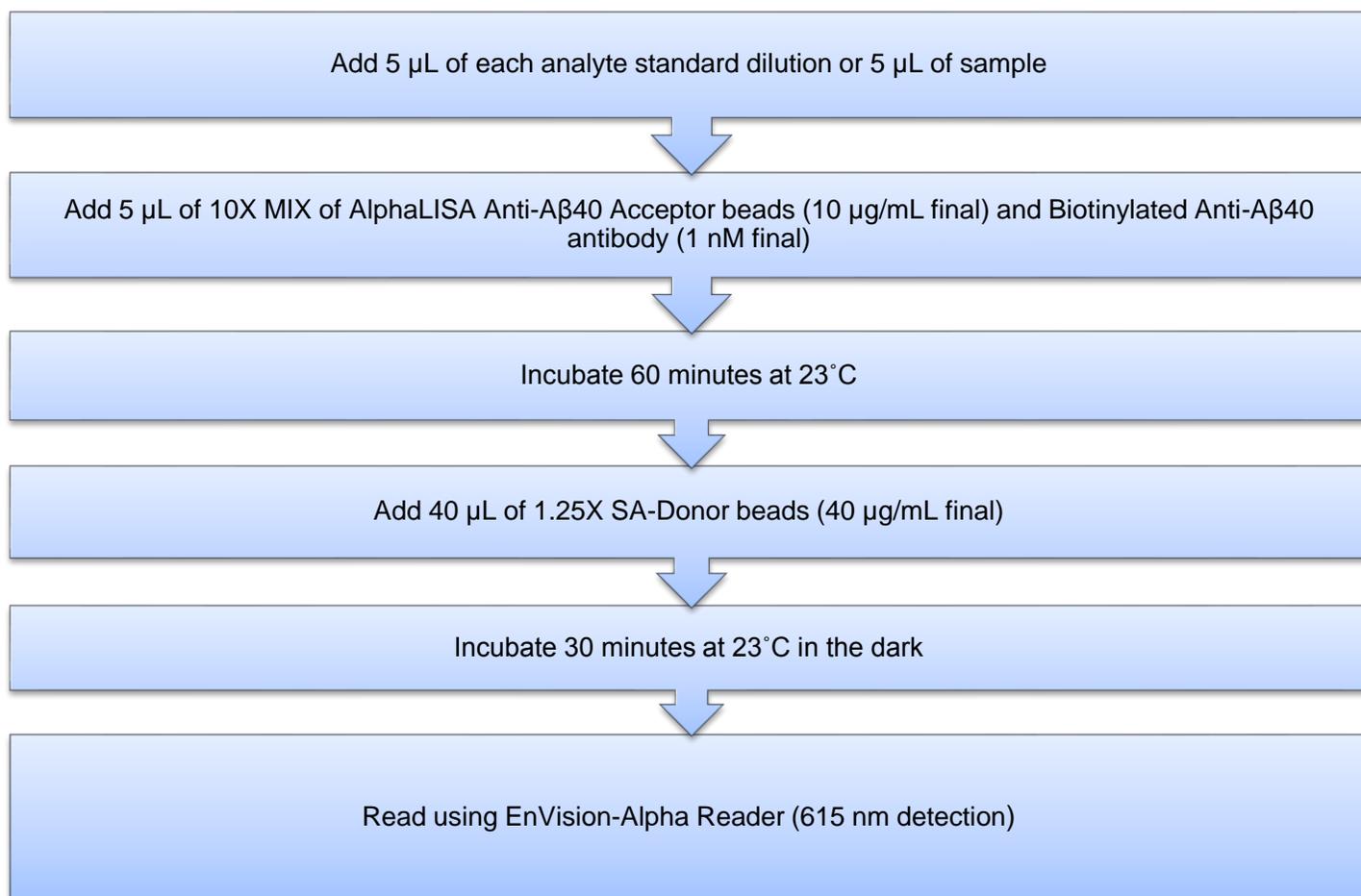
Tube	Vol. of A β 40 (μ L)	Vol. of diluent (μ L) *	[A β 40] in standard curve	
			(g/mL in 5 μ L)	(pg/mL in 5 μ L)
A	10 μ L of provided A β 40	190	1.00E-06	1 000 000
B	60 μ L of tube A	140	3.00E-07	300 000
C	60 μ L of tube B	120	1.00E-07	100 000
D	60 μ L of tube C	140	3.00E-08	30 000
E	60 μ L of tube D	120	1.00E-08	10 000
F	60 μ L of tube E	140	3.00E-09	3 000
G	60 μ L of tube F	120	1.00E-09	1 000
H	60 μ L of tube G	140	3.00E-10	300
I	60 μ L of tube H	120	1.00E-10	100
J	60 μ L of tube I	140	3.00E-11	30
K	60 μ L of tube J	120	1.00E-11	10
L	60 μ L of tube K	140	3.00E-12	3
M ** (background)	0	100	0	0
N ** (background)	0	100	0	0
O ** (background)	0	100	0	0
P ** (background)	0	100	0	0

* Dilute standards in diluent (e.g. 1X AlphaLISA Immunoassay Buffer).
At low concentrations of analyte, a significant amount of analyte can bind to the vial. Therefore, load the analyte

standard dilutions in the assay microplate within 60 minutes of preparation.

** Four background points in triplicate (12 wells) are used when LDL is calculated. If LDL does not need to be calculated, one background point in triplicate can be used (3 wells).

- 3) Preparation of 10X Anti-A β 40 AlphaLISA Acceptor beads (100 μ g/mL) and biotinylated Anti A β 40 Antibody (10 nM):
 - a. Add 50 μ L of 5 mg/mL **AlphaLISA Anti-A β 40 Acceptor beads** and 50 μ L of 500nM **Biotinylated Anti-A β 40 Antibody** to 2400 μ L of 1X AlphaLISA Immunoassay Buffer .
 - b. Prepare just before use.
- 4) Preparation of 1.25X Streptavidin (SA) Donor beads (50 μ g/mL):
 - a. Keep the beads under subdued laboratory lighting.
 - b. Add 200 μ L of 5 mg/mL SA-Donor beads to 19800 μ L of 1X AlphaLISA Immunoassay Buffer.
 - c. Prepare just before use.
- 5) In a white Optiplate (384 wells):



Read Settings: AlphaLISA signal is detected using an EnVision Multilabel Reader 2103 equipped with the Alpha option using the following settings: Total Measurement Time: 550 ms, Laser 680 nm Excitation Time: 180 ms, Mirror: D640as (Barcode# 444), Emission Filter: Wavelength 570nm, bandwidth: 100nm, Transmittance 75%, (Barcode# 244).

Data Analysis

- Calculate the average count value for the background wells.
- Generate a standard curve by plotting the AlphaLISA counts versus the concentration of analyte. A log scale can be used for either or both axes. No additional data transformation is required.
- Analyze data according to a nonlinear regression using the 4-parameter logistic equation (sigmoidal dose-response curve with variable slope) and a $1/Y^2$ data weighting (the values at maximal concentrations of analyte after the hook point should be removed for correct analysis).
- The LDL is calculated by interpolating the average background counts (12 wells without analyte) + 3 x standard deviation value (average background counts + (3xSD)) on the standard curve.
- The LLOQ as measured here is calculated by interpolating the average background counts (12 wells without analyte) + 10 x standard deviation value (average background counts + (10xSD)) on the standard curve. Alternatively, the true LLOQ can be determined by spiking known concentrations of analyte in the matrix and measuring the percent recovery, and then determining the minimal amount of spiked analyte that can be quantified within a given limit (usually +/- 20% or 30% of the real concentration).
- Read from the standard curve the concentration of analyte contained in the samples.
- If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.

Assay Performance Characteristics

AlphaLISA assay performance described below was determined using the 2 step protocol.

Assay Sensitivity:

The LDL and LLOQ were calculated as described above. The values correspond to the lowest concentration of analyte that can be detected in a volume of 5 μ L using the recommended assay conditions.

LDL (pg/mL)	LLOQ (pg/mL)	Buffer/Serum	# of experiments
42	89	AlphaLISA Immunoassay Buffer	6
77	134	Neurobasal A	6
99	201	DMEM/F12	6

* Note that LDL/ LLOQ can be decreased (i.e. sensitivity increased) by increasing the volume of analyte in the assay (e.g. use 10 μ L of analyte in a final assay volume of 50 μ L).

** Only the analytes were prepared in Cell Culture media. All of other components were prepared in Immunoassay Buffer.

Assay Precision:

The following assay precision data were calculated from the three independent assays using two different kit lots. In each lot, the analytes were prepared in AlphaLISA Immunoassay Buffer (IAB), Neurobasal A medium, or DMEM / F12 medium. Each assay consisted of one standard curve comprising 12 data points in triplicate and 12 background wells containing no analyte. The assays were performed in a 384-well format using AlphaLISA Immunoassay Buffer.

- Intra-assay precision:

The intra-assay precision was determined using 3 independent experiments for a total of 16 independent determinations in triplicate. CV% were calculated for each individual experiment then averaged. Shown is the average intra-experimental CV%.

A β 40	IAB	Neurobasal	DMEM/F12
CV%	5	8	8

- Inter-assay precision:

The inter-assay precision was determined using the data across 3 independent experiments with 16 measurements in triplicate. CV% was calculated by comparing the same measurement in each experiment. The CV% for all 16 measurements were then averaged. Shown is the inter-experimental CV%.

A β 40	IAB	Neurobasal	DMEM/F12
CV%	13	23	21

- Spike Recovery:

Three known concentrations of A β 40 were spiked into AlphaLISA Immunoassay Buffer (IAB), Neurobasal A medium, or DMEM/F12 medium. All samples, including non-spiked Immunoassay Buffers were measured in the assay. The average recovery was reported from 4 independent experiments each with 3 measurements in triplicate.

Spiked A β 40 (ng/mL)	% Recovery		
	IAB	Neurobasal	DMEM/F12
3	95	80	87
1	87	76	86
0.3	88	80	84

- Specificity for A β 40:

Cross-reactivity of the AlphaLISA A β 40 Kit was tested against the following proteins and was calculated at the A β 40 standard curve EC₅₀ point when performed in AlphaLISA Immunoassay Buffer. Both monomers and aggregates of A β 40 are detected by this assay kit.

Protein	% Cross-reactivity
Amyloid beta 1-38	0
Amyloid beta 1-42	0
Mouse/Rat Amyloid beta 1-40	0.2

Cerebrospinal Fluid Experiments

Pooled normal Cerebrospinal fluid (CSF) was utilized and AlphaLISA Immunoassay Buffer (IAB) was used as the diluent. A β 40 was detected in the normal Human CSF (data not shown). A β 40 is expected to be present at detectable levels in CSF from normal healthy subjects.

- Dilutional Linearity:

Dilutional linearity was determined by serial dilutions of Human CSF supplemented with 10 ng/mL of A β 40 then diluted with IAB.

Dilution Factor	% Recovery
1	43%
2	60%
4	82%
8	95%
16	100%
32	102%
64	101%
128	90%
256	100%

Troubleshooting Guide

You will find detailed recommendations for common situations you might encounter with your AlphaLISA Assay kit at:

http://www.perkinelmer.com/in/resources/technicalresources/applicationsupportknowledgebase/alphalisa-alphascreen-no-washassays/alpha_troubleshoot.xhtml

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

This product is not for resale or distribution except by authorized distributors.

This product is not for resale or distribution except by authorized distributors.

LIMITED WARRANTY: PerkinElmer BioSignal Inc. warrants that, at the time of shipment, the products sold by it are free from defects in material and workmanship and conform to specifications which accompany the product. PerkinElmer BioSignal Inc. makes no other warranty, express or implied with respect to the products, including any warranty of merchantability or fitness for any particular purpose. Notification of any breach of warranty must be made within 60 days of receipt unless otherwise provided in writing by PerkinElmer BioSignal Inc. No claim shall be honored if the customer fails to notify PerkinElmer BioSignal Inc. within the period specified. The sole and exclusive remedy of the customer for any liability of PerkinElmer BioSignal Inc. of any kind including liability based upon warranty (express or implied whether contained herein or elsewhere), strict liability contract or otherwise is limited to the replacement of the goods or the refunds of the invoice price of goods. PerkinElmer BioSignal Inc. shall not in any case be liable for special, incidental or consequential damages of any kind.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright© 2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.