

Human Amyloid beta 1-40 ELISA Kit ab193692

[3 References](#) [1 Image](#)

Overview

Product name Human Amyloid beta 1-40 ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.1 ng/ml

Range 0.1 ng/ml - 100 ng/ml

Recovery Sample specific recovery

Sample type	Average %	Range
Serum	101.3	77% - 125%
Plasma	97.39	68% - 136%
Cell culture media	104.3	71% - 145%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Predicted to work with: Mouse 

Product overview Abcam's Amyloid beta 1-40 Human ELISA Kit (ab193692) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of Human Amyloid beta 1-40 in serum, plasma, and cell culture supernatants. (UniprotID: P05067)

This assay employs an antibody specific for Human Amyloid beta 1-40 coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures Amyloid beta 1-40 present in the samples. The wells are washed and biotinylated anti-Human Amyloid beta 1-40 antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of Amyloid beta 1-40 bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

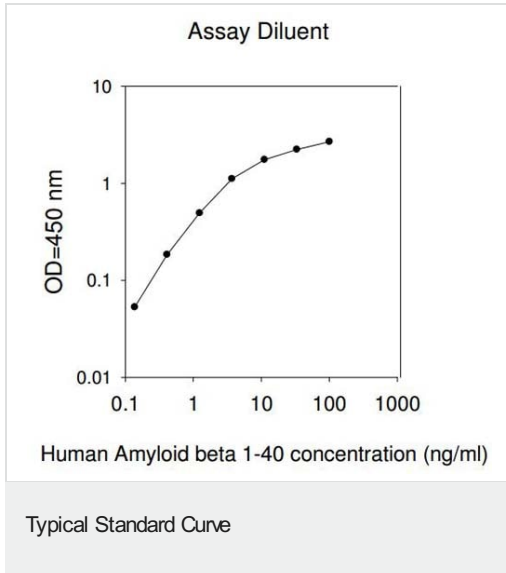
Platform Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
1100X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Biotinylated Human Amyloid beta detection antibody	2 vials
Human Amyloid beta Standards (lyophilized)	2 vials
Pre-coated Human amyloid beta 1-40 microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Images



Standard Curve using ab193692 Amyloid beta 1-40 Human ELISA Kit.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors