

AlphaLISA®

Human Amyloid β 1-40 (A β 1-40) Detection Kit**Product number:** AL202 C **Lot Number:** 3205725**Material provided:****Kit Format:** AL202C: 500 assay points AL202F: 5000 assay points

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

Analyte Format: AL202S: 0.3 μ g (one vial contains an amount of human Amyloid β 1-40 sufficient for performing 10 AlphaLISA standard curves)**Manufacturing date:** xx/xx/xx **Document version:** 1**Product Information**

Application: This kit is designed for the quantitative determination of human Amyloid β 1-40 in cerebrospinal fluid (CSF), buffered solution or cell culture medium using a homogeneous AlphaLISA assay (no wash steps). One antibody is specific to amino acids 17-24 (epitope VFFAE): mouse monoclonal antibody, clone number 4G8. The second antibody is specific to the C-terminus: mouse monoclonal antibody, clone number 11A50-B10.

Kit contents: The kit contains 5 components: AlphaLISA Acceptor beads coated with an Anti-Analyte Antibody, Streptavidin-coated Donor beads, Biotinylated Anti-Analyte Antibody, lyophilized analyte and 10X AlphaLISA Immunoassay Buffer.

Sensitivity: Lower Detection Limit (LDL): 87.9 pg/mL

Dynamic Range: 87.9 – 100 000 pg/mL

Storage: Store kit in the dark at +4°C. Store reconstituted analyte at -20°C.

Stability: This product is stable for at least 12 months from the manufacturing date when stored in its original packaging and the recommended storage conditions. Note: Once reconstituted, the human Amyloid β 1-40 analyte is stable for at least 45 days at -20°C

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. 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.

Maximum signal: 249560 counts

Minimum signal: 279 counts

EC₅₀: 11.100 ng/mL

LDL: 78.508 pg/mL

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

AL202-R Rev 01

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a modern, clean appearance.

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.