



Product Data Sheet

Product Name: β -Amyloid (1-42), mouse, rat

Catalog Number: AS-25381 (0.5mg) Lot Number: See label on vial
AS-25231 (1 mg)

Sequence: H-Asp-Ala-Glu-Phe-Gly-His-Asp-Ser-Gly-Phe-Glu-Val-Arg-His-Gln-Lys-
Leu-Val-Phe-Phe-Ala-Glu-Asp-Val-Gly-Ser-Asn-Lys-Gly-Ala-Ile-Ile-Gly-
Leu-Met-Val-Gly-Gly-Val-Val-Ile-Ala-OH (3-letter code)
DAEFGHDSGFVVRHQKLVFFAEDVGSNKGAIIGLMVGGVVIA
(1-letter code)

Molecular Weight: 4419.0

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Using DMSO, reconstitute by adding 50-60 μ l to 1 mg β -Amyloid (1-42). Dilute this stock solution with working buffer, such as PBS or other buffers.

Storage: β -Amyloid (1-42) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Authentic rodent A61718-05 was purchased from AnaSpec (San Jose, CA). A microtiter plate was precoated with BNT77 as the capture antibody. The captured A β was detected with horseradish peroxidase-labeled BA27 or BC05 for the quantitation of A β _{x-40} and A β _{x-42}, respectively. The brain tissue (about 120 mg) from each mouse was homogenized in 4 volumes of Tris-saline (50 mM Tris-HCl, pH 7.4, 150 mM NaCl) by a motor-driven homogenizer-
[Sawamura, N. et al. J. Biol. Chem. 275, 27901 \(2000\).](#)

Published Citations:

Sawamura, N. et al. *J. Biol. Chem.* **275**, 27901 (2000).

Related Product:

Name	Cat #	Size
DHL™ Fluorescent β -Amyloid (1-42) Sampler Kit	AS-72071	1 kit
SensoLyte® 520 β -Secretase Assay Kit	AS-71144	1 kit
SensoLyte® 520 TACE (α -Secretase) Activity Assay Kit *Fluorimetric*	AS-72085	1 kit

For Research Use Only