



Product Data Sheet

Product Name:	β-Amyloid (1-40)-Lys(Biotin-LC)	
Catalog Number:	AS-23518-01 (0.1 mg) AS-23517 (0.5 mg)	Lot Number: See label on vial
Sequence:	H-Asp-Ala-Glu-Phe-Arg-His-Asp-Ser-Gly-Tyr-Glu-Val-His-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-Lys(Biotin-LC)-OH (3-letter code) DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K (Biotin-LC) (1-letter code)	
Molecular Weight:	4797.5	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: β-Amyloid (1-40)-Lys(Biotin-LC) peptide is freely soluble in 1% NH₄OH.

Storage: β-Amyloid (1-40)-Lys(Biotin-LC) 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.

Description: This C-terminal biotinylated Aβ(1-40) contains a 6-carbon long chain (LC) to provide more accessibility for avidin attachment.

Related Products:

Name	Cat #	Size
β-Amyloid (1-40) Binding Peptide, Biotin-labeled Biotin-DWGKGGRWRLWPGASGKTEA	AS-62427	1 mg
β-Amyloid (1-40)-Lys(Biotin)-NH ₂ DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH ₂	AS-61483-01 AS-61483-05	0.1 mg 0.5 mg
Biotin-LC-β-Amyloid (1-40) Biotin-LC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	AS-24645-01 AS-24648 AS-24645	0.1 mg 0.5 mg 1 mg
f3-Amyloid (1-40)-Lys(LC-biotin)-NH ₂ , FAM-labeled 5-FAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(LC-BIOTIN)-NH ₂	AS-61962-01	0.1 mg
Biotin-f3-Amyloid (1-40) Biotin-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	AS-23512-01 AS-23512	0.1 mg 1 mg

For Research Use Only

