

Product Data Sheet

Product Name: GMP beta-Amyloid (1-42), human (Ammonium Salt)

Catalog Number: AS-GMP-20276-1 Lot Number: See label on vial

Size 1 mg NET peptide

Sequence H-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA-

(one-letter code) OH

Sequence NH₂ - Asp - Ala - Glu - Phe - Arg - His - Asp - Ser - Gly - Tyr - Glu - Val - (three-letter code) 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 – Ile - Ala – COOH

Molecular Weight: 4514.1 Da±0.2%

Storage: This peptide is shipped on dry ice. Upon receipt, store lyophilized peptide at

-15°C or lower. Reconstituted peptide can be aliquoted and stored at -15°C

or lower.

Description: This highly pure, well-characterized GMP beta-amyloid (1-42) peptide

has been produced under conditions compliant with 21 CFR part 820 and

ISO 13485.

Features & Benefits

Added assurance for your high profile projects		
Compliant with 21 CFR part 820 and ISO 13485		
Manufactured in controlled areas including ISO 7 cleanroom for downstream		
processing		
Lot to lot reproducibility, consistency, and traceability		

Applications*

Ideal for:

idodi ior.
In Vitro Diagnostics
QC Controls & Standards
Pre-Clinical animal studies

^{*}not to be used as an API in it's current form

Testing

Additional test results available on Certificate of Analysis

Attribute	Result
Peptide Purity	≥ 97%
Endotoxin	< 0.05 EU/mg
Bioburden	<1 CFU/mg
Solubility	Aqueous media (see CoA)
Monomer content	≥ 97%

References:

- 1. Wang X., et.al., Scientific Reports. (8): 4634 (2018)
- 2. Zhao Y., et.al., Neuroimage. (148):296-304 (2017)
- 3. Wang CY.,et.al., Alzheimer's Dement. (3):262-272 (2017)
- 4. Sun L., et al. Nanomedicine. (13):843 (2018)
- 5. Carneiro P., et al. Sensors & Actuators B:Chem. (239):157-65 (2017).
- 6. Wang JS., et al. J Am Soc Mass Spec. (4):786-795 (2018)
- 7. Leinbach A., et al. Clin Chem. 60(7):987-94 (2014).