



Product Data Sheet

Product Name:	GMP beta-Amyloid (1-42), human (Ammonium Salt)	
Catalog Number:	AS-GMP-20276-5	Lot Number: See label on vial
Size	5 mg NET peptide	
Sequence (one-letter code)	H- DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA-OH	
Sequence (three-letter code)	NH ₂ - 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 - 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:

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

*not to be used as an API

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. Nanmedicine.(**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).

Not to be used as an API