



## Product Data Sheet

<b>Product Name:</b>	[Gly <sup>22</sup> ]-β-Amyloid (1-40), Arctic Mutation	
<b>Catalog Number:</b>	AS-61262 (0.5 mg)	<b>Lot Number:</b> See label on vial
<b>Sequence:</b>	H-Asp-Ala-Glu-Phe-Arg-His-Asp-Ser-Gly-Tyr-Glu-Val-His-His-Gln-Lys-Leu-Val-Phe-Phe-Ala-Gly-Asp-Val-Gly-Ser-Asn-Lys-Gly-Ala-Ile-Ile-Gly-Leu-Met-Val-Gly-Gly-Val-Val-OH (3-letter code) DAEFRHDSGYEVHHQKLVFFAGDVGSNKGAIIGLMVGGVV (1-letter code)	
<b>Molecular Weight:</b>	4257.8	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** [Gly<sup>22</sup>]- β-Amyloid peptide is freely soluble in basic buffer.

**Storage:** [Gly<sup>22</sup>]-β-Amyloid 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.

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
[Gly22]- β-Amyloid (1-42), Arctic Mutation DAEFRHDSGYEVHHQKLVFFAGDVGSNKGAIIGLMVGGVVIA	AS-61967-05	0.5 mg
[Gly21]-β-Amyloid (1-40), A21G, Flemish Mutation DAEFRHDSGYEVHHQKLVFFGEDVGSNKGAIIGLMVGGVV	AS-62150	0.5 mg
[Asn7]-β-Amyloid (1-42), Tottori-Japanese Mutation DAEFRHNSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	AS-63324	0.5 mg
[Asn23]-β-Amyloid (1-40), Iowa Mutation DAEFRHDSGYEVHHQKLVFFAENVGSNKGAIIGLMVGGVV	AS-62145	0.5 mg
[Lys22]-β-Amyloid (1-40), Italian Mutation DAEFRHDSGYEVHHQKLVFFAKDVGSNKGAIIGLMVGGVV	AS-62147	0.5 mg

[Gln22, Asn23]- $\beta$ -Amyloid (1-40), E22Q/D23N Dutch/Iowa double mutation DAEFRHDSGYEVHHQKLVFFAQNVGSNKGAIIGLMVGGVV	AS-62146	0.5 mg
[Gln22]- $\beta$ -Amyloid (1-40), Dutch Mutation DAEFRHDSGYEVHHQKLVFFAQDVGSNKGAIIGLMVGGVV	AS-61261	0.5 mg

*For Research Use Only*