



## Product Data Sheet

---

<b>Product Name:</b>	Amyloid Bri (1-23)	
<b>Catalog Number:</b>	AS-65579 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	Glu-Ala-Ser-Asn-Cys-Phe-Ala-Ile-Arg-His-Phe-Glu-Asn-Lys-Phe-Ala-Val-Glu-Thr-Leu-Ile-Cys-Ser-OH  EASNCFAIRHFENKFAVETLICS-OH	
<b>Molecular Weight:</b>	2631.9	
<b>Peptide Purity:</b>	Peak Area by HPLC $\geq$ 95%	
<b>Storage:</b>	This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at $-20^{\circ}\text{C}$ or lower. Reconstituted peptide can be aliquoted and stored at $-20^{\circ}\text{C}$ or lower.	
<b>Description:</b>	Amyloid Bri (1-23) and Amyloid Bri (1-34) peptides are post-translational products of mutated <i>BRI2</i> gene. Amyloid Bri peptides are found as plaques in brains of patients with Familiar British Dementia (FBD) that shares many common features with Alzheimer's Disease.	
<b>Related Products:</b>	Amyloid Bri (1-34), Cat# AS-65580 Beta-Amyloid (1-40), Cat#AS-24236 Beta-Amyloid (1-42), Cat#AS-20276	
<b>References:</b>	<ol style="list-style-type: none"><li>1. Lelyveld V., et.al., <i>Spectroscopy</i> <b>(15)</b>, 129-139 (2001)</li><li>2. Cantlon A., et.al., <i>J Biol Chem</i> ,1-27 (2015)</li><li>3. Cantlon A., et.al., <i>J Parkin Dis Alzh Dis.</i> <b>2(1)</b>, 1-21 (2015)</li></ol>	

*For Research Use Only*