

## **Product Data Sheet**

Product Name: Amyloid Bri (1-34)

Catalog Number: AS-65580 (1 mg) Lot Number: See label on vial

Sequence: Glu-Ala-Ser-Asn-Cys-Phe-Ala-Ile-Arg-His-Phe-Glu-Asn-Lys-Phe-Ala-Val-

Glu-Thr-Leu-Ile-Cys-Ser-Arg-Thr-Val-Lys-Lys-Asn-Ile-Ile-Glu-Glu-Asn-OH

EASNCFAIRHFENKFAVETLICSRTVKKNIIEEN-OH

Molecular Weight: 3955.3

Peptide Purity: Peak Area by HPLC ≥95%

Storage: This 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: Amyloid Bri (1-23) and Amyloid Bri (1-34) peptides are post-translational

products of the mutated *BRI2* gene. Amyloid Bri peptides are found as plaques in the brains of patients with Familiar British Dementia (FBD), a

condition that shares common features with Alzheimer's Disease.

Related Products: Amyloid Bri (1-23), Cat# AS-65579

Beta-Amyloid (1-40), Cat#AS-24236 Beta-Amyloid (1-42), Cat#AS-20276

References: 1. Lelyveld V., et.al., Spectroscopy (15), 129-139 (2001)

2. Cantlon A., et.al., J Biol Chem ,1-27 (2015)

3. Cantlon A., et.al., J Parkin Dis Alzh Dis. 2(1), 1-21 (2015)

For Research Use Only