



Product Data Sheet

Product Name: [Glu1]-Fibrinopeptide B, Glufib

Catalog Number: AS-60501-1 (1 mg) **Lot Number:** See label on vial

Sequence: H-Glu-Gly-Val-Asn-Asp-Asn-Glu-Glu-Gly-Phe-Phe-Ser-Ala-Arg-OH (3-letter code)
EGVNDNEEGFFSAR (1-letter code)

Molecular Weight: 1570.6

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: [Glu1]-Fibrinopeptide is freely soluble in water.

Storage: [Glu1]-Fibrinopeptide 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: This peptide is derived from fibrinopeptide B amino acid residues 1-14. It is used as a mass spec (MS) standard in proteomic research.

Published Citations:

Fu, C. et al. *Mol. Cell. Proteomics* doi:10.1074/mcp.M800580-MCP200 (2009).

Related Products:

Name	Cat #	Size
Fibrinopeptide B, human Pyr-GVNDNEEGFFSAR	AS-60697	1 mg
Fibrinopeptide A, human ADSGEGDFLAEGGGVR	AS-20735	1 mg

For Research Use Only