

PRODUCT INFORMATION



[Glu¹] Fibrinopeptide B (trifluoroacetate salt)

Item No. 44042

Formal Name: (2S,5S,8S,11S,14S,20S,23S,26S,29S,32S,35S,41S)-41-amino-26,32-bis(2-amino-2-oxoethyl)-11,14-dibenzyl-20,23-bis(2-carboxyethyl)-29-(carboxymethyl)-2-(3-guanidinopropyl)-8-(hydroxymethyl)-35-isopropyl-5-methyl-4,7,10,13,16,19,22,25,28,31,34,37,40-tridecaoxo-3,6,9,12,15,18,21,24,27,30,33,36,39-tridecaazatetracontanedioic acid, trifluoroacetate salt

H—Glu—Gly—Val—Asn—Asp—Asn—Glu—Glu—Gly—Phe—

Phe—Ser—Ala—Arg—OH

• XCF₃COOH

Synonyms: GFB, GluFib

Peptide Sequence: EGVNDNEEGFFSAR-OH

MF: C₆₆H₉₅N₁₉O₂₆ • XCF₃COOH

FW: 1,570.6

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

[Glu¹] Fibrinopeptide B is supplied as a solid. A stock solution may be made by dissolving the [Glu¹] fibrinopeptide B in the solvent of choice, which should be purged with an inert gas. [Glu¹] Fibrinopeptide B is slightly soluble (0.1-1 mg/ml) in DMSO.

[Glu¹] Fibrinopeptide B is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

[Glu¹] Fibrinopeptide B is a synthetic peptide that corresponds to amino acids 1-14 of fibrinopeptide B with a glutamate instead of pyro-glutamine in position 1. It has been used as a model peptide and instrument control for peptide analysis by mass spectrometry (MS).^{1,2}

References

1. Kristensen, D.B., Imamura, K., Miyamoto, Y., *et al.* Mass spectrometric approaches for the characterization of proteins on a hybrid quadrupole time-of-flight (Q-TOF) mass spectrometer. *Electrophoresis* **21**(2), 430-439 (2000).
2. Davaliev, K., Kiprijanovska, S., Dimovski, A., *et al.* Comparative evaluation of two methods for LC-MS/MS proteomic analysis of formalin fixed and paraffin embedded tissues. *J. Proteomics* **235**, 104117 (2021).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/17/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM