

PRODUCT INFORMATION



FcIII Peptide (trifluoroacetate salt)

Item No. 44809

Formal Name: L- α -aspartyl-L-cysteinyl-L-alanyl-L-tryptophyl-L-histidyl-L-leucylglycyl-L- α -glutamyl-L-leucyl-L-valyl-L-tryptophyl-L-cysteinyl-L-threonine, cyclic (2 \rightarrow 12)-disulfide, trifluoroacetate salt

Peptide Sequence: D[CAWHLGELVWC]T-OH

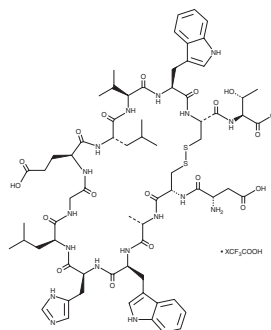
MF: C₆₉H₉₅N₁₇O₁₉S₂ • XCF₃COOH
FW: 1,530.7

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

FcIII peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the FcIII peptide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. FcIII peptide (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in DMSO.

Description

FcIII peptide is a synthetic, cyclic 13-amino acid peptide and fusion tag.¹⁻³ It inhibits binding of the protein A Z domain to IgG Fc with a K_i value of 25 nM.¹ FcIII peptide can be added to the N- or C-terminus of a protein and has been used as an epitope tag on recombinant proteins to facilitate their detection and purification.^{2,3} The inclusion of FcIII peptide on recombinant proteins also increases their plasma half-life in mice.² FcIII peptide has also been used as a building block in the synthesis of artificial human Fc receptors to facilitate IgG uptake in Jurkat cells, which do not intrinsically express Fc receptors.⁴

References

1. DeLano, W.L., Ultsch, M.H., de Vos, A.M., *et al.* Convergent solutions to binding at a protein-protein interface. *Science* **287**(5456), 1279-1283 (2000).
2. Sockolosky, J.T., Kivimäe, S., and Szoka, F.C. Fusion of a short peptide that binds immunoglobulin G to a recombinant protein substantially increases its plasma half-life in mice. *PLoS One* **9**(7), e102566 (2014).
3. Feng, S., Tian, E., Zhang, L., *et al.* Development of the Fc-III tagged protein expression system for protein purification and detection. *PLoS One* **7**(8), e44208 (2012).
4. Boonyarattanakalin, S., Martin, S.E., Sun, Q., *et al.* A synthetic mimic of human Fc receptors: Defined chemical modification of cell surfaces enables efficient endocytic uptake of human immunoglobulin-G. *J. Am. Chem. Soc.* **128**(35), 11463-11470 (2006).

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, 02/10/2026

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