

PRODUCT INFORMATION

HBP08 (trifluoroacetate salt)

Item No. 36961

Formal Name: glycyl-L-tyrosyl-L-histidyl-L-tyrosyl-L-
α-glutamyl-L-arginyl-L-tryptophyl-L-
isoleucyl-L-histidine, trifluoroacetate salt

Peptide Sequence: GYHYERWIIH-OH

MF: C₆₀H₇₇N₁₇O₁₄ • XCF₃COOH

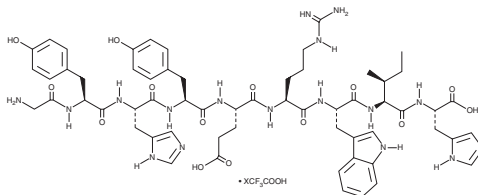
FW: 1,260.4

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

HBP08 (trifluoroacetate salt) is supplied as a solid. HBP08 (trifluoroacetate salt) is sparingly soluble (1-10 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

HBP08 is a peptide inhibitor of the protein-protein interaction between chemokine (C-X-C motif) ligand 12 (CXCL12) and high mobility group protein B1 (HMGB1).¹ It binds to HMGB1 (K_d = 0.8 μM) and inhibits HMGB1-induced inhibition of CXCL12-induced migration of 300-19 pre-B cells expressing human chemokine (C-X-C motif) receptor 4 (CXCR4).

Reference

1. Sgrignani, J., Cecchinato, V., Fassi, E.M.A., *et al.* Systematic development of peptide inhibitors targeting the CXCL12/HMGB1 interaction. *J. Med. Chem.* **64**(18), 13439-13450 (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, 04/04/2024

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