

PRODUCT INFORMATION



Interleukin-8 (54-72) (human) (scrambled) (trifluoroacetate salt)

Item No. 36979

Synonyms: C-X-C Motif Chemokine Ligand (54-72),
CXCL8 (54-72), IL-8 (54-72)

Peptide Sequence: Ac-KVREKNEKWFVEQRVALNS-NH₂ Ac-Lys-Val-Arg-Glu-Lys-Asn-Glu-Lys-Trp-Phe-
Val-Glu-Gln-Arg-Val-Ala-Leu-Asn-Ser-NH₂

MF: C₁₀₇H₁₇₃N₃₃O₃₀ • XCF₃COOH

FW: 2,401.8

Purity: ≥98% • XCF₃COOH

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

Interleukin-8 (IL-8) (54-72) (human) (scrambled) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the IL-8 (54-72) (human) (scrambled) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. IL-8 (54-72) (human) (scrambled) (trifluoroacetate salt) is soluble in the organic solvent DMSO. The solubility of IL-8 (54-72) (human) (scrambled) (trifluoroacetate salt) in DMSO is approximately 5 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of IL-8 (54-72) (human) (scrambled) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of IL-8 (54-72) (human) (scrambled) (trifluoroacetate salt) in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

IL-8 (54-72) (human) (scrambled) is a synthetic peptide with identical amino acid composition to IL-8 (54-72) (human) (Item No. 36978) with 15 out of 19 amino acids in a different position than the original sequence. It has been used as a negative control for IL-8 (54-72) (human) in neutrophil migration assays.¹

Reference

1. Martínez-Burgo, B., Cobb, S.L., Pohl, E., *et al.* A C-terminal CXCL8 peptide based on chemokine-glycosaminoglycan interactions reduces neutrophil adhesion and migration during inflammation. *Immunology* **157**(2), 173-184 (2019).

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, 10/18/2023

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