

PRODUCT INFORMATION



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

Item No. 36978

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

Peptide Sequence: Ac-KENWVQRVVEKFLKRAENS-NH₂ Ac-Lys-Glu-Asn-Trp-Val-Gln-Arg-Val-Val-Glu-

MF: C₁₀₇H₁₇₃N₃₃O₃₀ • XCF₃COOH Lys-Phe-Leu-Lys-Arg-Ala-Glu-Asn-Ser-NH₂

FW: 2,401.8 • XCF₃COOH

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

Interleukin-8 (IL-8) (54-72) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the IL-8 (54-72) (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. IL-8 (54-72) (human) (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of IL-8 (54-72) (human) (trifluoroacetate salt) in these solvents is approximately 10 and 1 mg/ml, respectively.

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) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of IL-8 (54-72) (human) (trifluoroacetate salt) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

IL-8 (54-72) is a peptide that corresponds to amino acids 54-72 of the mature human IL-8 sequence and 81-99 of the pre-processed sequence.¹ It inhibits flow-based adhesion of primary human neutrophils to human umbilical vein endothelial cells (HUVECs).

References

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, 10/19/2023

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