

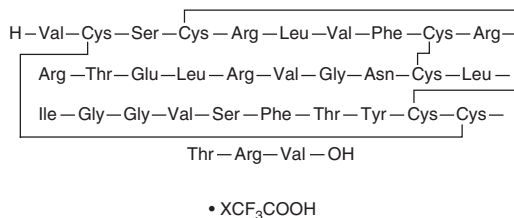
PRODUCT INFORMATION



Corticostatin (human) (trifluoroacetate salt)

Item No. 24888

Synonyms: DEFA4 Protein, α -Defensin 4, Defensin HNP-4, Human Neutrophil Peptide 4, Neutrophil Defensin 4
MF: $C_{157}H_{255}N_{49}O_{43}S_6 \cdot XCF_3COOH$
FW: 3,709.4
Purity: $\geq 95\%$
Supplied as: A lyophilized powder
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Corticostatin (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the corticostatin (human) (trifluoroacetate salt) in water. The solubility of corticostatin (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Corticostatin is a corticostatic peptide that has been found in human granulocytes.¹ It inhibits corticosterone production stimulated by adrenocorticotrophic hormone (ACTH) in rat adrenal cells ($IC_{50} = 0.7 \mu M$). Corticostatin also reduces mitogen-stimulated release of IFN- γ and IL-6 in human peripheral blood mononuclear cells (PBMCs).²

References

1. Singh, A., Bateman, A., Zhu, Q., *et al.* Structure of a novel human granulocyte peptide with anti-ACTH activity. *Biochem. Biophys. Res. Commun.* **155(1)**, 524-529 (1988).
2. Masera, R.G., Bateman, A., Muscettola, M., *et al.* Corticostatins/defensins inhibit in vitro NK activity and cytokine production by human peripheral blood mononuclear cells. *Regul. Pept.* **62(1)**, 13-21 (1996).

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, 11/18/2022

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