

PRODUCT INFORMATION



IDR 1002 (trifluoroacetate salt)

Item No. 36893

Formal Name: L-valyl-L-glutamyl-L-arginyl-L-tryptophyl-L-leucyl-L-isoleucyl-L-valyl-L-tryptophyl-L-arginyl-L-isoleucyl-L-arginyl-L-lysineamide, trifluoroacetate salt

Peptide Sequence: VQRWLIVWRIRK-NH₂

MF: C₇₉H₁₃₀N₂₆O₁₃ • XCF₃COOH

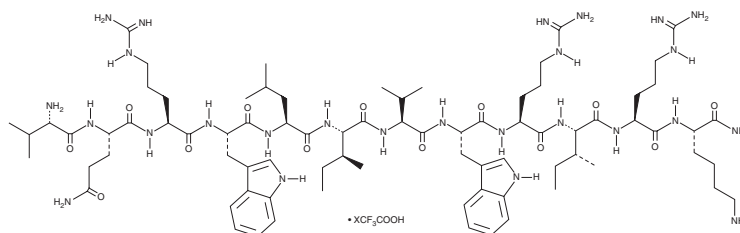
FW: 1,652.0

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

IDR 1002 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the IDR 1002 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. IDR 1002 (trifluoroacetate salt) is slightly soluble in the organic solvent DMSO.

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 IDR 1002 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. IDR 1002 (trifluoroacetate salt) is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

IDR 1002 is a synthetic cationic innate defense regulator peptide.¹ It increases the production of chemokine (C-C motif) ligand 2 (CCL2), CCL7, and chemokine (C-X-C motif) ligand 1 (CXCL1) in isolated human peripheral blood mononuclear cells (PBMCs) when used at a concentration of 100 µg/ml. IDR 1002 (200 µg/animal) increases the numbers of neutrophils and monocytes in mouse peritoneum and protects against *S. aureus* and *E. coli* infections in mice.

Reference

1. Nijnik, A., Madera, L., Ma, S., *et al.* Synthetic cationic peptide IDR-1002 provides protection against bacterial infections through chemokine induction and enhanced leukocyte recruitment. *J. Immunol.* **184**(5), 2539-2550 (2010).

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, 12/21/2022

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