

PRODUCT INFORMATION

Urumin (trifluoroacetate salt)

Item No. 24277

MF: $C_{129}H_{198}N_{42}O_{35}S_2 \cdot XCF_3COOH$ H-Ile-Pro-Leu-Arg-Gly-Ala-Phe-Ile-Asn-Gly-
 FW: 2,961.4 Arg-Trp-Asp-Ser-Gln-Cys-His-Arg-Phe-Ser-
 Purity: $\geq 90\%$ Asn-Gly-Ala-Ile-Ala-Cys-Ala-OH
 Supplied as: A lyophilized powder • XCF_3COOH
 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

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

Description

Urumin is a peptide originally isolated from *H. bahuvistara* (South Indian frog) skin that has antiviral activity.¹ It inhibits H1 hemagglutinin-containing human influenza A viruses by 60 to >90% and H3 hemagglutinin-containing viruses by less than 50%. Urumin specifically targets the H1 hemagglutinin stalk in an antiviral assay using a chimeric virus and in an ELISA assay and acts by actively destroying influenza virions. It reduces viral titers of drug-resistant and -sensitive strains of H1N1 virus in a plaque assay and in mice. Intranasal administration of urumin (20 μg) prior to a lethal dose of PR8 influenza increases survival, with 70% of urumin-treated mice surviving after 14 days compared with 20% of control animals.

Reference

1. Holthausen, D.J., Lee, S.H., Kumar, V.T.V., et al. An amphibian host defense peptide is virucidal for human H1 hemagglutinin-bearing influenza viruses. *Immunity* **46**(4), 587-595 (2017).

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, 10/21/2022

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