

PRODUCT INFORMATION



Acetyl Hexapeptide-38 (trifluoroacetate salt)

Item No. 27258

Formal Name: (S)-N-((2S,3R)-1-amino-3-hydroxy-1-oxobutan-2-yl)-5-guanidino-2-((2S,5S,8S,11S)-11-(hydroxymethyl)-2,5,8-triisopropyl-4,7,10,13-tetraoxo-3,6,9,1-tetraazatetradecanamido) pentanamide, trifluoroacetate

MF: $C_{30}H_{56}N_{10}O_9 \cdot XCF_3COOH$

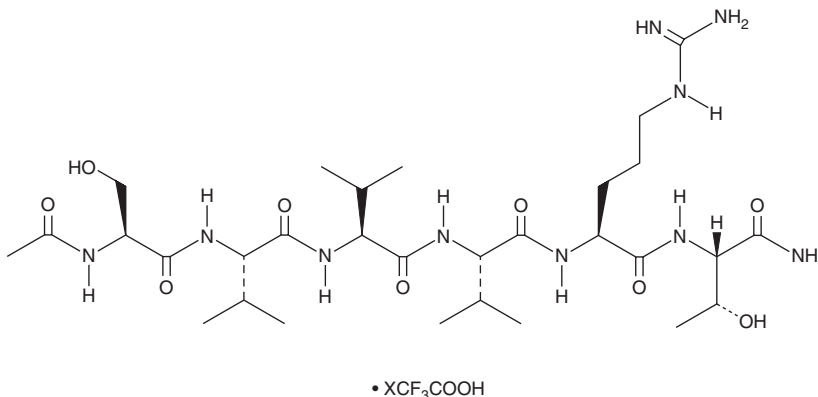
FW: 700.8

Purity: $\geq 98\%$

Supplied as: A solid

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

Acetyl hexapeptide-38 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the acetyl hexapeptide-38 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Acetyl hexapeptide-38 (trifluoroacetate salt) is slightly soluble in DMSO and dimethyl formamide.

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 acetyl hexapeptide-38 (trifluoroacetate salt) can be prepared by directly dissolving the acetyl hexapeptide-38 (trifluoroacetate salt) in aqueous buffers. The solubility of acetyl hexapeptide-38 (trifluoroacetate salt) in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Acetyl hexapeptide-38 is a hexapeptide that decreases IL-8 secretion in human keratinocytes when used alone or in combination with acetyl pentapeptide-46.¹ Formulations containing acetyl hexapeptide-38 have been used to diminish the appearance of skin aging.

Reference

1. Tittl, J. and Wolf, C. Compositions and methods for mitigating skin irritation and enhancing skin barrier function. Gogo Industries, Inc. U.S. 2016/0000858 A1 (2016).

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