

PRODUCT INFORMATION



HCKFWW (trifluoroacetate salt)

Item No. 39157

Formal Name: L-histidyl-L-cysteinyl-L-lysyl-L-phenylalanyl-L-tryptophyl-L-tryptophan, trifluoroacetate salt

Synonyms: His-Cys-Lys-Phe-Trp-Trp-OH, HW6

Peptide Sequence: HCKFWW-OH

MF: C₄₆H₅₅N₁₁O₇S • XCF₃COOH

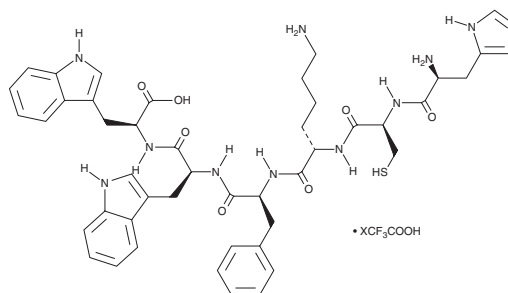
FW: 906.1

Purity: ≥98%

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

HCKFWW (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the HCKFWW (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

HCKFWW is a thiol-containing peptide and form of the HIV-1 integrase inhibitor HCKFWW-NH₂. It has been used as an analyte in the development of zirconium nanoparticles as probes and matrices for mass spectrometry applications.¹

Reference

1. Kailasa, S.K. and Wu, H.-F. Multifunctional ZrO₂ nanoparticles and ZrO₂-SiO₂ nanorods for improved MALDI-MS analysis of cyclodextrins, peptides, and phosphoproteins. *Anal. Bioanal. Chem.* **396**(3), 1115-1125 (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, 01/09/2024

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