

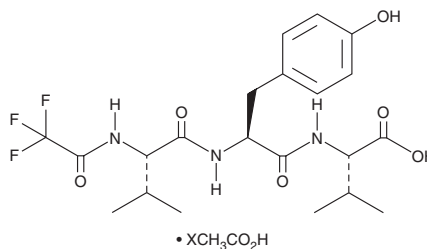
PRODUCT INFORMATION



Trifluoroacetyl Tripeptide-2 (acetate)

Item No. 35389

Formal Name:	N-(2,2,2-trifluoroacetyl)-L-valyl-L-tyrosyl-L-valine, acetate
Synonym:	F ₃ Ac-Val-Tyr-Val
MF:	C ₂₁ H ₂₈ F ₃ N ₃ O ₆ • XC ₂ H ₄ O ₂
FW:	475.5
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

Trifluoroacetyl tripeptide-2 (acetate) is supplied as a solid. A stock solution may be made by dissolving the trifluoroacetyl tripeptide-2 (acetate) in the solvent of choice, which should be purged with an inert gas. Trifluoroacetyl tripeptide-2 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of trifluoroacetyl tripeptide-2 (acetate) in these solvents is approximately 15, 2, and 10 mg/ml, respectively.

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 trifluoroacetyl tripeptide-2 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of trifluoroacetyl tripeptide-2 (acetate) in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Trifluoroacetyl tripeptide-2 is a peptide inhibitor of pancreatic and neutrophil elastases (K_i s = 0.00056 and 0.5 mM, respectively).¹

Reference

1. Dimicoli, J.L. and Bieth, J. On the inhibition of human leukocyte elastase by trifluoroacetyl tripeptides. *J. Biol. Chem.* **253**(10), 3459-3460 (1978).

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, 11/08/2022

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