

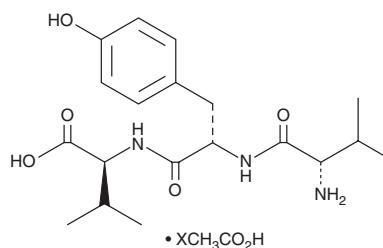
PRODUCT INFORMATION



Val-Tyr-Val-OH (acetate)

Item No. 43414

Formal Name: L-valyl-L-tyrosyl-L-valine, acetate
Synonyms: NSC 339922, Valine-Tyrosine-Valine
Peptide Sequence: VYV-OH
MF: $C_{19}H_{29}N_3O_5 \cdot XC_2H_4O_2$
FW: 379.5
Purity: $\geq 98\%$
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

Val-Tyr-Val-OH (acetate) is supplied as a solid. A stock solution may be made by dissolving the Val-Tyr-Val-OH (acetate) in the solvent of choice, which should be purged with an inert gas. Val-Tyr-Val-OH (acetate) is soluble (≥ 10 mg/ml) in 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 Val-Tyr-Val-OH (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Val-Tyr-Val-OH (acetate) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Val-Tyr-Val-OH is a tripeptide that has been used as a standard in HPLC, electrospray ionization mass spectroscopy (ESI-MS), or capillary electrophoresis (CE) analytical method development.¹⁻³

References

1. Uchiho, Y., Goto, Y., Kamahori, M., *et al.* New baseline correction method using near-infrared absorption of water in water/acetonitrile gradient high-performance liquid chromatography with far-ultraviolet absorbance detection. *Chromatographia* **80**, 329-333 (2017).
2. Wang, N.-H., Lee, W.-L., and Her, G.-R. Signal enhancement for peptide analysis in liquid chromatography-electrospray ionization mass spectrometry with trifluoroacetic acid containing mobile phase by postcolumn electrophoretic mobility control. *Anal. Chem.* **83**(16), 6163-6168 (2011).
3. Ferris, S.S., Lou, G., and Ewing, A.G. Scanning electrochemical detection in capillary electrophoresis. *J. Microcolumn. Sep.* **6**(3), 263-268 (1994).

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, 05/12/2025

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