

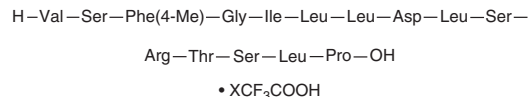
PRODUCT INFORMATION



VPM-p15 (trifluoroacetate salt)

Item No. 37470

Synonym: T1V/F3Phe(4-Me)
Peptide Sequence: VSXGILLDLRSTSLP-OH (X = 4-methylphenylalanine)
MF: C₇₅H₁₂₆N₁₈O₂₂ • XCF₃COOH
FW: 1,631.9
Purity: ≥95%
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

VPM-p15 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the VPM-p15 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. VPM-p15 (trifluoroacetate salt) is slightly soluble in 30% acetonitrile/water.

VPM-p15 (trifluoroacetate salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

VPM-p15 is a synthetic peptide agonist for the adhesion G protein-coupled receptor (GPCR) G₂ (ADGRG2), an orphan GPCR involved in male infertility.¹ It induces cAMP accumulation in HEK293 cells expressing human ADGRG2 (EC₅₀ = 1.41 μM).

Reference

1. Sun, Y., Zhang, D., Ma, M.L., *et al.* Optimization of a peptide ligand for the adhesion GPCR ADGRG2 provides a potent tool to explore receptor biology. *J. Biol. Chem.* **296**, 100174 (2021).

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, 02/15/2024

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