

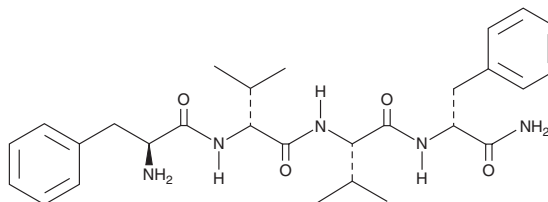
PRODUCT INFORMATION



Bilaid A1

Item No. 31243

CAS Registry No.: 2393866-02-7
Formal Name: L-phenylalanyl-D-valyl-L-valyl-D-phenylalaninamide
Synonyms: H-FVVF-NH₂, H-L-Phe-D-Val-L-Val-D-Phe-NH₂
MF: C₂₈H₃₉N₅O₄
FW: 509.6
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

Bilaid A1 is supplied as a solid. A stock solution may be made by dissolving the bilaid A1 in the solvent of choice, which should be purged with an inert gas. Bilaid A1 is soluble in DMSO and methanol.

Description

Bilaid A1 is a tetrapeptide agonist of the μ -opioid receptor ($K_i = 750$ nM in HEK293 cell membranes expressing the human receptor) and a derivative of bilaid A (Item No. 31242).¹ It inhibits forskolin-induced cAMP accumulation by 47% in HEK293 cells expressing the human μ -opioid receptor when used at a concentration of 10 μ M.

Reference

1. Dekan, Z., Sianati, S., Yousuf, A., *et al.* A tetrapeptide class of biased analgesics from an Australian fungus targets the μ -opioid receptor. *Proc. Natl. Acad. Sci. USA* **116**(44), 22353-22358 (2019).

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, 11/04/2022

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