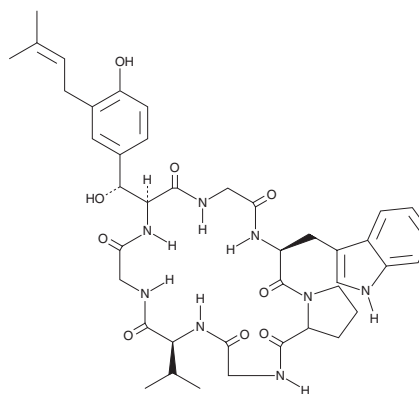


PRODUCT INFORMATION



WIN 66,306
Item No. 33161

CAS Registry No.: 151928-32-4
Formal Name: cyclo[glycyl-L-tryptophyl-L-prolylglycyl-L-valylglycyl-(βR)-β-hydroxy-3-(3-methyl-2-butenyl)-L-tyrosyl]
MF: C₄₁H₅₂N₈O₉
FW: 800.9
Purity: ≥95%
Supplied as: A residue
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Aspergillus nanangensis*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

WIN 66,306 is supplied as a residue. A stock solution may be made by dissolving the WIN 66,306 in the solvent of choice, which should be purged with an inert gas. WIN 66,306 is soluble in organic solvents such as methanol and DMSO.

Description

WIN 66,306 is a cyclic heptapeptide antagonist of neurokinin-1 (NK₁) and NK₂ receptors originally isolated from *A. flavipes*.¹ It binds to NK₁ and NK₂ receptors (K_is = 8 and 3.8 μM, respectively, for the human receptors) and inhibits contractions induced by substance P (Item No. 24035) in isolated guinea pig ileum in a concentration-dependent manner.

Reference

1. Barrow, C.J., Doleman, M.S., Bobko, M.A., *et al.* Structure determination, pharmacological evaluation, and structure-activity studies of a new cyclic peptide substance P antagonist containing the new amino acid 3-prenyl-β-hydroxytyrosine, isolated from *Aspergillus flavipes*. *J. Med. Chem.* **37(3)**, 356-363 (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, 08/24/2023

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