

PRODUCT INFORMATION



PG-106 (trifluoroacetate salt)

Item No. 41553

Formal Name: N-acetyl-L-norleucyl-L- α -aspartyl- β -alanyl-3-(2-naphthalenyl)-D-alanyl-L-arginyl-L-tryptophyl-L-lysynamide, (2 \rightarrow 7)-lactam, trifluoroacetate salt

Synonym: Ac-Nle-c[Asp- β Ala-DNal(2')-Arg-Trp-Lys]-NH₂

MF: C₅₁H₆₉N₁₃O₉ • XCF₃COOH

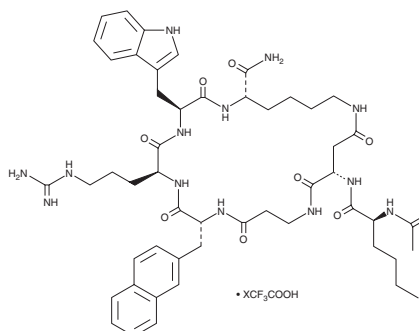
FW: 1,008.2

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

PG-106 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PG-106 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. PG-106 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in DMSO.

Description

PG-106 is a cyclic peptidomimetic antagonist of melanocortin receptor 3 (MC3R; IC₅₀ = 210 nM).¹ It is selective for MC3R over MC4R and MC5R at 10 μ M.

Reference

1. Grieco, P., Cai, M., Han, G., *et al.* Further structure-activity studies of lactam derivatives of MT-II and SHU-9119: Their activity and selectivity at human melanocortin receptors 3, 4, and 5. *Peptides* **28(6)**, 1191-1196 (2007).

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, 11/04/2025

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