

PRODUCT INFORMATION



GR94800

Item No. 38773

CAS Registry No.: 141636-65-9

Formal Name: N-benzoyl-L-alanyl-L-alanyl-D-tryptophyl-L-phenylalanyl-D-prolyl-L-prolyl-L-norleucinamide

Synonyms: Benzoyl-Ala-Ala-trp-Phe-pro-Pro-Nle-NH₂, GR98400

Peptide Sequence: X-AAwFpPZ-NH₂ (X = Benzoyl, Z = Norleucine)

MF: C₄₉H₆₁N₉O₈

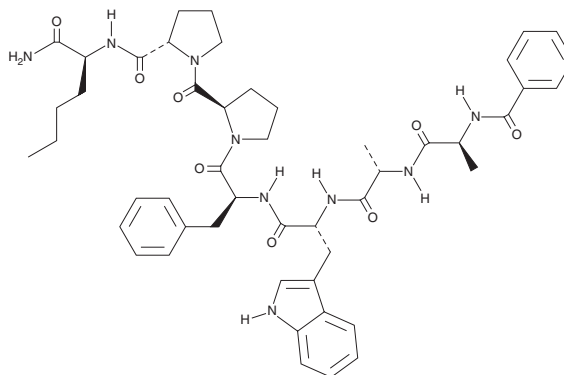
FW: 904.1

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

GR94800 is supplied as a solid. A stock solution may be made by dissolving the GR94800 in the solvent of choice, which should be purged with an inert gas. GR94800 is soluble in the organic solvent acetonitrile.

Description

GR94800 is a synthetic peptide neurokinin-2 (NK₂) receptor antagonist (K_i = 1 nM).¹ It is selective for NK₂ over NK₃ (K_i = 1,000 nM). GR94800 (0.1 and 1 μM) decreases balloon distension-induced submaximal and maximal atropine-resistant ascending excitatory reflex (AER) in isolated guinea pig ileum circular muscles.²

References

1. Antoniou, M. and Poulos, C. Analogues of the C-terminal fragments of neurokinins with modifications at their C-terminal methionyl residue. Structure-activity studies. *Int. J. Pept. Protein Res.* **43(4)**, 344-350 (1994).
2. Maggi, C.A., Patacchini, R., Bartho, L., *et al.* Tachykinin NK₁ and NK₂ receptor antagonists and atropine-resistant ascending excitatory reflex to the circular muscle of the guinea-pig ileum. *Br. J. Pharmacol.* **112(1)**, 161-168 (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, 12/20/2023

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