

PRODUCT INFORMATION



GR83074 (trifluoroacetate salt)

Item No. 44017

Formal Name: N²-[(1,1-dimethylethoxy)carbonyl]-L-arginyl-L-alanyl-D-tryptophyl-L-phenylalanyl-D-prolyl-L-prolyl-L-norleucinamide, trifluoroacetate salt

Synonym: Boc-Arg-Ala-D-Trp-Phe-D-Pro-Pro-Nle-NH₂

Peptide Sequence: Boc-RAwFpPX-NH₂ (X = norisoleucine)

MF: C₅₀H₇₂N₁₂O₉ • XCF₃COOH

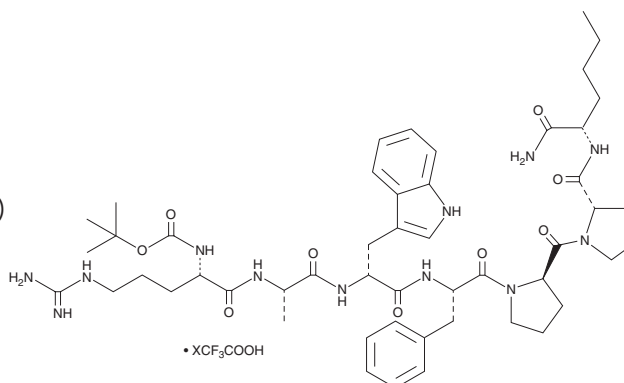
FW: 985.2

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

GR83074 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the GR83074 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. GR83074 (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

GR83074 (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

GR83074 is a peptide antagonist of the neurokinin-2 (NK₂) receptor.¹ It selectively inhibits neurokinin A-induced contractions in rat colon muscularis mucosae, which express NK₂ receptors, over substance P-induced contractions in guinea pig ileum longitudinal smooth muscle strips, which express NK₁ receptors (pK_bs = 8.2 and 5.8, respectively). GR83074 (0.3 μmol/kg) inhibits neurokinin A-induced bronchoconstriction in guinea pigs.

Reference

1. McElroy, A.B., Clegg, S.P., Deal, M.J., *et al.* Highly potent and selective heptapeptide antagonists of the neurokinin NK-2 receptor. *J. Med. Chem.* **35**(14), 2582-2591 (1992).

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, 10/14/2025

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