

PRODUCT INFORMATION



R 396 (trifluoroacetate salt)

Item No. 43083

Formal Name: N-acetyl-L-leucyl-L- α -aspartyl-L-glutamyl-L-tryptophyl-L-phenylalanyl-glycinamide, trifluoroacetate salt

Synonyms: Ac-Leu-Asp-Gln-Trp-Phe-Gly-NH₂

Peptide Sequence: Ac-LDQWFG-NH₂

MF: C₃₉H₅₁N₉O₁₀ • XCF₃COOH

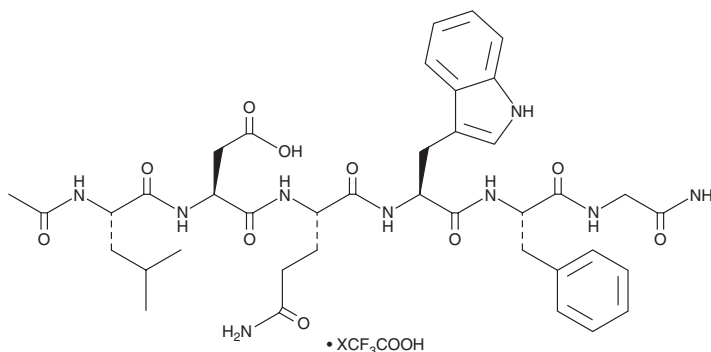
FW: 805.9

Purity: ≥90%

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

R 396 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the R 396 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. R 396 (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO.

Description

R 396 is a peptide antagonist of the neurokinin-2B (NK_{2B}) receptor (pA₂ = 7.63).¹ It is selective for the NK_{2B} receptor over NK₁, NK_{2A}, and NK₃ receptors (pA₂s = <5, 5.41, and <5, respectively). Intracerebroventricular administration of R 396 (650 pmol/animal) inhibits increases in mean arterial blood pressure, heart rate, face washing, and grooming induced by neurokinin A (NKA) but not substance P (Item No. 24035) in conscious rats.²

References

1. Rovero, P., Astolfi, M., Manzini, S., *et al.* Structure-activity relationship study of R396, an NK₂ tachykinin antagonist selective for the NK_{2B} receptor subtype. *Neuropeptides* **23(3)**, 143-145 (1992).
2. Tschöpe, C., Picard, P., Culman, J., *et al.* Use of selective antagonists to dissociate the central cardiovascular and behavioural effects of tachykinins on NK₁ and NK₂ receptors in the rat. *Br. J. Pharmacol.* **107(3)**, 750-755 (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, 06/30/2025

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