

PRODUCT INFORMATION



GR64349 (acetate)

Item No. 42915

Formal Name: (S)-4-(((S)-1-(((S)-1-(((S)-1-(((R)-1-((S)-1-(((S)-1-amino-4-(methylthio)-1-oxobutan-2-yl)amino)-4-methyl-1-oxopentan-2-yl)-2-oxopyrrolidin-3-yl)amino)-3-methyl-1-oxobutan-2-yl)amino)-1-oxo-3-phenylpropan-2-yl)amino)-3-hydroxy-1-oxopropan-2-yl)amino)-3-((S)-2,6-diaminohexanamido)-4-oxobutanoic acid

Synonyms: Lys-Asp-Ser-Phe-Val-[R-γ-Lac]-Leu-Met-NH₂
[Lys³,Gly⁸-R-Lac-Leu⁹]NKA (3-10),
[Lys³,Gly⁸-R-γ-lactam-Leu⁹]NKA (3-10),
[Lys³,Gly⁸-R-Lac-Leu⁹]Neurokinin A (3-10),

Peptide Sequence: KDSFV-[R-γ-Lac]-LM-NH₂

MF: C₄₂H₆₈N₁₀O₁₁S • C₂H₄O₂

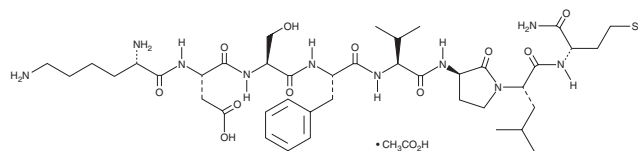
FW: 981.2

Purity: ≥98%

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

GR64349 (acetate) is supplied as a solid. A stock solution may be made by dissolving the GR64349 (acetate) in the solvent of choice, which should be purged with an inert gas. GR64349 (acetate) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

GR64349 (acetate) is slightly soluble (0.1-1 mg/ml) 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

GR64349 is a peptide agonist of the neurokinin-2 (NK₂) receptor.¹ It selectively induces contractions in isolated rat colon muscularis mucosae, which endogenously express NK₂ receptors, over isolated guinea pig ileum longitudinal smooth muscle and everted rat portal vein, which endogenously express NK₁ and NK₃ receptors, respectively (EC₅₀s = 3.7, 4,237, and 1,177 nM, respectively). GR64349 also induces contractions in isolated guinea pig trachea in a concentration-dependent manner.² *In vivo*, GR64349 (1.1 nmol/animal) induces contralateral rotations in rats when infused into the substantia nigra.³

References

1. Deal, M.J., Hagan, R.M., Ireland, S.J., *et al.* Conformationally constrained tachykinin analogues: Potent and highly selective neurokinin NK-2 receptor agonists. *J. Med. Chem.* **35(22)**, 4195-4204 (1992).
2. Beresford, I.J., Sheldrick, R.L., Ball, D.I., *et al.* GR159897, a potent non-peptide antagonist at tachykinin NK2 receptors. *Eur. J. Pharmacol.* **272(2-3)**, 241-248 (1995).
3. Elliott, P.J., Mason, G.S., Stephens-Smith, M., *et al.* Behavioural and biochemical responses following activation of midbrain dopamine pathways by receptor selective neurokinin agonists. *Neuropeptides* **19(2)**, 119-126 (1991).

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, 05/12/2025

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