

PRODUCT INFORMATION

Galanin (2-11) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt)

Item No. 34065

Formal Name: L-tryptophyl-L-threonyl-L-leucyl-L-asparaginyl-L-seryl-L-alanylglycyl-L-tyrosyl-L-leucyl-L-leucinamide, trifluoroacetate salt

Synonyms: AR-M1896, Gal(2-11)-NH₂, WTLNSAGYLL-NH₂

MF: C₅₄H₈₁N₁₃O₁₄ • XCF₃COOH

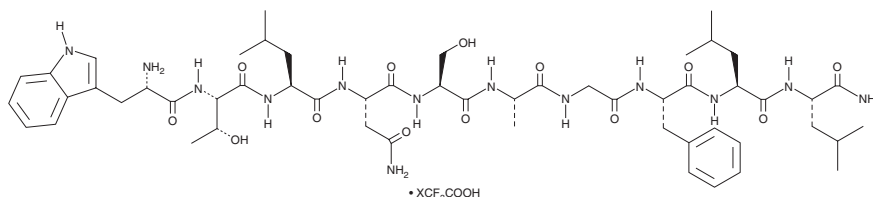
FW: 1,136.3

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

Galanin (2-11) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the galanin (2-11) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Galanin (2-11) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of galanin (2-11) amide (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in these solvents is approximately 5 mg/ml.

Description

Galanin (2-11) amide is a synthetic peptide fragment of the neuropeptide galanin and an agonist of the galanin-2 (GAL₂) receptor (EC₅₀ = 9.32 nM in a fluorescence imaging plate reader assay).¹ It selectively binds to the GAL₂ receptor (IC₅₀ = 1.76 nM for the rat receptor) over the GAL₁ receptor (IC₅₀ = 879 nM for the human receptor) but does also bind to the GAL₃ receptor (K_i = 271 nM for the rat receptor).^{1,2} Intracerebroventricular administration of galanin (2-11) amide (1 nmol/animal) decreases immobility in the forced swim test in rats.³ It also reduces the hind paw mechanical pain threshold and increases the hind paw cold sensitivity threshold in rats.¹

References

1. Liu, H.-X., Brumovsky, P., Schmidt, R., *et al.* Receptor subtype-specific pronociceptive and analgesic actions of galanin in the spinal cord: Selective actions via GalR1 and GalR2 receptors. *Proc. Natl. Acad. Sci. USA* **98(17)**, 9960-9964 (2001).
2. Lu, X., Lundström, L., and Bartfai, T. Galanin (2-11) binds to GalR3 in transfected cell lines: Limitations for pharmacological definition of receptor subtypes. *Neuropeptides* **39(3)**, 165-167 (2005).
3. Kuteeva, E., Wardi, T., Lundström, L., *et al.* Differential role of galanin receptors in the regulation of depression-like behavior and monoamine/stress-related genes at the cell body level. *Neuropsychopharmacology* **33(11)**, 2573-2585 (2008).

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.

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