

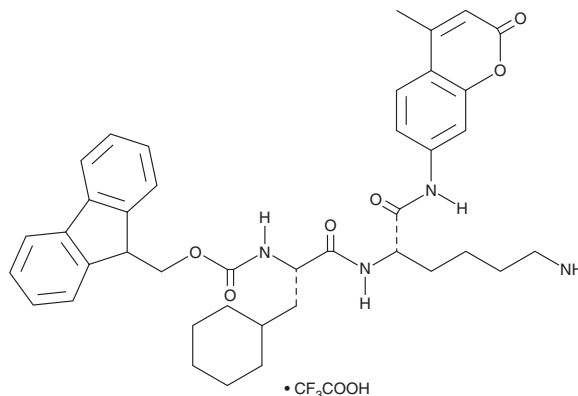
PRODUCT INFORMATION



Galnon (trifluoroacetate salt)

Item No. 29925

CAS Registry No.: 1217448-19-5
Formal Name: 3-cyclohexyl-N-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-alanyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-lysine, trifluoroacetate salt
Synonym: Fmoc-β-Cha-Lys-AMC
MF: C₄₀H₄₆N₄O₆ • CF₃COOH
FW: 792.9
Purity: ≥98%
Supplied as: A crystalline 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

Galnon (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the galnon (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Galnon (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of galnon (trifluoroacetate salt) in these solvents is approximately 10 mg/ml.

Galnon (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, galnon (trifluoroacetate salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Galnon (trifluoroacetate salt) has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Galnon is a galanin (GAL) receptor agonist ($K_D = 2.9 \mu\text{M}$ in Bowes cells expressing GAL₁ receptors).¹ It inhibits forskolin-induced adenylate cyclase activity in rat hippocampal membranes ($EC_{50} = 10 \mu\text{M}$). Galnon (2 mg/kg) reduces the maximal seizure score and increases the latency to first seizure in a mouse model of seizure induced by pentylenetetrazole (PTZ; Item No. 18682). It reduces cocaine-induced motor activity and frontal cortex, but not nucleus accumbens, dopamine overflow, as well as blocks cocaine-primed reinstatement of cocaine-seeking behavior in rats.² Galnon (1.5-5 mg/kg) reduces food intake in rats and mice.³

References

1. Saar, K., Mazarati, A.M., Mahlapuu, R., *et al.* Anticonvulsant activity of a nonpeptide galanin receptor agonist. *Proc. Natl. Acad. Sci. USA* **99**(10), 7136-7141 (2002).
2. Ogbonmwan, Y.E., Sciolino, N.R., Groves-Chapman, J.L., *et al.* The galanin receptor agonist, galnon, attenuates cocaine-induced reinstatement and dopamine overflow in the frontal cortex. *Addict. Biol.* **20**(4), 701-713 (2015).
3. Abramov, U., Florén, A., Echevarria, D.J., *et al.* Regulation of feeding by galnon. *Neuropeptides* **38**(1), 55-61 (2004).

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, 04/01/2020

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