

PRODUCT INFORMATION



Neuropeptide γ (trifluoroacetate salt)

Item No. 43052

Formal Name:	L- α -aspartyl-L-alanylglycyl-L-histidylglycyl-L-glutaminy-L-isoleucyl-L-seryl-L-histidyl-L-lysyl-L-arginyl-L-histidyl-L-lysyl-L-threonyl-L- α -aspartyl-L-seryl-L-phenylalanyl-L-valylglycyl-L-leucyl-L-methioninamide, trifluoroacetate salt	H-Asp-Ala-Gly-His-Gly-Gln-Ile-Ser-His-Lys-Arg-His-Lys-Thr-Asp-Ser-Phe-Val-Gly-Leu-Met-NH ₂
Synonyms:	γ -Neuropeptide, NPG, γ -PPT (72-92), γ -Preprotachykinin-1 (72-92)	
Peptide Sequence:	DAGHGQISHKRHKTDSEFVGLM-NH ₂	
MF:	C ₉₉ H ₁₅₈ N ₃₄ O ₂₉ S • XCF ₃ COOH	
FW:	2,320.6	
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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of neuropeptide γ (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Neuropeptide γ (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Neuropeptide γ is a peptide agonist of the neurokinin-2 (NK₂) receptor.¹ It induces phosphatidylinositol turnover in hamster urinary bladder preparations (EC₅₀ = 7.1 nM) and smooth muscle contractions in isolated hamster bladder (EC₅₀ = 112 nM). Neuropeptide γ selectively binds to NK₂ receptors over NK₁ and NK₃ receptors (IC₅₀s = 0.2, 3.16, and 50.1 nM, respectively). It induces salivation in rats in a dose-dependent manner.² Intrathecal administration of neuropeptide γ (0.078-78 nmol/animal) increases blood pressure and heart rate in conscious rats.³

References

1. van Giersbergen, P.L., Shatzer, S.A., Burcher, E., *et al.* Comparison of the effects of neuropeptide K and neuropeptide γ with neurokinin A at NK₂ receptors in the hamster urinary bladder. *Naunyn Schmiedebergs Arch. Pharmacol.* **345**(1), 51-56 (1992).
2. Takeda, Y. and Krause, J.E. γ -Preprotachykinin-(72-92)-peptide amide potentiates substance P-induced salivation. *Eur. J. Pharmacol.* **161**(2-3), 267-271 (1989).
3. Poulat, P., de Champlain, J., and Couture, R. Cardiovascular responses to intrathecal neuropeptide γ in conscious rats: Receptor characterization and mechanism of action. *Br. J. Pharmacol.* **117**(2), 250-257 (1996).

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, 06/13/2025

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