

PRODUCT INFORMATION



Neuromedin N (trifluoroacetate salt)

Item No. 24707

Formal Name:	L-lysyl-L-isoleucyl-L-prolyl-L-tyrosyl-L-isoleucyl-L-leucine, trifluoroacetate salt	
Synonym:	NMN	
MF:	$C_{38}H_{63}N_7O_8 \cdot XCF_3COOH$	H—Lys—Ile—Pro—Tyr—Ile—Leu—OH
FW:	745.9	• XCF_3COOH
Purity:	≥95%	
Supplied as:	A lyophilized powder	
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

Neuromedin N (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the neuromedin N (trifluoroacetate salt) in water. The solubility of neuromedin N (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Neuromedin N is a neurotensin-like hexapeptide that is synthesized from the same precursor as neurotensin.¹ It is highly expressed in the brain but is also expressed in the gastrointestinal system and adrenal glands. It competitively inhibits neurotensin binding to rat brain synaptic membranes and has increased potency in the presence of the peptidase inhibitor bestatin ($IC_{50}S = 16.7$ and 3 nM, respectively).² Neuromedin N (10 pM) increases the phagocytic function of isolated murine peritoneal macrophages and increases chemotaxis two-fold compared with control, an effect that is blocked by the calcium-channel blocker ryanodine.³ It induces hypotension in rats and contraction of isolated guinea pig ileum.⁴ Neuromedin N ($0.05 - 5$ µg, i.c.v.) also elicits hypothermia in rats in a dose- and time-dependent manner, which is potentiated by bestatin (Item No. 21217).⁵ Neuromedin N is more potent than neurotensin in stimulating spontaneous motor activity and increasing dopamine metabolism in several brain regions following injection into the rat ventral tegmental area but is less potent than neurotensin when injected into the nucleus accumbens.⁶

References

1. Kitabgi, P. Neurotensin and neuromedin N are differentially processed from a common precursor by prohormone convertases in tissues and cell lines. *Results Probl. Cell Differ.* **50**, 85-96 (2010).
2. Checler, F., Vincent, J.P., and Kitabgi, P. Neuromedin N: High affinity interaction with brain neurotensin receptors and rapid inactivation by brain synaptic peptidases. *Eur. J. Pharmacol.* **126**(3), 239-244 (1986).
3. De la Fuente, M., Garrido, J.J., Arahuetes, R.M., et al. Stimulation of phagocytic function in mouse macrophages by neurotensin and neuromedin N. *J. Neuroimmunol.* **42**(1), 97-104 (1993).
4. Minamino, N., Kangawa, K., and Matsuo, H. Neuromedin N: A novel neurotensin-like peptide identified in porcine spinal cord. *Biochem. Biophys. Res. Commun.* **122**(2), 542-549 (1984).
5. Dubuc, I., Nouel, D., Coquerel, A., et al. Hypothermic effect of neuromedin N in mice and its potentiation by peptidase inhibitors. *Eur. J. Pharmacol.* **151**(1), 117-121 (1988).
6. Kalivas, P.W., Richardson-Carlson, R., and Duffy, P. Neuromedin N mimics the actions of neurotensin in the ventral tegmental area but not in the nucleus accumbens. *J. Pharmacol. Exp. Ther.* **238**(3), 1126-1131 (1986).

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, 11/17/2022

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