

# PRODUCT INFORMATION



## Neuromedin S (human) (trifluoroacetate salt)

Item No. 24544

**Formal Name:** L-isoleucyl-L-leucyl-L-glutamyl-L-arginylglycyl-L-serylglycyl-L-threonyl-L-alanyl-L-alanyl-L-valyl-L- $\alpha$ -aspartyl-L-phenylalanyl-L-threonyl-L-lysyl-L-lysyl-L- $\alpha$ -aspartyl-L-histidyl-L-threonyl-L-alanyl-L-threonyl-L-tryptophylglycyl-L-arginyl-L-prolyl-L-phenylalanyl-L-phenylalanyl-L-leucyl-L-phenylalanyl-L-arginyl-L-prolyl-L-arginyl-L-aspartamide, trifluoroacetate salt

H-Ile-Leu-Gln-Arg-Gly-Ser-Gly-Thr-Ala-Ala-Val-

Asp-Phe-Thr-Lys-Lys-Asp-His-Thr-Ala-Thr-Trp-

Gly-Arg-Pro-Phe-Phe-Leu-Phe-Arg-Pro-Arg-Asn-NH<sub>2</sub>

**Synonym:** NMS

**MF:** C<sub>173</sub>H<sub>265</sub>N<sub>53</sub>O<sub>44</sub> • XCF<sub>3</sub>COOH

• XCF<sub>3</sub>COOH

**FW:** 3,791.3

**Purity:**  $\geq 95\%$

**Supplied as:** A lyophilized powder

**Storage:** -20°C

**Stability:**  $\geq 4$  years

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

### Laboratory Procedures

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

### Description

Neuromedin S is a neuropeptide expressed in brain.<sup>1</sup> *In vivo*, neuromedin S (44  $\mu$ g/day) reduces body weight and food intake in diet-induced obese (DIO) mice.

### Reference

1. Peier, A., Kosinski, J., Cox-York, K., et al. The antiobesity effects of centrally administered neuromedin U and neuromedin S are mediated predominantly by the neuromedin U receptor 2 (NMUR2). *Endocrinology* **150**(7), 3101-3109 (2009).

#### 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, 11/21/2022

#### 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](http://WWW.CAYMANCHEM.COM)