

PRODUCT INFORMATION

[Lys⁵,Tyr⁶,mLeu⁹,Nle¹⁰]-Neurokinin A (4-10) (trifluoroacetate salt)

Item No. 42740

Formal Name:	(3S,6S,9S,12S,18S,21S)-3-amino-6-(4-aminobutyl)-21-hydrazineyl-9-(4-hydroxybenzyl)-18-isobutyl-12-isopropyl-19-methyl-4,7,10,13,16,17,20-hepta-oxo-5,8,11,14,19-pentaazapentacosanoic acid, trifluoroacetate salt	H-Asp-Lys-Tyr-Val-Gly-mLeu-Nle-NH ₂ • XCF ₃ COOH
Synonyms:	Neurokinin A (4-10)-S5K/F6Y/L9mL/M10Nle, [Lys ⁵ ,Tyr ⁶ ,mLeu ⁹ ,Nle ¹⁰]-NKA (4-10)	
Peptide Sequence:	DKYVGmLNle-NH ₂	
MF:	C ₃₉ H ₆₅ N ₉ O ₁₀ • XCF ₃ COOH	
FW:	820.0	
Purity:	≥95%	
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

[Lys⁵,Tyr⁶,mLeu⁹,Nle¹⁰]-Neurokinin A (4-10) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the XX in the solvent of choice, which should be purged with an inert gas. [Lys⁵,Tyr⁶,mLeu⁹,Nle¹⁰]-Neurokinin A (4-10) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

[Lys⁵,Tyr⁶,mLeu⁹,Nle¹⁰]-Neurokinin A (4-10) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

[Lys⁵,Tyr⁶,mLeu⁹,Nle¹⁰]-Neurokinin A (4-10) is a derivative of neurokinin A (Item No. 24221) containing serine-to-lysine, phenylalanine-to-tyrosine, leucine-to-methyleucine, and methionine-to-norleucine substitutions at positions 5, 6, 9, and 10, respectively, and a peptide fragment of the neurokinin-2 (NK₂) receptor agonist EB1001.¹

Reference

1. Sass, F., Ma, T., Ekberg, J.H., *et al.* NK2R control of energy expenditure and feeding to treat metabolic diseases. *Nature* **635**(8040), 987-1000 (2024).

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, 03/11/2025

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