

PRODUCT INFORMATION



DKYVG(methyl-Leu)Methionine Sulfone-NH₂ (trifluoroacetate salt)

Item No. 42741

Formal Name: (3S,6S,12S,15S,18S,21S)-21-amino-18-(4-aminobutyl)-3-carbamoyl-15-(4-hydroxybenzyl)-6-isobutyl-12-isopropyl-7-methyl-1-(methylsulfonyl)-5,8,11,14,17,20-hexaoxo-4,7,10,13,16,19-hexaazatricosan-23-oic acid, trifluoroacetate salt

Synonyms: Asp-Lys-Tyr-Val-mLeu-Mox-NH₂, [Lys⁵,Tyr⁶,mLeu⁹,Mox¹⁰]-NKA (4-10), Neurokinin A (4-10)-S5K/F6Y/L9mL/M10Mox

Peptide Sequence: DKYVGmLX-NH₂ (X = methionine sulfone)

MF: C₃₈H₆₃N₉O₁₂S • XCF₃COOH

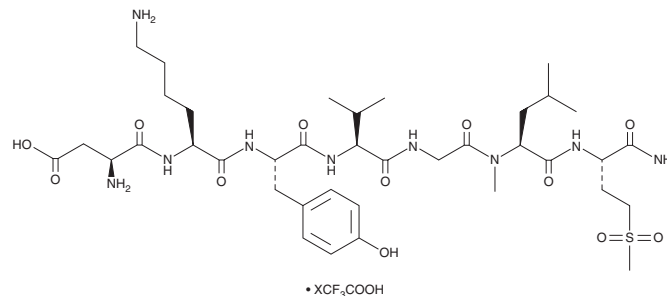
FW: 870.0

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

DKYVG(methyl-Leu)Methionine sulfone-NH₂ (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the DKYVG(methyl-Leu)methionine sulfone-NH₂ in the solvent of choice, which should be purged with an inert gas. DKYVG(methyl-Leu)Methionine sulfone-NH₂ is slightly soluble (0.1-1 mg/ml) in acetonitrile.

DKYVG(methyl-Leu)Methionine sulfone-NH₂ 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

DKYVG(methyl-Leu)Methionine sulfone-NH₂ is a derivative of neurokinin A (Item No. 24221) containing serine-to-lysine, phenylalanine-to-tyrosine, leucine-to-methylleucine, and methionine-to-methionine sulfone substitutions at positions 5, 6, 9, and 10, respectively, and a peptide fragment of the neurokinin-2 (NK₂) receptor agonist EB1002.¹

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, 06/24/2025

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