

PRODUCT INFORMATION



Tat-M2NX (trifluoroacetate salt)

Item No. 37447

Peptide Sequence: H-YGRKKRRQRRRGSREPGEMLPRLKRVL RQEFWV-OH H-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-Arg-Gly-Ser-Arg-Glu-Pro-Gly-Glu-Met-Leu-Pro-Arg-Lys-Leu-Lys-Arg-Val-Leu-Arg-Gln-Glu-Phe-Trp-Val-OH
MF: C₁₉₀H₃₂₃N₇₁O₄₅S • XCF₃COOH
FW: 4,354.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
• XCF₃COOH

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

Laboratory Procedures

Tat-M2NX (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Tat-M2NX (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Tat-M2NX is a peptide antagonist of transient receptor potential melastatin 2 (TRPM2).¹ It inhibits hydrogen peroxide-induced calcium influx in HEK293 cells expressing human TRPM2 when used at concentrations ranging from 25 to 100 μM. *In vivo*, Tat-M2NX (20 mg/kg) reduces brain infarct volume in male, but not female, mice in a model of stroke induced by middle cerebral artery occlusion (MCAO) when administered prior to or 3 hours-post occlusion.

Reference

1. Shimizu, T., Dietz, R.M., Cruz-Torres, I., *et al.* Extended therapeutic window of a novel peptide inhibitor of TRPM2 channels following focal cerebral ischemia. *Exp. Neurol.* **283(Pt A)**, 151-156 (2016).

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, 06/14/2023

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