

PRODUCT INFORMATION



Tat-QFNP12 (trifluoroacetate salt)

Item No. 38845

Formal Name: L-tyrosylglycyl-L-arginyl-L-lysyl-L-lysyl-L-arginyl-L-arginyl-L-glutaminyl-L-arginyl-L-arginyl-L-arginyl-L-valyl-L-lysyl-L-lysyl-L-glutamyl-L-alanyl-L-glutamyl-L-leucyl-L-aspartyl-L-lysyl-L-tyrosyl-L-leucyl-L-glutamic acid, trifluoroacetate salt

H-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-Arg-Val-Lys-Lys-Glu-Ala-Glu-Leu-Asp-Lys-Tyr-Leu-Glu-OH
• XCF₃COOH

Peptide Sequence: YGRKKRRQRRRVKKEAELDKYLE-OH

MF: C₁₃₀H₂₂₅N₄₇O₃₅ • XCF₃COOH

FW: 3,006.5

Purity: ≥90%

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

Tat-QFNP12 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Tat-QFNP12 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Tat-QFNP12 is a peptide containing a transcriptional transactivator (Tat) transmembrane domain and an inhibitor of the protein-protein interaction between N-Myc downstream regulated gene 2 (NDRG2) and the protein phosphatase Mg²⁺/Mn²⁺ dependent 1A (PPM1A).¹ It decreases the loss of blood-brain barrier endothelial tight junctions induced by increased levels of matrix metalloproteinase-9 (MMP-9), reduces cerebral edema, as well as increases spontaneous activity and symmetry of limb movement, in an endovascular puncture-induced mouse model of subarachnoid hemorrhage when administered at a dose of 20 mg/kg.

Reference

1. Feng, D., Zhou, J., Liu, H., *et al.* Astrocytic NDRG2-PPM1A interaction exacerbates blood-brain barrier disruption after subarachnoid hemorrhage. *Sci. Adv.* **8(39)**, eabq2423 (2022).

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, 08/08/2023

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