

PRODUCT INFORMATION



T100-Mut (trifluoroacetate salt)

Item No. 37464

Peptide Sequence: Myr-WKVRQRNKKVQQESQTALVVNQRLFA-

OH (where Myr = myristoyl)

MF: $C_{161}H_{271}N_{49}O_{42}S \cdot XCF_3COOH$

FW: 3,597.3

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

Myr-Trp-Lys-Val-Arg-Gln-Arg-Asn-Lys-Lys-Val-

Gln-Gln-Gln-Glu-Ser-Gln-Thr-Ala-Leu-Val-

Val-Asn-Gln-Arg-Cys-Leu-Phe-Ala-OH

• XCF_3COOH

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

Laboratory Procedures

T100-Mut (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the T100-Mut (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

T100-Mut is a cell-permeable peptide and transient receptor potential ankyrin 1 (TRPA1) inhibitor that contains the C-terminal sequence of Tmem100-3Q, a mutant form of the endogenous peptide Tmem100.¹ It reduces calcium flux induced by mustard oil, a TRPA1 agonist, in primary mouse dorsal root ganglia (DRG) when used at a concentration of 200 nM. Subcutaneous administration of T100-Mut (5 μ l of a 2 mM solution) into the hind paw reduces paw licking induced by mustard oil but not the TRP vanilloid 1 (TRPV1) agonist capsaicin (Item Nos. 92350 | 10010743) in mice. It also reduces mechanical hyperalgesia induced by complete Freund's adjuvant (CFA) or paclitaxel (Item No. 10461) in wild-type, but not TRPV1 knockout, mice.

Reference

1. Weng, H.J., Patel, K.N., Jeske, N.A., *et al.* Tmem100 is a regulator of TRPA1-TRPV1 complex and contributes to persistent pain. *Neuron* **85**(4), 833-846 (2015).

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, 07/11/2025

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