

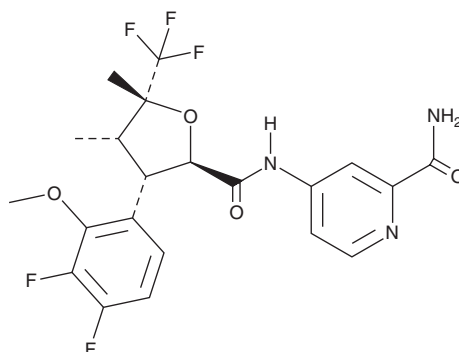
PRODUCT INFORMATION



Suzetrigine

Item No. 40799

CAS Registry No.: 2649467-58-1
Formal Name: 4-[[[(2R,3S,4S,5R)-3-(3,4-difluoro-2-methoxyphenyl)tetrahydro-4,5-dimethyl-5-(trifluoromethyl)-2-furanyl]carbonyl]amino]-2-pyridinecarboxamide
Synonym: VX-548
MF: C₂₁H₂₀F₅N₃O₄
FW: 473.4
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

Suzetrigine is supplied as a solid. A stock solution may be made by dissolving the suzetrigine in the solvent of choice, which should be purged with an inert gas. Suzetrigine is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

Suzetrigine is an inhibitor of voltage-gated sodium channel 1.8 (Na_v1.8; IC₅₀ = 0.7 nM for the human channel).¹ It is selective for Na_v1.8 over a panel of eight additional voltage-gated sodium channels (IC₅₀s = >10,000 nM for all).

Reference

1. Jones, J., Correll, D.J., Lechner, S.M., *et al.* Selective inhibition of Na_v1.8 with VX-548 for acute pain. *N. Engl. J. Med.* **389**(5), 393-405 (2023).

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, 05/24/2024

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