

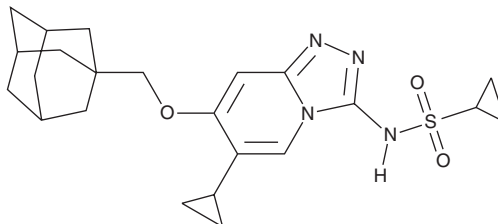
PRODUCT INFORMATION



GNE-131

Item No. 45386

CAS Registry No.: 1629063-81-5
Formal Name: N-[6-cyclopropyl-7-(tricyclo[3.3.1.1^{3,7}]-dec-1-ylmethoxy)-1,2,4-triazolo[4,3-a]pyridin-3-yl]-cyclopropanesulfonamide
MF: C₂₃H₃₀N₄O₃S
FW: 442.6
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

GNE-131 is supplied as a solid. A stock solution may be made by dissolving the GNE-131 in the solvent of choice, which should be purged with an inert gas. GNE-131 is soluble (≥10 mg/ml) in DMSO.

Description

GNE-131 is an inhibitor of voltage-gated sodium channel 1.7 (Na_v1.7; IC₅₀ = 0.003 μM).¹ It is selective for Na_v1.7 over Na_v1.1, Na_v1.2, Na_v1.5, and Na_v1.6 (IC₅₀s = 0.045, 0.007, 0.110, and 0.092 μM, respectively). *In vivo*, GNE-131 (30 mg/kg) reduces aconitine-induced nocifensive behavior in a transgenic mouse model of inherited erythromelalgia (IEM).

Reference

1. Focken, T., Chowdhury, S., Zenova, A., *et al.* Design of conformationally constrained acyl sulfonamide isosteres: Identification of N-([1,2,4]triazolo[4,3-a]pyridin-3-yl)methane-sulfonamides as potent and selective hNa_v1.7 inhibitors for the treatment of pain. *J. Med. Chem.* **61**(11), 4810-4831 (2018).

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, 03/30/2026

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