

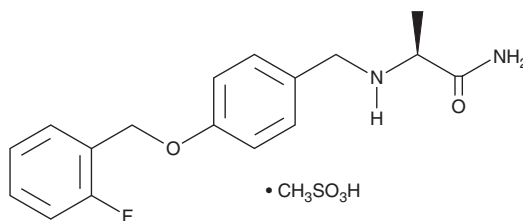
PRODUCT INFORMATION



Ralfinamide (mesylate)

Item No. 45954

CAS Registry No.: 202825-45-4
Formal Name: 2S-[[[4-[(2-fluorophenyl)methoxy]phenyl]methyl]amino]propanamide, methanesulfonate
Synonym: NW-1029
MF: C₁₇H₁₉FN₂O₂ • CH₃SO₃H
FW: 398.4
Purity: ≥98%
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

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

Description

Ralfinamide is an inhibitor of voltage-gated sodium channels (Na_v).¹ It induces tonic blocks of tetrodotoxin-resistant and -sensitive sodium currents in isolated rat dorsal root ganglion (DRG) neurons (IC₅₀s = 55 and 22 μM, respectively). Ralfinamide reduces mechanical allodynia in a rat model of arthritis induced by complete Freund's adjuvant (CFA) and in a rat model of neuropathic pain induced by chronic constriction injury (CCI).² It has anticonvulsant activity in models of seizures induced by bicuculline (Item No. 11727) or maximal electroshock (MES) in mice (ED₅₀s = 25.8 and 7 mg/kg, respectively).³

References

1. Stummann, T.C., Salvati, P., Fariello, R.G., *et al.* The anti-nociceptive agent ralfinamide inhibits tetrodotoxin-resistant and tetrodotoxin-sensitive Na⁺ currents in dorsal root ganglion neurons. *Eur. J. Pharmacol.* **510**(3), 197-208 (2005).
2. Veneroni, O., Maj, R., Calabresi, M., *et al.* Anti-allodynic effect of NW-1029, a novel Na⁺ channel blocker, in experimental animal models of inflammatory and neuropathic pain. *Pain* **102**(1-2), 17-25 (2003).
3. Pevarello, P., Bonsignori, A., Dostert, P., *et al.* Synthesis and anticonvulsant activity of a new class of 2-[(arylalkyl)amino]alkanamide derivatives. *J. Med. Chem.* **41**(4), 579-590 (1998).

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, 07/13/2026

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