

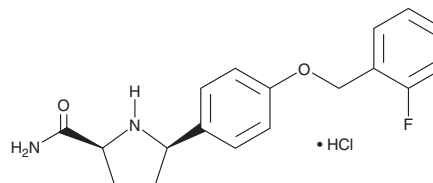
PRODUCT INFORMATION



Raxatrigine (hydrochloride)

Item No. 23490

CAS Registry No.: 934240-31-0
Formal Name: (2S)-5R-[4-[(2-fluorophenyl)methoxy]phenyl]-2-pyrrolidinecarboxamide, monohydrochloride
Synonyms: BIIB074, CNV1014802, GSK1014802A, Vixotrigine
MF: C₁₈H₁₉FN₂O₂ • HCl
FW: 350.8
Purity: ≥98%
UV/Vis.: λ_{max}: 229 nm
Supplied as: A crystalline 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

Raxatrigine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the raxatrigine (hydrochloride) in the solvent of choice. Raxatrigine (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of raxatrigine (hydrochloride) in these solvents is approximately 33 and 16 mg/ml, respectively.

Raxatrigine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, raxatrigine (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Raxatrigine (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Raxatrigine is an inhibitor of voltage-gated sodium channels (Na_v; IC₅₀s = 22, 13, 23, 10, 84, 15, 32, and 6 μM for Na_v1.1-Na_v1.8 channels in a FLIPR membrane potential assay using HEK293 cells).¹ Its activity is state-dependent at the Na_v1.7 channel, with a nine-fold increase in potency for the open/inactivated state (IC₅₀ = 6.3 μM) compared to the closed/resting state (IC₅₀ = 54 μM). Raxatrigine (30 mg/kg, i.p.) decreases spontaneous pain behavior in an OD1 mouse pain model.

Reference

1. Deuis, J.R., Wingerd, J.S., Winter, Z., *et al.* Analgesic effects of GpTx-1, PF-04856264 and CNV1014802 in a mouse model of Na_v1.7-mediated pain. *Toxins (Basel)* **8**(3), 78 (2016).

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, 12/22/2022

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