

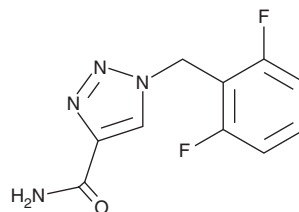
PRODUCT INFORMATION



Rufinamide

Item No. 18870

CAS Registry No.: 106308-44-5
Formal Name: 1-[(2,6-difluorophenyl)methyl]-1H-1,2,3-triazole-4-carboxamide
Synonyms: CGP 33101, RUF 331
MF: C₁₀H₈F₂N₄O
FW: 238.2
Purity: ≥98%
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

Rufinamide is supplied as a crystalline solid. A stock solution may be made by dissolving the rufinamide in the solvent of choice, which should be purged with an inert gas. Rufinamide is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of rufinamide in these solvents is approximately 10 and 2 mg/ml, respectively.

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

Description

Rufinamide is an anticonvulsant.¹ It inhibits the activation of voltage-gated sodium channel 1.1 (Na_v1.1) when used at a concentration of 100 μM.³ Rufinamide inhibits Na_v1.1, but not Na_v1.2, Na_v1.3, and Na_v1.6, opening and increases the action potential threshold in primary rat hippocampal neurons. It is an inhibitor of carbonic anhydrase VA (CAVA; K_i = 343.8 nM) that is selective for CAVA over CAI and CAII (K_is = >10,000 nM for both).² Rufinamide (100 μM) prolongs the preictal phase and reduces seizure-like event frequency in an *in vitro* model of epileptiform activity in rat hippocampal slices.⁴ It inhibits seizures induced by pentylenetetrazole (Item No. 18682) in a mouse model of epilepsy (ED₅₀ = 54 mg/kg, i.p.) and reduces kainic acid-induced neuronal cell death in the mouse hippocampal CA3 region when used at doses of 25, 50, and 100 mg/kg.^{5,6} Formulations containing rufinamide have been used in the treatment of seizures associated with Lennox-Gastaut Syndrome (LGS).

References

1. Wheless, J.W. and Vazquez, B. Rufinamide: A novel broad-spectrum antiepileptic drug. *Epilepsy Curr.* **10**(1), 1-6 (2010).
2. Costa, G., Carta, F., Ambrosio, F.A., *et al.* A computer-assisted discovery of novel potential anti-obesity compounds as selective carbonic anhydrase VA inhibitors. *Eur. J. Med. Chem.* **181**, 111565 (2019).
3. Gilchrist, J.J., Dutton, S., Diaz-Bustamante, M., *et al.* Na_v1.1 modulation by a novel triazole compound attenuates epileptic seizures in rodents. *ACS Chem. Biol.* **9**(5), 1204-1212 (2014).
4. Gáll, Z., Orbán-Kis, K. and Szilágyi, T. Differential effects of sodium channel blockers on *in vitro* induced epileptiform activities. *Arch. Pharm. Res.* **40**(1), 112-121 (2017).
5. White, H.S., Franklin, M.R., Kupferberg, H.J., *et al.* The anticonvulsant profile of rufinamide (CGP 33101) in rodent seizure models. *Epilepsia* **49**(7), 1213-1220 (2008).
6. Park, J.-A. and Lee, C.-H. Effect of Rufinamide on the kainic acid-induced excitotoxic neuronal death in the mouse hippocampus. *Arch. Pharm. Res.* **41**(7), 776-783 (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, 11/03/2022

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