

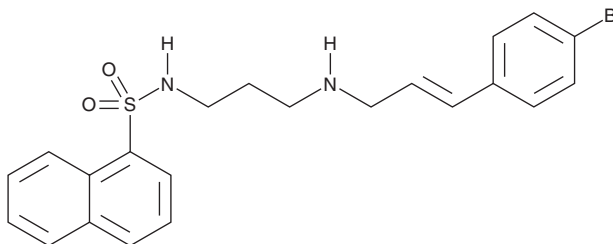
PRODUCT INFORMATION



Sulfonadyn-47

Item No. 39402

Formal Name: (E)-N-(3-((3-(4-bromophenyl)allyl)amino)propyl)naphthalene-1-sulfonamide
MF: C₂₂H₂₃BrN₂O₂S
FW: 459.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 261 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Sulfonadyn-47 is supplied as a solid. A stock solution may be made by dissolving the sulfonadyn-47 in the solvent of choice, which should be purged with an inert gas. Sulfonadyn-47 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of sulfonadyn-47 in these solvents is approximately 10 mg/ml.

Description

Sulfonadyn-47 is a dynamin inhibitor.¹ It inhibits dynamin GTPase activity in a cell-free assay (IC₅₀ = 3.5 μM). Sulfonadyn-47 inhibits clathrin-mediated endocytosis in U2OS cells (IC₅₀ = 27.3 μM), as well depolarization-stimulated uptake of FM4-64, a marker of synaptic vesicle endocytosis (SVE), in isolated rat brain synaptosomes (IC₅₀ = 12.3 μM). *In vivo*, sulfonadyn-47 (30 and 60 mg/kg) has anticonvulsant activity in the 6 Hz corneal kindling mouse model of seizures.

Reference

1. Odell, L.R., Jones, N.C., Chau, N., *et al.* The sulfonadyns: A class of aryl sulfonamides inhibiting dynamin I GTPase and clathrin mediated endocytosis are anti-seizure in animal models. *RSC Med. Chem.* **14**(8), 1492-1511 (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, 09/13/2023

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