

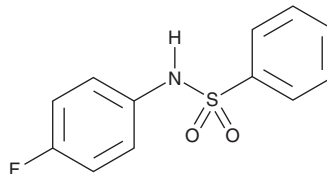
PRODUCT INFORMATION

ELN484228

Item No. 29398



CAS Registry No.: 312-63-0
Formal Name: N-(4-fluorophenyl)-benzenesulfonamide
Synonym: NSC 164389
MF: $C_{12}H_{10}FNO_2S$
FW: 251.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

ELN484228 is supplied as a solid. A stock solution may be made by dissolving the ELN484228 in the solvent of choice, which should be purged with an inert gas. ELN484228 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ELN484228 in these solvents is approximately 1 and 10 mg/ml, respectively.

Description

ELN484228 is an inhibitor of α -synuclein-induced toxicity.¹ It prevents inhibition of phagocytosis induced by overexpression of α -synuclein in H4 neuroglioma cells when used at concentrations of 30, 100, and 300 μM . ELN484228 (10 μM) protects dopaminergic neurons from cell death and neurite retraction induced by A53T mutant α -synuclein in a rat midbrain culture model of Parkinson's disease.

Reference

1. Toth, G., Gardai, S.J., Zago, W., *et al.* Targeting the intrinsically disordered structural ensemble of α -synuclein by small molecules as a potential therapeutic strategy for Parkinson's disease. *PLoS One* 9(2), e87133 (2014).

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, 11/23/2022

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