

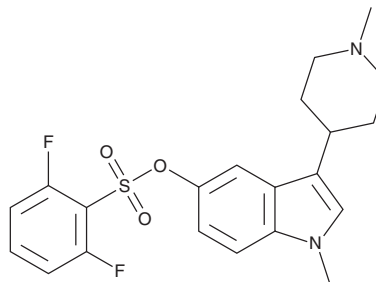
PRODUCT INFORMATION



SGS 518

Item No. 43670

CAS Registry No.: 445441-26-9
Formal Name: 2,6-difluoro-benzenesulfonic acid, 1-methyl-3-(1-methyl-4-piperidiny)-1H-indol-5-yl ester
Synonym: LY483518
MF: C₂₁H₂₂F₂N₂O₃S
FW: 420.5
Purity: ≥95%
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

SGS 518 is supplied as a solid. A stock solution may be made by dissolving the SGS 518 in the solvent of choice, which should be purged with an inert gas. SGS 518 is sparingly soluble (1-10 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Description

SGS 518 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT₆.¹

Reference

1. Upton, N., Chuang, T.T., Hunter, A.J., *et al.* 5-HT₆ receptor antagonists as novel cognitive enhancing agents for Alzheimer's disease. *Neurotherapeutics* **5**(3), 458-469 (2008).

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, 08/06/2025

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