

PRODUCT INFORMATION



GSM2

Item No. 41322

CAS Registry No.: 1093978-89-2
Formal Name: 6-fluoro-1,3,4,5-tetrahydro-3-[4-[3-methoxy-4-(4-methyl-1H-imidazol-1-yl)phenyl]-1H-1,2,3-triazol-1-yl]-1-(2,2,2-trifluoroethyl)-2H-1-benzazepin-2-one

Synonyms: γ -Secretase Modulator 2, γ -SM2

MF: $C_{25}H_{22}F_4N_6O_2$

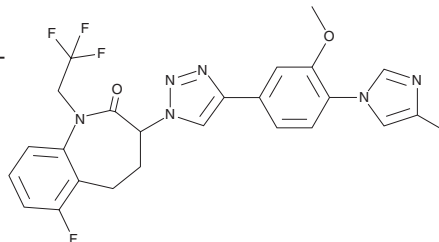
FW: 514.5

Purity: $\geq 98\%$

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

GSM2 is supplied as a solid. A stock solution may be made by dissolving the GSM2 in the solvent of choice, which should be purged with an inert gas. GSM2 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

GSM2 is an inhibitor of γ -secretase.¹ It inhibits the production of amyloid- β (1-42) (A β 42) in SH-SY5Y neuroblastoma cells overexpressing SPA4CT, which is a peptide containing the N-terminal signal peptide and C-terminal residues of amyloid precursor protein (APP).

Reference

1. Fischer, C., Munoz, B., Zultanski, S., *et al.* Triazole derivatives for treating Alzheimer's disease and related conditions. **WO2008/156580A1** (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, 09/10/2024

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