

PRODUCT INFORMATION

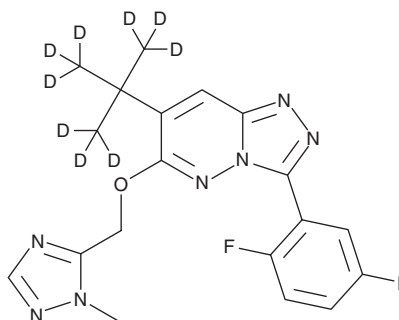


CTP354

Item No. 36618

CAS Registry No.: 1213669-91-0
Formal Name: 3-(2,5-difluorophenyl)-7-[1,1-di(methyl-d₃)ethyl-2,2,2-d₃]-6-[(1-methyl-1H-1,2,4-triazol-5-yl)methoxy]-1,2,4-triazolo[4,3-b]pyridazine
MF: C₁₉H₁₀D₉F₂N₇O
FW: 408.5

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
UV/Vis.: λ_{max}: 248 nm
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

CTP354 is supplied as a solid. A stock solution may be made by dissolving the CTP354 in the solvent of choice, which should be purged with an inert gas. CTP354 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CTP354 in ethanol is approximately 2 mg/ml and approximately 3 mg/ml in DMSO and DMF.

Description

CTP354 is a deuterated derivative of the GABA_A receptor modulator L-838,417.¹ It binds to the GABA_A receptor benzodiazepine binding site in non-α₁ subunit-containing GABA_A receptors. CTP354 has a longer duration of action than L-838,417 in a rat model of neuropathic pain induced by sciatic nerve ligation.

Reference

1. Harbeson, S.L. and Tung, R.D. Deuterium medicinal chemistry: A new approach to drug discovery and development. *Medchem News* **24(2)**, 8-22 (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, 12/20/2022

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