

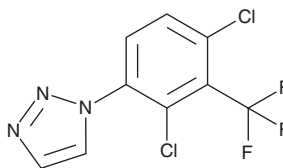
PRODUCT INFORMATION



ARUK3001185

Item No. 37113

CAS Registry No.: 2411969-39-4
Formal Name: 1-[2,4-dichloro-3-(trifluoromethyl)phenyl]-1H-1,2,3-triazole
Synonym: 1-Phenyl-1H-1,2,3-triazole 1
MF: C₉H₄Cl₂F₃N₃
FW: 282.1
Purity: ≥98%
UV/Vis.: λ_{max}: 288 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

ARUK3001185 is supplied as a crystalline solid. A stock solution may be made by dissolving the ARUK3001185 in the solvent of choice, which should be purged with an inert gas. ARUK3001185 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Description

ARUK3001185 is an inhibitor of palmitoleoyl-protein carboxylesterase NOTUM (IC₅₀ = 6.7 nM), an enzyme involved in the deacylation of Wnt.¹ It is selective for NOTUM over a panel of 39 serine hydrolases and 485 kinases, including casein kinase 1 (CK1) and glycogen synthase kinase 3 (GSK3), at 10 μM and a panel of 47 receptors, ion channels, transporters, and enzymes at 30 μM, but does inhibit the serotonin (5-HT) receptor subtype 5-HT_{3A}, α4β2 subunit-containing nicotinic acetylcholine receptors (nAChRs), and monoamine oxidase A (MAO-A; IC₅₀s = 4.5, 10, and 28 μM, respectively). ARUK3001185 prevents NOTUM-induced decreases in Wnt3a signaling in a cell-based reporter assay (EC₅₀ = 110 nM).

Reference

1. Willis, N.J., Mahy, W., Sipthorp, J., *et al.* Design of a potent, selective, and brain-penetrant inhibitor of Wnt-deactivating enzyme notum by optimization of a crystallographic fragment hit. *J. Med. Chem.* **65**(10), 7212-7230 (2122).

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, 01/17/2023

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