

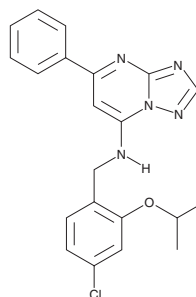
PRODUCT INFORMATION



AF64394

Item No. 36826

CAS Registry No.: 1637300-25-4
Formal Name: N-[[4-chloro-2-(1-methylethoxy)phenyl]methyl]-5-phenyl-[1,2,4]triazolo[1,5-a]pyrimidin-7-amine
MF: C₂₁H₂₀ClN₅O
FW: 393.9
Purity: ≥98%
UV/Vis.: λ_{max}: 253 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

AF64394 is supplied as a solid. A stock solution may be made by dissolving the AF64394 in the solvent of choice, which should be purged with an inert gas. AF64394 is soluble in DMSO.

Description

AF64394 is an inverse agonist of the orphan G protein-coupled receptor GPR3.¹ It selectively inhibits cAMP accumulation in HEK293 cells expressing GPR3 over GPR6 or GPR12 (IC₅₀s = 0.5, 7.9, and 12 μM, respectively).

Reference

1. Jensen, T., Elster, L., Nielsen, S.M., *et al.* The identification of GPR3 inverse agonist AF64394; The first small molecule inhibitor of GPR3 receptor function. *Bioorg. Med. Chem. Lett.* **24**(22), 5195-5198 (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, 03/20/2023

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