

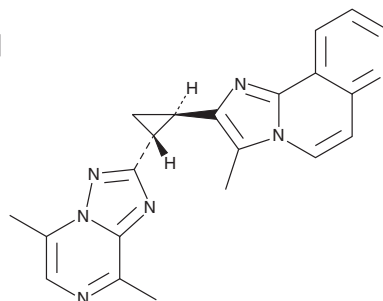
PRODUCT INFORMATION



SEP-0372814

Item No. 32836

CAS Registry No.: 1516895-53-6
Formal Name: *rel*-2-[(1R,2R)-2-(5,8-dimethyl[1,2,4]triazolo[1,5-a]pyrazin-2-yl)cyclopropyl]-3-methyl-imidazo[2,1-f][1,6]naphthyridine
MF: C₂₁H₁₉N₇
FW: 369.4
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 271 nm
Supplied as: A crystalline 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

SEP-0372814 is supplied as a crystalline solid. A stock solution may be made by dissolving the SEP-0372814 in the solvent of choice, which should be purged with an inert gas. SEP-0372814 is soluble in the organic solvent methanol at a concentration of approximately 1 mg/ml. SEP-0372814 is slightly soluble in chloroform and dimethyl formamide.

Description

SEP-0372814 is an inhibitor of phosphodiesterase 10A (PDE10A; IC₅₀ = ≤10 nM for the recombinant human enzyme).¹

Reference

1. Campbell, J.E., Jones, P.G., and Malcolm, S. Heteroaryl compounds and methods of use thereof. *Sunovion Pharmaceuticals*. US20150322090A1 (2015).

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, 12/09/2022

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