

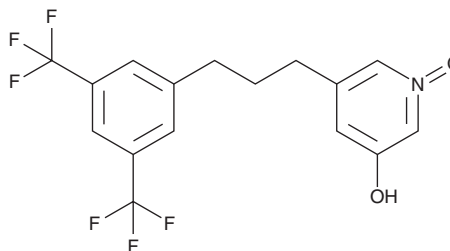
PRODUCT INFORMATION



OX04528

Item No. 42377

CAS Registry No.: 3028055-45-7
Formal Name: 5-[3-[3,5-bis(trifluoromethyl)phenyl]propyl]-3-pyridinol, 1-oxide
MF: C₁₆H₁₃F₆NO₂
FW: 365.3
Purity: ≥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

OX04528 is supplied as a solid. A stock solution may be made by dissolving the OX04528 in the solvent of choice, which should be purged with an inert gas. OX04528 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

OX04528 is a G protein-biased agonist of G protein-coupled receptor 84 (GPR84).¹ It inhibits forskolin-induced cAMP production in CHO cells expressing human GPR84 (EC₅₀ = 0.00598 nM) and does not induce β-arrestin recruitment when used at concentrations up to 80 μM.

Reference

1. Wang, P., Raja, A., Luscombe, V.B., *et al.* Development of highly potent, G-protein pathway biased, selective, and orally bioavailable GPR84 agonists. *J. Med. Chem.* **67**(1), 110-137 (2024).

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, 12/16/2024

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