

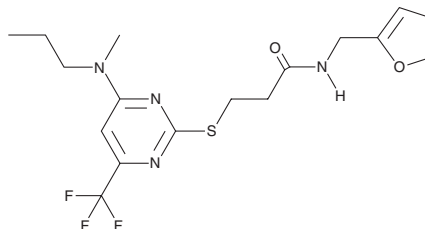
PRODUCT INFORMATION



SET2

Item No. 40360

CAS Registry No.: 2313525-20-9
Formal Name: N-(2-furanylmethyl)-3-[[4-(methylpropylamino)-6-(trifluoromethyl)-2-pyrimidinyl]thio]-propanamide
MF: C₁₇H₂₁F₃N₄O₂S
FW: 402.4
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

SET2 is supplied as a solid. A stock solution may be made by dissolving the SET2 in the solvent of choice, which should be purged with an inert gas. SET2 is soluble in (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

SET2 is an inhibitor of transient receptor potential vanilloid 2 (TRPV2; IC₅₀ = 0.46 μM).¹ It selectively inhibits currents induced by the TRP channel modulator 2-APB (Item No. 64970) in a patch-clamp assay using HEK293T cells overexpressing TRPV2 at 0.3 or 1 μM over HEK293T cells overexpressing TRPV1 or TRPV3 at 3 or 10 μM. SET2 (20 μM) reduces migration induced by lysophosphatidic acid (LPA) in PC-3M prostate cancer cells to a similar extent as the LPA receptor 1 (LPA1) antagonist AM966 (Item No. 22048). It also reduces LPA-induced increases in current in a patch-clamp assay using HEK293T cells when used at a concentration of 10 μM.

Reference

1. Chai, H., Cheng, X., Zhou, B., *et al.* Structure-based discovery of a subtype-selective inhibitor targeting a transient receptor potential vanilloid channel. *J. Med. Chem.* **62(3)**, 1373-1384 (2019).

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, 03/28/2024

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