

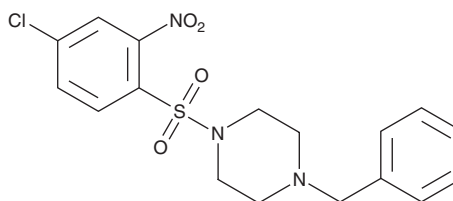
PRODUCT INFORMATION



RN-1747

Item No. 37573

CAS Registry No.: 1024448-59-6
Formal Name: 1-[(4-chloro-2-nitrophenyl)sulfonyl]-4-(phenylmethyl)piperazine
MF: C₁₇H₁₈ClN₃O₄S
FW: 395.9
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

RN-1747 is supplied as a solid. A stock solution may be made by dissolving the RN-1747 in the solvent of choice, which should be purged with an inert gas. RN-1747 is soluble in methanol and DMSO.

Description

RN-1747 is an agonist of transient receptor potential vanilloid 4 (TRPV4).¹ It selectively induces calcium influx in HEK293 cells expressing human, rat, or mouse TRPV4 (EC₅₀s = 0.77, 4.1, and 4 μM, respectively) over HEK293 cells expressing human TRPV1, TRPV3, or transient receptor potential melastatin 8 (TRPM8; EC₅₀s = >100, >30, and >30 μM, respectively). RN-1747 (10 μM) induces inward current in *Xenopus* oocytes expressing human TRPV4. Topical, but not intravenous, administration of RN-1747 induces hypothermia in rats.²

References

1. Vincent, F., Acevedo, A., Nguyen, M.T., *et al.* Identification and characterization of novel TRPV4 modulators. *Biochem. Biophys. Res. Commun.* **389**(3), 490-494 (2009).
2. Vizin, R.C., Scarpellini, C.S., Ishikawa, D.T., *et al.* TRPV4 activates autonomic and behavioural warmth-defence responses in Wistar rats. *Acta Physiol. (Oxf.)* **214**(2), 275-289 (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, 10/16/2023

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