

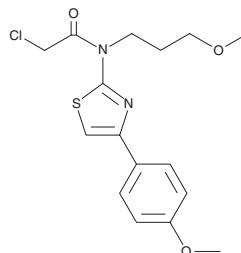
PRODUCT INFORMATION



JT010

Item No. 29732

CAS Registry No.: 917562-33-5
Formal Name: 2-chloro-N-[4-(4-methoxyphenyl)-2-thiazolyl]-N-(3-methoxypropyl)-acetamide
MF: C₁₆H₁₉ClN₂O₃S
FW: 354.9
Purity: ≥95%
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

JT010 is supplied as a solid. A stock solution may be made by dissolving the JT010 in the solvent of choice, which should be purged with an inert gas. JT010 is soluble in organic solvents such as ethanol and DMSO. The solubility of JT010 in these solvents is approximately 50 and 100 mM, respectively.

Description

JT010 is a transient receptor potential ankyrin 1 (TRPA1) agonist ($EC_{50} = 0.65$ nM in a calcium influx assay).¹ It is selective for TRPA1 over TRP vanilloid 1 (TRPV1), TRPV3, TRPV4, TRP canonical channel 5 (TRPC5), TRP melastatin 2 (TRPM2), and TRPM8 channels at 1 μ M.

Reference

1. Takaya, J., Mio, K., Shiraishi, T., *et al.* A potent and site-selective agonist of TRPA1. *J. Am. Chem. Soc.* **137**(50), 15859-15864 (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, 02/07/2024

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