

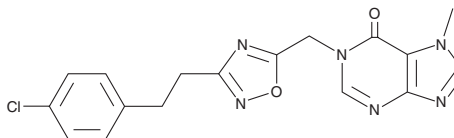
PRODUCT INFORMATION



AM0902

Item No. 37290

CAS Registry No.: 1883711-97-4
Formal Name: 1-[[3-[2-(4-chlorophenyl)ethyl]-1,2,4-oxadiazol-5-yl]methyl]-1,7-dihydro-7-methyl-6H-purin-6-one
MF: C₁₇H₁₅ClN₆O₂
FW: 370.8
Purity: ≥98%
UV/Vis.: λ_{max}: 214 nm
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

AM0902 is supplied as a solid. A stock solution may be made by dissolving the AM0902 in the solvent of choice, which should be purged with an inert gas. AM0902 is soluble in DMSO.

Description

AM0902 is a transient receptor potential ankyrin 1 (TRPA1) antagonist (IC₅₀ = 0.131 μM for the human receptor in a calcium mobilization assay).¹ It is selective for human TRPA1 over human transient receptor potential vanilloid 1 (TRPV1) and TRPV4, rat TRPV1, TRPV3, and transient receptor potential melastatin 8 (TRPM8), and a panel of 40 additional targets at 10 μM. *In vivo*, AM0902 (10 and 30 mg/kg) reduces the number of tail flinches induced by the TRPA1 agonist allyl isothiocyanate (AITC) in rats.

Reference

1. Schenkel, L.B., Olivieri, P.R., Boezio, A.A., *et al.* Optimization of a novel quinazolinone-based series of transient receptor potential A1 (TRPA1) antagonists demonstrating potent *in vivo* activity. *J. Med. Chem.* **59**(6), 2794-2809 (2016).

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, 02/07/2023

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