

PRODUCT INFORMATION



YQ-128

Item No. 36451

CAS Registry No.: 2454246-18-3

Formal Name: N-[(5-chloro-2-propoxyphenyl)methyl]-N-[2-[4-[(2-propyn-1-ylamino)sulfonyl]phenyl]ethyl]-3-thiopheneacetamide

MF: C₂₇H₂₉ClN₂O₄S₂

FW: 545.1

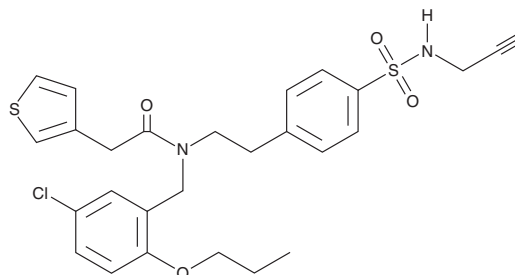
Purity: ≥98%

UV/Vis.: λ_{max}: 229 nm

Supplied as: A crystalline 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

YQ-128 is supplied as a crystalline solid. A stock solution may be made by dissolving the YQ-128 in the solvent of choice, which should be purged with an inert gas. YQ-128 is soluble in the organic solvent DMSO.

Description

YQ-128 is an inhibitor of the NOD-like receptor protein 3 (NLRP3) inflammasome (IC₅₀ = 0.3 μM).¹ It inhibits LPS- and ATP-induced IL-1β release from isolated mouse peritoneal macrophages (IC₅₀ = 1.59 μM). YQ-128 (10 mg/kg) inhibits LPS-induced IL-1β production in wild-type but not *Nlrp3*^{-/-} mice.

Reference

1. Jiang, Y., He, L., Green, J., *et al.* Discovery of second-generation NLRP3 inflammasome inhibitors: Design, synthesis, and biological characterization. *J. Med. Chem.* **62**(21), 9718-9731 (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, 12/19/2022

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