

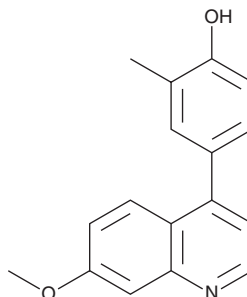
PRODUCT INFORMATION



CU-CPT9a

Item No. 28722

CAS Registry No.: 2165340-32-7
Formal Name: 4-(7-methoxy-4-quinolinyl)-2-methyl-phenol
MF: C₁₇H₁₅NO₂
FW: 265.3
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 316 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

CU-CPT9a is supplied as a crystalline solid. A stock solution may be made by dissolving the CU-CPT9a in the solvent of choice, which should be purged with an inert gas. CU-CPT9a is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CU-CPT9a in these solvents is approximately 5, 10, and 30 mg/ml, respectively.

CU-CPT9a is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CU-CPT9a should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CU-CPT9a has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CU-CPT9a is an antagonist of toll-like receptor 8 (TLR8).¹ It inhibits activation of NF-κB induced by the TLR8 agonist R-848 (Item No. 14806) in TLR8-overexpressing HEK-Blue cells (IC₅₀ = 0.5 nM). CU-CPT9a reverses R-848-induced increases in NF-κB p65, IRAK4, and TRAF3 protein levels in HEK-Blue cells.

Reference

1. Zhang, S., Hu, Z., Tanji, H., et al. Small-molecule inhibition of TLR8 through stabilization of its resting state. *Nat. Chem. Biol.* **14**(1), 58-64 (2018).

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, 09/03/2024

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