

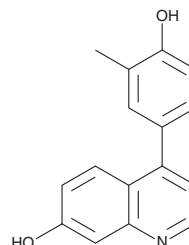
PRODUCT INFORMATION



CU-CPT9b

Item No. 28723

CAS Registry No.: 2162962-69-6
Formal Name: 4-(4-hydroxy-3-methylphenyl)-7-quinolinol
MF: C₁₆H₁₃NO₂
FW: 251.3
Purity: ≥98%
UV/Vis.: λ_{max}: 238 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-CPT9b is supplied as a crystalline solid. A stock solution may be made by dissolving the CU-CPT9b in the solvent of choice, which should be purged with an inert gas. CU-CPT9b is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CU-CPT9b in these solvents is approximately 30 mg/ml.

Description

CU-CPT9b is an antagonist of toll-like receptor 8 (TLR8; K_d = 21 nM).¹ It inhibits activation of NF-κB induced by the TLR8 agonist R-848 (Item No. 14806) in TLR8-overexpressing HEK-Blue cells with an IC₅₀ value of 0.7 nM.

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, 11/02/2022

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