

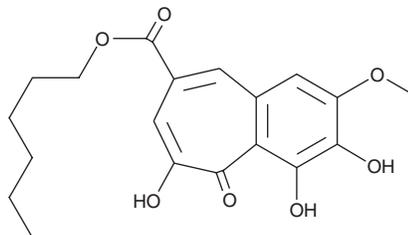
PRODUCT INFORMATION



CU CPT 22

Item No. 15244

CAS Registry No.: 1416324-85-0
Formal Name: 3,4,6-trihydroxy-2-methoxy-5-oxo-5H-benzocycloheptene-8-carboxylic acid, hexyl ester
MF: C₁₉H₂₂O₇
FW: 362.4
Purity: ≥98%
UV/Vis.: λ_{max}: 309, 403 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 CPT 22 is supplied as a crystalline solid. A stock solution may be made by dissolving the CU CPT 22 in the solvent of choice, which should be purged with an inert gas. CU CPT 22 is soluble in the organic solvent DMSO.

Description

CU CPT 22 is an antagonist of toll-like receptor 1 (TLR1) and TLR2 heterodimers (TLR1/TLR2; $K_i = 0.41 \mu\text{M}$).¹ It inhibits nitric oxide (NO) production induced by the TLR2 agonist Pam₃CSK₄ (Item No. 24126) in RAW 264.7 cells ($\text{IC}_{50} = 0.58 \mu\text{M}$). It is selective for TLR1/TLR2 over TLR2/TLR6 and TLR3, -4, and -7 homodimers, as well as 10 kinases at 0.5 μM . CU CPT 22 (8 μM) inhibits Pam₃CSK₄-induced release of TNF- α and IL-1 β in RAW 264.7 cells by 60 and 95%, respectively.

Reference

1. Cheng, K., Wang, X., Zhang, S., *et al.* Discovery of small-molecule inhibitors of the TLR1/TLR2 complex. *Angew Chem. Int. Ed. Engl.* **51(49)**, 12246-12249 (2012).

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, 11/28/2022

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