

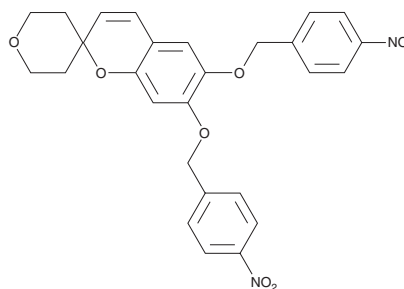
PRODUCT INFORMATION



CU-CPT17e

Item No. 39534

CAS Registry No.: 2109805-75-4
Formal Name: 2',3',5',6'-tetrahydro-6,7-bis[(4-nitrophenyl)methoxy]-spiro[2H-1-benzopyran-2,4'-[4H]pyran]
MF: C₂₇H₂₄N₂O₈
FW: 504.5
Purity: ≥95%
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

CU-CPT17e is supplied as a solid. A stock solution may be made by dissolving the CU-CPT17e in the solvent of choice, which should be purged with an inert gas. CU-CPT17e is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CU-CPT17e in these solvents is approximately 2 mg/ml. CU-CPT17e is slightly soluble in ethanol.

Description

CU-CPT17e is an agonist of toll-like receptor 3 (TLR3), TLR8, and TLR9.¹ It selectively induces NF-κB activation in HEK293 cells expressing recombinant human TLR3, TLR8, or TLR9 (EC₅₀s = 4.8, 13.45, and 5.66 μM, respectively) over HEK293 cells expressing recombinant human TLR2, TLR4, TLR5, or TLR7 (EC₅₀s = >100 μM for all). CU-CPT17e (5-40 μM) induces TNF-α, IL-6, and IL-8 production in THP-1 cells. It also inhibits the growth of HeLa cells (IC₅₀ = 2.71 μM).

Reference

1. Zhang, L., Dewan, V., and Yin, H. Discovery of small molecules as multi-toll-like receptor agonists with proinflammatory and anticancer activities. *J. Med. Chem.* **60**(12), 5029-5044 (2017).

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, 10/26/2023

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