

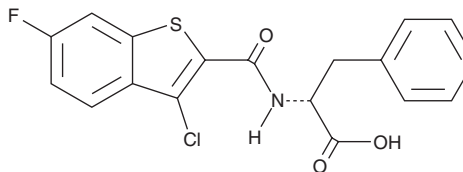
PRODUCT INFORMATION



CU-CPT4a

Item No. 30951

CAS Registry No.: 1279713-77-7
Formal Name: N-[(3-chloro-6-fluorobenzo[b]thien-2-yl)carbonyl]-D-phenylalanine
MF: C₁₈H₁₃ClFNO₃S
FW: 377.8
Purity: ≥98%
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-CPT4a is supplied as a solid. A stock solution may be made by dissolving the CU-CPT4a in the solvent of choice, which should be purged with an inert gas. CU-CPT4a is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

CU-CPT4a is a toll-like receptor 3 (TLR3) antagonist ($IC_{50} = 3.44 \mu\text{M}$).¹ It is selective for TLR3 over TLR4, TLR2/6, TLR1/2, and TLR7 at 27 μM . CU-CPT4a inhibits dsRNA binding to TLR3 with a K_i value of 2.96 μM in a fluorescence anisotropy assay. It decreases poly(I:C)-induced production of TNF- α and IL-1 β in RAW 264.7 cells when used at a concentration of 27 μM .

Reference

1. Cheng, K., Wang, X., and Yin, H. Small-molecule inhibitors of the TLR3/dsRNA complex. *J. Am. Chem. Soc.* **133**(11), 3764-3767 (2011).

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, 12/07/2022

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