

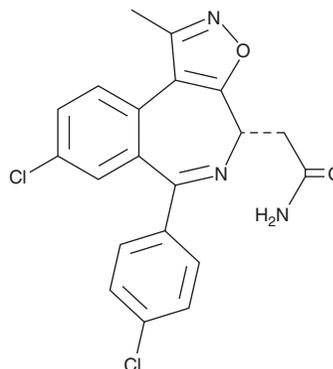
PRODUCT INFORMATION



CPI-268456

Item No. 20973

CAS Registry No.: 1380087-86-4
Formal Name: (4S)-8-chloro-6-(4-chlorophenyl)-1-methyl-4H-isoxazolo[5,4-d][2]benzazepine-4-acetamide
MF: C₂₀H₁₅Cl₂N₃O₂
FW: 400.3
Purity: ≥98%
UV/Vis.: λ_{max}: 256 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

CPI-268456 is supplied as a crystalline solid. A stock solution may be made by dissolving the CPI-268456 in the solvent of choice, which should be purged with an inert gas. CPI-268456 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CPI-268456 in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

CPI-268456 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CPI-268456 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CPI-268456 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CPI-268456 is a ligand of bromodomain-containing protein 4 (BRD4).¹ It binds to BRD4 (IC₅₀ = <0.5 μM in a cell-free assay) and inhibits LPS-induced IL-6 secretion in THP-1 cells (IC₅₀ = <0.5 μM).

Reference

1. Albrecht, B.K., Audia, J.E., Côté, A., *et al.* Bromodomain inhibitors and uses thereof. *Constellation Pharmaceuticals, Inc. US9522920B2* (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, 12/15/2022

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