

PRODUCT INFORMATION

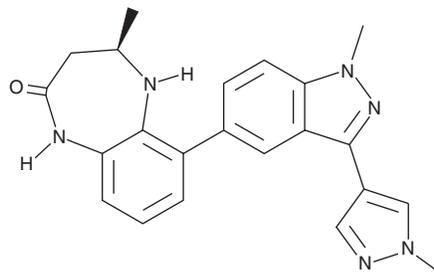


CPI-637

Item No. 23500

CAS Registry No.: 1884712-47-3
Formal Name: (4R)-1,3,4,5-tetrahydro-4-methyl-6-[1-methyl-3-(1-methyl-1H-pyrazol-4-yl)-1H-indazol-5-yl]-2H-1,5-benzodiazepin-2-one

MF: C₂₂H₂₂N₆O
FW: 386.5
Purity: ≥98%
UV/Vis.: λ_{max}: 201, 229, 320 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-637 is supplied as a crystalline solid. A stock solution may be made by dissolving the CPI-637 in the solvent of choice. CPI-637 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of CPI-637 in these solvents is approximately 0.1, 0.5, and 0.2 mg/ml, respectively.

CPI-637 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CPI-637 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CPI-637 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CPI-637 is an inhibitor of cyclic-AMP response element binding protein (CBP) and adenoviral E1A binding protein of 300 kDa (EP300), a homologous pair of bromodomain-containing proteins (IC₅₀s = 30 and 51 nM in a time-resolved-FRET (TR-FRET) assay).¹ It is selective for CBP and EP300 over the BET family of bromodomains (IC₅₀s = >10 μM), however, CPI-637 is active at bromodomain BRD9 (IC₅₀ = 0.73 μM). CPI-637 also inhibits MYC expression (EC₅₀ = 0.6 μM) in AMO-1 cells.

Reference

1. Taylor, A.M., Côté, A., Hewitt, M.C., *et al.* Fragment-based discovery of a selective and cell-active benzodiazepinone CBP/EP300 bromodomain inhibitor (CPI-637). *ACS Med. Chem. Lett.* **7**(5), 531-536 (2016).

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, 12/22/2022

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