

PRODUCT INFORMATION



LH1307

Item No. 28728

CAS Registry No.: 2375720-38-8

Formal Name: N,N'-((((((2,2'-dimethyl-[1,1'-biphenyl]-3,3'-diyl)bis(methylene))bis(oxy))bis(2-((5-cyanopyridin-3-yl)methoxy)-5-methyl-4,1-phenylene))bis(methylene))bis(azanediyl))bis(ethane-2,1-diyl)) diacetamide

MF: C₅₄H₅₈N₈O₆

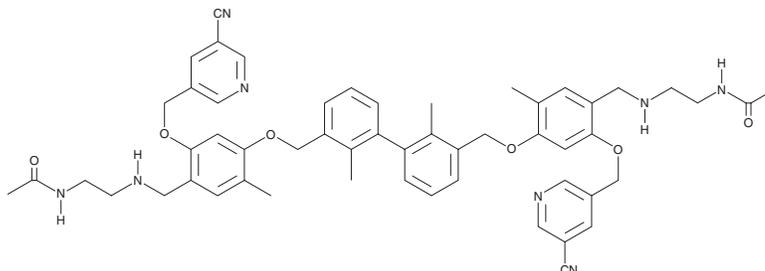
FW: 915.1

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

LH1307 is supplied as a crystalline solid. A stock solution may be made by dissolving the LH1307 in the solvent of choice, which should be purged with an inert gas. LH1307 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of LH1307 in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

LH1307 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LH1307 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LH1307 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LH1307 is an inhibitor of the interaction between programmed cell death 1 (PD-1) and its ligand PD-L1 that has an IC₅₀ value of 3 nM in a homologous time-resolved fluorescence (HTRF) assay.¹ It increases the activation of Jurkat cells expressing PD-1 in co-culture with U2OS or CHO cells expressing PD-L1 (EC₅₀s = 79 and 763 nM, respectively, in reporter assays).

Reference

1. Basu, S., Yang, J., Xu, B., *et al.* Design, synthesis, evaluation, and structural studies of C₂-symmetric small molecule inhibitors of programmed cell death-1/programmed death-ligand 1 protein-protein interaction. *J. Med. Chem.* **62**(15), 7250-7263 (2019).

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, 12/23/2019

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