

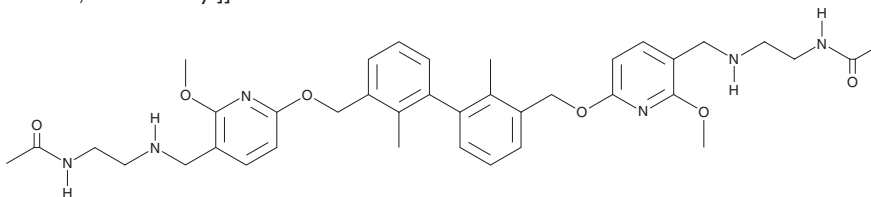
PRODUCT INFORMATION



LH1306

Item No. 28727

CAS Registry No.: 2182653-84-3
Formal Name: N,N'-[(2,2'-dimethyl[1,1'-biphenyl]-3,3'-diyl)bis[methyleneoxy(6-methoxy-2,5-pyridinediyl)methyleneimino-2,1-ethanediyl]]bis-acetamide
MF: C₃₈H₄₈N₆O₆
FW: 684.8
Purity: ≥95%
UV/Vis.: λ_{max}: 283 nm
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

LH1306 is supplied as a crystalline solid. A stock solution may be made by dissolving the LH1306 in the solvent of choice, which should be purged with an inert gas. LH1306 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LH1306 in these solvents is approximately 0.25 and 1 mg/ml, respectively.

Description

LH1306 is an inhibitor of the interaction between programmed cell death 1 (PD-1) and its ligand PD-L1 that has an IC₅₀ value of 25 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 = 334 and 4,214 nM, respectively, in reporter assays).

References

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, 10/21/2019

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