

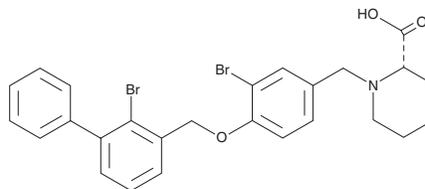
PRODUCT INFORMATION



CAY10774

Item No. 32569

CAS Registry No.: 2714327-65-6
Formal Name: (2S)-1-[[3-bromo-4-[(2-bromo[1,1'-biphenyl]-3-yl)methoxy]phenyl]methyl]-2-piperidinecarboxylic acid
MF: C₂₆H₂₅Br₂NO₃
FW: 559.3
Purity: ≥95%
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

CAY10774 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10774 in the solvent of choice, which should be purged with an inert gas. CAY10774 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10774 in these solvents is approximately 1 mg/ml.

CAY10774 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10774 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10774 has a solubility of approximately 0.33 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

CAY10774 is an inhibitor of the protein-protein interaction between programmed cell death 1 (PD-1) and its ligand PD-L1 (IC₅₀ = 15 nM in a homologous time-resolved fret (HTRF) assay).¹ It increases the activation of Jurkat cells expressing PD-1 in co-culture with artificial antigen-presenting cells (aAPCs) expressing PD-L1 (EC₅₀ = 6.6 μM in a reporter assay). CAY10774 increases surface expression of PD-1 on primary human CD4⁺ and CD8⁺ T cells co-cultured with aAPCs.

Reference

1. Konieczny, M., Musielak, B., Kocik, J., *et al.* Di-bromo-based small-molecule inhibitors of the PD-1/PD-L1 immune checkpoint. *J. Med. Chem.* **63**(19), 11271-11285 (2020).

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, 03/21/2024

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