

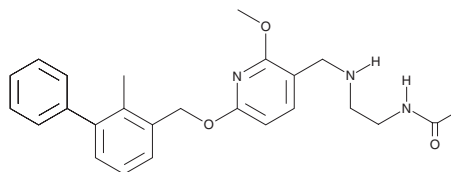
PRODUCT INFORMATION



PD-1/PD-L1 Inhibitor 2

Item No. 19132

CAS Registry No.: 1675203-84-5
Formal Name: N-[2-[[[2-methoxy-6-[(2-methyl[1,1'-biphenyl]-3-yl)methoxy]-3-pyridinyl)methyl]amino]ethyl]-acetamide
Synonym: Programmed Cell Death 1/Programmed Cell Death-Ligand 1 Inhibitor 2
MF: C₂₅H₂₉N₃O₃
FW: 419.5
Purity: ≥98%
UV/Vis.: λ_{max}: 281 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

PD-1/PD-L1 inhibitor 2 is supplied as a crystalline solid. A stock solution may be made by dissolving the PD-1/PD-L1 inhibitor 2 in the solvent of choice, which should be purged with an inert gas. PD-1/PD-L1 inhibitor 2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PD-1/PD-L1 inhibitor 2 in these solvents is approximately 5, 2, and 10 mg/ml, respectively.

PD-1/PD-L1 inhibitor 2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PD-1/PD-L1 inhibitor 2 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PD-1/PD-L1 inhibitor 2 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The programmed death-1/programmed death-ligand 1 (PD-1/PD-L1) interaction plays a dominant role in the suppression of T cell responses, especially in a tumor microenvironment, protecting tumor cells from lysis.¹ PD-1/PD-L1 inhibitor 2 is reported to prevent the interaction of PD-L1 with PD-1 with an IC₅₀ value of 18 nM.²

References

1. Pardoll, D.M. The blockade of immune checkpoints in cancer immunotherapy. *Nat. Rev. Cancer* **12**(4), 252-264 (2012).
2. Chupak, L.S. and Zheng, X. Compounds useful as immunomodulators. *Bristol-Myers Squibb WO2015/034820A1* (2015)

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, 09/30/2022

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