

PRODUCT INFORMATION



PD-1/PD-L1 Inhibitor N11

Item No. 39663

Formal Name: 2-methyl-2-(((4-(4-phenylindolin-1-yl)pyrido[3,2-d]pyrimidin-7-yl)methyl)amino)propan-1-ol

Synonyms: PD-1/PD-L1 Interaction Inhibitor, Programmed Cell Death 1/Programmed Cell Death-Ligand 1 Inhibitor N11

MF: C₂₆H₂₇N₅O

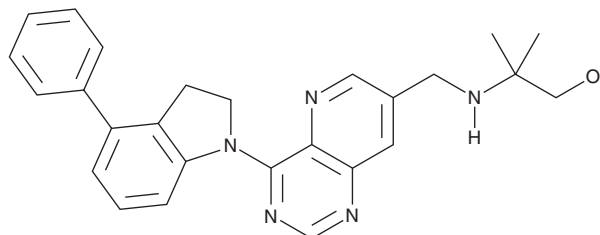
FW: 425.5

Purity: ≥95%

Supplied as: A 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

PD-1/PD-L1 inhibitor N11 is supplied as a solid. A stock solution may be made by dissolving the PD-1/PD-L1 inhibitor N11 in the solvent of choice, which should be purged with an inert gas. PD-1/PD-L1 inhibitor N11 is soluble in acetonitrile and DMSO.

Description

PD-1/PD-L1 inhibitor N11 is an inhibitor of the protein-protein interaction between programmed cell death protein 1 (PD-1) and PD-1 ligand (PD-L1; IC₅₀ = 6.3 nM).¹ It increases T cell proliferation in a PD-L1-overexpressing CT26 colon cancer cell and T cell co-culture when used at a concentration of 20 nM. PD-1/PD-L1 inhibitor N11 (20 and 80 nM) promotes T cell infiltration into an *in vitro* PD-L1-overexpressing murine Lewis lung carcinoma tumor matrix. *In vivo*, PD-1/PD-L1 inhibitor N11 (20 mg/kg) reduces tumor growth and promotes CD8⁺ and CD4⁺ T cell tumor infiltration in a PD-L1-overexpressing murine Lewis lung carcinoma model.

Reference

- Meng, Y., Wang, Y., Zhang, S., *et al.* Novel 4-arylindolines containing a pyrido[3,2-d]pyrimidine moiety as the programmed cell death-1/programmed cell death-ligand 1 interaction inhibitors for tumor immunotherapy. *J. Med. Chem.* **66**(17), 11815-11830 (2023).

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, 01/04/2024

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