

PRODUCT INFORMATION



ARB-272572

Item No. 42547

CAS Registry No.: 2368182-63-0
Formal Name: N,N'-(2,2'-dimethyl[1,1'-

MF: C₃₂H₃₆N₆O₄

FW: 568.7

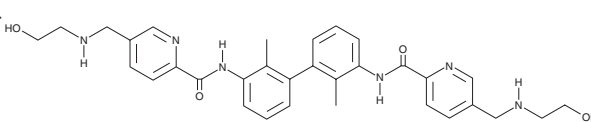
Purity: ≥98%

Supplied as: A 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

ARB-272572 is supplied as a solid. A stock solution may be made by dissolving the ARB-272572 in the solvent of choice, which should be purged with an inert gas. ARB-272572 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

ARB-272572 is an inhibitor of the protein-protein interaction between programmed cell death protein 1 (PD-1) and PD-1 ligand (PD-L1; IC₅₀ = 400 pM in a homogenous time-resolved fluorescence (HTRF) assay).¹ It also inhibits the PD-1/PD-L1 interaction in a cell-based nuclear factor of activated T cells (NFAT) reporter assay (IC₅₀ = 17 nM) and decreases cell surface levels of PD-L1 in primary human dendritic cells and B cells. ARB-272572 reduces cytomegalovirus antigen-induced IFN-γ production in a cytomegalovirus (CMV) recall assay using peripheral blood mononuclear cells (PBMCs) isolated from CMV seropositive patients (IC₅₀ = 3 nM). In the presence of hepatitis B virus (HBV) peptides, it induces T cell proliferation and IFN-γ production in PBMCs isolated from patients with chronic HBV infection when used at a concentration of 500 nM. *In vivo*, ARB-272572 (10 mg/kg) reduces tumor volume and PD-L1 levels in CD45⁺ tumor cells in a humanized MC-38 murine model of colon cancer.

Reference

1. Park, J.-J., Thi, E.P., Carpio, V.H., *et al.* Checkpoint inhibition through small molecule-induced internalization of programmed death-ligand 1. *Nat. Commun.* **12**(1), 1222 (2021).

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, 02/21/2025

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