

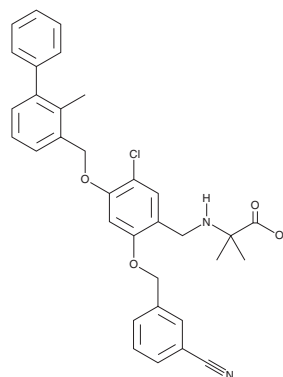
# PRODUCT INFORMATION



## NP19

Item No. 31791

**CAS Registry No.:** 2377916-66-8  
**Formal Name:** N-[[5-chloro-2-[(3-cyanophenyl)methoxy]-4-[[2-methyl[1,1'-biphenyl]-3-yl)methoxy]phenyl]methyl]-2-methyl-alanine  
**MF:** C<sub>33</sub>H<sub>31</sub>ClN<sub>2</sub>O<sub>4</sub>  
**FW:** 555.1  
**Purity:** ≥98%  
**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

NP19 is supplied as a crystalline solid. A stock solution may be made by dissolving the NP19 in the solvent of choice, which should be purged with an inert gas. NP19 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of NP19 in these solvents is approximately 30 mg/ml.

### Description

NP19 is an inhibitor of the protein-protein interaction between programmed cell death 1 (PD-1) and its ligand PD-L1 (IC<sub>50</sub> = 12.5 nM in a homogenous time-resolved fluorescence (HTRF) assay).<sup>1</sup> It increases IFN-γ production by isolated human peripheral blood mononuclear cells (PBMCs) in co-culture with anti-CD3 single-chain variable fragment- and PD-L1-expressing Hep3B hepatoma cells (EC<sub>50</sub> = ~3 μM). NP19 (25, 50, and 100 mg/kg) reduces tumor growth in a B16/F10 murine melanoma model and an H22 murine hepatocellular carcinoma model.

### Reference

1. Cheng, B., Ren, Y., Niu, X., *et al.* Discovery of novel resorcinol dibenzyl ethers targeting the programmed cell death-1/programmed cell death-ligand 1 interaction as potential anticancer agents. *J. Med. Chem.* **63**(15), 8338-8358 (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