

# PRODUCT INFORMATION



## BMS-1001 (hydrochloride)

Item No. 27891

**CAS Registry No.:** 2113650-04-5  
**Formal Name:** N-[[2-[[3-(3-cyanophenyl)methoxy]-4-[[3-(2,3-dihydro-1,4-benzodioxin-6-yl)-2-methylphenyl]methoxy]-5-methylphenyl]methyl]-D-serine, monohydrochloride

**MF:** C<sub>35</sub>H<sub>34</sub>N<sub>2</sub>O<sub>7</sub> • HCl

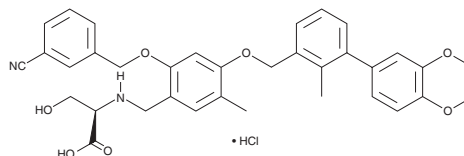
**FW:** 631.1

**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

BMS-1001 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS-1001 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BMS-1001 (hydrochloride) is soluble in DMSO.

### Description

BMS-1001 is a small molecule inhibitor of the interaction between programmed death protein 1 (PD-1) and its ligand programmed cell death ligand 1 (PD-L1) with an IC<sub>50</sub> value of 2.25 nM in a homogeneous time-resolved fluorescence (HTRF) binding assay.<sup>1</sup>

### Reference

1. Guzik, K., Zak, K.M., Grudnik, P., *et al.* Small-molecule inhibitors of the programmed cell death-1/programmed death-ligand 1 (PD-1/PD-L1) interaction via transiently induced protein states and dimerization of PD-L1. *J. Med. Chem.* **60**(13), 5857-5867 (2017).

#### 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, 12/19/2022

#### 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