

PRODUCT INFORMATION



BMS 1166 (hydrochloride)

Item No. 27774

CAS Registry No.: 2113650-05-6
Formal Name: (4R)-1-[[5-chloro-2-[(3-cyanophenyl)methoxy]-4-[[3-(2,3-dihydro-1,4-benzodioxin-6-yl)-2-methylphenyl]methoxy]phenyl]methyl]-4-hydroxy-D-proline, monohydrochloride

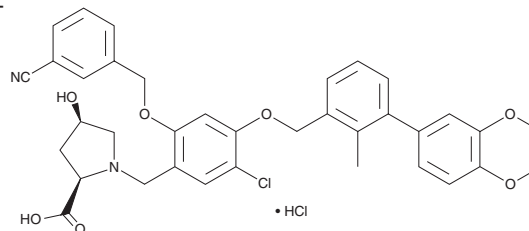
MF: C₃₆H₃₃ClN₂O₇ • HCl
FW: 677.6

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

BMS 1166 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the BMS 1166 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BMS 1166 (hydrochloride) is soluble in the organic solvent DMSO.

Description

BMS 1166 is an inhibitor of the protein-protein interaction between programmed cell death 1 (PD-1) and its ligand PD-L1 that has an IC₅₀ value of 1.4 nM in a homologous time-resolved fluorescence (HTRF) assay.¹ It increases the activation of Jurkat cells expressing PD-1 in co-culture with CHO cells expressing PD-L1 (EC₅₀ = 276 nM in a reporter assay).²

References

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).
2. Skalniak, L., Zak, K.M., Guzik, K., *et al.* Small-molecule inhibitors of PD-1/PD-L1 immune checkpoint alleviate the PD-L1-induced exhaustion of T-cells. *Oncotarget* **8**(42), 72167-72181 (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