

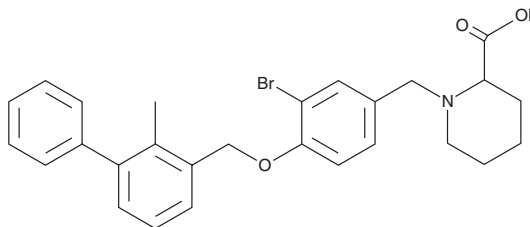
PRODUCT INFORMATION



BMS-8

Item No. 22355

CAS Registry No.: 1675201-90-7
Formal Name: 1-[[3-bromo-4-[(2-methyl[1,1'-biphenyl]-3-yl)methoxy]phenyl]methyl]-2-piperidinecarboxylic acid
MF: C₂₇H₂₈BrNO₃
FW: 494.4
Purity: ≥98%
UV/Vis.: λ_{max}: 234 nm
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-8 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS-8 in the solvent of choice, which should be purged with an inert gas. BMS-8 is slightly soluble in organic solvent methanol.

Description

BMS-8 is a small molecule inhibitor of the protein-protein interaction between programmed death protein 1 (PD-1) and its ligand programmed cell death ligand 1 (PD-L1) with an IC₅₀ value of 146 nM in a homogeneous time-resolved fluorescence (HTRF) binding assay.¹

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