

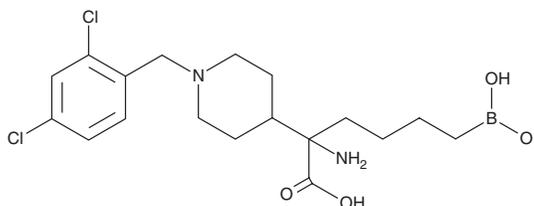
PRODUCT INFORMATION



CB-1158

Item No. 27297

CAS Registry No.: 1345810-21-0
Formal Name: α -amino- α -(4-boronobutyl)-1-
[(2,4-dichlorophenyl)methyl]-4-
piperidineacetic acid
Synonyms: INCB 01158, Numidargistat
MF: C₁₈H₂₇BCl₂N₂O₄
FW: 417.1
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CB-1158 is supplied as a solid. A stock solution may be made by dissolving the CB-1158 in the solvent of choice, which should be purged with an inert gas. CB-1158 is soluble in the organic solvent DMSO.

Description

CB-1158 is an inhibitor of arginase I and II (IC₅₀s = 0.086 and 0.296 μ M for the human enzymes, respectively).¹ It is selective for arginase I and II over neuronal nitric oxide synthase (nNOS), endothelial NOS (eNOS), and inducible NOS (iNOS; IC₅₀s = >50 μ M for all). CB-1158 reverses inhibition of T cell proliferation induced by granulocytes *in vitro* (IC₅₀ = 0.204 μ M). It reduces tumor growth in a CT26 murine colon cancer model when administered at a dose of 100 mg/kg twice per day.

Reference

1. Steggerda, S.M., Bennett, M.K., Chen, J., *et al.* Inhibition of arginase by CB-1158 blocks myeloid cell-mediated immune suppression in the tumor microenvironment. *J. Immunother. Cancer* **5**(1), 101 (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, 12/22/2022

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