

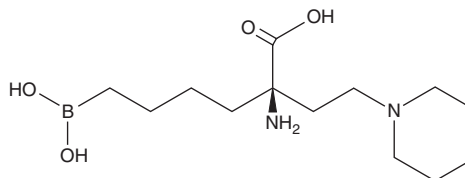
PRODUCT INFORMATION



Arginase Inhibitor 1

Item No. 28418

CAS Registry No.: 1345808-25-4
Formal Name: (αR)-α-amino-α-(4-boronobutyl)-
1-piperidinebutanoic acid
MF: C₁₃H₂₇BN₂O₄
FW: 286.2
Purity: ≥90%
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

Arginase inhibitor 1 is supplied as a solid. A stock solution may be made by dissolving the arginase inhibitor 1 in the solvent of choice, which should be purged with an inert gas. Arginase inhibitor 1 is soluble in organic solvents such as DMSO. It is also soluble in water. The solubility of arginase inhibitor 1 in DMSO and water is approximately 50 and 30 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Arginase inhibitor 1 is an inhibitor of arginase I and II (IC₅₀s = 223 and 509 nM, respectively, for the recombinant human enzymes).¹ It decreases arginase I activity in CHOK cells expressing human arginase I (IC₅₀ = 8 μM). Arginase inhibitor 1 (100 mg/kg, i.v.) reduces infarct size in a rat model of myocardial ischemia and reperfusion injury induced by left-main coronary artery occlusion.

Reference

1. Van Zandt, M.C., Whitehouse, D.L., Golebiowski, A., *et al.* Discovery of (R)-2-amino-6-borono-2-(2-(piperidin-1-yl)ethyl)hexanoic acid and congeners as highly potent inhibitors of human arginases I and II for treatment of myocardial reperfusion injury. *J. Med. Chem.* **56(6)**, 2568-2580 (2013).

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, 12/08/2022

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