

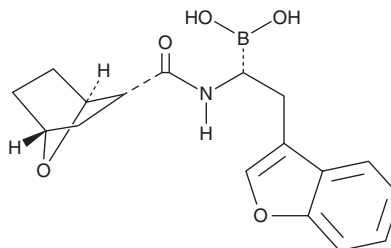
PRODUCT INFORMATION



M-3258

Item No. 38160

CAS Registry No.: 2285330-15-4
Formal Name: B-[(1R)-2-(3-benzofuranyl)-1-[[[(1S,2R,4R)-7-oxabicyclo[2.2.1]hept-2-ylcarbonyl]amino]ethyl]-boronic acid
MF: C₁₇H₂₀BNO₅
FW: 329.2
Purity: ≥98%
UV/Vis.: λ_{max}: 248 nm
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

M-3258 is supplied as a solid. A stock solution may be made by dissolving the M-3258 in the solvent of choice, which should be purged with an inert gas. M-3258 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

M-3258 is an inhibitor of the immunoproteasome subunit large multifunctional peptidase 7 (LMP7; IC₅₀ = 4.1 nM).¹ It is selective for LMP7 over the 20S proteasomal catalytic subunits β1, -β2, and -β5 (IC₅₀s = >10,000, 2,519, and >10,000 nM, respectively), as well as LMP2 and proteasome subunit beta type 10 (PSMB10; IC₅₀s = >10,000 nM for both). It increases ubiquitinated protein levels (EC₅₀ = 1,980 nM) and induces apoptosis in MM.1S multiple myeloma cells (EC₅₀ = 420 nM). M-3258 (10 mg/kg per day) reduces tumor volume in U266B1, OPM-2, and RPMI-8226 mouse xenograft models of multiple myeloma.

Reference

1. Sanderson, M.P., Friese-Hamim, M., Walter-Bausch, G., *et al.* M3258 is a selective inhibitor of the immunoproteasome subunit LMP7 (β5i) delivering efficacy in multiple myeloma models. *Mol. Cancer Ther.* 20(8), 1378-1387 (2021).

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, 07/26/2023

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