

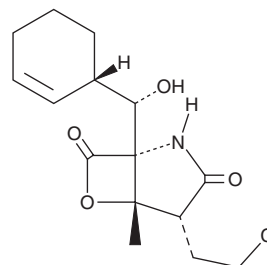
PRODUCT INFORMATION



Salinosporamide A

Item No. 10007311

CAS Registry No.: 437742-34-2
Formal Name: (1R,4R,5S)-4-(2-chloroethyl)-1-[(S)-(1S)-2-cyclohexen-1-ylhydroxymethyl]-5-methyl-6-oxa-2-azabicyclo[3.2.0]heptane-3,7-dione
Synonyms: Marizomib, ML-858, NPI-0052
MF: C₁₅H₂₀ClNO₄
FW: 313.8
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥1 year



Special Conditions: Keep under inert gas. Protect from light. Protect from moisture and oxygen.

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

Laboratory Procedures

Salinosporamide A is supplied as a solid. A stock solution may be made by dissolving the salinosporamide A in the solvent of choice. Salinosporamide A is soluble in the organic solvent DMSO.

Description

Salinosporamide A is a natural proteasome inhibitor.¹ It is a β -lactone- γ -lactam that irreversibly inhibits all three protease activities within the proteasome (IC_{50} s = 3.5, 28, and 430 nM for β 5 (chymotrypsin), β 2 (macropain), and β 1 (C5) subunits, respectively), both *in vitro* and *in vivo*.^{2,3} Salinosporamide A has potential applications in cancer therapy, particularly in combination with other chemotherapeutics.^{1,2}

References

1. Kubiczikova, L., Pour, L., Sedlarikova, L., *et al.* Proteasome inhibitors - molecular basis and current perspectives in multiple myeloma. *J. Cell. Mol. Med.* **18(6)**, 947-961 (2014).
2. Chauhan, D., Catley, L., Li, G., *et al.* A novel orally active proteasome inhibitor induces apoptosis in multiple myeloma cells with mechanisms distinct from bortezomib. *Cancer Cell* **8(5)**, 407-419 (2005).
3. Kale, A.J. and Moore, B.S. Molecular mechanisms of acquired proteasome inhibitor resistance. *J. Med. Chem.* **55(23)**, 10317-10327 (2012).

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, 09/27/2017

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