

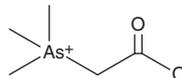
PRODUCT INFORMATION



Arsenobetaine

Item No. 34515

CAS Registry No.: 64436-13-1
Formal Name: (carboxymethyl)trimethyl-arsonium, inner salt
MF: C₅H₁₁AsO₂
FW: 178.1
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Arsenobetaine is supplied as a solid. A stock solution may be made by dissolving the arsenobetaine in water. We do not recommend storing the aqueous solution for more than one day.

Description

Arsenobetaine is an organoarsenical and a compatible solute that has been found in various marine animals, such as lobsters and crabs, as well as terrestrial organisms, including earthworms and lichens.^{1,2} It is protective against *B. subtilis* cell death induced by high osmolarity or extreme temperatures when used at a concentration of 1 mM.² Arsenobetaine is non-toxic to mice (LD₅₀ = 10 g/kg).^{1,2}

References

1. Popowich, A., Zhang, Q., and Le, X. Arsenobetaine: The ongoing mystery. *Natl. Sci. Rev.* **3(4)**, 451-458 (2016).
2. Hoffmann, T., Warmbold, B., Smits, S.H.J., *et al.* Arsenobetaine: An ecophysiological important organoarsenical confers cytoprotection against osmotic stress and growth temperature extremes. *Environ. Microbiol.* **20(1)**, 305-323 (2018).

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, 11/14/2022

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