

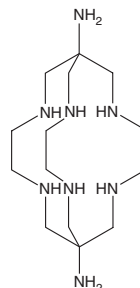
PRODUCT INFORMATION



DiAmSar

Item No. 43826

CAS Registry No.: 91002-72-1
Formal Name: 3,6,10,13,16,19-hexaazabicyclo[6.6.6]
eicosane-1,8-diamine
Synonyms: Diaminosarcophagine, (NH₂)₂-Sar
MF: C₁₄H₃₄N₈
FW: 314.5
Purity: ≥95%
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

DiAmSar is supplied as a solid. A stock solution may be made by dissolving the DiAmSar in the solvent of choice, which should be purged with an inert gas. DiAmSar is soluble (≥10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of DiAmSar can be prepared by directly dissolving the solid in aqueous buffers. DiAmSar is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

DiAmSar is a metal chelator.^{1,2} It has been conjugated to various tumor-targeting ligands and complexed with the copper radioisotope ⁶⁴Cu for PET imaging *in vivo*.

References

1. Wei, L., Ye, Y., Wadas, T.J., *et al.* ⁶⁴Cu-labeled CB-TE2A and diamsar-conjugated RGD peptide analogs for targeting angiogenesis: comparison of their biological activity. *Nucl. Med. Biol.* **36(3)**, 277-285 (2009).
2. Huang, C.-W., Li, Z., Cai, H., *et al.* Biological stability evaluation of the α2β1 receptor imaging agents: Diamsar and DOTA conjugated DGEA peptide. *Bioconjug. Chem.* **22(2)**, 256-263 (2011).

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, 08/28/2025

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