

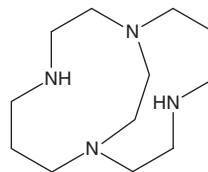
PRODUCT INFORMATION



CB-Cyclam

Item No. 43825

CAS Registry No.: 130701-19-8
Formal Name: 1,4,8,11-tetraazabicyclo[6.6.2]hexadecane
Synonym: Cross-bridged Cyclam
MF: C₁₂H₂₆N₄
FW: 226.4
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

CB-cyclam is supplied as a solid. A stock solution may be made by dissolving the CB-cyclam in the solvent of choice, which should be purged with an inert gas. CB-cyclam is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in 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 CB-cyclam can be prepared by directly dissolving the solid powder in aqueous buffers. CB-cyclam is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

CB-cyclam is a metal chelator.¹ It has also been used in the synthesis of bifunctional chelators for copper radioisotopes.²

References

1. Lelong, E., Suh, J.-M., Kim, G., *et al.* Complexation of C-functionalized cyclams with copper(II) and zinc(II): Similarities and changes when compared to parent cyclam analogues. *Inorg. Chem.* **60**(15), 10857-10872 (2021).
2. Urbanovský, P., David, T., Hlinová, V., *et al.* Cross-bridged cyclam derivatives with bis(phosphinate) and phosphinate-phosphonate pendant arms (cb-BPC) as chelators for copper radioisotopes. *Org. Biomol. Chem.* **23**(3), 699-720 (2025).

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, 08/29/2025

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