

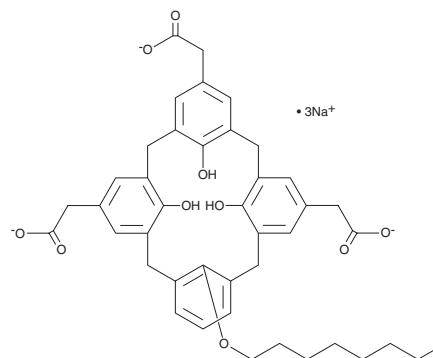
PRODUCT INFORMATION



CALX8

Item No. 24777

CAS Registry No.: 1872296-37-1
Formal Name: 26,27,28-trihydroxy-25-(octyloxy)-pentacyclo[19.3.1.1^{3,7}.1^{9,13}.1^{15,19}]octacos-1(25),3,5,7(28),9,11,13(27),15,17,19(26),21,23-dodecaene-5,11,17-triacetic acid, trisodium salt
MF: C₄₂H₄₃O₁₀ • 3Na
FW: 776.8
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

CALX8 is supplied as a powder. A stock solution may be made by dissolving the CALX8 in the solvent of choice, which should be purged with an inert gas. CALX8 is soluble in organic solvents such as methanol and DMSO. CALX8 is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

CALX8 is a calixarene detergent that has been used to solubilize membrane proteins.¹ It has a critical micelle concentration (CMC) of 2.2 mM.

Image



Membrane protein solubilization from Sf9 membranes.
The target was extracted from Sf9 membranes by using CALX8 reagent at 10-fold the critical micelle concentration (CMC). After solubilization, samples were centrifuged at 100,000 g. Proteins from pellets (P) and supernatants (S) were separated on a 4-15% Tris-glycine SDS-PAGE, transferred to PVDF membrane and immunodetected with a specific antibody. T = total, P = pellet, S = supernatant.

Reference

1. Desuzinges Mandon, E., Agez, M., Pellegrin, R., et al. Novel systematic detergent screening method for membrane proteins solubilization. *Anal. Biochem.* **517**, 40-49 (2017).

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, 07/21/2025

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