

# PRODUCT INFORMATION



## CALXCHOL

Item No. 24783

**CAS Registry No.:** 2121543-66-4  
**Formal Name:** 25-[[1-(3 $\alpha$ )-cholest-5-en-3-yl-1H-1,2,3-triazol-4-yl]methoxy]-26,27,28-trihydroxy-pentacyclo[19.3.1.1<sup>3,7</sup>.1<sup>9,13</sup>.1<sup>15,19</sup>]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:** C<sub>64</sub>H<sub>74</sub>N<sub>3</sub>O<sub>10</sub> • 3Na

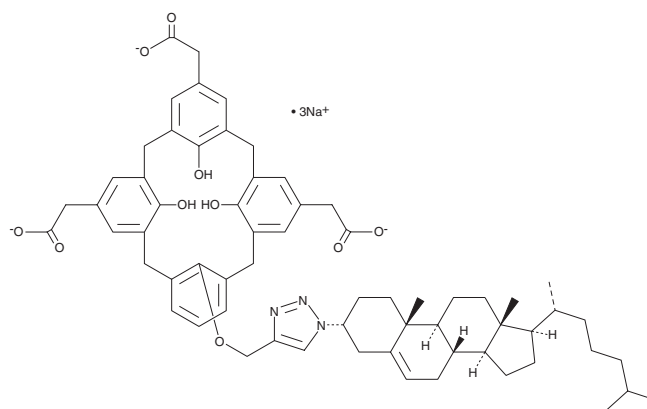
**FW:** 1,114.3

**Purity:** ≥85%

**Supplied as:** A powder

**Storage:** -20°C

**Stability:** ≥4 years



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

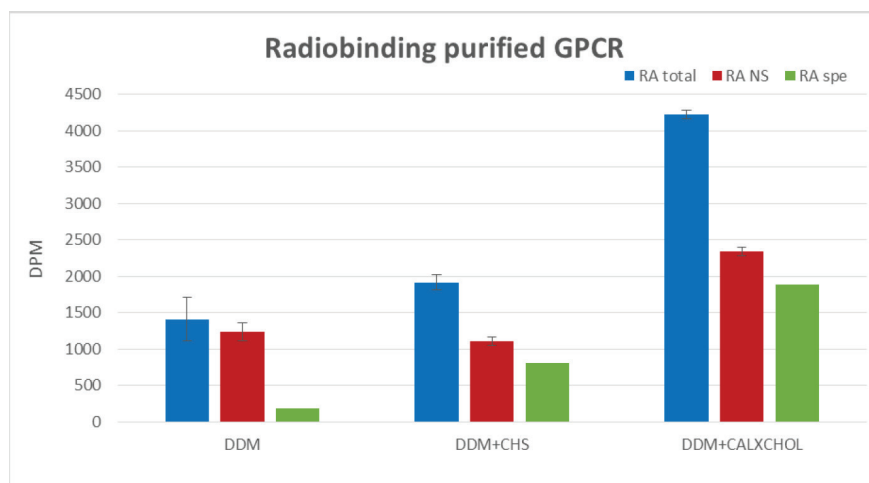
### Laboratory Procedures

CALXCHOL is supplied as a powder. A stock solution may be made by dissolving the CALXCHOL in water (up to 4%). We do not recommend storing the aqueous solution for more than one day.

### Description

CALXCHOL is a calixarene detergent that can be used to solubilize and stabilize membrane proteins.

### Image



Binding of radioligand on GPCR protein, purified in reference detergent (DDM) with or without CHS or CALXCHOL. Purified protein was incubated with radioligand in absence (total, blue bars) or presence (non-specific signal, red bars) of an excess of cold ligand. After filtration on GF/C membranes and washing, scintillation agent was added and radioactivity was detected using a Microbeta2. Specific radioactivity (green bars) corresponds to (total signal) - (non-specific signal).

#### 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, 12/19/2022

#### 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