



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

---

A Geno Technology, Inc. (USA) brand name

# CALXCHOL

(Cat. # 786-1521, 786-1522 & 786-1523)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1521	CALXCHOL	10 mg
786-1522	CALXCHOL	50 mg
786-1523	CALXCHOL	100mg

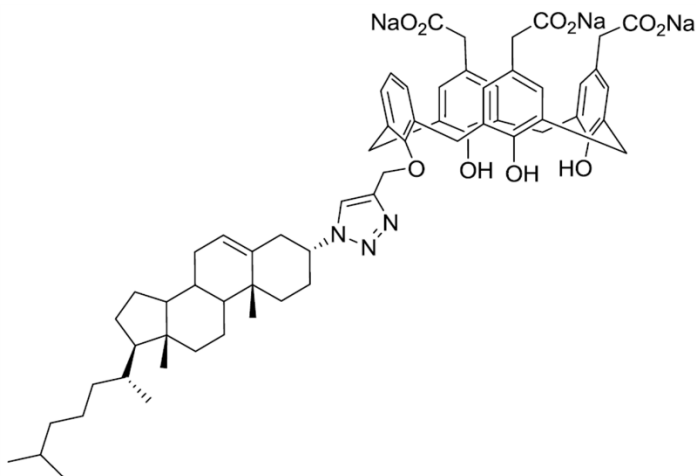
## REAGENT INFORMATION

- Compound Name: CALXCHOL
- IUPAC Name: 5,11,17-tris[(carboxy)methyl]-25-monomethoxytriazolo-(3 $\alpha$ -cholesteryl)-26,27,28-trihydroxycalix[4]arene
- Application: Membrane Protein Extraction
- Surfactant type: Calixarene
- Molecular Formula: C<sub>64</sub>H<sub>74</sub>N<sub>3</sub>Na<sub>3</sub>O<sub>10</sub>
- Molecular Weight: 1114.2 g/mol
- Percent composition: C (64.94%), H (5.58%), N (8.88%), O (20.60%)
- Physical state: Beige powder
- Purity (HPLC, 214nm): 85%
- Retention Time (RP<sub>18</sub> HPLC): t<sub>R</sub> = 22.9 min

## STORAGE AND HANDLING

- Storage conditions: Store in <-20°C freezer for up to one year
- Solubility: Soluble in water (up to 4%)

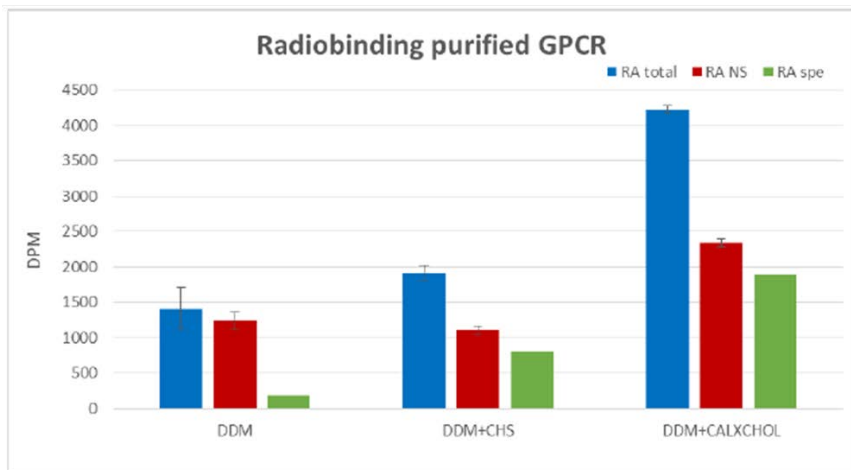
## STRUCTURE



## BIOCHEMICAL VALIDATION DATA

### 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).



## REFERENCES

1. Proprietary patent: "Method for the synthesis of steroidal calixarene compounds and their use for extraction, solubilization and/or stabilization of native and functional membrane proteins"; WO 2017134087 A1 ; Publication date: August 10, 2017.

Last saved: 7/17/2018 EM



[www.GBiosciences.com](http://www.GBiosciences.com)