



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

---

A Geno Technology, Inc. (USA) brand name

# DDLAC

(Cat. # 786-1527, 786-1528 & 786-1529)



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

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1527	DDLAC	250 mg
786-1528	DDLAC	500 mg
786-1529	DDLAC	1 g

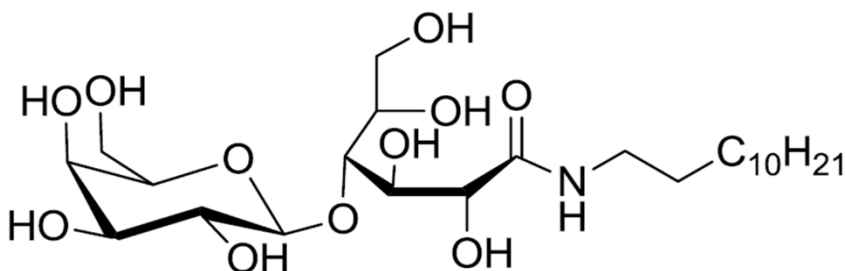
## REAGENT INFORMATION

- Compound Name: DDLAC
- IUPAC Name: N-dodecylamine lactobionamide
- Application: Membrane Protein Extraction
- Molecular Formula:  $C_{24}H_{47}NO_{11}$
- Molecular Weight: 525.6 g/mol
- Percent composition: C, 54.84; H, 9.01; N, 2.66; O, 33.48
- Physical state: White powder
- Purity (HPLC, 214nm): <95%
- Retention Time (RP<sub>18</sub> HPLC):  $t_R = 14.9$  min
- Critical Micelle Concentration: 0.25mM

## STORAGE AND HANDLING

- Storage conditions: Store in <-20°C freezer for up to one year
- Solubility: Soluble in water (1.25mM), methanol and DMSO.

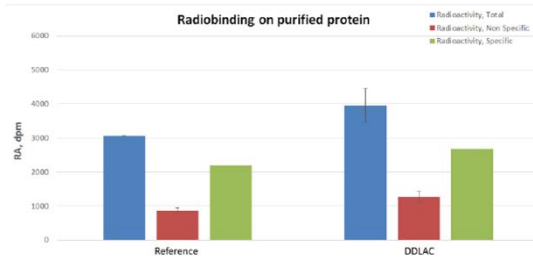
## STRUCTURE



## BIOCHEMICAL VALIDATION DATA

### Binding of radioligand on GPCR protein, purified in reference detergent with or without addition of DDLAC as an additive.

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. Lebaupain et al., Langmuir, 2006, 22 (21), pp 8881–8890

Last saved: 7/17/2018 EM



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