



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

---

A Geno Technology, Inc. (USA) brand name

# DLAC

(Cat. # 786-1509, 786-1510 & 786-1511)



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

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1509	DLAC	250 mg
786-1510	DLAC	500 mg
786-1511	DLAC	1 g

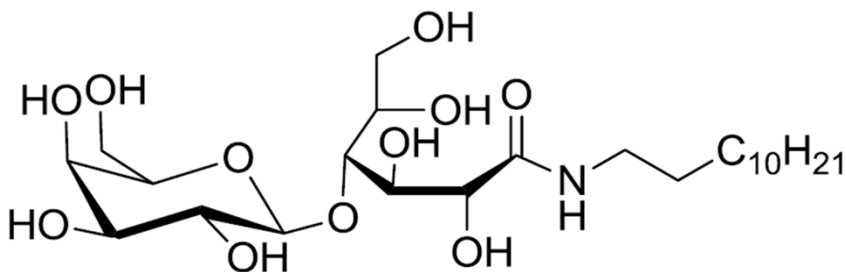
## REAGENT INFORMATION

- Compound Name: DLAC
- IUPAC Name: N-decylamine lactobionamide
- Application: Membrane Protein Extraction
- Molecular Formula:  $C_{22}H_{43}NO_{11}$
- Molecular Weight: 497.6 g/mol
- Percent composition: C, 53.1; H, 8.71; N, 2.81; O, 35.37
- Physical state: White powder
- Purity (HPLC, 214nm): 97%
- Retention Time (RP<sub>18</sub> HPLC):  $t_R = 10.6$  min
- Critical Micelle Concentration: 1.3mM

## STORAGE AND HANDLING

- Storage conditions: Store in  $<-20^{\circ}C$  freezer for up to one year
- Solubility: Soluble in water (65mM), methanol and DMSO.

## STRUCTURE

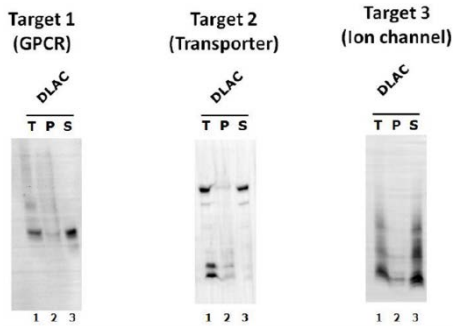


## BIOCHEMICAL VALIDATION DATA

### Membrane proteins solubilization.

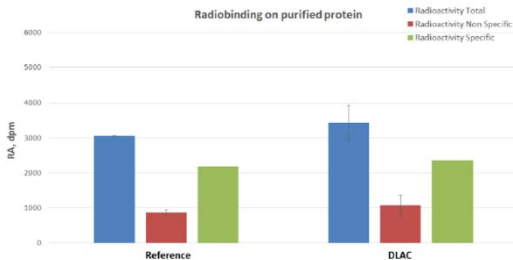
The 3 targets were extracted from Sf9 membranes (GPCR), E. coli membranes (Transporter) and mammalian membranes (ion channel) by using DLAC reagent at 10-fold the critical micelle concentration (cmc). After solubilization, samples were centrifuged at 100000g. Proteins from pellets (P) and supernatants (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.



### Binding of radioligand on GPCR protein, purified in reference detergent or in DLAC.

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)