

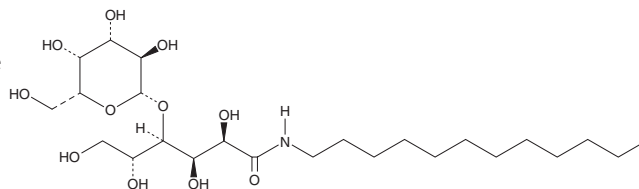
# PRODUCT INFORMATION



## DDLAC

Item No. 24785

**CAS Registry No.:** 69313-68-4  
**Formal Name:** N-dodecyl-4-O-β-D-galactopyranosyl-D-gluconamide  
**MF:** C<sub>22</sub>H<sub>47</sub>NO<sub>11</sub>  
**FW:** 525.6  
**Purity:** ≥95%  
**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.

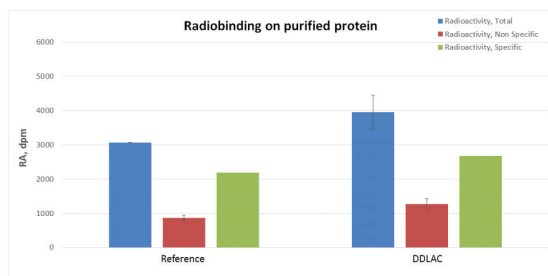
## Laboratory Procedures

DDLAC is supplied as a powder. A stock solution may be made by dissolving the DDLAC in the solvent of choice, which should be purged with an inert gas. DDLAC is soluble in organic solvents such as methanol and DMSO. DDLAC is also soluble in water at a concentration of approximately 1.25 mM. We do not recommend storing the aqueous solution for more than one day.

## Description

DDLAC is a detergent synthesized from lactobionic acid.<sup>1,2</sup> It can be used to stabilize membrane proteins and has a critical micelle concentration (CMC) of 0.25 mM.

## Image



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. El Ghoul, M., Escoula, B., Rico, I., et al. *J. Fluorine Chem.* **59(1)**, 107-112 (1992).
2. Lebaupain, F., Salvay, A.G., Oliver, B.B., et al. *Langmuir* **22(21)**, 8881-8890 (2006).

### 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