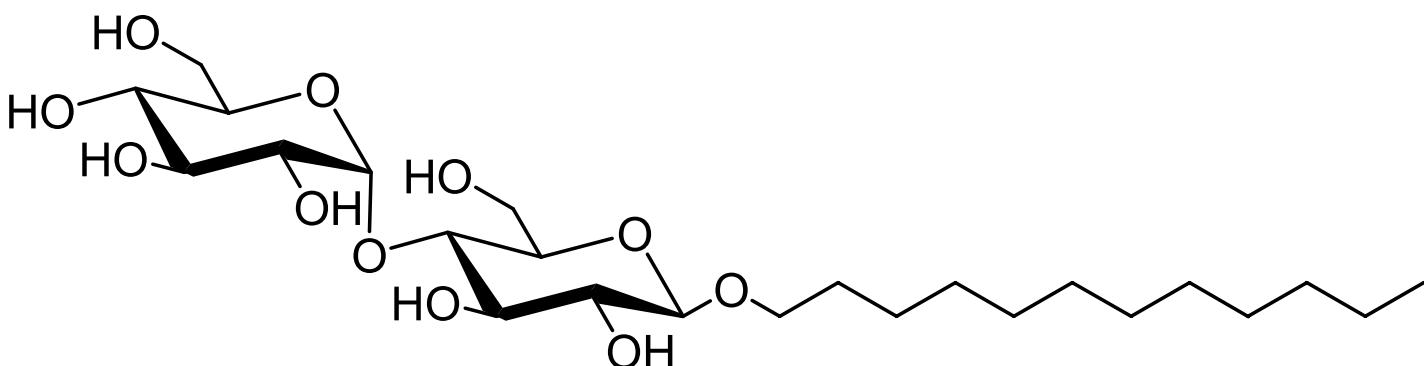


# TECHNICAL DATA SHEET

## n-dodecyl- $\beta$ -D-maltopyranoside

Catalog Number	850520	Physical state	Powder
Purity	> 99%; contains <2% $\alpha$ isomer	Transition temp.	No data
CAS	69227-93-6	CMC	~0.2 mM
Synonyms	n-dodecyl- $\beta$ -D-maltoside; DDM; lauryl maltoside; dodecyl 4-O- $\alpha$ -D-glucopyranosyl- $\beta$ -D-glucopyranoside	pK <sub>a</sub>	No data
Molec. Formula	C <sub>24</sub> H <sub>46</sub> O <sub>11</sub>	TLC mobile phase	C:M*, 80:20, v/v
MW	510.615	Exact Mass	510.304
Percent composition	C 56.45% H 9.08% O 34.47%		
Stability	Store in <-20°C freezer for up to one year		
Solubility	Soluble in ethanol, methanol, water and chloroform		
Web link	<a href="#">850520</a>		

\*C, chloroform; M, methanol



### Description:

Non-ionizing detergents, such as n-dodecyl- $\beta$ -D-maltopyranoside, help solubilize membrane proteins (Reisinger and Eickacker, 2008). To determine the structure of a membrane protein, it must keep its native structure when extracted. DDM can preserve the activity of such diverse membrane proteins as Bacillus multidrug resistance ATP protein (Matar-Merheb et al, 2011), bacteriorhodopsin (Sasaki et al, 2011) and a G-protein coupled receptor (O'Malley et al, 2011) among others. Also, DDM is used as a reference for surfactants (Hovers et al, 2011).

### References:

- Reisinger V, Eichacker LA (2008) Isolation of membrane protein complexes by blue native electrophoresis. Methods Mol Biol. 424: 423-31
- Matar-Merheb et al (2011) Structuring detergents for extracting and stabilizing functional membrane proteins. PLoS One. 6(3): e18036
- Sasaki et al (2011) Sensitive detection of protein-lipid interaction change on bacteriorhodopsin using dodecyl  $\beta$ -D-maltoside. Biochemistry. 50(12): 2283-90
- O'Malley et al (2011) The morphology and composition of cholesterol-rich micellar nanostructures determine transmembrane protein (GPCR) activity. Biophys J. 100(2): L11-3
- Hovers et al (2011) A class of mild surfactants that keep integral membrane proteins water-soluble for functional studies and crystallization. Mol Membr Biol. 28(3): 171-81

### Related products: [Detergents](#)

**MSDS:** Available at [www.avantilipids.com](http://www.avantilipids.com) for Product Number 850520

Avanti Polar Lipids, Inc., 700 Industrial Park Drive, Alabaster, AL 35007-9105, U.S.A.

Phone •(800) 227-0651 •(205) 663-2494 •Fax (800) 229-1004 •(205) 663-0756

Email •[orders@avantilipids.com](mailto:orders@avantilipids.com) •Technical Questions: [technical@avantilipids.com](mailto:technical@avantilipids.com)

•Inquiries: [info@avantilipids.com](mailto:info@avantilipids.com) •Analytical: [analytical@avantilipids.com](mailto:analytical@avantilipids.com)

