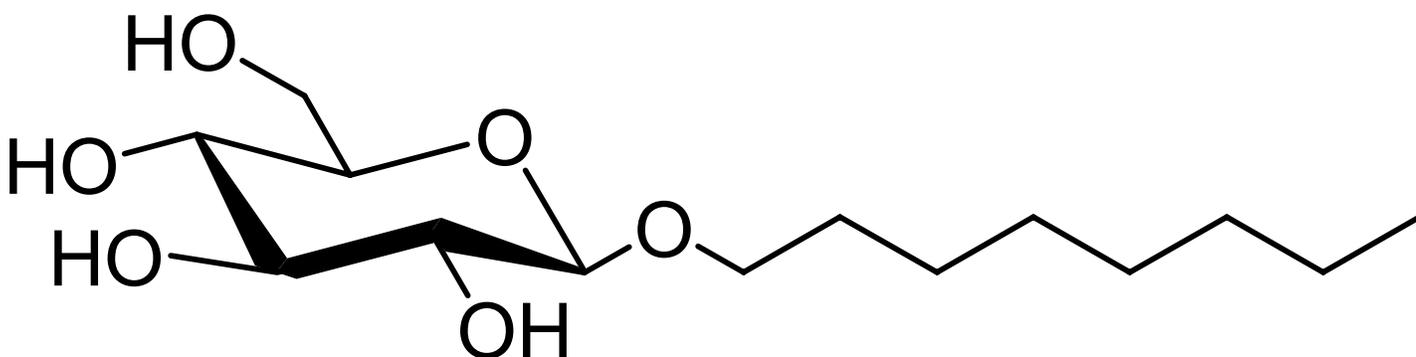


# TECHNICAL DATA SHEET

## n-octyl- $\beta$ -D-glucoside

Catalog Number	850511	Physical state	Powder
Purity	> 99%; contains <2% $\alpha$ isomer	Transition temp.	No data
CAS	29836-26-8	CMC	~20 mM
Synonyms	octyl glucoside; OG; OGP	pK <sub>a</sub>	No data
Molec. Formula	C <sub>14</sub> H <sub>28</sub> O <sub>6</sub>	TLC mobile phase	C:M*, 80:20, v/v
MW	292.369	Exact Mass	292.189
Percent composition	C 57.51% H 9.65% O 32.83%		
Stability	Store in <-20°C freezer for up to one year		
Solubility	Soluble in ethanol, methanol, water and chloroform		
Web link	<a href="#">850511</a>		

\*C, chloroform; M, methanol



### Description:

Non-ionizing detergents help extract, solubilize and denature membrane proteins (Stroud, 2011; Antharavally et al, 2011). To determine the structure of a membrane protein, it must keep its native structure when extracted. Many researchers use combinations of non-ionizing detergents to isolate membrane proteins in vitro, but these detergents must be removed to analyze the protein (Antharavally et al, 2011). Octyl glucoside has versatile uses including: reconstituting an ion channel (Raja and Vales, 2011), solubilizing an enzyme (Wu et al, 2011) and incorporating a protein into liposomes (Madani, Peralvarez-Marín and Graslund, 2011). Also, non-ionizing detergent complexes are used in protein crystallization studies (Sennoga et al, 2003).

### References:

- Stroud RM (2011) New tools in membrane protein determination. *F1000 Biol Rep*.3:8
- Antharavally BS et al (2011) Efficient removal of detergents from proteins and peptides in a spin column format. *Anal Biochem*. 421(1): 39-44
- Raja M, Vales E. Improved technique for reconstituting incredibly high and soluble amounts of tetrameric K<sup>+</sup> channel in natural membranes. *J Membr Biol*. 241(3): 141-4
- Wu LC et al (2011). Purification, identification and cloning of lysoplasmalogenase, the enzyme that catalyzes hydrolysis of the vinyl ether bond of lysoplasmalogen. *J Biol Chem*. 286(28):24916-30
- Madani F, Peralvarez-Marín A, Graslund A. Liposome model systems to study the endosomal escape of cell-penetrating peptides: transport across phospholipid membranes induced by a proton gradient. *J Drug Deliv*. 2011:897592
- Sennoga C et al (2003) Membrane-crystallization in cubo: temperature-dependent phase behavior of monoolein-detergent mixtures. *Acta Crystallogr D Biol Crystallogr*. 59(Pt 2): 239-46

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

**MSDS:** Available at [www.avantilipids.com](http://www.avantilipids.com) for Product Number 850511