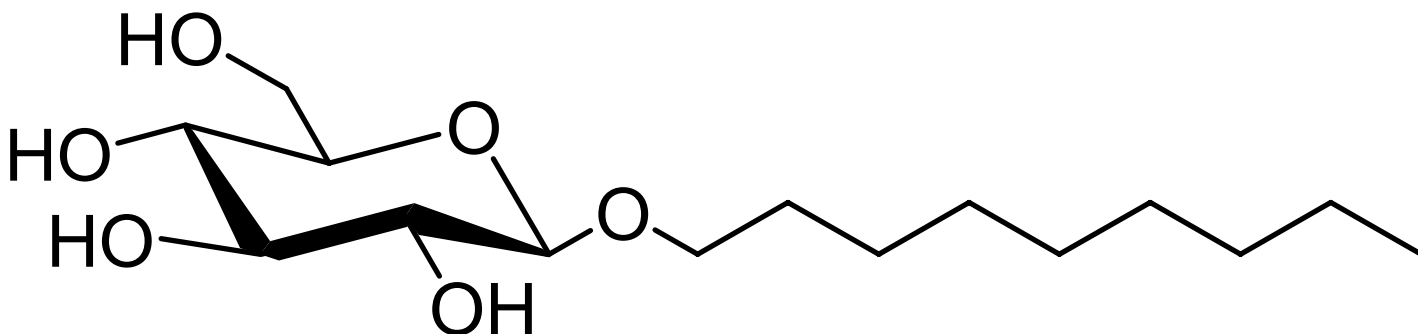


TECHNICAL DATA SHEET

n-nonyl- β -D-glucoside

Catalog Number	850510	Physical state	Powder
Purity	> 99%; contains <2% α isomer	Transition temp.	No data
CAS	69984-73-2	CMC	6.5 mM
Synonyms	n-nonyl- β -D-glucoside; nonyl glucoside; NG; NGP; n-nonyl- β -D-glucopyranoside	pK_a	No data
Molec. Formula	C ₁₅ H ₃₀ O ₆	TLC mobile phase	C:M*, 80:20, v/v
MW	306.395	Exact Mass	306.204
Percent composition	C 58.80% H 9.87% O 31.33%		
Stability	Store in <-20°C freezer for up to one year		
Solubility	Soluble in ethanol, methanol, water and chloroform		
Web link	850510		

*C, chloroform; M, methanol



Description:

Non-ionizing detergents, such as n-nonyl- β -D-glucoside, help solubilize membrane proteins (Reisinger and Eichacker, 2008). 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. Different proteins show distinct accessibility patterns to lipids like nonyl glucoside (Opekarova et al, 2005), suggesting these detergents may play a role in stabilizing membrane proteins also. Finally, non-ionizing detergent complexes are used in protein crystallization studies (Sennoga et al, 2003)

References:

- Reisinger V, Eichacker LA (2008) Isolation of membrane protein complexes by blue native electrophoresis. *Methods Mol Biol.* 424: 423-31
- Opekarova M, Malinska K, Novakova L, Tanner W (2005) Differential effect of phosphatidylethanolamine depletion on raft proteins: further evidence for diversity of rafts in *Saccharomyces cerevisiae*. *Biochim Biophys Acta.* 1711(1): 87-95
- 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 for Product Number 850510