

# n-Dodecyl- $\beta$ -D-maltoside (high purity)

Non-ionic detergent free of  $\alpha$ -isomer (HPLC)

Non-ionic detergent similar to n-octyl- $\beta$ -D-glucopyranoside (Prod. No. ALX-500-001). However, increased alkyl chain length has been shown to increase thermal stability of lipid-free rhodopsin in solution. Applicable to membrane-protein solubilization studies.

Citations: 8

[View Online »](#)

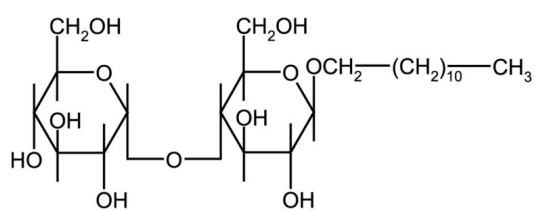
## Ordering Information

[Order Online »](#)

ALX-500-007-G001	1g
ALX-500-007-G005	5g
ALX-500-007-G025	25g

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Handling** Protect from light and moisture. Hygroscopic.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** n-Dodecyl-β-D-maltopyranoside

**Appearance** White crystal powder.

**CAS** 69227-93-6

**Formula**  $C_{24}H_{46}O_{11}$

**MW** 510.6

**Purity** ≥99% (HPLC)(free of α-isomer)

**Solubility** 100mg/ml soluble in water and 100% ethanol.

**Last modified:** May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)