

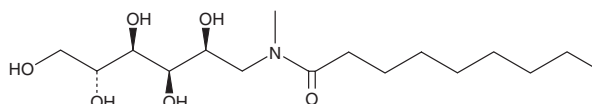
PRODUCT INFORMATION



MEGA-9

Item No. 25702

CAS Registry No.: 85261-19-4
Formal Name: 1-deoxy-1-[methyl(1-oxononyl)amino]-D-glucitol
Synonym: Nonanoyl-N-Methylglucamide
MF: C₁₆H₃₃NO₆
FW: 335.4
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MEGA-9 is supplied as a crystalline solid. A stock solution may be made by dissolving the MEGA-9 in the solvent of choice, which should be purged with an inert gas. MEGA-9 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MEGA-9 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

MEGA-9 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MEGA-9 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MEGA-9 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MEGA-9 is a nonionic detergent that can be used to solubilize membrane proteins.¹ It has a critical micelle concentration (CMC) of 20 mM under no-salt conditions and CMCs ranging from 2 to 17.1 mM under high and low salt conditions for a variety of salts.² It has been used to solubilize the melibiose transport carrier from *E. coli* membranes and reconstitute it into liposomes.³

References

1. Hildreth, J.E.K. N-D-Gluco-N-methylalkanamide compounds, a new class of non-ionic detergents for membrane biochemistry. *Biochem. J.* **207**(2), 363-366 (1982).
2. Miyagishi, S., Okada, K., and Asakawa, T. Salt effect on critical micelle concentrations of nonionic surfactants, N-acyl-N-methylglucamides (MEGA-n). *J. Colloid. Interface Sci.* **238**(1), 91-95 (2001).
3. Hanatani, M., Nishifuji, K., Futai, M., *et al.* Solubilization and reconstitution of membrane proteins of *Escherichia coli* using alkanoyl-N-methylglucamides. *J. Biochem.* **95**(5), 1349-1353 (1984).

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, 10/28/2022

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