

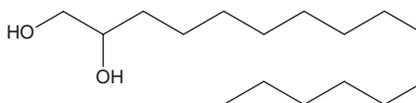
PRODUCT INFORMATION



1,2-Hexadecanediol

Item No. 17445

CAS Registry No.: 6920-24-7
Formal Name: hexadecane-1,2-diol
Synonyms: Cetyl glycol, NSC 71525
MF: C₁₆H₃₄O₂
FW: 258.4
Purity: ≥98%
Supplied as: A 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

1,2-Hexadecanediol is supplied as a solid. A stock solution may be made by dissolving the 1,2-hexadecanediol in the solvent of choice, which should be purged with an inert gas. 1,2-Hexadecanediol is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

1,2-Hexadecanediol is a hydroxyl-substituted hexadecane.^{1,2} It has been used with iron (III) acetylacetonate in the synthesis of monodisperse iron-, cobalt-, or manganese-containing lipid nanoparticles (LNPs) or cobalt-iron nanocrystals.^{1,2} Formulations containing 1,2-hexadecanediol have been used as hair and skin conditioners.

References

1. Sun, S., Zeng, H., Robinson, D.B., *et al.* Monodisperse MFe₂O₄ (M = Fe, Co, Mn) nanoparticles. *J. Am. Chem. Soc.* **126(1)**, 273-279 (2004).
2. Shemer, G., Tirosh, E., Livneh, T., *et al.* Tuning a colloidal synthesis to control Co²⁺ doping in ferrite nanocrystals. *J. Phys. Chem.* **111(39)**, 14334-14338 (2007).

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, 05/09/2024

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