

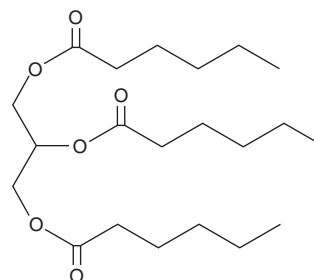
PRODUCT INFORMATION



1,2,3-Trihexanoyl Glycerol

Item No. 27281

CAS Registry No.: 621-70-5
Formal Name: hexanoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glycerol trihexanoate, NSC 406885, Tricaproin, Trihexanoin
MF: C₂₁H₃₈O₆
FW: 386.5
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,2,3-Trihexanoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-trihexanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Trihexanoyl glycerol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1,2,3-trihexanoyl glycerol in these solvents is approximately 30 mg/ml.

Description

1,2,3-Trihexanoyl glycerol is a triacylglycerol that contains hexanoic acid at the *sn*-1, *sn*-2, and *sn*-3 positions. Dietary administration of 1,2,3-trihexanoyl glycerol reduces serum cholesterol levels in rats compared to corn oil-fed control animals.¹ It also decreases plasma cholesterol levels and glutamyl transpeptidase activity, increases plasma LDL, HDL, and protein levels, and reduces diarrhea rate in weaned piglets when administered as 0.5% of the diet.²

References

1. Clifford, A.J., Smith, L.M., Creveling, R.K., *et al.* Effects of dietary triglycerides on serum and liver lipids and sterol excretion of rats. *J. Nutr.* **116(6)**, 944-956 (1986).
2. Wu, T., Li, K., Yi, D., *et al.* Dietary supplementation with trihexanoin enhances intestinal function of weaned piglets. *Int. J. Mol. Sci.* **19(10)**, E3277 (2018).

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, 01/21/2026

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