

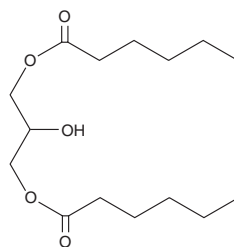
PRODUCT INFORMATION



Dihexanoin

Item No. 27293

CAS Registry No.: 33774-66-2
Formal Name: hexanoic acid, diester with
1,2,3-propanetriol
Synonym: Dicaproin
MF: $C_{15}H_{28}O_5$
FW: 288.4
Purity: $\geq 98\%$
Supplied as: A liquid
Storage: -20°C
Stability: ≥ 4 years



Unspecified isomer

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

Laboratory Procedures

Dihexanoin is supplied as a liquid. A stock solution may be made by dissolving the dihexanoin in the solvent of choice, which should be purged with an inert gas. Dihexanoin is soluble in organic solvents such as ethyl acetate and chloroform.

Description

Dihexanoin is a diacylglycerol that contains the saturated short-chain fatty acid hexanoic acid at two positions.

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