

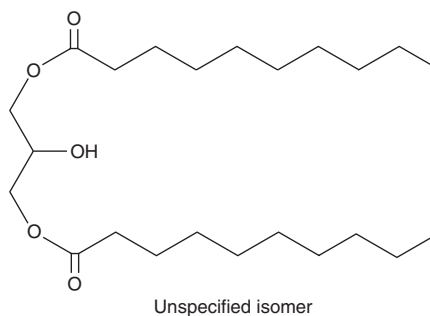
PRODUCT INFORMATION



Didecanoin

Item No. 26992

CAS Registry No.: 53988-07-1
Formal Name: decanoic acid, diester with 1,2,3-propanetriol
Synonym: Dicaprin
MF: C₂₃H₄₄O₅
FW: 400.6
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Didecanoin is supplied as a liquid. A stock solution may be made by dissolving the didecanoin in the solvent of choice, which should be purged with an inert gas. Didecanoin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of didecanoin in these solvents is approximately 30, 5, and 20 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of didecanoin can be prepared by directly dissolving the liquid in aqueous buffers. The solubility of didecanoin in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Didecanoin is a diacylglycerol that contains the saturated medium-chain fatty acid decanoic acid (Item No. 20838) 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, 01/16/2020

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