

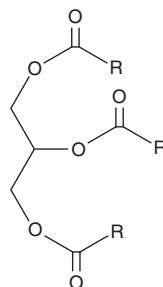
PRODUCT INFORMATION



Caprylic/Capric Triglyceride

Item No. 45757

CAS Registry No.: 65381-09-1
Formal Name: decanoic acid, ester with
1,2,3-propanetriol octanoate
Synonyms: Octanoic/Decanoic Triacylglycerol,
Octanoyl/Decanoyl Triacylglycerol
MF: $C_{29}H_{54}O_6$ (for TG(10:0/8:0/8:0))
FW: 498.7
Purity: $\geq 98\%$
Supplied as: A neat oil
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Caprylic/capric triglyceride is supplied as a neat oil. A stock solution may be made by dissolving the caprylic/capric triglyceride in the solvent of choice, which should be purged with an inert gas. Caprylic/capric triglyceride is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Caprylic/capric triglyceride is a medium-chain triacylglycerol (MCT) that primarily contains caprylic acid (octanoic acid; Item No. 33674) and capric acid (decanoic acid; Item No. 20838) fatty acyl chains. MCTs containing caprylic and capric fatty acyl chains are components of coconut oil, palm kernel oil, and milk fat.¹ Caprylic and capric acid-containing MCTs have been used in the preparation of lipid nanoparticles (LNPs) and drug delivery systems for use *in vitro*.^{2,3} This product contains triacylglycerol molecular species with C8:0 and C10:0 acyl chains, including a dominant species containing capric acid in the *sn*-1 position and caprylic acid in the *sn*-2 and *sn*-3 positions. Formulations containing caprylic/capric triglyceride have been used in creams, lotions, and cosmetics.

References

1. Jadhav, H.B. and Annapure, U.S. Triglycerides of medium-chain fatty acids: A concise review. *J. Food Sci. Technol.* **60**(8), 2143-2152 (2022).
2. Azhar Shekoufeh Bahari, L. and Hamishehkar, H. The impact of variables on particle size of solid lipid nanoparticles and nanostructured lipid carriers; a comparative literature review. *Adv. Pharm. Bull.* **6**(2), 143-151 (2016).
3. Zulfakar, M.H., Pubadi, H., Ibrahim, S.I., et al. Medium-chain triacylglycerols (MCTs) and their fractions in drug delivery systems: A systematic review. *J. Oleo. Sci.* **73**(3), 293-310 (2024).

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, 06/29/2026

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