

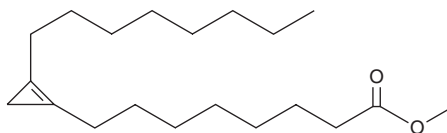
PRODUCT INFORMATION



Sterculic Acid methyl ester

Item No. 26736

CAS Registry No.: 3220-60-8
Formal Name: 2-octyl-1-cyclopropene-1-octanoic acid, methyl ester
Synonyms: Methyl Sterculate, SFE 20:2
MF: C₂₀H₃₆O₂
FW: 308.5
Purity: ≥95%
Supplied as: A solution in hexane
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Sterculic Acid methyl ester is supplied as a solution in hexane. To change the solvent, simply evaporate the hexane under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol, ethyl ether, and chloroform purged with an inert gas can be used.

Description

Sterculic acid methyl ester is an ester form of sterculic acid (Item No. 26735), which is an inhibitor of Δ^9 desaturase.¹ Sterculic acid methyl ester (0.75 mM) inhibits the growth of, and is toxic to, the bacteria *R. opacus*.² It decreases the fatty acid content, increases the ratio of palmitate to other fatty acids, and decreases the levels of stearate and oleate in *R. opacus* when used at concentrations of 0.25 or 0.5 mM. Sterculic acid methyl ester (50 ppm) has a synergistic effect on increased tumor growth induced by aflatoxin Q₁ in rainbow trout.

References

1. Bichi, E., Toral, P.G., Hervás, G., *et al.* Inhibition of Δ^9 -desaturase activity with sterculic acid: Effect on the endogenous synthesis of *cis*-9 18:1 and *cis*-9, *trans*-11 18:2 in dairy sheep. *J. Dairy Sci.* **95**(9), 5242-5252 (2012).
2. Wältermann, M. and Steinbüchel, A. In vitro effects of sterculic acid on lipid biosynthesis in *Rhodococcus opacus* strain PD630 and isolation of mutants defective in fatty acid desaturation. *FEMS Microbiol. Lett.* **190**(1), 45-50 (2000).

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, 03/04/2024

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