

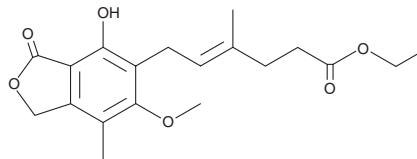
PRODUCT INFORMATION



Ethyl Mycophenolate

Item No. 35065

CAS Registry No.: 32483-51-5
Formal Name: (4E)-6-(1,3-dihydro-4-hydroxy-6-methoxy-7-methyl-3-oxo-5-isobenzofuranyl)-4-methyl-4-hexenoic acid, ethyl ester
MF: C₁₉H₂₄O₆
FW: 348.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Ethyl mycophenolate is supplied as a solid. A stock solution may be made by dissolving the ethyl mycophenolate in the solvent of choice, which should be purged with an inert gas. Ethyl mycophenolate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ethyl mycophenolate in DMSO and DMF is approximately 16 and 20 mg/ml, respectively. Ethyl mycophenolate is slightly soluble in ethanol.

Description

Ethyl mycophenolate is a potential impurity found in commercial preparations of the mycophenolic acid prodrug mycophenolate mofetil.¹

Reference

1. Lu, Y.T., Qi, W.Z., Chen, Y.F., *et al.* Separation and identification of process-related substances and degradation products in mycophenolate mofetil by liquid chromatography coupled with quadrupole-time of flight mass spectrometry. *J. Sep. Sci.* **42(16)**, 2650-2659 (2019).

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, 10/10/2022

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