

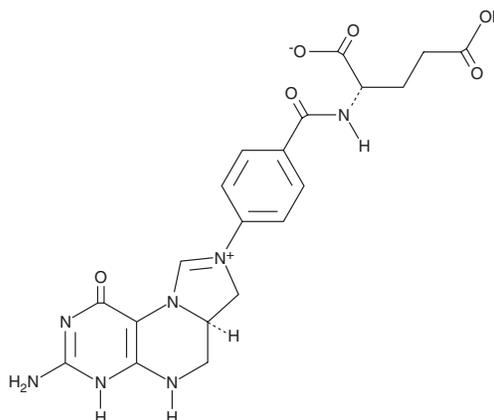
PRODUCT INFORMATION



(6R,S)-5,10-Methenyl-5,6,7,8-tetrahydrofolic Acid

Item No. 31333

CAS Registry No.: 7444-29-3
Formal Name: (6aR)-3-amino-8-[4-[[[(1S)-1,3-dicarboxypropyl]amino]carbonyl]phenyl]-1,2,5,6,6a,7-hexahydro-1-oxo-imidazo[1,5-f]pteridinium, inner salt
Synonyms: Anhydroleucovorin, 6R-N⁵-N¹⁰-methylene-5,6,7,8-tetrahydrofolate
MF: C₂₀H₂₁N₇O₆
FW: 455.4
Purity: ≥95%
UV/Vis.: λ_{max}: 370 nm
Supplied as: A crystalline 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

(6R,S)-5,10-Methenyl-5,6,7,8-tetrahydrofolic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the (6R,S)-5,10-methenyl-5,6,7,8-tetrahydrofolic acid in the solvent of choice, which should be purged with an inert gas. (6R,S)-5,10-Methenyl-5,6,7,8-tetrahydrofolic acid is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

(6R,S)-5,10-Methenyl-5,6,7,8-tetrahydrofolic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (6R,S)-5,10-methenyl-5,6,7,8-tetrahydrofolic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (6R,S)-5,10-Methenyl-5,6,7,8-tetrahydrofolic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(6R,S)-5,10-Methenyl-5,6,7,8-tetrahydrofolic acid is an intermediate in the biosynthesis of thymidylate.¹

Reference

1. Karunaratne, K., Luedtke, N., Quinn, D.M., *et al.* Flavin-dependent thymidylate synthase: N5 of flavin as a Methylene carrier. *Arch. Biochem. Biophys.* **632**, 11-19 (2017).

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, 12/13/2022

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