

PRODUCT INFORMATION



(6R,S)-5,6,7,8-Tetrahydrofolic Acid (hydrochloride)

Item No. 20526

CAS Registry No.: 150731-85-4
Formal Name: N-[4-[[[(2-amino-1,4,5,6,7,8-hexahydro-4-oxo-6-pteridiny]methyl)amino]benzoyl]-L-glutamic acid, trihydrochloride

Synonyms: Tetrahydrofolate, THFA

MF: C₁₉H₂₃N₇O₆ • 3HCl

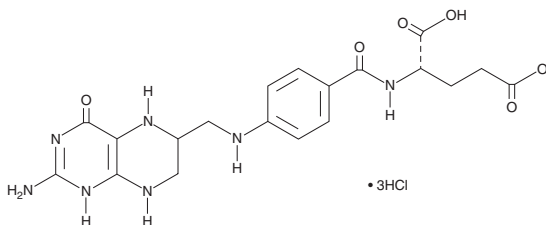
FW: 554.8

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

(6R,S)-5,6,7,8-Tetrahydrofolic acid (THFA) (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the THFA (hydrochloride) in the solvent of choice. THFA (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of THFA (hydrochloride) in these solvents is approximately 20 and 10 mg/ml, respectively.

THFA (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, THFA (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. THFA (hydrochloride) 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

THFA, the reduced form of folic acid, serves as a cofactor in methyltransferase reactions and is the major one-carbon carrier in one carbon metabolism.¹ Its metabolites participate in the synthesis of thymidine for incorporation into DNA or the synthesis of purines, as well as in the formation of methionine, which can be converted to the ubiquitous methyl donor, S-adenosylmethionine (Item No. 13956).¹ THFA has also been used to study the activation of metabolite-binding riboswitches.²

References

1. Ragsdale, S.W. Catalysis of methyl group transfers involving tetrahydrofolate and B12. *Vitam. Horm.* **79**, 293-324 (2008).
2. Breaker, R.R. Prospects for riboswitch discovery and analysis. *Mol. Cell* **43(6)**, 867-879 (2011).

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, 11/14/2016

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