

PRODUCT INFORMATION



L-Folinic Acid (calcium salt hydrate)

Item No. 23833

Formal Name: N-[4-[[[(6S)-2-amino-5-formyl-3,4,5,6,7,8-hexahydro-4-oxo-6-pteridiny]]methyl]amino]benzoyl]-L-glutamic acid, monocalcium salt, hydrate

Synonyms: L-Folinate, 5-Formyltetrahydrofolate

MF: C₂₀H₂₁N₇O₇ • Ca [XH₂O]

FW: 511.5

Purity: ≥98%

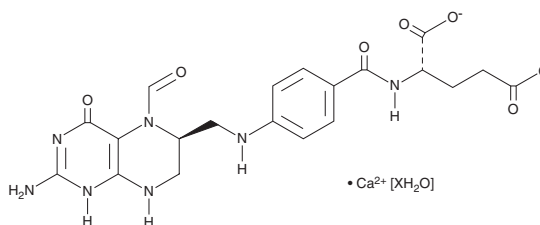
UV/Vis.: λ_{max}: 221, 287 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Synthetic



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

Laboratory Procedures

L-Folinic acid (calcium salt hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the L-folinic acid (calcium salt hydrate) in water. The solubility of L-folinic acid (calcium salt hydrate) in water is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Folinic acid is a reduced folic acid and isomer of DL-folinic acid (Item No. 20383).¹ Formulations containing DL-folinic acid, a racemic mixture of D-folinic acid and L-folinic acid, that have been used to prevent toxicity following therapy with high-dose methotrexate, to enhance the antitumor activity of 5-fluorouracil, and as a folate supplement.^{2,3} Clinical studies demonstrate that formulations containing L-folinic acid have similar efficacy and tolerability as those containing D,L-folinic acid and have been used interchangeably for modification of antitumor activity of 5-fluorouracil, while D-folinic acid is inactive.¹

References

1. Koor, P.A., Karim, S.M., and Marshall, J.L. Is levoleucovorin an alternative to racemic leucovorin? A literature review. *Clin. Colorectal Cancer* **8**(4), 200-206 (2009).
2. Danenberg, P.V., Gustavsson, B., Johnston, P., et al. Folates as adjuvants to anticancer agents: Chemical rationale and mechanism of action. *Crit. Rev. Oncol. Hematol.* **106**, 118-131 (2016).
3. Frye, R.E., Rossignol, D., Casanova, M.F., et al. A review of traditional and novel treatments for seizures in autism spectrum disorder: Findings from a systematic review and expert panel. *Front. Public Health* **1**, 31 (2013).

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, 10/27/2023

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