

PRODUCT INFORMATION



DL-Folinic Acid (calcium salt)

Item No. 20383

CAS Registry No.: 1492-18-8
Formal Name: N-[4-[[[(2-amino-5-formyl-1,4,5,6,7,8-hexahydro-4-oxo-6-pteridiny]methyl]amino]benzoyl]-L-glutamic acid, monocalcium salt

Synonyms: DL-Folate, NSC 3590, 5-Formyltetrahydrofolate

MF: C₂₀H₂₁N₇O₇ • Ca

FW: 511.5

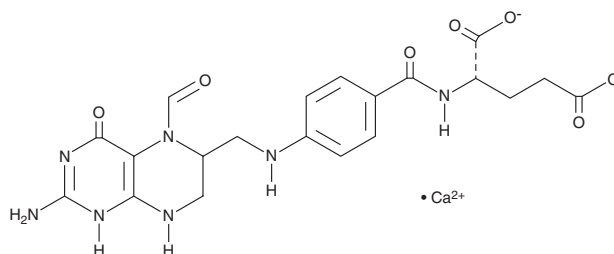
Purity: ≥98%

UV/Vis.: λ_{max}: 220, 287 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

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

Description

DL-Folinic acid is a formylated form of tetrahydrofolic acid (Item No. 18263), the reduced form of folic acid, a B vitamin. DL-Folinic acid is used to prevent toxicity following therapy with high-dose methotrexate (Item No. 13960), to enhance the antitumor activity of 5-fluorouracil (Item No. 14416), and as a folate supplement.^{1,2}

References

1. 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).
2. 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, 05/03/2022

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