

PRODUCT INFORMATION

N-(4-Aminobenzoyl)-L-glutamic Acid (hydrate)

Item No. 33919

Synonyms: pABG, *p*-Aminobenzoyl-glutamic Acid, PABA-Glu

MF: C₁₂H₁₄N₂O₅ • XH₂O

FW: 266.3

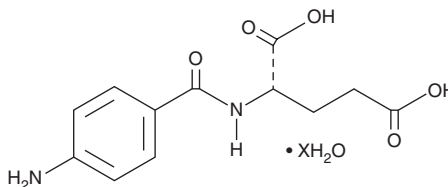
Purity: ≥98%

UV/Vis.: λ_{max}: 287 nm

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

N-(4-Aminobenzoyl)-L-glutamic acid (hydrate) is supplied as a solid. A stock solution may be made by dissolving the N-(4-aminobenzoyl)-L-glutamic acid (hydrate) in the solvent of choice, which should be purged with an inert gas. N-(4-Aminobenzoyl)-L-glutamic acid (hydrate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of N-(4-aminobenzoyl)-L-glutamic acid (hydrate) in these solvents is approximately 1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of N-(4-aminobenzoyl)-L-glutamic acid (hydrate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N-(4-aminobenzoyl)-L-glutamic acid (hydrate) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-(4-Aminobenzoyl)-L-glutamic acid is an oxidation product and a metabolite of tetrahydrofolate.¹ Levels of N-(4-aminobenzoyl)-L-glutamic acid are increased in the plasma of mice fed a high folic acid-containing diet.²

References

1. Reed, L.S. and Archer, M.C. Oxidation of tetrahydrofolic acid by air. *J. Agric. Food Chem.* **28**(4), 801-805 (1980).
2. Burton, M.A., Antoun, E., Penailillo, R.S., *et al.* Folic acid induces intake-related changes in the mammary tissue transcriptome of C57BL/6 mice. *Nutrients* **12**(9), 2821 (2020).

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, 10/27/2022

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