

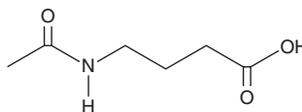
PRODUCT INFORMATION



4-Acetamidobutyric Acid

Item No. 29881

CAS Registry No.: 3025-96-5
Formal Name: 4-(acetylamino)-butanoic acid
Synonyms: 4-Acetamidobutanoic Acid,
N-acetyl-4-Aminobutanoic Acid,
N-acetyl-γ-Aminobutyric Acid,
NSC 27423
MF: C₆H₁₁NO₃
FW: 145.2
Purity: ≥95%
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

4-Acetamidobutyric acid is supplied as a solid. A stock solution may be made by dissolving the 4-acetamidobutyric acid in the solvent of choice, which should be purged with an inert gas. 4-Acetamidobutyric acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-acetamidobutyric acid in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

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 4-acetamidobutyric acid can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 4-acetamidobutyric acid in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-Acetamidobutyric acid is a GABA derivative.¹ Urinary levels of 4-acetamidobutyric acid are elevated in patients with stage 5 chronic kidney disease compared to patients with stage 3 kidney disease and those without kidney disease.

Reference

1. Boelaert, J., t'Kindt, R., Schepers, E., *et al.* State-of-the-art non-targeted metabolomics in the study of chronic kidney disease. *Metabolomics* **10(3)**, 425-442 (2014).

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, 12/16/2022

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