

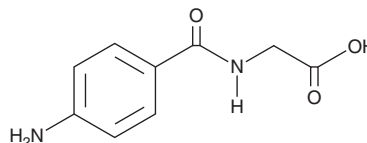
PRODUCT INFORMATION



4-Aminohippuric Acid

Item No. 23726

CAS Registry No.: 61-78-9
Formal Name: N-(4-aminobenzoyl)-glycine
Synonyms: *p*-Aminohippuric Acid, *para*-Aminohippuric Acid, NSC 7550, NSC 13064
MF: C₉H₁₀N₂O₃
FW: 194.2
Purity: ≥98%
UV/Vis.: λ_{max}: 283 nm
Supplied as: A solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

4-Aminohippuric acid is supplied as a solid. A stock solution may be made by dissolving the 4-aminohippuric acid in the solvent of choice. 4-Aminohippuric acid is slightly soluble in chloroform, methanol, and DMSO.

Description

4-Aminohippuric acid is a substrate for various renal transporters that has been used in the study of anion uptake in the kidney. It is a substrate for organic anion transporter 1 (OAT1; K_m = 14.3 μM in *X. laevis* oocytes), the human inorganic phosphate transporter (NPT1; K_m = 2.66 mM in HEK293 cells), and apical multidrug resistance protein (MDR2; K_m = 880 μM in HEK293 cells).¹⁻³ 4-Aminohippuric acid is a glycine amide form of 4-aminobenzoic acid (Item No. 18659). Formulations containing 4-aminohippuric acid have been used as markers to determine renal plasma flow.

References

1. Sekine, T., Watanabe, N., Hosoyamada, M., *et al.* Expression cloning and characterization of a novel multispecific organic anion transporter. *J. Biol. Chem.* **272**(30), 18526-18529 (1997).
2. Uchino, H., Tamai, I., Yamashita, K., *et al.* *p*-aminohippuric acid transport at renal apical membrane mediated by human inorganic phosphate transporter NPT1. *Biochem. Biophys. Res. Commun.* **270**(1), 254-259 (2000).
3. Leier, I., Hummel-Eisenbeiss, J., Cui, Y., *et al.* ATP-dependent *para*-aminohippurate transport by apical multidrug resistance protein MRP2. *Kidney Int.* **57**(4), 1636-1642 (2000).

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, 11/28/2022

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