

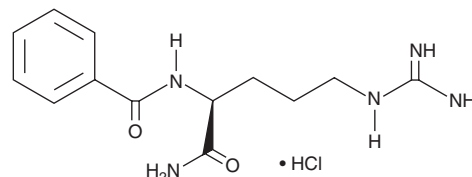
PRODUCT INFORMATION



N- α -Benzoyl-L-Argininamide (hydrochloride)

Item No. 35015

CAS Registry No.: 4299-03-0
Formal Name: (S)-N-[1-(aminocarbonyl)-4-[(aminoiminomethyl)amino]butyl]-benzamide, monohydrochloride
Synonyms: BAA, Bz-Arg-NH₂, Benzoylarginamide, NSC 343719
MF: C₁₃H₁₉N₅O₂ • HCl
FW: 313.8
Purity: ≥98%
UV/Vis.: λ_{max} : 227 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- α -Benzoyl-L-argininamide (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the N- α -benzoyl-L-argininamide (hydrochloride) in the solvent of choice, which should be purged with an inert gas. N- α -Benzoyl-L-argininamide (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N- α -benzoyl-L-argininamide (hydrochloride) in these solvents is approximately 1, 2, and 3 mg/ml, respectively.

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- α -benzoyl-L-argininamide (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N- α -benzoyl-L-argininamide (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N- α -Benzoyl-L-argininamide is a nonpeptide substrate of proteases and deiminases.¹⁻³ It has been used as a substrate for the quantification of papain, carboxypeptidase B, protein arginine deiminase 2 (PAD2), and PAD4 activities.

References

- Whitaker, J.R. and Bender, M.L. Kinetics of papain-catalyzed hydrolysis of α -N-Benzoyl-L-arginine ethyl ester and α -N-Benzoyl-L-argininamide. *J. Am. Chem. Soc.* **87**(12), 2728-2737 (1965).
- Lones, M., Chatterjee, R., Singh, H., et al. Lysosomal carboxypeptidase B from rabbit lung purification and characterization. *Arch. Biochem. Biophys.* **221**(1), 64-78 (1983).
- Nakayama-Hamada, M., Suzuki, A., Kubota, K., et al. Comparison of enzymatic properties between hPADI2 and hPADI4. *Biochem. Biophys. Res. Commun.* **327**(1), 192-200 (2005).

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, 12/09/2022

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