

PRODUCT INFORMATION



Na-Benzoyl-L-Arginine-pNA (hydrochloride)

Item No. 36331

CAS Registry No.: 21653-40-7

Formal Name: N-[(1S)-4-[(aminoiminomethyl)amino]-1-[[[4-(4-nitrophenyl)amino]carbonyl]butyl]-benzamide, monohydrochloride

Synonyms: Na-Benzoyl-L-Arginine-4-Nitroanilide, Na-Benzoyl-L-Arginine-p-Nitroanilide

MF: $C_{19}H_{22}N_6O_4 \cdot HCl$

FW: 434.9

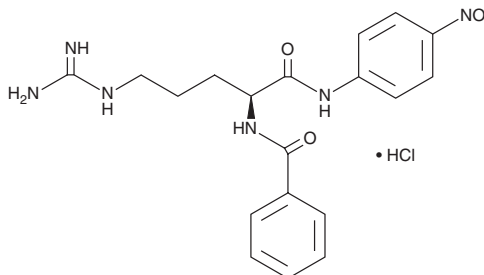
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 228, 314 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Na-Benzoyl-L-arginine-pNA (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the Na-benzoyl-L-arginine-pNA (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Na-Benzoyl-L-arginine-pNA (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Na-benzoyl-L-arginine-pNA (hydrochloride) in these solvents is approximately 2 mg/ml.

Description

Na-Benzoyl-L-arginine-pNA is a colorimetric substrate for trypsin and papain.¹ Enzymatic cleavage of Na-benzoyl-L-arginine-pNA releases p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of trypsin or papain activity.

Reference

1. Nishi, N., Tokura, S., and Noguchi, J. The synthesis of benzoyl-L-arginine-p-nitroanilide. *B. Chem. Soc. Jpn.* **43**(9), 2900-2907 (1970).

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, 11/30/2022

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