

PRODUCT INFORMATION

Cbz-Gly-Pro-Arg-pNA (acetate)

Item No. 42990

Formal Name: N-[(phenylmethoxy)carbonyl]glycyl-L-prolyl-

Synonyms: N-(4-nitrophenyl)-L-argininamide, acetate

Cbz-GPR-p-nitroanilide, Cbz-GPR-pNA, Protease Chromogenic Substrate

MF: $C_{27}H_{34}N_8O_7 \cdot XC_2H_4O_2$

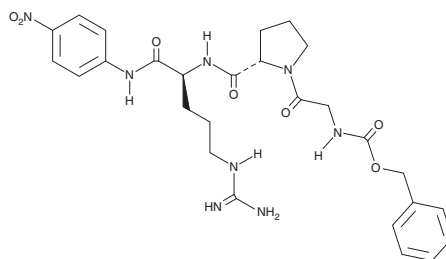
FW: 582.6

Purity: $\geq 98\%$

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

Cbz-Gly-Pro-Arg-pNA is supplied as a solid. A stock solution may be made by dissolving the Cbz-Gly-Pro-Arg-pNA in the solvent of choice, which should be purged with an inert gas. Cbz-Gly-Pro-Arg-pNA is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Cbz-Gly-Pro-Arg-pNA is a colorimetric substrate for proteases, including thrombin and mast cell tryptase.^{1,2} Proteases preferentially bind to and cleave the Gly-Pro-Arg (GPR) peptide sequence to release p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of protease activity.

References

1. Hogg, P.J. and Jackson, C.M. Formation of a ternary complex between thrombin, fibrin monomer, and heparin influences the action of thrombin on its substrates. *J. Biol. Chem.* **265**(1), 248-255 (1990).
2. Harvima, I.T., Harvima, R.J., Eloranta, T.O., *et al.* The allosteric effect of salt on human mast cell tryptase. *Biochim. Biophys. Acta* **956**(2), 133-139 (1988).

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, 04/04/2025

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