

PRODUCT INFORMATION

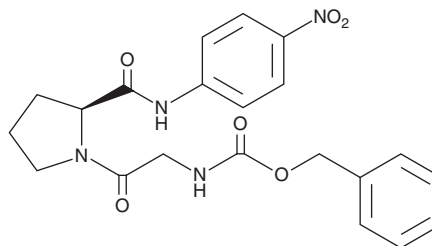


Z-Gly-Pro-pNA

Item No. 36135

CAS Registry No.: 65022-15-3
Formal Name: N-[(phenylmethoxy)carbonyl]glycyl-L-(4-nitrophenyl)-L-prolinamide
Synonyms: Carbobenzoxy-glycyl-L-prolyl-p-Nitroanilide, N-CBZ-Glycyl-L-proline 4-Nitroanilide, Z-Gly-Pro-4-Nitroanilide

MF: C₂₁H₂₂N₄O₆
FW: 426.4
Purity: ≥98%
UV/Vis.: λ_{max}: 314 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

Z-Gly-Pro-pNA is supplied as a solid. A stock solution may be made by dissolving the Z-Gly-Pro-pNA in the solvent of choice, which should be purged with an inert gas. Z-Gly-Pro-pNA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Z-Gly-Pro-pNA in ethanol is approximately 12 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 Z-Gly-Pro-pNA can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Z-Gly-Pro-pNA in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Z-Gly-Pro-pNA is a colorimetric substrate for prolyl endopeptidase.¹ Enzymatic cleavage of Z-Gly-Pro-pNA releases p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nM as a measure of prolyl endopeptidase activity.

Reference

1. Kalwant, S. and Porter, A.G. Purification and characterization of human brain prolyl endopeptidase. *Biochem. J.* **276**, 237-244 (1991).

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, 12/19/2022

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