PRODUCT INFORMATION



Z-D-Arg-Gly-Arg-pNA (hydrochloride)

Item No. 27583

CAS Registry No.: 113711-77-6

Formal Name: N²-[(phenylmethoxy)carbonyl]-D-

arginylglycyl-N-(4-nitrophenyl)-L-

argininamide, dihydrochloride

Synonyms: Benzyl-D-Arg-Gly-Arg-pNA,

> Benzyl-D-Arg-Gly-Arg-p-nitroanilide, Factor Xa Chromogenic Substrate,

S-2765

MF: C₂₈H₃₉N₁₁O₇ • 2HCl

FW: 714.6 **Purity:** ≥98% UV/Vis.: λ_{max} : 315 nm Supplied as: A solid -20°C Storage: Stability: ≥4 years

2HCl

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

Laboratory Procedures

Z-D-Arg-Gly-Arg-pNA (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the Z-D-Arg-Gly-Arg-pNA (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Z-D-Arg-Gly-Arg-pNA (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Z-D-Arg-Gly-Arg-pNA (hydrochloride) in these solvents is approximately 14 mg/ml. Z-D-Arg-Gly-Arg-pNA (hydrochloride) is slightly soluble in ethanol.

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-D-Arg-Gly-Arg-pNA (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Z-D-Arg-Gly-Arg-pNA (hydrochloride) in PBS, pH 7.2, is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Z-D-Arg-Gly-Arg-pNA is a colorimetric substrate for Factor Xa.1,2 Factor Xa preferentially binds to and cleaves the Arg-Gly-Arg (RGR) peptide sequence to release p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of Factor Xa activity.

References

- 1. Gustavsson, S.I. and Arielly, S. New Peptide Derivatives. KabiVitrum AB 4,797,472 (1989).
- 2. Joo, H.-S., Park, G.-C., Kim, K.-M., et al. Novel alkaline protease from the polychaeta, Periserrula leucophryna: Purification and characterization. Proc. Biochem. 36(8-9), 893-900 (2001).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information Buyer agrees to purchase the mater can be found on our website.

Copyright Cayman Chemical Company, 10/04/2022

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