

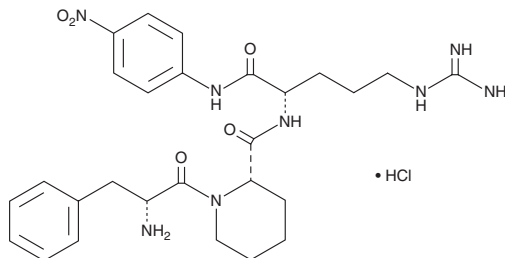
PRODUCT INFORMATION



S-2238 (hydrochloride)

Item No. 34944

CAS Registry No.: 160192-34-7
Formal Name: D-phenylalanyl-(2S)-2-piperidinecarbonyl-N-(4-nitrophenyl)-L-argininamide, monohydrochloride
Synonyms: H-D-Phe-Pip-Arg-pNA,
D-Phe-Piperidinecarbonyl-Arg-p-nitroanilide
MF: C₂₇H₃₆N₈O₅ • HCl
FW: 589.1
Purity: ≥98%
UV/Vis.: λ_{max}: 313 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

S-2238 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the S-2238 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. S-2238 (hydrochloride) is soluble in DMSO.

Description

S-2238 is a chromogenic substrate for thrombin.^{1,2} Upon enzymatic cleavage, *p*-nitroanilide (*p*NA) is released which can be quantified by colorimetric detection at 405 nm as a measure of thrombin activity.

References

1. Abildgaard, U., Lie, M., and Ødegård, O.R. Antithrombin (heparin cofactor) assay with “new” chromogenic substrates (S-2238 and Chromozym TH). *Thromb. Res.* **11(4)**, 549-553 (1977).
2. Ohno, H., Kosaki, G., Kambayashi, J., *et al.* FOY: [ethyl P-(6-guanidinohexanoyloxy) benzoate] methanesulfonate as a serine proteinase inhibitor. I. Inhibition of thrombin and factor Xa in vitro. *Thromb. Res.* **19(4-5)**, 579-588 (1980).

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, 12/02/2022

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