

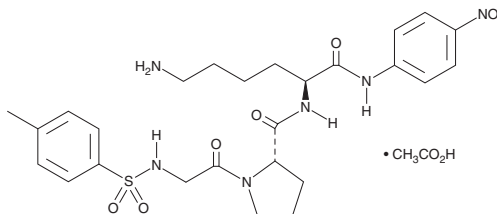
PRODUCT INFORMATION



N-p-Tosyl-Gly-Pro-Lys-pNA (acetate)

Item No. 31544

CAS Registry No.: 88793-79-7
Formal Name: N-[(4-methylphenyl)sulfonyl]glycyl-L-prolyl-N-(4-nitrophenyl)-L-lysine, monoacetate
Synonym: N-p-Tosyl-Gly-Pro-Lys-p-nitroanilide
MF: $C_{26}H_{34}N_6O_7S \cdot C_2H_4O_2$
FW: 634.7
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 226 nm
Supplied as: A crystalline 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

N-p-Tosyl-Gly-Pro-Lys-pNA (acetate) is supplied as a crystalline solid. A stock solution may be made by dissolving the N-p-Tosyl-Gly-Pro-Lys-pNA (acetate) in the solvent of choice, which should be purged with an inert gas. N-p-Tosyl-Gly-Pro-Lys-pNA (acetate) is soluble in the organic solvent ethanol at a concentration of approximately 50 mg/ml.

Description

N-p-Tosyl-Gly-Pro-Lys-pNA is a colorimetric substrate for plasmin.¹⁻³ Plasmin binds and hydrolyzes N-p-tosyl-Gly-Pro-Lys-pNA to release p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of plasmin activity.

References

1. Jespersen, J., Gram, J., and Astrup, T. The autodigestion of human plasmin follows a bimolecular mode of reaction subject to product inhibition. *Thromb. Res.* **41(3)**, 395-404 (1986).
2. Kim, S.H. and Choi, N.S. Electrophoretic analysis of protease inhibitors in fibrin zymography. *Anal. Biochem.* **270(1)**, 179-181 (1999).
3. Chen, T. and Rael, E.D. Purification of M5, a fibrinolytic proteinase from *Crotalus molossus molossus* venom that attacks complement. *Int. J. Biochem. Cell Biol.* **29(5)**, 789-799 (1997).

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, 12/08/2022

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