

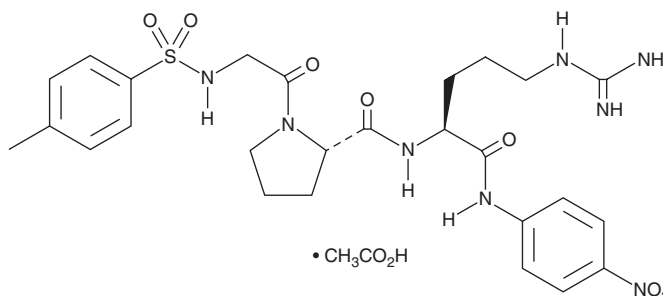
# PRODUCT INFORMATION



## N-(p-Tosyl)-GPR-pNA (acetate)

Item No. 27568

**CAS Registry No.:** 86890-95-1  
**Formal Name:** N-[(4-methylphenyl)sulfonyl]glycyl-L-prolyl-N-(4-nitrophenyl)-L-argininamide, monoacetate  
**Synonyms:** N-(p-Tosyl)-Gly-Pro-Arg-p-Nitroanilide, Tos-Gly-Pro-Arg-pNA  
**MF:** C<sub>26</sub>H<sub>34</sub>N<sub>8</sub>O<sub>7</sub>S • C<sub>2</sub>H<sub>4</sub>O<sub>2</sub>  
**FW:** 662.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 225, 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

N-(p-Tosyl)-GPR-pNA (acetate) is supplied as a solid. A stock solution may be made by dissolving the N-(p-tosyl)-GPR-pNA (acetate) in the solvent of choice, which should be purged with an inert gas. N-(p-Tosyl)-GPR-pNA (acetate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of N-(p-tosyl)-GPR-pNA (acetate) in these solvents is approximately 3 and 20 mg/ml, respectively.

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 N-(p-tosyl)-GPR-pNA (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N-(p-tosyl)-GPR-pNA (acetate) in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

N-(p-Tosyl)-GPR-pNA is a colorimetric substrate for thrombin.<sup>1</sup> Thrombin preferentially binds to and cleaves 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 thrombin activity.

### Reference

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).

#### 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, 10/07/2022

#### 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