

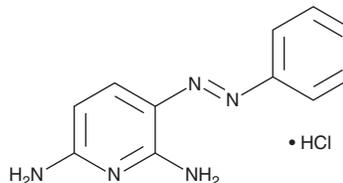
# PRODUCT INFORMATION



## Phenazopyridine (hydrochloride)

Item No. 29683

**CAS Registry No.:** 136-40-3  
**Formal Name:** 3-(2-phenyldiazenyl)-2,6-pyridinediamine, monohydrochloride  
**MF:** C<sub>11</sub>H<sub>11</sub>N<sub>5</sub> • HCl  
**FW:** 249.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 238, 394 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

Phenazopyridine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the phenazopyridine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Phenazopyridine (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of phenazopyridine (hydrochloride) in these solvents is approximately 1 mg/ml.

Phenazopyridine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, phenazopyridine (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Phenazopyridine (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Phenazopyridine is an azo dye with analgesic properties.<sup>1</sup> It decreases the bladder distension-induced firing rate of afferent bladder A $\gamma$ -fibers, but not C-fibers, in anesthetized rats when administered at a doses of 0.3, 1, and 3 mg/kg.<sup>2</sup> Formulations containing phenazopyridine have been used as analgesics in the treatment of urinary tract infections.

### References

1. Iqbal, J., Gupta, A., and Husain, A. Photochemistry of phenazopyridine hydrochloride. *Pharmazie* **61(9)**, 747-750 (2006).
2. Aizawa, N. and Wyndaele, J.J. Effects of phenazopyridine on rat bladder primary afferent activity, and comparison with lidocaine and acetaminophen. *Neurol. Urodyn.* **29(8)**, 1445-1450 (2010).

#### 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, 10/21/2022

#### 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