

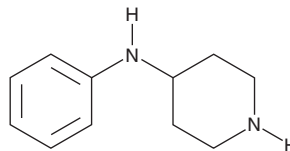
# PRODUCT INFORMATION



## N-Phenylpiperidin-4-amine

Item No. 30898

**CAS Registry No.:** 23056-29-3  
**Formal Name:** N-phenyl-4-piperidinamine  
**Synonyms:** 4-Anilinopiperidine, NSC 80678  
**MF:** C<sub>11</sub>H<sub>16</sub>N<sub>2</sub>  
**FW:** 176.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 248 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-Phenylpiperidin-4-amine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-phenylpiperidin-4-amine in the solvent of choice, which should be purged with an inert gas. N-Phenylpiperidin-4-amine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N-phenylpiperidin-4-amine in these solvents is approximately 30 mg/ml.

### Description

N-Phenylpiperidin-4-amine is a building block.<sup>1</sup> It has been used in the synthesis of *M. tuberculosis* decaprenylphosphoryl-β-D-ribose 2'-epimerase (DprE1) inhibitors.

### Reference

1. Borthwick, J.A., Alemparte, C., Wall, I., et al. *Mycobacterium tuberculosis* decaprenylphosphoryl-β-D-ribose oxidase inhibitors: Expedient reconstruction of suboptimal hits into a series with potent in vivo activity. *J. Med. Chem.* **63**(5), 2557-2576 (2020).

#### 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, 12/01/2022

#### 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](http://WWW.CAYMANCHEM.COM)