

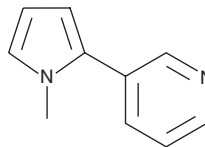
# PRODUCT INFORMATION



## β-Nicotyrine

Item No. 16333

**CAS Registry No.:** 487-19-4  
**Formal Name:** 3-(1-methyl-1H-pyrrol-2-yl)-pyridine  
**Synonyms:** NSC 127943, NSC 407276  
**MF:** C<sub>10</sub>H<sub>10</sub>N<sub>2</sub>  
**FW:** 158.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 288 nm  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

β-Nicotyrine is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of β-nicotyrine in these solvents is approximately 30 and 50 mg/ml, respectively.

β-Nicotyrine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of β-nicotyrine should be diluted with the aqueous buffer of choice. The solubility of β-nicotyrine in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

β-Nicotyrine is an alkaloid that has been found in tobacco.<sup>1,2</sup> It reduces the number of dsDNA breaks induced by the tobacco-specific carcinogen nitrosamine (4-(methylnitrosamino)-1-(3-pyridyl)-1-butanone; Item No. 16414) in a comet assay using HepaRG hepatocellular carcinoma cells when used at concentrations of 5 and 50 μM.<sup>2</sup>

### References

1. Burns, D.T. Rapid determination of certain alkaloids, other than nicotine, in tobacco. *J. Chromatogr.* **133(2)**, 379-381 (1977).
2. Ordonez, P., Sierra, A.B., Camacho, O.M., *et al.* Nicotine, cotinine, and β-nicotyrine inhibit NNK-induced DNA-strand break in the hepatic cell line HepaRG. *Toxicol. In Vitro* **28(7)**, 1329-1337 (2014).

#### 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, 12/21/2023

#### 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