

PRODUCT INFORMATION

NSC 19005

Item No. 32898

CAS Registry No.: 5452-87-9

Formal Name: N-methyl-N-[2-(2-pyridinyl)ethyl]-
2-pyridineethanamine

Synonym: Betahistine Impurity C

MF: C₁₅H₁₉N₃

FW: 241.3

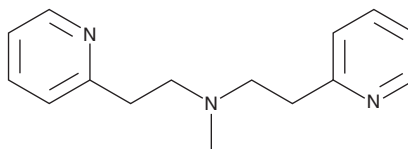
Purity: ≥95%

UV/Vis.: λ_{max}: 261 nm

Supplied as: A neat oil

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

NSC 19005 is supplied as a neat oil. A stock solution may be made by dissolving the NSC 19005 in the solvent of choice, which should be purged with an inert gas. NSC 19005 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NSC 19005 in these solvents is approximately 30 mg/ml.

NSC 19005 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NSC 19005 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. NSC 19005 has a solubility of approximately 0.12 mg/ml in a 1:7 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NSC 19005 is a potential impurity found in commercial preparations of betahistine.¹ It is a degradation product formed by exposure to oxidative, thermal, or photolytic stress.

Reference

1. de Mello, M.B.M., de Oliveira, A.F., de Oliveira, C.L., et al. Characterization and *in silico* mutagenic assessment of a new betahistine degradation impurity. *J. Braz. Chem. Soc.* **30(7)**, 1415-1424 (2019).

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, 12/09/2022

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