

# PRODUCT INFORMATION



## 1-methoxy-5-Methylphenazinium (methyl sulfate)

Item No. 21258

CAS Registry No.: 65162-13-2

Formal Name: 1-methoxy-5-methyl-phenazinium, methyl sulfate

Synonym: 1-Methoxyphenazine methosulfate, 1-methoxy PMS

MF:  $C_{14}H_{13}N_2O \cdot CH_3SO_4$

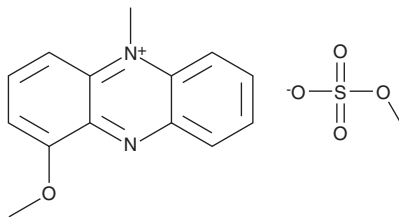
FW: 336.4

Purity:  $\geq 98\%$

Supplied as: A crystalline solid

Storage:  $-20^\circ C$

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



### Laboratory Procedures

1-methoxy-5-Methylphenazinium (methyl sulfate) (1-methoxy PMS) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-methoxy PMS in the solvent of choice. 1-methoxy PMS is soluble in water. The solubility of 1-methoxy PMS in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

1-methoxy PMS is a photochemically stable mediator of electron transport between NADPH and various electron acceptors such as tetrazolium dyes.<sup>1</sup> It can be used to visualize NAD-linked dehydrogenase activities.

### Reference

1. Hisada, R., and Yagi, T. 1-Methoxy-5-methylphenazinium methyl sulfate. A photochemically stable electron mediator between NADH and various electron acceptors. *J. Biochem.* **82(5)**, 1469-1473 (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, 02/14/2017

#### 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