

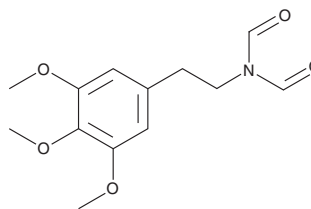
# PRODUCT INFORMATION



## N,N-Diformylmescaline

Item No. 9003607

**Formal Name:** N-formyl-N-(3,4,5-trimethoxyphenethyl) formamide  
**MF:** C<sub>13</sub>H<sub>17</sub>NO<sub>5</sub>  
**FW:** 267.3  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Laboratory Procedures

N,N-Diformylmescaline is supplied as a crystalline solid. A stock solution may be made by dissolving the N,N-diformylmescaline in the solvent of choice, which should be purged with an inert gas. N,N-Diformylmescaline is soluble in organic solvents such as acetonitrile, DMSO, and dimethyl formamide (DMF). The solubility of N,N-diformylmescaline in acetonitrile and DMF is approximately 30 mg/ml and approximately 10 mg/ml in DMSO.

### Description

N,N-Diformylmescaline (Item No. 9003607) is an analytical reference standard categorized as a phenethylamine. It has been found in seized samples suspected of containing novel psychoactive substances (NPS).<sup>1</sup> This product is intended for research and forensic applications.

### Reference

1. Gallagher, R., McLaughlin, M.G., Blakey, K., *et al.* N,N-Diformylmescaline: A novel analogue of mescaline detected in Queensland. *Drug Test. Anal.* **15(2)**, 204-212 (2022).

#### 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, 07/07/2023

#### 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