

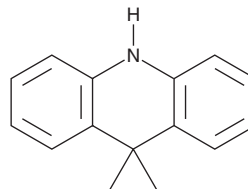
PRODUCT INFORMATION



9,9-Dimethyl-9,10-dihydroacridine

Item No. 45217

CAS Registry No.: 6267-02-3
Formal Name: 9,10-dihydro-9,9-dimethyl-acridine
Synonyms: 9,9-Dimethylcarbazine, DM-AD, DMAC, NSC 36671
MF: C₁₅H₁₅N
FW: 209.3
Purity: ≥98%
Supplied as: A 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

9,9-Dimethyl-9,10-dihydroacridine is supplied as a solid. A stock solution may be made by dissolving the 9,9-Dimethyl-9,10-dihydroacridine in the solvent of choice, which should be purged with an inert gas. 9,9-Dimethyl-9,10-dihydroacridine is soluble (≥10 mg/ml) in ethanol.

Description

9,9-Dimethyl-9,10-dihydroacridine is a synthetic intermediate and an electron donor.¹ It has been used in the synthesis of thermally activated delayed fluorescence (TADF) materials for use in organic light-emitting diodes (OLEDs).

Reference

1. Madushani, B., Mamada, M., Goushi, K., *et al.* Multiple donor-acceptor design for highly luminescent and stable thermally activated delayed fluorescence emitters. *Sci. Rep.* **13**(1), 7644 (2023).

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, 02/27/2026

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