

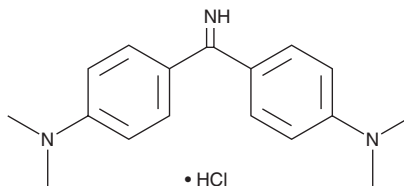
PRODUCT INFORMATION



Auramine O

Item No. 35787

CAS Registry No.: 2465-27-2
Formal Name: 4,4'-carbonimidoylbis[N,N-dimethyl-benzenamine, monohydrochloride
Synonyms: AO, AuO
MF: $C_{17}H_{21}N_3 \cdot HCl$
FW: 303.8
UV/Vis.: λ_{max} : 204, 307, 430 nm
Abs./Em. Max: 432/550 nm
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

Auramine O is supplied as a solid. A stock solution may be made by dissolving the auramine O in the solvent of choice, which should be purged with an inert gas. Auramine O is soluble (≥ 10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Auramine O is a diphenylmethane dye.¹ It displays absorbance/emission maxima of 432/550 nm, respectively. Auramine O is an acid-fast dye that has been used in the staining of various bacteria, including *M. tuberculosis*.² It induces autophagy in, and promotes migration of, A549 lung cancer cells.³ Formulations containing auramine O were previously used as food additives.

References

1. Sabnis, R.W. Handbook of biological dyes and stains: Synthesis and industrial applications. (2010).
2. Ryan, G.J., Shapiro, H.M., and Lenaerts, A.J. Improving acid-fast fluorescent staining for the detection of mycobacteria using a new nucleic acid staining approach. *Tuberculosis (Edinb)* **94**(5), 511-518 (2014).
3. Tung, J.-C., Huang, W.-C., Yang, J.-C., et al. Auramine O, an incense smoke ingredient, promotes lung cancer malignancy. *Environ. Toxicol.* **32**(11), 2379-2391 (2017).

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, 10/29/2025

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