

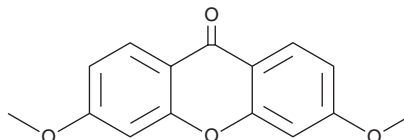
PRODUCT INFORMATION



3,6-Dimethoxy-9H-xanthen-9-one

Item No. 34806

CAS Registry No.: 15007-07-5
Synonym: 3,6-Dimethoxyxanthone
MF: $C_{15}H_{12}O_4$
FW: 256.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 241, 307 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

3,6-Dimethoxy-9H-xanthen-9-one is supplied as a solid. A stock solution may be made by dissolving the 3,6-dimethoxy-9H-xanthen-9-one in the solvent of choice, which should be purged with an inert gas. 3,6-Dimethoxy-9H-xanthen-9-one is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 3,6-dimethoxy-9H-xanthen-9-one these solvents is approximately 2 and 1 mg/ml, respectively.

Description

3,6-Dimethoxy-9H-xanthen-9-one is a synthetic intermediate and a derivative of xanthone (Item No. 34098).¹⁻³ It has been used in the synthesis of fluorescein derivatives, as well as compounds with anticancer or acetylcholinesterase inhibitory activities.

References

1. Šebej, P., Wintner, J., Müller, P., *et al.* Fluorescein analogues as photoremovable protecting groups absorbing at ~ 520 nm. *J. Org. Chem.* **78**(5), 1833-1843 (2013).
2. Liu, C., Zhang, M., Zhang, Z., *et al.* Synthesis and anticancer potential of novel xanthone derivatives with 3,6-substituted chains. *Bioorg. Med. Chem.* **24**(18), 4263-4271 (2016).
3. Alawi, M.S., Awad, T.A., Mohamed, M.A., *et al.* Insights into the molecular basis of acetylcholinesterase inhibition by xanthenes: An integrative *in silico* and *in vitro* approach. *Mol. Simul.* **46**(4), 253-261 (2020).

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, 09/22/2021

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