

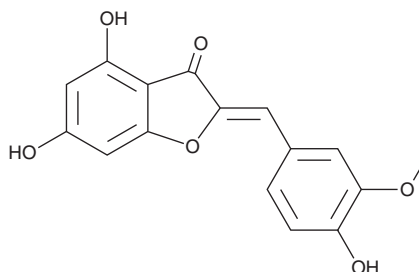
PRODUCT INFORMATION



4,6,4'-Trihydroxy-3'-methoxyaurone

Item No. 37937

CAS Registry No.: 873787-99-6
Formal Name: 4,6-dihydroxy-2Z-[(4-hydroxy-3-methoxyphenyl)methylene]-3(2H)-benzofuranone
MF: C₁₆H₁₂O₆
FW: 300.3
Purity: ≥95%
UV/Vis.: λ_{max}: 395 nm
Supplied as: A crystalline 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

4,6,4'-Trihydroxy-3'-methoxyaurone is supplied as a crystalline solid. A stock solution may be made by dissolving the 4,6,4'-trihydroxy-3'-methoxyaurone in the solvent of choice, which should be purged with an inert gas. 4,6,4'-Trihydroxy-3'-methoxyaurone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 4,6,4'-trihydroxy-3'-methoxyaurone in these solvents is approximately 30 mg/ml. 4,6,4'-Trihydroxy-3'-methoxyaurone is slightly soluble in ethanol.

4,6,4'-Trihydroxy-3'-methoxyaurone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4,6,4'-trihydroxy-3'-methoxyaurone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 4,6,4'-Trihydroxy-3'-methoxyaurone has a solubility of approximately 0.50 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4,6,4'-Trihydroxy-3'-methoxyaurone is an herbicide.¹ It inhibits the growth of rape root and barnyardgrass seedlings when used at concentrations of 100 and 1,000 µg/ml.

Reference

1. Zhang, M., Xu, X.-H., Cui, Y., *et al.* Synthesis and herbicidal potential of substituted aurones. *Pest Manag. Sci.* **68(11)**, 1512-1522 (2012).

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, 11/17/2022

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