

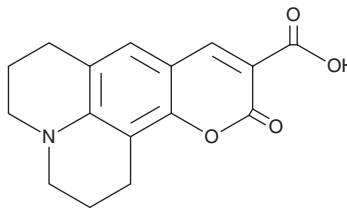
PRODUCT INFORMATION



Coumarin 343

Item No. 41970

CAS Registry No.: 55804-65-4
Formal Name: 2,3,6,7-tetrahydro-11-oxo-1H,5H,11H-[1]benzopyrano[6,7,8-ij]quinolizine-10-carboxylic acid
MF: C₁₆H₁₅NO₄
FW: 285.3
Purity: ≥95%
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

Coumarin 343 is supplied as a solid. A stock solution may be made by dissolving the coumarin 343 in the solvent of choice, which should be purged with an inert gas. Coumarin 343 is slightly soluble (0.1-1 mg/ml) in chloroform and DMSO.

Description

Coumarin 343 is a coumarin laser dye.¹ It displays absorption/emission maxima of 446/519 nm, respectively.

Reference

1. Reynolds, G.A. and Drexhage, K.H. New coumarin dyes with rigidized structure for flashlamp-pumped dye lasers. *Opt. Commun.* **13(3)**, 222-225 (1975).

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, 10/17/2024

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