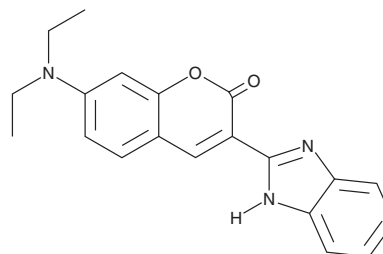


PRODUCT INFORMATION



Coumarin 7 Item No. 35762

CAS Registry No.: 27425-55-4
Formal Name: 3-(1H-benzimidazol-2-yl)-7-(diethylamino)-2H-1-benzopyran-2-one
Synonyms: C.I. 551200, NSC 303254
MF: C₂₀H₁₉N₃O₂
FW: 333.4
Purity: ≥98%
UV/Vis.: λ_{max}: 250, 435 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

Coumarin 7 is supplied as a solid. A stock solution may be made by dissolving the coumarin 7 in the solvent of choice, which should be purged with an inert gas. Coumarin 7 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of coumarin 7 in these solvents is approximately 1 mg/ml.

Description

Coumarin 7 is a coumarin laser dye.^{1,2}

References

1. Mishra, V.R. and Sekar, N. Photostability of coumarin laser dyes - a mechanistic study using global and local reactivity descriptors. *J. Fluoresc.* **27(3)**, 1101-1108 (2017).
2. Yenagi, J.V., Gorbali, M.R., Savildatti, M.I., *et al.* Study of photophysical parameters and measurement of laser characteristics of coumarin 7 and acriflavine. *Spectrosc. Lett.* **25(1)**, 63-72 (1992).

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, 12/13/2022

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