

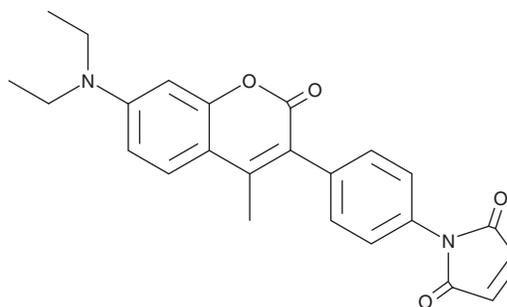
# PRODUCT INFORMATION



## 7-Diethylamino-3-(4-maleimidophenyl)-4-methylcoumarin

Item No. 18209

**CAS Registry No.:** 76877-33-3  
**Formal Name:** 1-[4-[7-(diethylamino)-4-methyl-2-oxo-2H-1-benzopyran-3-yl]phenyl]-1H-pyrrole-2,5-dione  
**MF:** C<sub>24</sub>H<sub>22</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 402.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 210, 378 nm  
**Ex./Em. Max:** 387/468 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

7-Diethylamino-3-(4-maleimidophenyl)-4-methylcoumarin is supplied as a solid. A stock solution may be made by dissolving the 7-diethylamino-3-(4-maleimidophenyl)-4-methylcoumarin in the solvent of choice, which should be purged with an inert gas. 7-Diethylamino-3-(4-maleimidophenyl)-4-methylcoumarin is slightly soluble in DMSO and methanol.

### Description

7-Diethylamino-3-(4-maleimidophenyl)-4-methylcoumarin is a thiol-reactive fluorescent probe.<sup>1</sup> It displays excitation/emission maxima of 387/468 nm, respectively, and fluorescence intensity increases when bound to cysteine residues.<sup>2</sup>

### References

1. Sippel, T.O. New fluorochromes for thiols: Maleimide and iodoacetamide derivatives of a 3-phenylcoumarin fluorophore. *J. Histochem. Cytochem.* **29(2)**, 314-316 (1981).
2. Chen, R.F. and Scott, C.H. Atlas of fluorescence spectra and lifetimes of dyes attached to protein. *Anal. Lett.* **18(4)**, 393-421 (1985).

#### 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, 12/07/2022

#### 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