

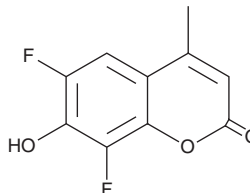
PRODUCT INFORMATION



6,8-Difluoro-7-hydroxy-4-methylcoumarin

Item No. 36017

CAS Registry No.: 215868-23-8
Formal Name: 6,8-difluoro-7-hydroxy-4-methyl-2H-1-benzopyran-2-one
MF: C₁₀H₆F₂O₃
FW: 212.2
Purity: ≥98%
Abs./Em. Max: 358/455 nm
UV/Vis.: λ_{max}: 320 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

6,8-Difluoro-7-hydroxy-4-methylcoumarin is supplied as a solid. A stock solution may be made by dissolving the 6,8-difluoro-7-hydroxy-4-methylcoumarin in the solvent of choice, which should be purged with an inert gas. 6,8-Difluoro-7-hydroxy-4-methylcoumarin is soluble in methanol, DMSO, and dimethyl formamide.

Description

6,8-Difluoro-7-hydroxy-4-methylcoumarin is a fluorophore.¹ It has been conjugated to streptavidin or secondary antibodies for flow cytometry and used to label bacteria for multi-color fluorescence *in situ* hybridization (FISH).^{1,2} 6,8-Difluoro-7-hydroxy-4-methylcoumarin has also been coupled to peptides to improve electrospray ionization and MS sensitivity in MALDI-TOF MS.³ It displays absorption/emission maxima of 358/455 nm, respectively.⁴

References

1. Takada, T., Matsumoto, K., and Nomoto, K. Development of multi-color FISH method for analysis of seven *Bifidobacterium* species in human feces. *J. Microbiol. Methods* **58(3)**, 413-421 (2004).
2. Telford, W.G. Analysis of UV-excited fluorochromes by flow cytometry using near-ultraviolet laser diodes. *Cytometry A*. **61(1)**, 9-17 (2004).
3. Pashkova, A., Moskovets, E., and Karger, B.L. Coumarin tags for improved analysis of peptides by MALDI-TOF MS and MS/MS. 1. Enhancement in MALDI MS signal intensities. *Anal. Chem.* **76(15)**, 4550-4557 (2004).
4. Sun, W.C., Gee, K.R., and Haugland, R.P. Synthesis of novel fluorinated coumarins: Excellent UV-light excitable fluorescent dyes. *Bioorg. Med. Chem. Lett.* **8(22)**, 3107-3110 (1998).

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, 12/15/2022

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