

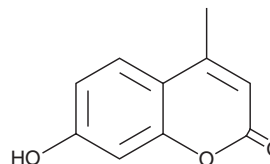
PRODUCT INFORMATION



4-Methylumbelliferone

Item No. 29603

CAS Registry No.: 90-33-5
Formal Name: 7-hydroxy-4-methyl-2H-1-benzopyran-2-one
Synonyms: 4-MU, Hymecromone, 4-Methyl-7-hydroxycoumarin, 7-Hydroxy-4-methylcoumarin
MF: C₁₀H₈O₃
FW: 176.2
Purity: ≥98%
Ex./Em. Max: 320 and 360 nm at low and high pH, respectively/
445 and 455 nm at low and high pH, respectively
UV/Vis.: λ_{max}: 218, 325 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

4-Methylumbelliferone is supplied as a solid. A stock solution may be made by dissolving the 4-methylumbelliferone in the solvent of choice, which should be purged with an inert gas. 4-Methylumbelliferone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-methylumbelliferone in ethanol is approximately 0.25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

4-Methylumbelliferone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-methylumbelliferone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 4-Methylumbelliferone has a solubility of approximately 0.5 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-Methylumbelliferone (4-MU) is a synthetic coumarin with diverse biological activities.¹⁻³ It inhibits hyaluronic acid synthesis in PC3-ML and DU145 cells (IC₅₀s = ~0.4 mM for both).¹ 4-MU inhibits the growth of PC3-ML, DU145, C4-2B, LNCaP, and LAPC-4 prostate cancer cells (IC₅₀s = 0.2-0.4 mM) and invasion and chemotactic motility of PC3-ML and DU145 cells *in vitro*. It reduces tumor growth and microvessel density in a PC3-ML mouse xenograft model when administered at doses of 225 and 450 mg/kg twice per day. 4-MU is active against *H. pylori* and *V. cholerae in vitro* (MICs = 59 and 100-200 µg/ml, respectively).² Derivatives and conjugates of 4-MU have been used as fluorogenic substrates to measure enzyme activity.⁴⁻⁶ Upon enzymatic cleavage, 4-MU is released and its fluorescence can be used to quantify enzyme activity. 4-MU fluorescence is pH-dependent with excitation maxima of 320 and 360 nm at low (1.97-6.72) and high pH (7.12-10.3), respectively, and an emission maximum ranging from 445 to 455 nm, increasing as pH decreases.³

References

1. Lokeshwar, V.B., Lopez, L.E., Munoz, D., *et al. Cancer Res.* **70(7)**, 2613-2623 (2010).
2. Kawase, M., Varu, B., Shah, A., *et al. Arzneimittelforschung* **51(1)**, 67-71 (2001).
3. Zhi, H., Wang, J., Wang, S., *et al. J. Spectrosc.* **1(1)**, 147128 (2013).
4. Karpova, E.A., Voznyi, Y.V., Keulemans, J.L.M., *et al. J. Inherit. Metab. Dis.* **19(3)**, 278-285 (1996).
5. Kelly, S. and Bakhru-Kishore, R. *Clinica Chimica Acta* **97(2-3)**, 239-242 (1979).
6. Wendeler, M., and Sandhoff, K. *Glycoconj. J.* **26(8)**, 945-952 (2009).

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