

MUG

MUG (4-MU-glcUA; 4-Methylumbelliferyl- β -D-glucuronic acid dihydrate), also known as MUGlcU, is a fluorogenic β -glucuronidase substrate that releases the blue fluorescent 4-methyl-7-hydroxycoumarin on enzymatic hydrolysis. The substrate is commonly used for identifying *E. coli* contamination and for detecting marker GUS gene expression in plants with high sensitivity.



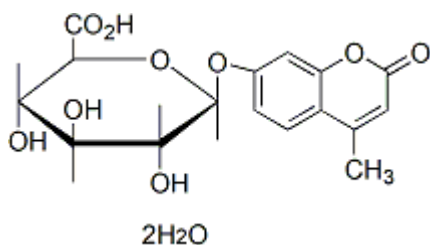
Product attributes

CAS number	6160-80-1
Excitation/Emission	360/450 nm (end product)

Product Description

MUG (4-MU-glcUA; 4-Methylumbelliferyl- β -D-glucuronic acid dihydrate), also known as MUGlcU, is a fluorogenic β -glucuronidase substrate that releases the blue fluorescent 4-methyl-7-hydroxycoumarin on enzymatic hydrolysis. The substrate is commonly used for identifying *E. coli* contamination and for detecting marker GUS gene expression in plants with high sensitivity.

- $\lambda_{Ex}/\lambda_{Em}$ = 360/450 nm (end product)
- White solid soluble in DMSO
- Store desiccated at $\leq 4^{\circ}\text{C}$ and protect from light
- $\text{C}_{16}\text{H}_{16}\text{O}_9 \cdot 2\text{H}_2\text{O}$
- MW: 388.3
- [6160-80-1]



References

1. BioTechniques 8, 39 (1990).
2. Plant Mol. Biol. Rep. 5, 387 (1987).

This datasheet was generated on January 2, 2026 at 09:53:14 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mug-muglcu-4-mu-glcua-4-methylumbelliferyl-b-d-glucuronic-acid-dihydrate/>