

## MUG

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



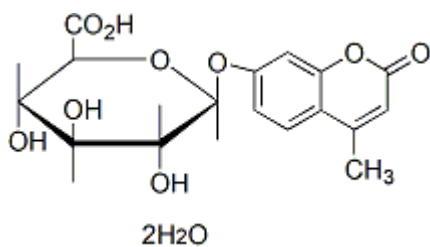
### Product attributes

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

## Product Description

MUG (4-MU-glcUA; 4-Methylumbelliferyl- $\beta$ -D-glucuronic acid dihydrate), also known as MUGlcU, is a fluorogenic  $\beta$ -glucuronidase substrate that releases the blue fluorescent 4-methyl-7-hydroxycoumarin 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).