46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: MUG (MUGlcU; 4-MU-glcUA; 4-methylumbelliferyl-β-D-glucuronic acid

dihydrate)

CATALOG # 10020

MOLECULAR $C_{16}H_{16}O_9 \cdot 2H_2O$ **INFORMATION:** MWt: 388.3

[6160-80-1]

OH OH OH OH OH

PROPERTIES:

Color & Form White solid

Soluble in water at 100mg/mL

Absorption/Emission 360/450 nm

Extinction Coefficient

Extinction

STORAGE AND HANDLING:

Store at 4°C. For long-term storage, store at -20°C. MUG at 100 mg/100 ml of H₂O plus 2 drops of Triton X-100, which is sterile filtered, is stable for 6 months when

stored refrigerated.

APPLICATION: MUG is a fluorogenic β-glucuronidase substrate that releases the blue fluorescent 4-

methyl-7-hydroxycoumarin ($\lambda_{ex} \setminus \lambda_{em} = 360 \setminus 450$ nm) 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.

MUG is also used for identifying E. coli contamination in drinking water and for rapid

bacterial identification in blood cultures.

Ref: BioTechnique 8, 39(1990); Plant Mol. Biol. Rep. 5, 387(1987)

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on

clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.