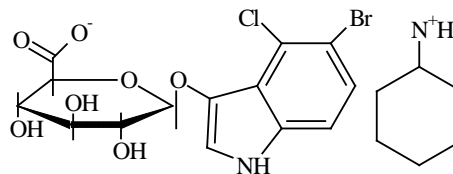


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **X-GlcU, CHA** (5-bromo-4-chloro-3-indolyl- β -D-glucuronide, cyclohexylammonium salt)

CATALOG NUMBER: 10016

MOLECULAR INFORMATION: $C_{20}H_{26}BrClN_2O_7$
Mwt: 521.8
[18656-96-7]

**PROPERTIES:**

Color & Form White solid
Purity $\geq 99\%$ by HPLC
Solubility Soluble in water and DMF (2%, DMF)

STORAGE AND HANDLING: Store desiccated at $-20\text{ }^{\circ}\text{C}$ upon receipt. Protect from light, especially in solution.

APPLICATION: Chromogenic β -glucuronidase substrate widely used for monitoring *E. coli* contamination in food (such as meat, dairy products and shellfish, etc) and water and for the detection of GUS gene expression in plants. X-GlcU also has clinical applications in the assessment of urinary tract infection by detecting the presence of *E. coli*. The substrate yields a dark blue precipitate after enzymatic hydrolysis.

Ref: *J. Food Protect.* **53**(6), 508(1990); *Appl. Environ. Microbiol.* **54**, 1874(1988); *J. Clin. Microbiol.* **27**(4), 778(1989); *Plant cell Reports* **13**(1), 17(1993); *J. Am. Soc. Horticult. Sci.* **115**(4), 686(1990).

TOXICITY: Unknown

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.