

X-GlcU, CHA

Alternate name: BCIG. X-GlcU (5-Bromo-4-chloro-3-indolyl- β -D-glucuronide, cyclohexylammonium salt) is a 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.



Product attributes

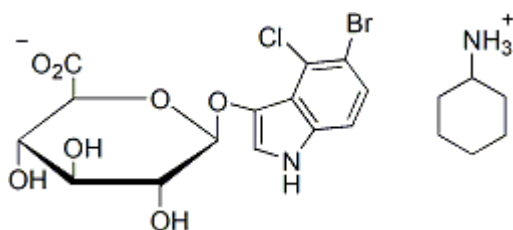
CAS number	18656-96-7
------------	------------

Product Description

Alternate name: BCIG

X-GlcU (5-Bromo-4-chloro-3-indolyl- β -D-glucuronide, cyclohexylammonium salt) is a 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 (λ_{max} 615 nm) precipitate after enzymatic hydrolysis.

- White solid soluble in water or DMSO
- Store desiccated at -20 °C and protect from light
- $20\text{H}_{26}\text{BrClN}_2\text{O}_7$
- MW: 521.8
- [18656-96-7]



References

1. J Food Protect 53(6), 508 (1990).
2. Appl Environ Microbiol 54, 1874 (1988).
3. J Clin Microbiol 27(4), 778 (1989).
4. Plant Cell Reports 13(1), 17 (1993).
5. J Am Soc Horticul Sci 115(4), 686 (1990).

This datasheet was generated on January 18, 2026 at 02:22:07 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/x-glcu-cha-also-called-bcig-or-5-bromo-4-chloro-3-indolyl-b-dglucuronidecyclohexylammonium-salt/>