X-GIcU, Na

X-GlcU, Na (5-Bromo-4-chloro-3-indolyl-β-D-glucuronide, sodium salt), also known as X-GlcA, Na, is similar to X-GlcU, CHA, but has better water solubility.



Product attributes

Call us: 800-304-5357

CAS number

12954-41-9

Product Description

X-GlcU, Na (5-Bromo-4-chloro-3-indolyl-β-D-glucuronide, sodium salt), also known as X-GlcA, Na, is similar to X-GlcU, CHA

(<https://biotium.com/product/x-glcu-cha-also-called-bcig-or-5-bromo-4-chloro-3-indolyl-b-dglucuronidecyclohexylammonium-salt/">10016), but has better water solubility.

- White solid soluble in water or DMSO
- Store desiccated at 4°C and protect from light
- C₁₄H₁₂BrClNNaO₇
- MW: 444.6
- [12954-41-9]

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 Horticult Sci 115(4), 686 (1990).

This datasheet was generated on November 29, 2025 at 11:23:43 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/x-glcu-na-x-glca-na-5-bromo-4-chloro-3-indolyl-b-d-glucuronide-sodium-salt/