Rose-β-D-GlcU,CHA

Rose-β-D-GlcU

(6-Chloro-3-indolyl-β-D-glucuronide,cyclohexylammonium salt), also known as Salmon-β-D-GlcUA, is similar to X-GlcU, but the product of the substrate is a rose-colored ($\lambda_{max} \sim 540$ nm) precipitate, which is easier to detect against the background of plant cells.

nictum 680

Call us : 800-304-5357

Product attributes

CAS number

138182-20-4

Product Description

Rose- β -D-GlcU (6-Chloro-3-indolyl- β -D-glucuronide,cyclohexylammonium salt), also known as Salmon- β -D-GlcUA, is similar to X-GlcU, but the product of the substrate is a rose-colored ($\lambda_{max} \sim 540$ nm) precipitate, which is easier to detect against the background of plant cells.

- White solid soluble in H2O or DMSO
- Store desiccated at 4°C and protect from light
- C₂₀H₂₇CIN₂O₇
- MW: 442.9
- [138182-20-4]

$$\begin{array}{c} -O_2C \\ OH \\ OH \\ OH \end{array}$$

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 11:24:48 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rose-b-d-glcua-6-chloro-3-indolyl-b-dglucuronidecyclohexylammonium-salt/