

X-β-D-cellobioside

X-β-D-cellobioside (5-Bromo-4-chloro-3-indolyl-beta-D-cellobioside) is a chromogenic substrate for β-cellobiosidase. Enzymatic cleavage of the substrate generates a dark blue-colored precipitate.



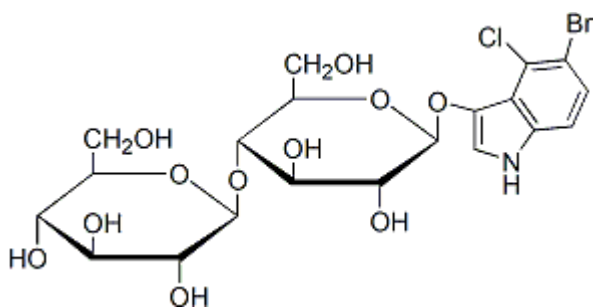
Product attributes

CAS number	177966-52-8
------------	-------------

Product Description

X-β-D-cellobioside (5-Bromo-4-chloro-3-indolyl-beta-D-cellobioside) is a chromogenic substrate for β-cellobiosidase. Enzymatic cleavage of the substrate generates a dark blue-colored precipitate.

- White solid soluble in DMSO
- Store at 4 °C and protect from light
- C₂₀H₂₅BrClNO₁₁
- MW: 570.78
- [177966-52-8]



This datasheet was generated on June 15, 2026 at 12:49:33 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/x-b-d-cellobioside-5-bromo-4-chloro-3-indolyl-beta-d-cellobioside/>