

Red- β -D-Gal

Red- β -D-Gal (5-Bromo-6-chloro-3-indolyl- β -D-galactopyranoside) is a chromogenic substrate for β -galactosidase. Similar to X-gal, but generates a red or magenta ($\lambda_{\text{max}} \sim 565 \text{ nm}$) colored precipitate on enzyme hydrolysis.



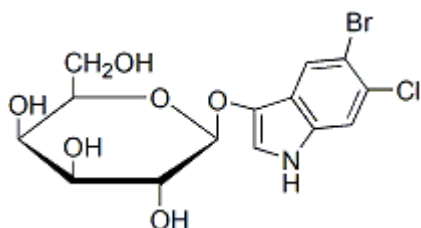
Product attributes

CAS number	93863-88-8
------------	------------

Product Description

Red- β -D-Gal (5-Bromo-6-chloro-3-indolyl- β -D-galactopyranoside) is a chromogenic substrate for β -galactosidase. Similar to X-gal, but generates a red or magenta ($\lambda_{\text{max}} \sim 565 \text{ nm}$) colored precipitate on enzyme hydrolysis.

- White solid soluble in DMSO
- Store at 4 °C and protect from light
- $\text{C}_{14}\text{H}_{15}\text{BrClNO}_6$
- MW: 408.6
- [93863-88-8]



References

1. Molecular Cloning: A Laboratory Manual, Cold Spring Harbor Laboratory, B. 14, p 186 (1989).

This datasheet was generated on January 8, 2026 at 12:58:06 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/red-b-d-gal-5-bromo-6-chloro-3-indolyl-b-d-galactopyranoside/>