## 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 ( $\lambda_{max} \sim 565$  nm) colored precipitate on enzyme hydrolysis.



## **Product attributes**

CAS number

93863-88-8

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

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

- White solid soluble in DMSO
- Store at 4°C and protect from light
- C<sub>14</sub>H<sub>15</sub>BrClNO<sub>6</sub>
- 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 November 8, 2025 at 09:40:45 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/red-b-d-gal-5-bromo-6-chloro-3-indolyl-b-d-galactopyranoside/">https://biotium.com/product/red-b-d-gal-5-bromo-6-chloro-3-indolyl-b-d-galactopyranoside/</a>