

## Purple- $\beta$ -D-Gal

Purple- $\beta$ -D-Gal (5-Iodo-3-indolyl- $\beta$ -D-galactopyranoside) is a chromogenic substrate for  $\beta$ -galactosidase. Similar to X-gal, but generates a purple ( $\lambda_{\max}$  ~575 nm) colored precipitate on enzyme hydrolysis.



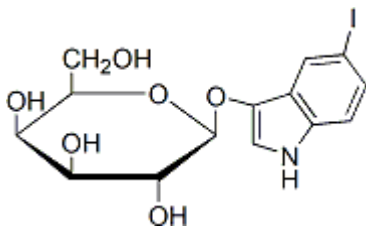
### Product attributes

CAS number	36473-36-6
------------	------------

## Product Description

Purple- $\beta$ -D-Gal (5-Iodo-3-indolyl- $\beta$ -D-galactopyranoside) is a chromogenic substrate for  $\beta$ -galactosidase. Similar to X-gal, but generates a purple ( $\lambda_{\max}$  ~575 nm) colored precipitate on enzyme hydrolysis.

- Off-white solid soluble in water and DMSO.
- Store at 4 °C and protect from light
- C<sub>14</sub>H<sub>16</sub>INO<sub>6</sub>



## References

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

This datasheet was generated on June 20, 2026 at 08:01:11 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/purple-b-d-gal-5-iodo-3-indolyl-b-d-galactopyranoside/>