

Purple- β -D-Gal

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



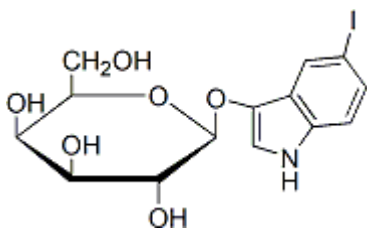
Product attributes

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

Product Description

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

- Off-white solid soluble in water and DMSO.
- Store at 4 °C and protect from light
- C₁₄H₁₆INO₆



References

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

This datasheet was generated on January 21, 2026 at 01:33:12 AM. 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/>