Glowing products for science

Purple-β-D-Gal

Purple-β-D-Gal (5-lodo-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

Call us: 800-304-5357

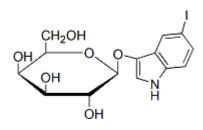
CAS number

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 November 23, 2025 at 03:59:23 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/