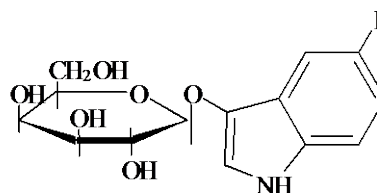


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Purple- $\beta$ -D-Gal (5-iodo-3-indolyl- $\beta$ -D-galactopyranoside)**CATALOG NUMBER:** 10014**MOLECULAR INFORMATION:** C<sub>14</sub>H<sub>16</sub>INO<sub>6</sub>  
Mwt: 421.6**PROPERTIES:**

<b>Color &amp; Form</b>	Off-white solid
<b>Purity</b>	≥ 99% by TLC
<b>Solubility</b>	Soluble in DMF or DMSO

**STORAGE AND HANDLING:** Store desiccated at ≤ 4 °C upon receipt. Protect from light, especially in solution.**APPLICATION:** Chromogenic substrate for  $\beta$ -galactosidase. Similar to X-gal, but generates a purple colored precipitate ( $\lambda_{\text{abs}} \sim 575 \text{ nm}$ ) on enzyme hydrolysis. Protocol for using the material is similar to that for X-gal.

Ref: 1) Molecular Cloning: A laboratory Manual, Cold Spring Harbor Laboratory, B. 14, p186(1989); 2) *Biotechniques*, **27**, 438(1999); 3) *Oncogen*, **10**, 2323(1995); 4) *Devel. Biol.*, **161**, 77(1994); 5) *J. Histochem. Cytochem.*, **42**, 1299(1994); 6) *BioTechniques*, **15**, 974(1993); 7) *BioTechniques*, **15**, 292(1993).

**TOXICITY:** Unknown

<b>FIRST AID:</b>	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

<b>Disclaimer:</b> Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.
---