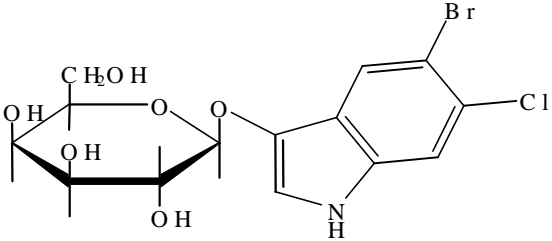


5-bromo-6-chloro-3-indolyl- β -D-galactopyranoside

Instruction Manual

| | |
|---------------------------------------|---|
| Catalog Number | PK-CA707-10012 |
| Description | Chromogenic substrate for β -galactosidase. Similar to X-gal, but generates a red or magenta ($\lambda_{\text{max}} \sim 565 \text{ nm}$) colored precipitate on enzyme hydrolysis. |
| Quantity | 100 mg |
| Excitation / Emission Maxima | NA |
| Molecular Structure |  |
| Molecular Weight / Molecular Formula | 408.6 Da; $\text{C}_{13}\text{H}_{15}\text{BrClNO}_6$ |
| Purity | $\geq 98\%$ by TLC |
| Appearance / Formulation / Solubility | White solid; soluble in DMSO and MeOH. |
| Storage & Stability | Store desiccated at -20°C upon receipt. Protect from light, especially when in solution. |
| Applications | β -galactosidase assays (e.g. lacZ reporter gene assays) |
| References | <ol style="list-style-type: none"> 1) <i>Molecular Cloning: A laboratory Manual</i>, Cold Spring Harbor Laboratory, B. 14, p186(1989) 2) <i>Biotechniques</i>, 27, 438(1999) 3) <i>Oncogen</i>, 10, 2323(1995) 4) <i>Devel. Biol.</i>, 161, 77(1994) 5) <i>J. Histochem. Cytochem.</i>, 42, 1299(1994) 6) <i>BioTechniques</i>, 15, 974(1993) 7) <i>BioTechniques</i>, 15, 292(1993) |
| Caution | 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. |

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.