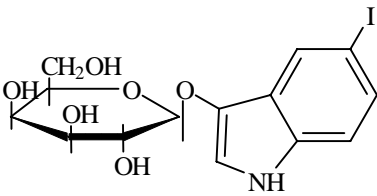


5-iodo-3-indolyl- β -D-galactopyranoside

Instruction Manual

Catalog Number	PK-CA707-10014
Description	Chromogenic substrate for β -galactosidase. Similar to X-gal, but generates a purple (λ_{max} ~ 575 nm) colored precipitate on enzyme hydrolysis.
Quantity	25 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	421.6 Da; C ₁₄ H ₁₆ ClNO ₆
Purity	≥ 99% by TLC
Appearance / Formulation / Solubility	Off-white solid; soluble in DMF, DMSO and water.
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.