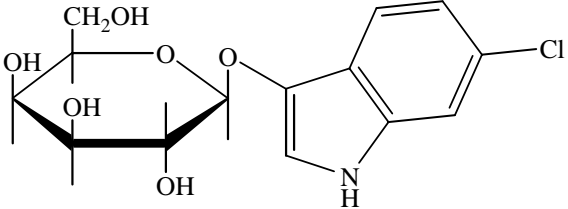


6-chloro-3-indoxyl-β-D-galactopyranoside

Instruction Manual

Catalog Number	PK-CA707-10013
Description	Rose-β-D-Gal is a chromogenic substrate for β-galactosidase. The product is similar to X-gal, but generates an intense pink colored precipitate ($\lambda_{max} \sim 540 \text{ nm}$) on enzymatic hydrolysis. Protocol for using the material is similar to that for X-gal (see references listed below).
Quantity	100 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	329.74 Da; C ₁₄ H ₁₆ ClNO ₆
Purity	≥ 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) 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)
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.