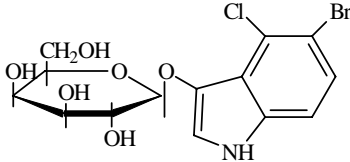


5-bromo-4-chloro-3-indolyl- β -galactopyranoside

Instruction Manual

Catalog Number	PK-CA707-10011
Description	X-gal is the most widely used chromogenic substrate for β -galactosidase. Enzyme hydrolysis of the substrate yields a dark blue colored precipitate (λ_{max} ~ 615 nm).
Quantity	100 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	408.6 Da; C ₁₃ H ₁₅ BrClNO ₆
Purity	≥ 99% by HPLC
Appearance / Formulation / Solubility	White solid; soluble in DMSO or DMF.
Storage & Stability	Store desiccated at -20°C. Hygroscopic! Keep tightly closed. Protect from light.
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.