

Chromogenic beta-Galactosidase Substrate Sampler Kit



Instruction Manual

Catalog Number	PK-CA707-10024
Description	For researchers wishing to investigate the suitability of PromoKine's various chromogenic β -galactosidase substrates, we offer this sampler kit, which contains 5 mg each of the following five different substrates: X-Gal (# PK-CA707-10011), Red- β -D-Gal (# PK-CA707-10012), Rose- β -D-Gal (# PK-CA707-10013), Purple- β -D-Gal (# PK-CA707-10014) and Green- β -D-Gal (# PK-CA707-10015).
Quantity	5 x 5 mg
Excitation / Emission Maxima	NA
Molecular Structure	See data sheets of the individual β -galactosidase substrates.
Molecular Weight / Molecular Formula	See data sheets of the individual β -galactosidase substrates.
Purity	See data sheets of the individual β -galactosidase substrates.
Appearance / Formulation / Solubility	Solids; soluble in DMSO, DMF and/or water. See data sheets of the individual β -galactosidase substrates
Storage & Stability	Store at -20°C. Protect from light. See data sheets of the individual β -galactosidase substrates.
Applications	β -galactosidase assays (e.g. lacZ reporter gene assays)
References	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	See data sheets of the individual β -galactosidase substrates.

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