



X-Gal
Catalog Number 170-3455
(5-bromo-4-chloro-3-indolyl- β -D-galactopyranoside)

Contents	250 mg
Storage	Desiccated below 0 °C. Protect from light.
Use	<p>X-Gal is used to indicate the presence of functional β-galactosidase. It is a substrate of the enzyme and forms a blue precipitate when cleaved. For convenience, Bio-Rad provides X-Gal in an oversized bottle. The total 250 mg can be dissolved directly in the bottle with 12.5 ml of dimethylformamide to give a final concentration of 2%. The solution can be stored in the original bottle at 4 °C for up to 1 year.</p> <p>This product is intended for research use only.</p>
Quality assurance	<p>X-Gal has been tested functionally with <i>E. coli</i> MV1190. Blue colonies form only in the presence of X-gal and IPTG (isopropyl-β-D-thiogalactopyranoside), an inducer of β-galactosidase.</p>