## Chromogenic β-Galactosidase Substrate Sampler

For researchers wishing to investigate the suitability of our various chromogenic  $\beta$ -galactosidase substrates, we offer this sampler kit, which contains 5 mg each of the following five different substrates: X-Gal (10011), Red- $\beta$ -D-Gal (10012), Rose- $\beta$ -D-Gal (10013), Purple- $\beta$ -D-Gal (10014) and Green- $\beta$ -D-Gal (10015).



## **Product attributes**

Call us: 800-304-5357

Email: btinfo@biotium.com

## **Product Description**

For researchers wishing to investigate the suitability of our various chromogenic  $\beta$ -galactosidase substrates, we offer this sampler kit, which contains 5 mg each of the following five different substrates: X-Gal (10011), Red- $\beta$ -D-Gal (10012), Rose- $\beta$ -D-Gal (10013), Purple- $\beta$ -D-Gal (10014) and Green- $\beta$ -D-Gal (10015).

This datasheet was generated on November 8, 2025 at 07:53:30 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/chromogenic-b-galactosidase-substrate-sampler-kit/">https://biotium.com/product/chromogenic-b-galactosidase-substrate-sampler-kit/</a>