



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## Chromogenic $\beta$ -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

### 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 June 4, 2026 at 06:53:39 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/chromogenic-b-galactosidase-substrate-sampler-kit/>