



Glowing products for science

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

Chromogenic β -Glucuronidase Substrate Sampler

For researchers wishing to investigate the suitability of our various chromogenic β -glucuronidase substrates, we offer this sampler kit, which contains 5 mg each of the following three different substrates: X-GlcU, CHA, Red- β -D-GlcU, CHA, and Rose- β -D-GlcU, CHA.



Product attributes

Product Description

For researchers wishing to investigate the suitability of our various chromogenic β -glucuronidase substrates, we offer this sampler kit, which contains 5 mg each of the following three different substrates: X-GlcU, CHA ([10016](#)), Red- β -D-GlcU, CHA ([10018](#)), and Rose- β -D-GlcU, CHA ([10019](#)).

This datasheet was generated on June 4, 2026 at 07:01:58 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/chromogenic-b-glucuronidase-substrate-sampler-kit/>