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

Call us: 800-304-5357

Email: btinfo@biotium.com

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 November 4, 2025 at 06:47:39 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/chromogenic-b-glucuronidase-substrate-sampler-kit/