



Glowing products for science

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

Chromogenic Phosphatase Substrate Sampler

The sampler kit contains 10 mg each of BCIP Na (10001), BCIP Pink (10006), BCIP Red (10004) and 250 mg of NBT (10008). NBT forms a co-precipitant and enhances the BCIP reaction.



Product attributes

Product Description

The sampler kit contains 10 mg each of BCIP Na ([10001](#)), BCIP Pink ([10006](#)), BCIP Red ([10004](#)) and 250 mg of NBT ([10008](#)). NBT forms a co-precipitant and enhances the BCIP reaction.

This datasheet was generated on January 8, 2026 at 10:20:12 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/chromogenic-phosphatase-substrate-sampler-kit/>