Glowing products for science

Calcium Calibration Buffer Kit

This kit provides buffers for generating a range of accurate calcium concentrations for the calibration of fluorescent Ca^{2+} indicators.



Product attributes

Storage Conditions

Store at 2 to 8 °C

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

This kit provides calibration buffers and a protocol for producing a range of accurate calcium concentrations for the calibration of fluorescent Ca²⁺ indicators.

This datasheet was generated on November 6, 2025 at 02:26:58 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/calcium-calibration-buffer-kit-zero-and-10-mm-caegta-2x50-ml/