



Glowing products for science

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

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

Product Description

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

This datasheet was generated on June 3, 2026 at 02:57:30 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/>