



Glowing products for science

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

MCB Glutathione Detection Kit

This product is discontinued and no longer available.



Product attributes

Product Description

This product is discontinued and no longer available.

This datasheet was generated on June 24, 2026 at 09:11:22 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mcb-glutathione-detection-kit/>