



Glowing products for science

Precast GelRed® Agarose Gels, 1% Agarose/TAE

This item is discontinued.



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

Product attributes

Assay type/options	DNA/RNA gel staining
--------------------	----------------------

Product Description

This item is discontinued.

References

Download a list of selected [References for GelRed® and GelGreen®](#).

This datasheet was generated on June 13, 2026 at 02:05:13 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/precast-gelred-agarose-gels-1-agarose-tae/>