



Glowing products for science

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

## Gel-Bright™ LED Gel Illuminator

This item is discontinued.



### Product attributes

## Product Description

This item is discontinued.

## References

1. PLoS ONE 14(4): e0208966, [DOI: 10.1371/journal.pone.0208966](https://doi.org/10.1371/journal.pone.0208966)

This datasheet was generated on January 18, 2026 at 06:25:52 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/gel-bright-led-gel-illuminator/>