



Glowing products for science

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

CF®640R-UTP

CF®640R-UTP can be enzymatically incorporated into RNA probes for use in many biological applications.



Product attributes

Product Description

CF®640R-UTP can be enzymatically incorporated into RNA probes for use in many biological applications. Far-red fluorescent CF®640R is spectrally similar to Alexa Fluor® 647, but is much more photostable.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.

References

Download a list of [CF® dye references](#).

This datasheet was generated on January 9, 2026 at 07:15:37 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf640r-utp-lyophilized-powder/>