

CF®640R-UTP

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



Product attributes

Call us: 800-304-5357

Email: btinfo@biotium.com

Product Description

 ${\sf CF} \\ @ 640 \\ R-{\sf UTP} \ can \ be \ enzymatically \ incorporated \ into \ RNA \ probes \ for \ use \ in \ many \ biological$ 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 November 17, 2025 at 07:47:18 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cf640r-utp-lyophilized-powder/