



Glowing products for science

## CF®555-ddCTP

This product is discontinued.



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

### Product attributes

Excitation/Emission	555/565 nm
Colors	Red
Molecular weight	1383

## Product Description

This product is discontinued.

## References

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

This datasheet was generated on January 18, 2026 at 08:58:14 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cf555-ddctp-lyophilized-powder/>