

## CF®555-ddCTP

This product is discontinued.



## Call us : 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 November 22, 2025 at 01:50:29 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf555-ddctp-lyophilized-powder/">https://biotium.com/product/cf555-ddctp-lyophilized-powder/</a>