

## CF® Dye cAMP

Fluorescently labeled cAMP analogs that can be used to probe cAMP receptors.



### Product attributes

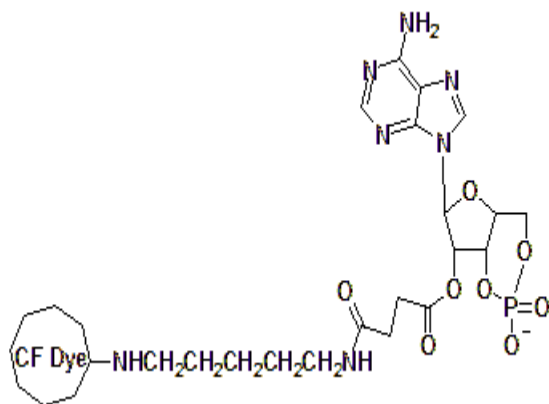
<b>Colors</b>	Green, Far-red
<b>Cell permeability</b>	Membrane impermeant
<b>Storage Conditions</b>	Store at -10 to -35 °C, Protect from light, Desiccate
<b>Reconstitution</b>	Do not reconstitute in buffer as it may cause the material to hydrolyze. Soluble in water, DMSO, or DMF. The stock solution is stable for at least 6 months when stored at -10 to -35 °C and protected from light.

## Product Description

CF®488A-cAMP and CF®640R-cAMP are fluorescently labeled cAMP analogs that can be used to probe cAMP receptors. Single molecules of fluorescently labeled cAMP have been imaged as the probe binds to the receptor on the surface of *Dictyostelium* cells (1).

- Superior CF® Dyes are bright, photostable, and water-soluble
- Available with green or far-red labels
- Soluble in H<sub>2</sub>O, DMF, or DMSO

Also see our [Biotin-cAMP](#) analogs.



### Superior CF® Dyes

Biotium's next-generation CF® dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about [CF® Dyes](#).

### CF® Dye cAMP

Product	Ex/Em	MW (g/mol)	Size	Catalog No.	Dye Features
<a href="#">CF®488A cAMP</a>	490/516 nm	1000	100 ug	<a href="#">00036</a>	<a href="#">CF®488A Features</a>
<a href="#">CF®640R cAMP</a>	642/663 nm	1285	100 ug	<a href="#">00037</a>	<a href="#">CF®640R Features</a>

CF is a registered trademark of Biotium, Inc. Alexa Fluor, Texas Red, and DyLight are registered trademarks of Thermo Fisher Scientific.

## References

1. Science (2001), <https://doi.org/10.1126/science.1063951>

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

This datasheet was generated on December 21, 2024 at 07:44:51 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-camp/>