

CF® Dye cAMP

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



Product attributes

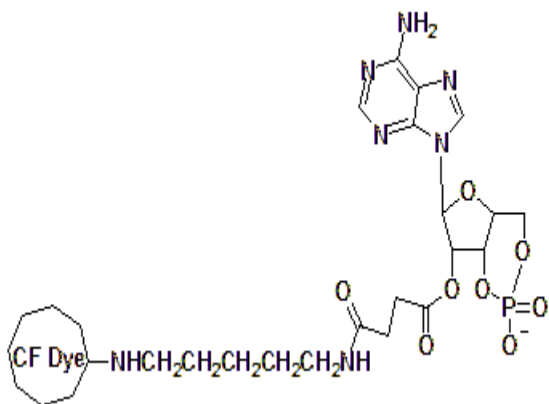
Colors	Green, Far-red
Cell permeability	Membrane impermeant
Storage Conditions	Store at -10 to -35 °C, Protect from light, Desiccate
Reconstitution	Do not reconstitute in buffer as it may cause the material to hydrolyze. Soluble in water, DMSO, or DMF. The stock solution is stable for at least 6 months 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 H₂O, 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
CF®488A cAMP	490/516 nm	1000	100 ug	00036	CF®488A Features
CF®640R cAMP	642/663 nm	1285	100 ug	00037	CF®640R Features

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 March 29, 2025 at 01:35:36 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-camp/>