Cyanine 644-cAMP

A fluorescently labeled cAMP analog that can be used to probe cAMP receptors.



Product attributes

| For live or fixed cells | For live/intact cells |
|-------------------------|-----------------------|
| Colors | Far-red |
| Excitation/Emission | 644/670 nm |
| Molecular weight | 1043.11 |

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Cyanine 644-cAMP is a fluorescently labeled cAMP analog 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).

- Dark Blue solid soluble in H2O, DMF, or DMSO
- Store desiccated at -20°C and protect from light

References

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

This datasheet was generated on November 17, 2025 at 04:10:22 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cyanine-644-camp/