Biotin-cAMP

A biotin-conjugated cAMP analog that can be used to probe cAMP receptors.



Product attributes

For live or fixed cells	For live/intact cells
Conjugation	Biotin
Molecular weight	826
Storage Conditions	Store at -10 to -35 °C. Protect from light. Desiccate

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

Product Description

Cyclic adenosine monophosphate is an important second messenger in many biological processes. This Biotin-cAMP, diisopropylethylammonium salt is a biotin-conjugated cAMP analog that can be used to probe cAMP receptors.

- Cyclic AMP conjugated with a biotin label
- Biotinylated for avidin or streptavidin reactivity and signal amplification
- White lyophilized solid soluble in water

Also see our fluorescently labeled CF® Dye cAMP analogs.

CF is a registered trademark of Biotium, Inc.

References

1. PLoS One, 12, 9, (2017), DOI: 10.1371/journal.pone.0184843

This datasheet was generated on November 4, 2025 at 06:26:57 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-camp-diisopropylethylammonium-salt/