

Biotin-cAMP

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



Product attributes

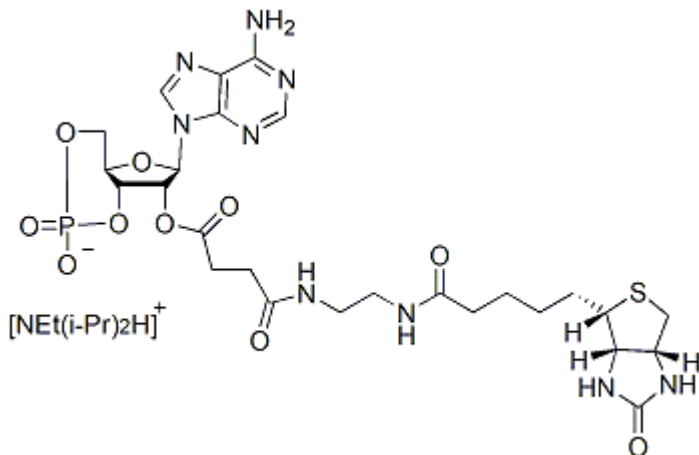
For live or fixed cells	For live/intact cells
Conjugation	Biotin
Molecular weight of antigen	826
Storage Conditions	Store at -10 to -35 °C, Protect from light, Desiccate

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](https://doi.org/10.1371/journal.pone.0184843)

This datasheet was generated on June 4, 2026 at 12:19:52 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-camp-diisopropylethylammonium-salt/>