

Biotin-cGMP, Diisopropylethylammonium Salt

A biotinylated cyclic nucleotide that is an important second messenger in many biological processes and has been used as a tracer in immunoassays.



Product attributes

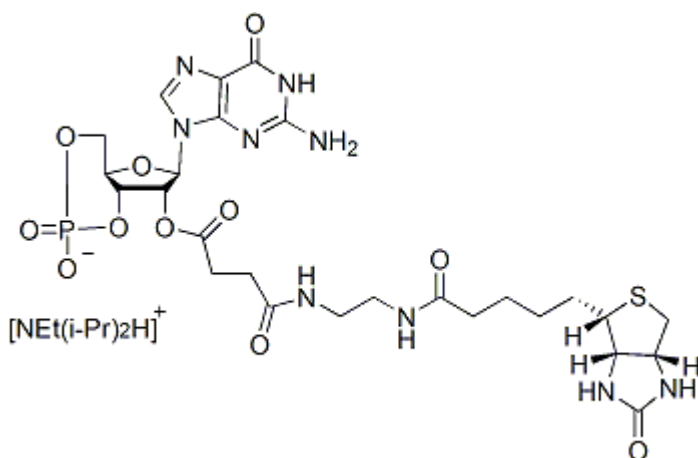
For live or fixed cells	For live/intact cells
Conjugation	Biotin
Molecular weight	832
Storage Conditions	Store at -10 to -35 °C, Protect from light, Desiccate

Product Description

Cyclic guanosine monophosphate (cGMP) is a cyclic nucleotide that is an important second messenger in many biological processes. Biotin-cGMP has been used as a tracer in immunoassays for the cGMP measurement in plasma or urine samples.

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

Chemical Structure:



References

1. PLoS One, 12, 9, (2017), [DOI: 10.1371/journal.pone.0184843](https://doi.org/10.1371/journal.pone.0184843)
2. J Immunoassay 16 (1), 37 (1995), [DOI: 10.1080/15321819508013548](https://doi.org/10.1080/15321819508013548)

This datasheet was generated on January 23, 2026 at 11:05:10 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-cgmp-diisopropylethylammonium-salt/>