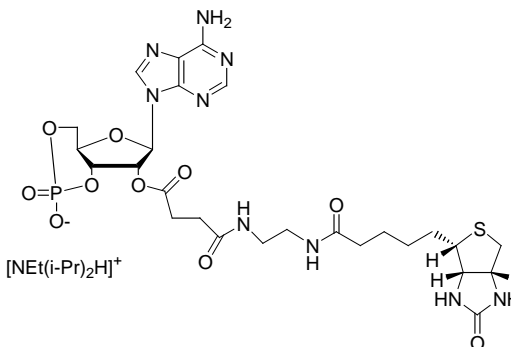


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Biotin-cAMP, diisopropylethylammonium salt**CATALOG #:** 00020/00020-1**MOLECULAR INFORMATION:** C₃₄H₅₅N₁₀O₁₀PS
Mwt: 826**PROPERTIES:**

Color & Form White solid
Purity 98% by HPLC
Solubility Soluble in H₂O, DMSO, and DMF. **Do not reconstitute in buffer as it may cause the material to hydrolyze.**

STORAGE AND HANDLING:

Stored desiccated at -20°C. The shelf life for solid is at least 12 months if stored desiccated at -20°C, and 6 months for solution if stored at -20°C.

APPLICATION:

Biotin-cAMP is a cAMP analog and can be used to probe cAMP receptors. In addition, we are in the process of developing fluorescently labeled cAMP and cGMP. Single molecule of fluorescently labeled cAMP has recently been imaged as the probe binds to the receptor on the surface of *Dictyostelium* cells (M. Ueda, et al. *Science* **294**, 864(2001)).

TOXICITY:

Unknown

FIRST AID:

Potentially harmful; avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.