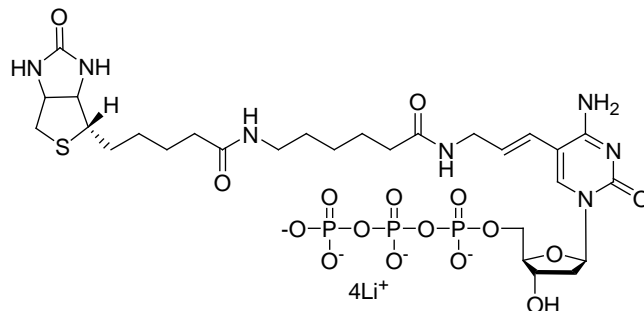


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Biotin-11-dCTP, 1 mM in pH 7.5 Tris-HCl buffer**CATALOG #** 40036**MOLECULAR INFORMATION:** C₂₈H₄₂N₇O₁₆P₃SLi₄
MWt: 885.4**PROPERTIES:**

Color & Form Colorless liquid
Purity ≥ 95% by TLC
Solubility N/A

STORAGE AND HANDLING:

Stored desiccated at -20 °C. This product is shipped on blue ice.

APPLICATION:

Biotin-11-dCTP can be enzymatically incorporated into DNA via nick translation, random priming, or 3'-end terminal labeling. The number "11" is the number of atoms in the linker between biotin and dCTP.

Please also see our streptavidin conjugates labeled with our exceptionally bright and photostable CF™ dyes. There are more than 20 CF dyes available with wavelengths spanning the entire visible and near-IR spectra.

Ref: 1) *Cell*: 73(2), 361 (1993); 2) *Anal Biochem*: 194(1),45(1991); 3) *Anal Biochem*: 169(2),376 (1988)**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.