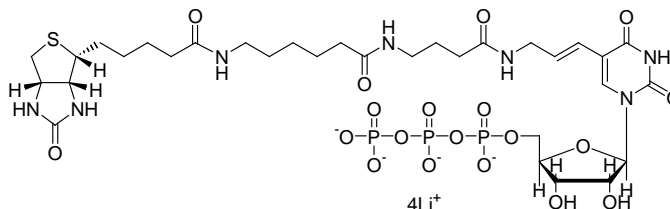


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Biotin-16-UTP, 10 mM in pH 7.5 Tris-HCl buffer**CATALOG #** 40023**MOLECULAR INFORMATION:** C₃₂H₄₈N₇O₁₉P₃SLi₄
MWt: 987.51**PROPERTIES:****Color & Form**
PurityColorless liquid
≥ 95% by TLC**STORAGE AND HANDLING:**

Stored at -20 °C. Decomposition of about 5% may occur within six months.

APPLICATION:

Biotin-16-UTP can be enzymatically incorporated into RNA. The number “16” is the number of atoms in the linker between biotin and UTP. The length of the linker affects the incorporation efficiency of the biotin-UTP probe into RNA using RNA polymerases, and it also affects biotin/avidin or biotin/streptavidin. In general, the shorter the linker, the more efficiently the biotin-UTP is incorporated into RNA by RNA polymerases. On the other hand, the longer the linker, the better biotin can interact with avidin or streptavidin.

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.

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.