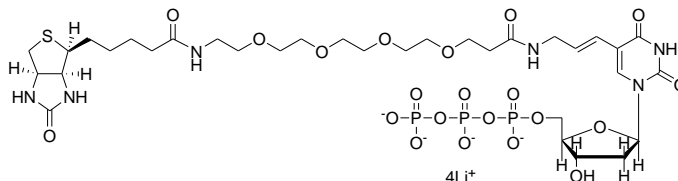


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Biotin-20-dUTP, lyophilized powder

CATALOG # 40030-1

MOLECULAR INFORMATION: $C_{33}H_{51}N_6O_{21}P_3SLi_4$
MWt: 1020.54



PROPERTIES:

Color & Form

Solid

Purity

≥ 95% by TLC

Contains lyophilized TE buffer

STORAGE AND HANDLING:

Store at -20 °C

APPLICATION:

This product is similar to biotin-11-dUTP (#40029) and biotin-16-dUTP (#40022) except that it has a longer and water-soluble spacer group, which should facilitate interaction between the biotin group and avidin or streptavidin. Note: for PCR applications, Taq polymerase should be used with dUTP conjugates, because dUTP inhibits archaeal polymerases such as Pfu and Vent.

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); 4) *Anal Biochem*. **211**,164(1993); 5) *PNAS* **99**,596(2002).

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.