

Product Information Sheet

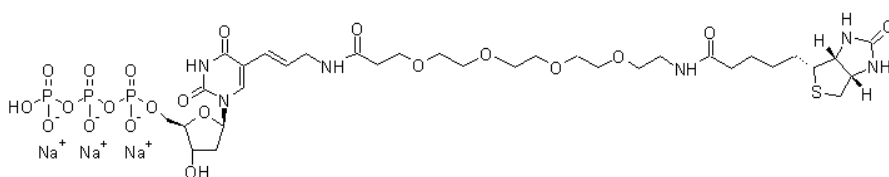
Ordering Information

Product Number:	17018
Product Name:	Biotin-20-dUTP *1 mM in Tris Buffer (pH 7.5)*
Unit Size:	25 nmoles
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 1062.75

Chemical Structure:



Soluble in:	Water
Excitation Wavelength:	NA
Emission Wavelength:	NA

Application Notes

The biotin-modified deoxyuridine 5'-triphosphates are widely used for a variety of non-radioactive DNA labeling reactions including nick translation, random prime labeling, cDNA labeling and 3'-end labeling. The biotinylated probes have been shown to hybridize to homologous nucleic acid at the same rate and to the same extent as non-biotinylated probes. The hybridized biotinylated DNA probes can be detected by avidin and streptavidin. Biotin-20-dUTP can be enzymatically incorporated into DNA via nick-translation, random priming, 3'-end terminal labeling or in the process of PCR. Biotin-20-dUTP is used to produce biotinylated DNA probes in a variety of hybridization applications including Southern blots, Northern blots, dot blots, fixed cells and tissues.