

Biotin-16-dUTP *1 mM in TE Buffer (pH 7.5)* *CAS 136632-31-0*

Catalog Number: 17017

Unit Size: 25 nmoles

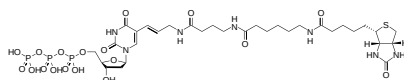
Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	6 months upon receiving

Chemical Properties

Appearance	Liquid colorless
Molecular Weight	947.78
Soluble In	Water

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

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-16-dUTP can be enzymatically incorporated into DNA via nick-translation, random priming, 3'-end terminal labeling or in the process of PCR. The number '16' is the number of carbon atoms in the backbone of the linker between dUTP and biotin. The longer the linker is, the more effective interaction of biotin with avidin occurs. On the other hand, the shorter the linker is, the more effective incorporation of dUTP into DNA. Biotin-16-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.