

Digoxigenin-11-dUTP *1 mM solution in TE Buffer*

Catalog Number: 17012

Unit Size: 25 nmoles

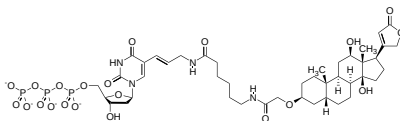
Product Details

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

Chemical Properties

Appearance	Liquid colorless
Molecular Weight	1066.92
Soluble In	Water

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Digoxigenin-11-dUTP can be enzymatically incorporated into DNA/cDNA as substitute for its natural counterpart dTTP. The resulting digoxigenin-labeled DNA/cDNA probes are subsequently detected using digoxigenin-antibodies conjugated with horseradish peroxidase (HRP), alkaline phosphatase (AP) or a fluorescent dye. The 11-atom linker is attached to the C5 position of uridine to minimize the interference of DIG on the substrate incorporation by PCR polymerase. Typically 1:2 Digoxigenin-11-dUTP/dTTP ratio for PCR and Nick Translation.