

Bio-16-UTP [Biotin-16-UTP] *1 mM*

Catalog Number: 17110, 17111

Unit Size: 25 nmoles, 100 nmoles

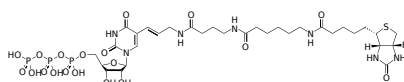
Product Details

| | |
|--------------------|--|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date | 12 months upon receiving |

Chemical Properties

| | |
|------------------|------------------|
| Appearance | Liquid colorless |
| Molecular Weight | 963.78 |
| Soluble In | Water |

Chemical Structure



Spectral Properties

| | |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength | N/A |

Applications

Bio-16-UTP is biotinylated UTP that can be used as a substrate in in vitro transcription reactions with a variety of RNA polymerases, including T7 RNA polymerase, SP6 RNA polymerase, or T3 RNA polymerase. Biotinylated RNA probes may be used in place of radioactively labeled RNA probes in many applications, including in situ hybridization and chromosome mapping. Bio-16-UTP can be detected using a variety of methods based on streptavidin binding. Biotin-bound streptavidin can be detected directly, using fluorescently labeled streptavidin, or indirectly, using a streptavidin-conjugated enzyme assay or appropriately labeled or conjugated anti-streptavidin antibody.