

ddGTP
[2',3'-Dideoxyguanosine-5'-triphosphate]
10 mM in ddH₂O

Catalog number: 17210
Unit size: 1 umole

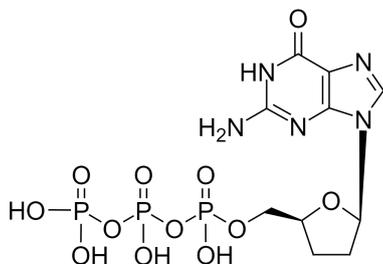
Product Details

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

Chemical Properties

| | |
|------------------|------------------|
| Appearance | Colorless Liquid |
| Molecular Weight | 491.18 |
| Soluble In | Water |

Chemical Structure



Applications

Sanger sequencing, also known as the chain termination method, is a technique for DNA sequencing based upon the selective incorporation of chain-terminating dideoxynucleotides (ddNTPs) by DNA polymerase. It was developed by Frederick Sanger and colleagues in 1977. Although the newer NGS technologies are becoming common in clinical research labs due to their higher throughput capabilities and lower costs per sample, Sanger sequencing with 99.99% accuracy is still the “gold standard” for clinical research sequencing. dd-GTP is one of the four critical ddNTP components for performing Sanger sequencing.