

## ReadiUse™ ddNTP Terminator Mix \*10 mM\*

Catalog number: 17205, 17206 Unit size: 100 nmoles, 10 umoles

**Product Details** 

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

**Chemical Properties** 

Appearance Liquid

Soluble In Water

## **Applications**

ReadiUse™ ddNTP Terminator Mix consists of 4 equimolar 2',3'-ddNTPs, including ddATP, ddGTP, ddTTP and ddCTP. They are provided in 100 mM solution and can be readily added to a sequencing solution or other biological samples. All the 2',3'-ddNTPs are highly purified.