



Oligo(dT)₂₀ Primer

Cat. No. 18418-020

Conc. 50 μ M

Size: 50 μ l

Store at -20°C

Description

Oligo(dT)₂₀ primer is a string of 20 deoxythymidylic acid residues that hybridizes to the poly(A) tail of mRNA and can be used as a primer for first-strand cDNA synthesis with reverse transcriptase. The primer is supplied in distilled water at a concentration of 50 μ M. The primer is phosphorylated at the 5' end.

Oligo(dT)₂₀ primer is recommended for use with SuperScript™ III RNase H⁻ Reverse Transcriptase (Cat. no. 18080-044), ThermoScript™ RNase H⁻ Reverse Transcriptase (Cat. no. 12236-014), or Cloned AMV Reverse Transcriptase (Cat. no. 12328-019).

We recommend using 1 μ l of oligo(dT)₂₀ primer per 20 μ l reaction volume.

Storage

Store at -20°C in a non-frost-free freezer.

Quality Control

This product is qualified in a first-strand cDNA reaction.

Part no. 18418020.pps

Rev. date.: 01/20/03

This product is distributed for laboratory research only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.