



Oligo(dT)₁₂₋₁₈ Primer

Cat. No. 18418-012

Conc.: 0.5 µg/µl

Size: 25 µg

Store at -20°C

Description

Oligo(dT)₁₂₋₁₈ Primer hybridizes to the poly(A) tail of mRNA and is suitable for use as a primer for first stand cDNA synthesis with reverse transcriptase. The primer is supplied in DEPC-treated water at a concentration of 0.5 µg/µl. The primer is phosphorylated at the 5' end.

Component

Oligo(dT)₁₂₋₁₈ Primer

Recommended Amount

We recommend using 0.5 µg of oligo(dT)₁₂₋₁₈ primer per µg of poly(A)⁺ RNA. For total RNA, we recommend 0.5 µg of primer for ≤ 5 µg of RNA.

Part no. 18418012.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.

Functional Assay Conditions

1 μ l oligo(dT)₁₂₋₁₈ primer (0.5 μ g/ μ l)

1.0 μ g poly(A)⁺ RNA

Sterile distilled water

Heat mixture to 70°C for 10 min, and quick chill on ice.

4 μ l 5X first strand buffer [250 mM Tris-HCl (pH 8.3), 375 mM KCl, 15 mM MgCl₂]

2 μ l 0.1 M DTT

1 μ l dNTPs (10 mM each dNTP)

1 μ l [α -³²P]dCTP (1 μ Ci/ μ l)

200 units SuperScriptTM RT (Cat. No. 18053-017) or M-MLV RT (Cat. No. 28025-013/28025-021)

Total volume: 20 μ l

Incubate at 37°C for 1 h.

Quality Control

This product is qualified in a first strand cDNA reaction.