

Anchored Oligo (dT)₂₀ Primer

Oligo (dT) Primers

Primers and Oligonucleotides

	Cat. No.	Size	Conc.
	PM-305S	200 µl	100 µM
	PM-305L	1 ml	100 µM

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C, avoid frequent thawing and freezing

Description

Anchored Oligo (dT) Primer hybridizes with highest efficiency to the poly(A) tail of mRNA and is recommended as primer for first stand cDNA synthesis, cDNA labeling or RT-PCR applications.

Sequence

5'-TTT TTT TTT TTT TTT TTT TTV N-3'

V: dA, dC, or dG

N: dA, dC, dG or dT

Molecular weight

6639

Package size and conversion

Anchored Oligo (dT) ₂₀ Primer			
PM-305S	200 µl / 100 µM	20 nmol	133 µg
PM-305L	1 ml / 100 µM	100 nmol	664 µg

Anchored Oligo (dT)₂₀ Primer (white cap)

100 µM in aqueous solution