

**Oligo (dT)₂₀**

Cat. No.	Amount
PM-304S	200 µl (20 nmol, 120 µg)
PM-304L	5 x 200 µl (100 nmol, 602 µg)

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 6022 g/mol

Purity: by HPLC

Form: clear aqueous solution

Concentration: 100 µM

Description:

Oligo (dT) Primer hybridizes to the poly(A) tail of mRNA and is used as primer for first stand cDNA synthesis with reverse transcriptases.

Sequence:

5'- TTT TTT TTT TTT TTT TTT TT -3'