

T7-Terminator

Sequencing Primer

Cat.-No.	Size	Conc.
PM-203	200 µl	10 µM

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

T7-Terminator, 2 nmol
HPLC purified, supplied as 10 µM aqueous solution

Sequence

5'-GCT AGT TAT TGC TCA GCG G-3'
19-mer

Description

The primer can be used for sequencing plasmids containing T7-Terminator sequences.