

M13/pUC forward

Sequencing Primer

Cat.-No.	Size	Conc.
PM-206	200 µl	10 µM

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

M13/pUC forward, 17-mer (-20), 2 nmol
HPLC purified, supplied as 10 µM aqueous solution

Sequence

5'-GTA AAA CGA CGG CCA GT-3'
17-mer

Description

The primer can be used for sequencing plasmids containing M13 or pUC forward sequences.