

M13/pUC reverse

Sequencing Primer

Cat.-No.	Size	Conc.
PM-207	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 reverse, 19-mer (-24), 2 nmol
HPLC purified, supplied as 10 µM aqueous solution

Sequence
5'-GGA AAC AGC TAT GAC CAT G-3'
19-mer

Description

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