

T3-Promoter

Sequencing Primer

Cat.-No.	Size	Conc.
PM-201	200 µl	10 µM

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

T3-Promoter, 2 nmol
HPLC purified, supplied as 10 µM aqueous solution

Sequence

5'-ATT AAC CCT CAC TAA AG-3'
17-mer

Description

The primer can be used for sequencing plasmids containing T3 RNA polymerase promoter sequences. It is complementary to the consensus sequence of the eleven class III promoter sequences for T3 RNA polymerase [1].

Note that primers cannot be used for certain plasmids that contain truncated promoters.

[1] McGraw, N. J. et al., Nucl. Acids Res. (1985) 13, 6753-6766