

# T7-Promoter

## Sequencing Primer

---

Cat.-No.	Size	Conc.
PM-202	200 µl	10 µM

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

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

### Sequence

5'-TAA TAC GAG TCA CTA TAG GGA-3'  
21-mer

### Description

The primer can be used for sequencing plasmids containing T7 RNA polymerase promoter sequences. It is complementary to conserved sequence in all of the class III promoters for T7 RNA polymerase [1].

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

[1] Dunn J. J. et al., J. Mol. Biol. (1983) 166, 477-535