

## FAQs for DNA End Modification: DNA A-Tailing

### 1. What enzymes can be used for A-Tailing?

Taq DNA Polymerase and Klenow exo-, are good choices for this application.

### 2. Does the DNA need to be purified after the PCR amplification and prior to the tailing reaction?

If the amplification has been carried out with a high fidelity polymerase, it is essential that the DNA be purified prior to the tailing reaction. The presence of any active high fidelity polymerase will result in the removal of any non-templated nucleotides added.