

## Properties of Thermostable DNA Polymerases

### Properties of Thermostable DNA Polymerases.

Organism	<i>Taq*</i> ( <i>Thermus aquaticus</i> )	<i>Tfl</i> ( <i>Thermus flavus</i> )	<i>Tth</i> ( <i>Thermus thermophilus</i> )	<i>Tli</i> ( <i>Thermococcus litoralis</i> )	<i>Pfu**</i> ( <i>Pyrococcus furiosus</i> )
MW	80kDa	94kDa	94kDa	90kDa	92kDa
Extension Temperature	74°C	74°C	74°C	74°C	75°C
5'→3' Exonuclease Activity	Yes	Yes	Yes	No	No
3'→5' Exonuclease Activity	No	No	No	Yes	Yes
Reverse Transcriptase Activity	Weak	Yes	Yes	No	N/A
Predominant PCR Product Ends	3'-A	3'-A	3'-A	70% Blunt; 30% Single-base overhangs	Blunt

\* GoTaq® DNA Polymerase and PCR Master Mix contain *Taq* DNA polymerase.

\*\* Not available in North America.

N/A = Not available.

### Reference

Newton, C.R. and Graham, A. (1994) In: *PCR, BIOS Scientific Publishers, Ltd., Oxford, UK.* 13.