

# Taq DNA Polymerase (1000 U) with 10 mM dNTP Mix (0.5 mL)

Version: 201807

**Cat.No.E00101**

**Size: 1PK**

	volume	quantity
Taq DNA Polymerase	0.2mL	1
10×Taq buffer	1.5mL	2
10 mM dNTP	0.5mL	1
15mM MgCl <sub>2</sub>	1.5mL	1

### Description

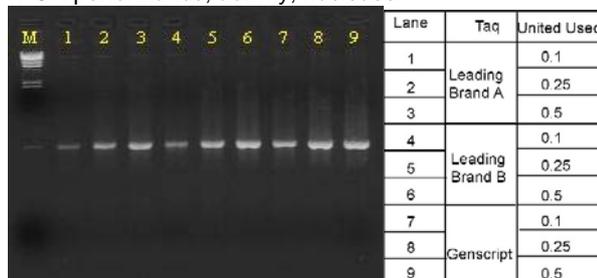
The Product contain 1000 U Taq DNA Polymerase (Cat. No. E00007) and 0.5 mL of 10 mM dNTP (Cat. No. D0056). The Taq DNA Polymerase is supplied with 10X Taq buffer (Cat. No. B0005) containing 15 mM magnesium chloride.

### Storage Buffer and Concentration

Supplied in 5 units/μL in 20 mM Tris HCl (pH 8.0), 0.1 mM EDTA, 1 mM DTT, 0.1% Triton X-100 and 50% glycerol.

### QC Tests

PCR performance, activity, nuclease.



\* The PCR process is covered by US. Patents Nos. 4683195 and 4683202, issued to Cetus and owned by Hoffman-La Roche Inc. GenScript does not encourage or support the unauthorized use of the PCR process. Use of this product is recommended for persons who either have a license to perform PCR or are not required to obtain a license. Sale of this product is restricted to regions or countries where native Taq DNA polymerase patents have been invalidated.