



Phi29 DNA Polymerase

Bacillus subtilis phage phi29, recombinant, *E. coli*

Cat. No.	Amount
PCR-381S	200 units
PCR-381L	1.000 units

Unit Definition: One unit is defined as the amount of the enzyme required to catalyze the incorporation of 10 nmol of dNTP into an acid-insoluble form in 30 minutes at 74 °C.

For *in vitro* use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 6 months

Form: liquid

Concentration: 10 units/μl

Applications:

- Rolling-circle amplification (RCA)
- Multiple displacement amplification (MDA)
- Whole genome amplification (WGA)
- Preparation of DNA template for sequencing

Description:

Phi29 DNA polymerase is a recombinant protein purified from *E. coli* cloned the gene encoding the DNA polymerase from *Phi29* phage. *Phi29* DNA polymerase is the replicative polymerase from the *Bacillus subtilis* phage *Phi29* and possesses the highest processivity and strand-displacement activity among the known DNA polymerase. *Phi29* DNA polymerase contains a 3'→5' exonuclease activity that enables proofreading capability.

Content:

Phi29 DNA Polymerase (brown cap)

10 units/μl *Phi29* DNA Polymerase in storage buffer (25 mM NaH₂PO₄ pH 7.0, 150 mM NaCl, 125 mM Imidazole, 50 % Glycerol, 2.5 mM 2-Mercaptoethanol)

Reaction Buffer (purple cap)

10x conc. complete reaction buffer containing 500 mM Tris-HCl pH 7.5, 100 mM MgCl₂, 100 mM (NH₂)₄SO₄ and 40 mM DTT

component	PCR-381S	PCR-381L
<i>Phi29</i> DNA Poly-merase	20 μl	100 μl
Reaction Buffer	100 μl	500 μl



Phi29 DNA Polymerase

Bacillus subtilis phage phi29, recombinant, *E. coli*

Recommended Assay Set-Up

component	Cat. No.	stock conc.	final conc.	1 assay @ 20 µl
PCR-grade Water	PCR-258			fill up to 20 µl
Reaction Buffer	PCR-381 (purple cap)	10x	1x	2 µl
dNTP Mix	NU-1006	10 mM	125 µM	0.25 µl
Random Hexamers	PM-301	100 µM	5 µM	1 µl
Target DNA				1 µl
Phi29 DNA Polymerase	PCR-381 (brown cap)	10 units/µl	5-10 units/assay	5-10 units

Incubation:

Incubate at 30 °C.

Inactivation:

Heat the mixture to 65 °C for 15 min.

Related Products:

Direct WGA Kit, #PCR-382

Phi29 WGA Kit, #APP-004