

	Package contents	Catalog number EP150B005	Size 5000 units	Kit contents
	Storage conditions	<ul style="list-style-type: none"> Store all contents at -20°C. 		
	Required materials	<ul style="list-style-type: none"> Template: cDNA, gDNA, λDNA Forward and reverse gene-specific primers Invitrogen™ 10 mM dNTP mix (Cat. no. 18427-088) 100 μM primer probe Water, nuclease-free Invitrogen™ ROX Reference Dye (Cat. no. 12223-012) 0.2 or 0.5-mL nuclease-free microcentrifuge tubes 		
	Timing	Varies depending on amplicon length		
	Product description	<ul style="list-style-type: none"> Low Glycerol LibertyTaq™ DNA Polymerase is a recombinant <i>Taq</i> polymerase complexed with a proprietary hot-start component that blocks polymerase activity at ambient temperatures. Activity is restored past 57°C, providing an automatic “hot start” and offering increased sensitivity, specificity, and yield, while allowing reaction assembly at room temperature. This enzyme has a non-template-dependent, terminal transferase activity that adds a single deoxyadenosine (A) to the 3' ends of PCR products. Like standard <i>Taq</i>, it has both 5' to 3' polymerase and 5' to 3' exonuclease activity, but lacks 3' to 5' exonuclease activity. Low Glycerol LibertyTaq™ DNA Polymerase is supplied at a concentration of 50 U/μL with minimal glycerol content, rendering it feasible for lyophilization applications. The Low Glycerol LibertyTaq™ DNA Polymerase is designed for qPCR applications. 		
	Important guidelines	Click here for important qPCR guidelines.		
	Online resources	For further information, contact outlicensing@lifetech.com .		

Enzyme characteristics

Hot-start:	Proprietary
Fidelity vs. <i>Taq</i>:	1X
Format:	Separate components

qPCR setup

Use the measurements below to prepare your qPCR experiment, or enter your own parameters in the column provided.

Component	25- μL rxn	Custom	Final conc. in 25- μL rxn
Water, nuclease-free	to 25 μL	to μL	—
20X Liberty PCR Buffer (– MgCl_2)	1.25 μL	μL	1X
50 mM MgCl_2	0.75 μL	μL	1.5 mM
10 mM dNTP mix	0.5 μL	μL	0.2 mM each
10 μM forward primer	0.75 μL	μL	0.3 μM
10 μM reverse primer	0.75 μL	μL	0.3 μM
100 μM primer probe	0.05 μL	μL	0.2 μM
30 μM ROX Reference Dye	0.025 μL	μL	30 nM
Template DNA	varies	μL	≤ 250 ng/rxn
Low Glycerol LibertyTaq™ DNA Polymerase (10 U/ μL) ¹	0.1 μL	μL	1 U/rxn

¹ Prior to use, dilute the Low Glycerol LibertyTaq™ DNA Polymerase to at least 10 U/ μL with Low Glycerol LibertyTaq™ Diluent included in the kit.

qPCR protocol

See pages 2 and 3 for instructions to prepare and run your qPCR experiment.


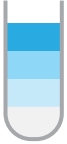

Optimization strategies


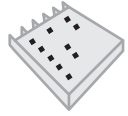
Click here for guidelines to optimize your qPCR experiment.

Purchaser notification

Click here for Limited Warranty, Disclaimer, and Licensing information.

The example qPCR procedure below shows appropriate volumes for a single **25- μ L** reaction. For multiple reactions, prepare a master mix of components common to all reactions to minimize pipetting error, and then dispense appropriate volumes into each 0.2–0.5 mL qPCR reaction tube or well of a MicroAmp™ EnduraPlate™ Optical 96- or 384-well plate prior to adding template DNA and primers. For 384-well plates, we recommend a maximum reaction volume of 10 μ L per well.

Steps	Action	Procedure details																																
1 	Thaw reagents	<p>Thaw, mix, and briefly centrifuge each component before use.</p> <p>Prior to use, dilute the Low Glycerol LibertyTaq™ DNA Polymerase (50 U/μL) to at least 10 U/μL in Low Glycerol LibertyTaq™ Diluent included in the kit.</p> <p>Note: After dilution, the Low Glycerol LibertyTaq™ DNA Polymerase (at 10 U/μL) can be stored at 4°C for use within 1–2 days. Do not store the diluted enzyme at –20°C.</p>																																
2 	Prepare qPCR master mix	<p>Add the following components to each qPCR reaction tube.</p> <p>Note: Consider the volumes for all components listed in steps 2 and 3 to determine the correct amount of water required to reach your final reaction volume.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>25-μL rxn</th> <th>Custom</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>Water, nuclease-free</td> <td>to 25 μL</td> <td>μL</td> <td>—</td> </tr> <tr> <td>20X Liberty PCR Buffer (– MgCl₂)</td> <td>1.25 μL</td> <td>μL</td> <td>1X</td> </tr> <tr> <td>50 mM MgCl₂</td> <td>0.75 μL</td> <td>μL</td> <td>1.5 mM</td> </tr> <tr> <td>10 mM dNTP mix</td> <td>0.5 μL</td> <td>μL</td> <td>0.2 mM each</td> </tr> <tr> <td>100 μM primer probe</td> <td>0.05 μL</td> <td>μL</td> <td>0.2 μM</td> </tr> <tr> <td>30 μM ROX Reference Dye</td> <td>0.025 μL</td> <td>μL</td> <td>30 nM</td> </tr> <tr> <td>Low Glycerol LibertyTaq™ DNA Polymerase (10 U/μL)</td> <td>0.1 μL</td> <td>μL</td> <td>1 U/rxn</td> </tr> </tbody> </table> <p>Mix and then briefly centrifuge the components.</p>	Component	25- μ L rxn	Custom	Final conc.	Water, nuclease-free	to 25 μ L	μ L	—	20X Liberty PCR Buffer (– MgCl ₂)	1.25 μ L	μ L	1X	50 mM MgCl ₂	0.75 μ L	μ L	1.5 mM	10 mM dNTP mix	0.5 μ L	μ L	0.2 mM each	100 μ M primer probe	0.05 μ L	μ L	0.2 μ M	30 μ M ROX Reference Dye	0.025 μ L	μ L	30 nM	Low Glycerol LibertyTaq™ DNA Polymerase (10 U/ μ L)	0.1 μ L	μ L	1 U/rxn
Component	25- μ L rxn	Custom	Final conc.																															
Water, nuclease-free	to 25 μ L	μ L	—																															
20X Liberty PCR Buffer (– MgCl ₂)	1.25 μ L	μ L	1X																															
50 mM MgCl ₂	0.75 μ L	μ L	1.5 mM																															
10 mM dNTP mix	0.5 μ L	μ L	0.2 mM each																															
100 μ M primer probe	0.05 μ L	μ L	0.2 μ M																															
30 μ M ROX Reference Dye	0.025 μ L	μ L	30 nM																															
Low Glycerol LibertyTaq™ DNA Polymerase (10 U/ μ L)	0.1 μ L	μ L	1 U/rxn																															
3 	Add template DNA and primers	<p>Add your template DNA and primers to each tube for a final reaction volume of 25 μL.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>25-μL rxn</th> <th>Custom</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>10 μM forward gene-specific primer</td> <td>0.75 μL</td> <td>μL</td> <td>0.3 μM</td> </tr> <tr> <td>10 μM reverse gene-specific primer</td> <td>0.75 μL</td> <td>μL</td> <td>0.3 μM</td> </tr> <tr> <td>Template DNA</td> <td>varies</td> <td>μL</td> <td>\leq250 ng/rxn (human gDNA)¹</td> </tr> </tbody> </table> <p>¹ See “Important guidelines”, page 1.</p> <p>Cap each tube, mix, and then briefly centrifuge the contents.</p>	Component	25- μ L rxn	Custom	Final conc.	10 μ M forward gene-specific primer	0.75 μ L	μ L	0.3 μ M	10 μ M reverse gene-specific primer	0.75 μ L	μ L	0.3 μ M	Template DNA	varies	μ L	\leq 250 ng/rxn (human gDNA) ¹																
Component	25- μ L rxn	Custom	Final conc.																															
10 μ M forward gene-specific primer	0.75 μ L	μ L	0.3 μ M																															
10 μ M reverse gene-specific primer	0.75 μ L	μ L	0.3 μ M																															
Template DNA	varies	μ L	\leq 250 ng/rxn (human gDNA) ¹																															

Steps	Action	Procedure details																											
4	 <p>Incubate reactions in a thermal cycler</p>	<p>qPCR cycling parameters (Two-step cycling protocol):</p> <table border="1"> <thead> <tr> <th>Step</th> <th>Temperature</th> <th>Time</th> </tr> </thead> <tbody> <tr> <td>Initial denaturation</td> <td>95°C</td> <td>5 minutes</td> </tr> <tr> <td rowspan="2">40 qPCR cycles</td> <td>Denature</td> <td>95°C</td> </tr> <tr> <td>Anneal/Extend¹</td> <td>~60°C (depending on primer T_m)</td> <td>1 minute</td> </tr> </tbody> </table> <p>¹ Data acquisition should be performed during the annealing/extension step for probe-based assays.</p> <p>qPCR cycling parameters (Three-step cycling protocol):</p> <table border="1"> <thead> <tr> <th>Step</th> <th>Temperature</th> <th>Time</th> </tr> </thead> <tbody> <tr> <td>Initial denaturation</td> <td>95°C</td> <td>5 minutes</td> </tr> <tr> <td rowspan="3">40 qPCR cycles</td> <td>Denature</td> <td>95°C</td> </tr> <tr> <td>Anneal</td> <td>~60°C (depending on primer T_m)</td> <td>30 seconds</td> </tr> <tr> <td>Extend¹</td> <td>72°C</td> <td>30 seconds</td> </tr> </tbody> </table> <p>¹ Data acquisition should be performed during the extension step for probe-based assays.</p>	Step	Temperature	Time	Initial denaturation	95°C	5 minutes	40 qPCR cycles	Denature	95°C	Anneal/Extend ¹	~60°C (depending on primer T _m)	1 minute	Step	Temperature	Time	Initial denaturation	95°C	5 minutes	40 qPCR cycles	Denature	95°C	Anneal	~60°C (depending on primer T _m)	30 seconds	Extend ¹	72°C	30 seconds
Step	Temperature	Time																											
Initial denaturation	95°C	5 minutes																											
40 qPCR cycles	Denature	95°C																											
	Anneal/Extend ¹	~60°C (depending on primer T _m)	1 minute																										
Step	Temperature	Time																											
Initial denaturation	95°C	5 minutes																											
40 qPCR cycles	Denature	95°C																											
	Anneal	~60°C (depending on primer T _m)	30 seconds																										
	Extend ¹	72°C	30 seconds																										
5	 <p>Analyze results</p>	<ul style="list-style-type: none"> Analyze results following your real-time instrument manufacturer's guidelines. You can store your samples overnight at 2–8°C, or at –20°C for longer storage. 																											

References

SantaLucia, Jr, J. (1998) A unified view of polymer, dumbbell and oligonucleotide DNA nearest-neighbor thermodynamics. Proc. Natl Acad. Sci., 95, 1460–1465.
 SantaLucia, Jr, J. and Hicks, D. (2004) The thermodynamics of DNA structural motifs. Annu. Rev. Biophys. Biomol. Struct., 33, 415–440.