

## NovaTaq™ Hot Start DNA Polymerase Kits

### Table of Contents

About the Kits.....	2
Description	2
Components	2
Storage	2
Standard Protocols.....	3
Standard PCR with NovaTaq Hot Start DNA Polymerase	3
Using NovaTaq™ Hot Start Master Mix	3

© 2011 EMD Chemicals, Inc., an affiliate of Merck KGaA, Darmstadt, Germany. All rights reserved. Perfectly Blunt® and Novagen® are registered trademarks of Merck KGaA, Darmstadt, Germany. AccepTOr™ and NovaTaq™ are trademarks of Merck KGaA, Darmstadt, Germany. TWEEN® is a registered trademark of ICI Americas, Inc.

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 6,127,155, 5,677,152 (claims 1 to 23 only), 5,773,258 (claims 1 and 6 only), and claims outside the US corresponding to expired US Patent 5,079,352. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California, 94404, USA

#### USA and Canada

Tel (800) 628-8470  
bioscienceshelp@  
emdchemicals.com

#### Europe

**France**  
Freephone  
0800 126 461

**Germany**  
Freecall  
0800 100 3496

**Ireland**  
Toll Free  
1800 409 445

**United Kingdom**  
Freephone  
0800 622 935

**All other  
European Countries**  
+44 115 943 0840

#### All Other Countries

**Contact Your Local Distributor**  
www.merck4biosciences.com  
bioscienceshelp@  
emdchemicals.com

techservice@merckbio.eu

www.merck4biosciences.com

FOR RESEARCH USE ONLY. NOT FOR HUMAN OR DIAGNOSTIC USE.

## About the Kits

NovaTaq Hot Start DNA Polymerase	250 U	71091-3
	1250 U	71091-4
NovaTaq Hot Start Master Mix Kit	200 rxn	71676-3
	1000 rxn	71676-4

## Description

NovaTaq Hot Start DNA Polymerase is a chemically modified form of *Taq* DNA Polymerase that is inactive at room temperature. The polymerase is a premium quality recombinant form of *Thermus aquaticus* DNA polymerase. The enzyme possesses 5'→3' DNA polymerase activity and lacks 3'→5' exonuclease activity. The enzyme preparation is > 95% homogenous by SDS-PAGE and lacks detectable RNase and endonuclease activities. The enzyme provides improved specificity when compared to standard *Taq* polymerase and can eliminate the presence of nonspecific amplification such as primer-dimers and misprimed products. The enzyme must be activated by heat treatment (7–10 min at 95°C), after which thermal cycling can proceed. NovaTaq Hot Start DNA Polymerase generates PCR products with 3'-dA overhangs, suitable for cloning with Novagen® Perfectly Blunt®, AccepTor™, and LIC Vector Kits.

NovaTaq Hot Start Master Mix Kit is a ready-to-use 2X mixture of NovaTaq Hot Start DNA Polymerase, ultrapure deoxynucleotides, and reaction buffer with MgCl<sub>2</sub>. The Master Mix simplifies the assembly of PCR reactions and offers advantages of time savings, consistency, and minimal risk of contamination. In addition to the Master Mix, the kit includes PCR Grade Water and MgCl<sub>2</sub> for increasing Mg<sup>2+</sup> concentrations. Simply add the NovaTaq Hot Start Master Mix to an equal volume containing DNA template, primers, and, if desired, additional MgCl<sub>2</sub>. The final diluted reaction contains 1.25 U of NovaTaq Hot Start DNA Polymerase per 50 µl. The two NovaTaq Hot Start Master Mix Kit sizes provide sufficient components for 200 or 1000 standard 50 µl reactions.

**NovaTaq Hot Start Unit definition:** One unit is defined as the amount of enzyme that will catalyze the incorporation of 10 nmol of dNTP into acid-insoluble form in 30 min at 72°C in a reaction containing 25 mM TAPS (tris-[hydroxymethyl]-methyl-amino-propane-sulfonic acid, sodium salt), pH 9.3 at 25°C, 50 mM KCl, 2 mM MgCl<sub>2</sub>, 1 mM β-mercaptoethanol, 0.2 mM dATP, dGTP, and dTTP, 0.1 µM [α-<sup>32</sup>P] dCTP, and 12.5 µg activated salmon sperm DNA in a 50 µl volume.

## Components

### NovaTaq™ Hot Start DNA Polymerase

- 1 x or 5 x 250 U NovaTaq Hot Start DNA Polymerase (5 U/µl)
- 1 x or 5 x 1.5 ml 10X NovaTaq Hot Start Buffer (10X = 670 mM Tris-HCl, 160 mM (NH<sub>4</sub>)<sub>2</sub>SO<sub>4</sub>, 0.1% TWEEN®-20)
- 1 x or 5 x 1.5 ml 25 mM MgCl<sub>2</sub>

### NovaTaq Hot Start Master Mix Kit

- 4 x or 20 x 1.25 ml NovaTaq Hot Start Master Mix (includes 250 U NovaTaq Hot Start DNA Polymerase)
- 3 x or 11 x 2 ml PCR Grade Water
- 1 x or 3 x 1.5 ml 25 mM MgCl<sub>2</sub>

## Storage

Store all components in a constant-temperature freezer at –20°C.

## Standard Protocols

This standard PCR protocol will provide satisfactory amplification in many cases. Include a negative control reaction lacking a template, and a positive control reaction using a template known to amplify with the primers. The optimal concentrations of enzyme, MgCl<sub>2</sub>, template, and primers are determined empirically.

### Standard PCR with NovaTaq Hot Start DNA Polymerase

- For each 50 µl reaction, assemble the following in a 0.5 ml PCR tube just prior to use:

X µl	PCR Grade Water (to final volume of 50 µl)
1 µl	dNTP Mix (10 mM each dATP, dCTP, dGTP, dTTP)
1 µl	5' primer, ~5 pmol/µl
1 µl	3' primer, ~5 pmol/µl
5 µl	10X NovaTaq Hot Start Buffer
3–5 µl	25 mM MgCl <sub>2</sub>
0.25 µl	NovaTaq Hot Start DNA Polymerase (1.25 U)
1 µl	DNA template (typically 10–100 ng)
50 µl	total volume
  - Mix gently. If necessary, centrifuge briefly. Add 2 drops mineral oil (if appropriate). Cap tubes. Place tubes in thermal cycler.
  - Activate the NovaTaq Hot Start DNA Polymerase with a 7–10 min 95°C incubation followed by the following cycling program (or other optimized cycling program):

Denature	30 s at 94°C
Anneal	30 s at the proper annealing temperature
Extend	1 min at 72°C

Repeat for 30 cycles  
Final extension for 10 min at 72°C
- Note: Any cycling program can be used with NovaTaq Hot Start DNA Polymerase with the addition of the polymerase activation step for 7–10 min at 95°C.*
- To remove oil overlay and inactivate polymerase, add 100 µl chloroform, mix 30 s, and centrifuge for 1 min. The top aqueous phase (which may appear cloudy) contains the DNA products. Transfer aqueous phase to fresh tube. Load and run a 1% agarose gel containing 0.5 µg/ml ethidium bromide. Visualize bands under UV illumination.

### Using NovaTaq™ Hot Start Master Mix

The NovaTaq Hot Start Master Mix is a 2X mixture of NovaTaq Hot Start DNA Polymerase, dNTPs and NovaTaq Buffer with MgCl<sub>2</sub>. When using NovaTaq Hot Start Master Mix, assemble the components of each 50 µl reaction as follows:

25 µl	NovaTaq Hot Start Master Mix
1 µl	5' primer, ~5 pmol/µl
1 µl	3' primer, ~5 pmol/µl
X µl	PCR Grade Water (to final volume of 50 µl)
1 µl	DNA template (typically 10 ng)
Y µl	25 mM MgCl <sub>2</sub> (if desired, see note below)
50 µl	total volume

Proceed with the standard cycling conditions as outlined in the previous section.

*Note: The reaction conditions above will contain a final concentration of 1.5 mM MgCl<sub>2</sub>. For some applications, increased MgCl<sub>2</sub> concentrations may be optimal. Table 1 below shows the volume of 25 mM MgCl<sub>2</sub> to be added to each 50 µl reaction to achieve the desired final concentration.*

Final MgCl <sub>2</sub> Concentration	Volume of 25 mM MgCl <sub>2</sub> to add per 50 µl reaction
1.5 mM	0 µl
2.0 mM	1 µl
2.5 mM	2 µl