

GeneAmp[®] Gold RNA PCR Reagent Kit

Protocol



© Copyright 2001, 2010 Applied Biosystems

For Research Use Only. Not for use in diagnostic procedures.

All rights reserved.

Printed in the U.S.A.

Notice to Purchaser: Limited License

Applied Biosystems and MicroAmp are registered trademarks of Applied Biosystems or its subsidiaries in the U.S. and certain other countries. MultiScribe is a trademark of Applied Biosystems or its subsidiaries in the U.S. and certain other countries.

GeneAmp and AmpliTaq Gold are registered trademarks of Roche Molecular Systems, Inc.

All other trademarks are the sole property of their respective owners.

Applied Biosystems is committed to providing the world's leading technology and information for life scientists.

Contents

Introduction	1
RNA-PCR Kit Features	1
RT-PCR Schematic	4
Materials and Equipment	5
GeneAmp Gold RNA PCR Reagent Kit	5
Kit Components	5
Control RNA and Primers	7
Storage and Stability	7
Enzyme Characteristics	8
Materials Required But Not Supplied	9
Safety Information	10
Document User Attention Words	10
Ordering Kits and Reagents	10
About MSDSs	10
Ordering MSDSs	10
Safety	11
Preventing Contamination	12
Introduction	12
General PCR Practices	12
Reagent Optimization Guidelines	13
General Guidelines	13
Preparing RT and PCR Reactions	13
Reaction Tubes	13
Magnesium Chloride Concentration	14
Primer Guidelines	14
RNA Template	15
Primers for RT	16

Cycle Optimization Guidelines	17
RT Thermal Cycling Conditions	17
PCR Amplification Conditions	17
Protocol for One-Step RNA PCR	19
Overview	19
Preparation of Reagents	19
Reaction Mix Preparation	19
One-Step Reaction Mix Preparation	20
Thermal Cycling Parameters for One-Step RNA PCR	21
Protocol for Two-Step RNA PCR	22
Overview	22
Preparation of Reagents	22
Reaction Mix Preparation	22
Two-Step Reaction Mix Preparation	22
RT Reaction Mix Preparation	23
RT Thermal Cycling Parameters for Two-Step RNA PCR	24
PCR Reaction Mix Preparation	25
PCR Thermal Cycling Parameters for Two-Step RNA PCR	27
Appendix A. Additional Primer and PCR Product Information	28
Additional Primers and PCR Products from pAW109	28
pAW109 Transcribed RNA	29
Performance Characteristics	29
Appendix B. References	30
Appendix C. Technical Support	32
Contacting Technical Support	32
To Contact Technical Support by E-mail	32
Hours for Telephone Technical Support	32
To Contact Technical Support by Telephone or Fax	33
To Reach Technical Support Through the Internet	36
To Obtain Documents on Demand	37

Introduction

RNA-PCR Kit Features

The GeneAmp® Gold RNA PCR Reagent Kit (P/N 4308206) is designed for the reverse transcription (RT) and polymerase chain reaction (PCR) amplification of a specific target RNA from total RNA, messenger RNA (mRNA), or viral RNA. The kit has the following features:

- ◆ Incorporation of the “Hot Start” technique by using AmpliTaq Gold® DNA Polymerase, which results in an increase in sensitivity, specificity, and yield of PCR products
- ◆ The capability to perform a unique, coupled one-step RNA PCR amplification, with reverse transcription followed by amplification in a single tube with a single buffer
- ◆ The capability to perform a more traditional two-step RNA PCR amplification, with reverse transcription followed by a separate PCR amplification step

The kit includes all the essential reagents to perform the reverse transcription of RNA to complementary DNA (cDNA) and the subsequent cDNA amplification using a single reaction tube and a single reaction buffer. This one-step process combines the versatility of MultiScribe™ Reverse Transcriptase, derived from Murine Leukemia Virus (MuLV), with the unique properties of AmpliTaq Gold® DNA Polymerase.

The provided RT-PCR Reaction Buffer has been optimized for both efficient reverse transcription by MultiScribe Reverse Transcriptase and efficient PCR amplification by AmpliTaq Gold DNA Polymerase, enabling it to be used in either the coupled one-step procedure or the traditional two-step procedure.

Hot Start Technique

The Hot Start technique (Faloona *et al*, 1990; Chou *et al*, 1992) is a simple modification of the original PCR reaction whereby the amplification reaction is initiated at a temperature above the optimal primer annealing temperature. In conventional PCR amplification, active reaction components are exposed to suboptimal annealing temperatures resulting in unintended priming and subsequent formation of non-specific products.

Because the unintended annealing occurs at the beginning of the PCR reaction, the non-specific products are efficiently amplified throughout the remaining PCR cycles. This unintended amplification results in poor yield of the desired product reducing the sensitivity of the experiment both by decreasing the desired amplification signal and by obscuring the signal with a high background.

Conversely, when the Hot Start PCR method is used, primers bind only to their specific target and the polymerase activity is directed to only the desired target. This results in increased sensitivity and yield, and decreased non-specific background amplification.

AmpliTaq Gold DNA Polymerase is a chemically modified form of *Thermus aquaticus* (Taq) DNA polymerase that requires high temperature activation. The modified enzyme is provided in an inactive state and remains so throughout the reverse transcription step. When the sample is heated to 95 °C, the modifier is permanently released, regenerating the active enzyme and initiating PCR amplification of the cDNA generated in the RT reaction. This thermal activation or “Hot Start” property of AmpliTaq Gold DNA Polymerase simplifies the procedure and provides increased sensitivity, specificity, and yield over conventional PCR techniques.

One-Step RNA-PCR

The coupled one-step RNA-PCR procedure eliminates the requirement for multiple manipulations and reagent additions following the reverse transcription reaction. This one-step procedure reduces the risk of contamination.

The one-step reaction can utilize any of these primers for the reverse transcription reaction: a sequence-specific primer, random hexamers, or oligo d(T)₁₆. In the one-step procedure, use sequence-specific primers (downstream or reverse PCR primers) to achieve the highest specificity.

Two-Step RNA-PCR

The two-step RNA-PCR procedure, in which an initial reverse transcription reaction is followed by an independent PCR amplification, is useful when detecting multiple transcripts from a single cDNA reaction, or when storing a portion of the cDNA for later use.

The two-step reaction can utilize any of these primers for the reverse transcription reaction: a sequence-specific primer, random hexamers, or oligo d(T)₁₆. In the two-step procedure, use either random hexamers

or oligo d(T)₁₆ as the cDNA synthesis primers to achieve the highest specificity in an RNA PCR amplification.

The table below compares the one-step and two-step features.

Comparison of RNA-PCR methods

Method	Primers for cDNA synthesis	Features
One-step	Sequence-specific primer, oligo d(T) ₁₆ or random hexamers	<ul style="list-style-type: none"> ◆ Requires single Reaction Mix ◆ Reduces risk of contamination
Two-step	Sequence-specific primer, oligo d(T) ₁₆ or random hexamers	<ul style="list-style-type: none"> ◆ Requires two Reaction Mixes ◆ cDNA can be stored or used for multiple amplifications

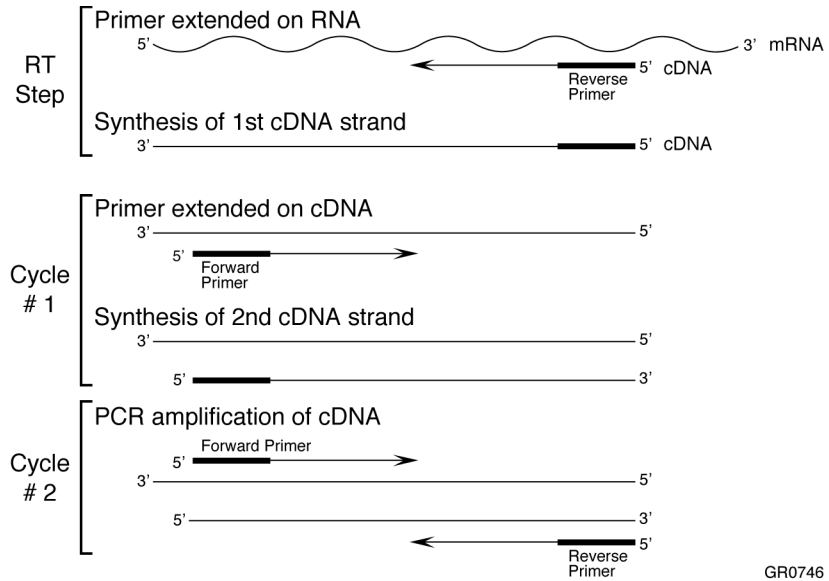
GeneAmp RNA PCR Control Kit

As a positive control, the kit contains a pAW109 complementary RNA (cRNA) template transcribed from the plasmid pAW109. The plasmid pAW109 contains an insert of a synthetic linear array of primer sequences for multiple target genes constructed such that “upstream” primers are followed by complementary sequences to their “downstream” primers in the same order.

The insert in pAW109 is identical to the insert previously described for the plasmid pAW108 (Wang *et al*, 1989) and is depicted in “A schematic diagram of the general arrangement of the RNA transcribed from plasmid pAW109.” on page 29.

The primers included in this kit flank an IL-1 α site and can be used to amplify a 308-bp sequence within the target. There is sufficient positive control pAW109 RNA and primers for at least 30 control RNA PCR amplifications.

RT-PCR Schematic As shown in the figure below, one-step RNA-PCR uses a single buffer that enables RT and PCR amplification to occur without interruption.



Schematic representation of RNA-PCR using the GeneAmp Gold RNA PCR Reagent Kit.

Materials and Equipment

GeneAmp Gold RNA PCR Reagent Kit The GeneAmp Gold RNA PCR Reagent Kit (P/N 4308206) is the complete kit providing materials to perform the reverse transcription of RNA and subsequent PCR amplification of samples plus control material (RNA and primers).

Kit Components The GeneAmp Gold RNA PCR Reagent Kit is comprised of two module kits that can be purchased separately:

Kit	P/N	Contents
GeneAmp Gold RNA PCR Reagent Kit	4308206	◆ GeneAmp® Gold RNA PCR Core Kit ◆ GeneAmp® Gold RNA PCR Control Kit ◆ Protocol (P/N 4310253)
GeneAmp Gold RNA PCR Core Kit	4312765	◆ RNA PCR Reagents ◆ Protocol (P/N 4310253)
GeneAmp Gold RNA PCR Control Kit	4308238	◆ pAW109 Control Reagents ◆ Protocol (P/N 4310253)

The contents of the modules are described in the tables on pages 6–7.

The GeneAmp Gold RNA PCR Reagent Kit and Core Kit each have a set of reagents sufficient to perform 100 RNA PCR amplifications of 50 μ L each. The GeneAmp Gold RNA PCR Control Kit (included in the Reagent Kit) has sufficient template and primers to perform at least 30 control reactions.

The following reagents are supplied with the GeneAmp Gold RNA PCR Core Kit:

The GeneAmp RNA PCR Core Kit reagents (P/N 4312765)

Reagent	Quantity (μL)	Description
MultiScribe™ Reverse Transcriptase	100	One tube containing 50 U/μL, MultiScribe (Murine Leukemia Virus) Reverse Transcriptase in DEPC water with 20 mM Tris-HCl, pH 7.5, 0.1 mM EDTA, 1 mM DTT, 0.01% Nonidet P40, 50% (v/v) glycerol.
RNase Inhibitor	100	One tube containing 20 U/μL RNase Inhibitor in DEPC water with 20 mM HEPES-KOH, pH 7.6, 50 mM KCL, 8 mM DTT, 50% (v/v) glycerol.
AmpliTaq Gold DNA Polymerase	50	One tube containing 5 U/μL of AmpliTaq Gold DNA Polymerase in 100 mM KCl, 20 mM Tris-HC, pH 9.0, 0.1 mM EDTA, 1 mM DTT, 0.5% Tween® 20, and 50% (v/v) glycerol.
dNTP Blend	1000	One tube containing 10 mM Deoxyribonucleotide triphosphates (2.5 mM each dATP, dCTP, dGTP, and dTTP) dissolved in DI water, titrated to pH 7.0 with NaOH.
100 mM DTT	500	One tube containing 100 mM Dithiothreitol in RNase free water.
5X RT-PCR Buffer	1500	One tube containing 150 mM Tris-HCl and 100 mM KCl, pH 8.3.
MgCl ₂ solution	1500	One tube containing 25 mM MgCl ₂ .
Random Hexamers	100	One tube containing 50 μM random primers for reverse transcription of RNA in 10 mM Tris-HCl, pH 8.3.
Oligo d(T) ₁₆	100	One tube containing 50 μM poly dT tailed primer for reverse transcription of RNA, in 10 mM Tris-HCL, pH 8.3. The sequence is: 5'-TTTTTTTTTTTTTTTT-3'.

The following reagents are supplied with the GeneAmp RNA PCR Control Kit.

The GeneAmp RNA PCR Control Kit reagents (P/N 4308238)

Reagent	Quantity (μL)	Description
pAW109 Control RNA	50	One tube containing 10 ⁶ copies/μL of RNA transcribed from pAW109 in 30 μg/mL <i>E. coli</i> rRNA, dissolved in 10 mM Tris-HCl, pH 7.0, 1 mM EDTA, and 100 mM NaCl.
Primer DM151	50	One tube containing 15 μM “upstream” IL-1α primer for amplification of pAW109 RNA, supplied in 10 mM Tris-HCl, pH 8.3. The sequence is: 5′-GTCTCTGAATCAGAAATCCTTCTATC-3′.
Primer DM152	50	One tube containing 15 μM “downstream” IL-1α primer for reverse transcription and amplification of pAW109 RNA, supplied in 10 mM Tris-HCl, pH 8.3. The sequence is: 5′-CATGTCAAATTTCACTGCTTCATCC-3′.

Control RNA and Primers The DM151 and DM152 primer pair amplifies a 308-bp product from the control pAW109 provided in the kit. The primer set amplifies a 421-bp product from authentic IL-1α mRNA. Refer to “Additional Primer and PCR Product Information” on page 28 for more information.

Storage and Stability Upon receipt, store the GeneAmp Gold RNA PCR Kit at -15 to -25 °C in a constant-temperature freezer. If stored under the recommended conditions, the product will maintain performance through the control date printed on the label.

Enzyme AmpliTaq Gold DNA Polymerase

Characteristics AmpliTaq Gold DNA Polymerase is a chemically modified form of AmpliTaq DNA Polymerase. This thermostable 94 kDa protein is encoded by a modified form of a *Thermus aquaticus* DNA polymerase gene and expressed in an *Escherichia coli* host (Lawyer, *et. al*, 1989).

Characteristics of AmpliTaq Gold DNA Polymerase

Item	Description
Unit Definition	The enzyme is provided at 5 Units/ μ L. One unit of enzyme will incorporate 10 nmoles of dNTPs into acid insoluble material per 30 min in a 10 min incubation at 74 °C under conditions listed under Analysis Conditions. Note The enzyme is shipped in an inactive form. Before analyzing, the enzyme is activated by heating for three hours at 80 °C.
Analysis Conditions	The analysis conditions are in the presence of 25 mM TAPS (tris-hydroxymethyl-methyl-amino-propanesulfonic) pH 9.3, 50 mM KCl, 2 mM MgCl ₂ , 1 mM β -mercaptoethanol, 200 μ M each of dATP, dGTP, dTTP, 100 μ M (α - ³² P)-dCTP (0.05–0.1 Ci/mmol), salmon sperm DNA activated by a modification of methods in Richardson <i>et al</i> , 1966; mixed in a final volume of 50 μ L and incubated at 74 °C for 10 min.
Storage Buffer	The storage buffer contains 20 mM Tris-HCl, pH 9.0 (room temperature), 100 mM KCl, 0.1 mM EDTA, 1.0 mM DTT (dithiothreitol), 0.5% Tween® 20, 50% (v/v) glycerol.
Storage Conditions	Store at –20 °C. If stored under proper conditions, the enzyme will remain active through the control date on the label.
Associated Activities	Non-specific endonuclease and exonuclease activities were not detectable after a 1 hour incubation of 600 ng of supercoiled pBR322 (dam ⁻ dcm ⁻) or 600 ng of <i>Msp</i> I-digested pBR322 DNA. The incubations were performed at 74 °C in the presence of 8 units of AmpliTaq Gold DNA polymerase. The enzyme has a fork-like structure-dependent polymerization enhanced 5' to 3' nuclease activity. The enzyme lacks a 3' to 5' exonuclease activity.

MultiScribe Reverse Transcriptase

MultiScribe Reverse Transcriptase, a recombinant form of the Murine Leukemia Virus (MuLV), is an RNA-dependent DNA polymerase that uses single-stranded RNA as a template in the presence of a primer to synthesize a complementary DNA strand. This enzyme is optimal for first strand cDNA synthesis. It is used in this kit for its robust reverse

transcription capability and its compatibility with random hexamers, oligo (dT), and gene specific primers.

Characteristics of MultiScribe Reverse Transcriptase

Item	Description
MultiScribe Reverse Transcriptase	The enzyme is provided at 50 Units/ μ L. One unit of enzyme will incorporate 1 nmol of dTTP into acid-precipitable material in 10 min at 37 °C using poly rA•oligo d(T) ₁₂₋₁₈ as a template primer. The ratio of oligo d(T) to poly rA is 3:1.
Source	The enzyme is purified from <i>E.coli</i> expressing the <i>pol</i> gene of MuLV on a plasmid.
Storage Buffer	The buffer contains 20 mM Tris–HCl, pH 7.5, 0.1 mM EDTA, 1 mM DTT, 0.01% (w/v) Nonidet® P40, 50% (v/v) glycerol, in diethylpyrocarbonate (DEPC) treated deionized (DI) water.

Materials Required But Not Supplied

The items listed below are required in addition to the reagents supplied in the GeneAmp Gold RNA PCR Reagent Kit.

User-supplied reagents

Equipment Item	Source
Centrifuge	Major laboratory suppliers (MLS)
Disposable gloves	MLS
MicroAmp Disposables	Applied Biosystems
Microcentrifuge	MLS
Pipettors, positive-displacement or air-displacement	MLS
Pipette tips, with filter plugs	MLS
Vortexer	MLS
RNase-free water	MLS
GeneAmp PCR Instrument System	Applied Biosystems

Safety Information

Document User Attention Words Four user words appear in the text of all Applied Biosystems user documentation. Each word implies a particular level of observation or action as follows.

Note This word is used to call attention to information.

IMPORTANT This word calls attention to information that is necessary for correct operation of the instrument.

CAUTION This word informs the user that damage to the instrument could occur if the user does not comply with the information. It also indicates a potentially hazardous situation that could result in minor or moderate injury to the user.

! WARNING ! This word informs the user that serious physical injury to the user or other persons could occur if these required precautions are not taken.

Ordering Kits and Reagents To order additional kits and reagents, please contact Applied Biosystems at one of the regional offices listed in this protocol. Have the part number of the kit or reagent you are ordering available when ordering.

About MSDSs Material and Safety Data Sheets (MSDSs) provide information about physical characteristics, health hazards, safety precautions, first aid, spill clean-up, and disposal procedures. Read the MSDSs before handling reagents or interacting with the instrument.

Ordering MSDSs Material Safety Data Sheets (MSDSs) for hazardous chemicals manufactured by Applied Biosystems will accompany your first shipment. To receive additional copies of MSDSs at no extra cost, call Applied Biosystems at (800) 327-3002.

! WARNING ! Some of the chemicals used in this manual may be hazardous. Follow the safety precautions stated in the specific chemical warnings that appear throughout the protocol.

! WARNING ! Some of the chemicals referred to in this manual may not have been provided with your kit. If the chemicals are not provided, they are not manufactured or sold by Applied Biosystems. Please obtain the Material Safety Data Sheets from their manufacturers.

! WARNING ! Some of the chemicals in this kit are sensitizers. While sensitizers are not hazardous in nature, some individuals may be sensitive to individual compounds.

Safety ! WARNING ! CHEMICAL HAZARD. Certain chemicals in this kit may be hazardous and therefore require special handling. Do not store, handle, or work with any chemicals or hazardous materials unless you have received appropriate safety training and have read and understood the Material Safety Data Sheet. Comply with all federal, state, and local laws related to chemical storage, handling, and disposal.

Preventing Contamination

Introduction Due to the high throughput and repetitive nature of PCR, special laboratory practices are necessary in order to avoid false positive amplifications (Kwok and Higuchi, 1989). This is because of the capability for single DNA molecule amplification provided by the PCR process (Saiki *et al*, 1985; Mullis *et al*, 1987).

General PCR Practices Please follow these recommended procedures:

- ◆ Wear a clean lab coat (not previously worn while handling amplified PCR products or used during sample preparation) and clean gloves when preparing samples for PCR amplification.
- ◆ Change gloves whenever you suspect that they are contaminated.
- ◆ Maintain separate areas and dedicated equipment and supplies for sample preparation, for PCR setup, and for PCR amplification and analysis of PCR products.
- ◆ Never bring amplified PCR products into the PCR setup area.
- ◆ Open and close all sample tubes carefully. Try not to splash or spray PCR samples.
- ◆ Keep reactions and components capped as much as possible.
- ◆ Clean lab benches and equipment periodically with 10% bleach solution.

Reagent Optimization Guidelines

General Guidelines These guidelines have been developed from the experience obtained at Applied Biosystems in the design of RT and PCR reactions.

Preparing RT and PCR Reactions Please follow these guidelines:

- ◆ Gently mix (avoid generating bubbles), Multiscribe Reverse Transcriptase, RNase inhibitor, AmpliTaq Gold DNA Polymerase and other thawed reagents, prior to pipetting.
- ◆ Spin the reagents down before pipetting.
- ◆ Pipette the enzymes carefully and slowly as they are in a viscous, 50%-glycerol solution.
- ◆ Prepare Master Mixes containing water, buffers, dNTPs, magnesium chloride, and enzyme, for both reverse transcription and PCR amplification.
- ◆ Use any combination of water, primer, and experimental volumes as long as the final combined volume is 50 μ L.
- ◆ Refrain from adding additional EDTA, citrate, or other chelators with the primers or experimental sample. If the final concentration of chelator exceeds 0.11 mM, additional $MgCl_2$ may need to be added.

Reaction Tubes The reaction tube guidelines are listed below.

- ◆ Use MicroAmp[®] 96 well Reaction Plates or MicroAmp[®] Reaction Tubes for your GeneAmp[®] Applied Biosystems tubes provide the best heat transfer because they fit uniformly into the wells of GeneAmp PCR Systems.
 - ◆ Use sterile reaction tubes and pipette tips to avoid nuclease contaminated plastic.
-

Magnesium Chloride Concentration Too little or too much magnesium chloride can reduce reverse transcriptase and amplification efficiency or result in nonspecific products. The optimal magnesium chloride concentration must be determined empirically. The guidelines for determining the optimum magnesium chloride concentration for each primer set are listed below.

- ◆ Test concentrations of magnesium chloride in the range of 1.25–5 mM in 0.5-mM increments.
- ◆ Raise the magnesium chloride concentration in the reaction mix proportionately if the samples contain EDTA, citrate, or other chelators.
- ◆ Adjust the magnesium chloride concentration in parallel with significant changes in the concentration (higher or lower) of sample RNA or dNTPs. This will keep the free magnesium ion concentration constant.

Primer Guidelines Guidelines for designing and using primers are listed below.

- ◆ Sequence-specific primers for RNA and cDNA should be 15–30 bases in length.
 - ◆ The % G +C should be near 50%.
 - ◆ Purify primers by gel electrophoresis or HPLC ion-exchange chromatography.
 - ◆ Determine the optimum primer concentration empirically by testing concentrations ranging 0.1–1.0 μ M primer. In general concentrations ranging 0.15–0.5 μ M will work for most amplifications.
 - Too low a concentration will result in little or no PCR product
 - Too high a concentration will result in amplification of non-target sequences.
 - ◆ Primer sequences should not complement within themselves or to each other, particularly at the 3' ends.
-
-

RNA Template This kit is designed to reverse transcribe and amplify RNA targets from mRNA or total RNA. The detection sensitivity of this kit is 500 copies of RNA.

A cDNA synthesis of up to 5 kb in length can be achieved with a subsequent amplification of up to 4 kb in length.

The guidelines for RT/PCR amplification of an RNA template are listed below.

- ◆ Start with at least 500 copies, but less than 1 μg total of the RNA template, to ensure a signal after 43 cycles. The cycle number can be reduced if you start with more than 500 copies. For example, 10,000 copies of RNA typically can be amplified with 35 cycles.
 - ◆ Amplify the positive control, pAW109, using a range of 500–1,000,000 copies diluted in the presence of 0.1 $\mu\text{g}/\mu\text{L}$ of carrier RNA. Always make fresh working solutions from stock preparations of pAW109.
-
-

Primers for RT The following primers can be used for cDNA synthesis:

- ◆ Random hexamers
- ◆ Oligo d(T)₁₆
- ◆ Sequence-specific reverse primers

The choice of primers for reverse transcription is best made after experimentally evaluating all three priming systems. For short RNA sequences containing no hairpin loops, any of the three priming systems should work equally well. For longer RNA transcripts or sequences containing hairpin loops, consider the following guidelines:

Primers for two-step RT-PCR

Primers	Selection Guidelines
Random hexamers	Try first for use with long reverse transcripts or reverse transcripts containing hairpin loops Use to transcribe all RNA (rRNA, mRNA, and tRNA)
Sequence-specific reverse primer	Use to reverse transcribe RNA-containing complementary sequences only
Oligo d(T) ₁₆	Use to reverse transcribe only eukaryotic mRNAs and retroviruses with poly-A tails Avoid long mRNA transcripts or amplicons greater than two kilobases upstream

Cycle Optimization Guidelines

RT Thermal Cycling Conditions The reverse transcriptase incubation condition guidelines are listed below.

- ◆ Incubate reactions for 12 minutes at 42 °C when reverse transcribing nonstructured RNA that is less than 2 kb in length.
- ◆ Raise the incubation temperature to 45 °C for 12 minutes for highly-structured RNA.
- ◆ If trying to achieve long transcripts, keep the temperature at 42 °C and increase the incubation time up to 20 minutes.

Note The half-life of MultiScribe reverse transcriptase is 10 minutes at 42 °C.

PCR Amplification Conditions **Anneal-Extend Temperature**

The selection of 60 °C for a combined anneal-extend temperature is optimal for amplification of the pAW109 cDNA. Optimal conditions for other primer-template pairs may require raising or lowering the annealing temperature in the range of 37–65 °C, followed by an extension temperature of 70–72 °C. In general, higher annealing temperatures result in a more specific product (Saiki *et al*, 1988 and Rychlik *et al*, 1990). The following are recommended guidelines.

- ◆ Determine the optimal annealing temperature empirically by testing increments of 5 °C or less until reaching the maximum in specificity and product.
- ◆ Amplification extension temperatures are typically 70–72 °C. When the anneal temperature is above 60 °C, it may be optimal to use two temperatures rather than three temperatures. In this case, complete the extension at the anneal temperature. The extension at lower temperatures may require time optimization for such combined anneal-extend steps.

Note Computer programs designed to calculate primer melting temperatures (T_m) such as Primer Express Applications-Based Primer Design Software (P/N 402809) can assist you in narrowing the suggested range of annealing temperatures for empirical determination.

Anneal-Extend Time

The length of the target sequence will affect the required extension time. AmpliTaq Gold DNA Polymerase has an extension rate of 2000–4000 bases per minute at 70–72 °C.

Note Polymerase activity is significant between 55–85 °C (Innis *et al*, 1988).

Typically, for products < 1 kb in length, the annealing times are 15–30 seconds and the extension times are 30–60 seconds. Using a combined anneal-extend step, 1 minute is sufficient for both steps.

As the amount of DNA increases in later cycles, the number of AmpliTaq Gold DNA Polymerase molecules may become limiting for the extension time allotted. To compensate for this, we recommend the following guidelines:

- ◆ Increase the extend times in later cycles to maintain the efficiency of amplification.
 - ◆ Use the AUTO program for the GeneAmp PCR Systems 2400, 9600, and 9700.
-

Protocol for One-Step RNA PCR

Overview Note The following procedure is optimized for one-step RNA PCR using the GeneAmp Gold RNA PCR Reagent Kit pAW109 control reagents to generate the 308 bp amplification product.

To design RNA PCR experiments for your samples and primers, refer to “Reagent Optimization Guidelines” on page 13 and “Cycle Optimization Guidelines” on page 17.

Preparation of Reagents The kit is stored at –15 to –25 °C. Prior to use, thaw all reagents except the enzymes. When the reagents are thawed, keep them on ice. Keep the enzymes (AmpliTaq Gold DNA Polymerase, MultiScribe Reverse Transcriptase, and RNase Inhibitor) in a freezer until immediately prior to use. Mix kit components by vortexing (or gentle mixing for enzymes), and using a microcentrifuge, briefly spin down the tube contents. When finished with the kit return it to the –15 to –25 °C freezer.

Reaction Mix Preparation Preparing a Reaction Mix of RNA PCR components is recommended in order to increase the precision of the results. The use of a Reaction Mix reduces the number of reagent transfers and minimizes volume loss due to pipetting.

General steps for Reaction Mix preparation

Step	Action
1	Prepare Reaction Mix by combining all the nonenzymatic components listed in the appropriate table.
2	Mix the components by pipetting up and down.
3	Vortex briefly.
4	Add the enzymatic components listed for the appropriate Reaction Mix.
5	Mix the components by inverting the microcentrifuge tube or by vortexing gently.

**One-Step
Reaction Mix
Preparation**

The ingredients of a 50- μ L reaction, one-step reaction mix are listed in the table below. To make the one-step Reaction Mix, follow the instructions described in “Reaction Mix Preparation” on page 19.

IMPORTANT MultiScribe Reverse Transcriptase and RNase Inhibitor are sensitive to air oxidation. Keep these reagents at -20°C until just prior to use.

One-Step Reaction Mix Components

Component	Volume/Tube (μ L)	Final Concentration
pAW109 RNA ^a	1.0	10^6 copies
RNase-free water	26.7	–
5X RT-PCR Buffer	10.0	1X
25 mM Magnesium Chloride	3.5	1.75 mM
10 mM dNTP Blend	4.0	0.8 mM (200 μ M of each dNTP)
RNase Inhibitor (20 Units/ μ L)	0.5	10 Units/50 μ L
100 mM DTT	2.5	5.0 mM
15 μ M Gene Specific Upstream Primer DM151	0.5	0.15 μ M
15 μ M Gene Specific Downstream Primer DM152	0.5	0.15 μ M
AmpliTaq Gold DNA Polymerase (5.0 Units/ μ L)	0.5	2.5 Units/50 μ L
MultiScribe Reverse Transcriptase (50 Units/ μ L)	0.3	15 Units/50 μ L
Total	50.0	

a. The sensitivity of the Control Reagent Kit reagents will allow successful amplification with as little as 1,000 copies of pAW109 RNA (refer to “Performance Characteristics” on page 29). When testing the kit at low copy number, make dilutions in the presence of 0.1 μ g/ μ L of carrier (such as *E. coli* ribosomal or transfer) RNA.

Thermal Cycling Parameters for One-Step RNA PCR

The thermal cycling parameters are optimized for the GeneAmp Gold RNA PCR Reagent Kit pAW109 control reagents to generate a 308-bp amplification product. The parameters described here are only for GeneAmp PCR Systems 2400, 9600, and 9700. See thermal cycler manuals for details on operation.

Some thermal cyclers may require longer (~60 sec) denaturation times. Refer to the manufacturers operating manuals.

Note All GeneAmp PCR Systems display the “Calculated Sample Temperature.”

Cycle parameters for one-step RNA PCR with pAW109 control reagents

Step	Reverse Transcription ^a	AmpliTaq Gold Activation	PCR CYCLE (43 cycles) ^b		PCR (Final Step)
	HOLD	HOLD	Denature	Anneal/Extend	HOLD
Time	12 min	10 min	20 sec	60 sec	7 min
Temperature	42 °C	95 °C	94 °C	62 °C	72 °C

a. When using random hexamers or oligo d(T)₁₆ for the RT priming, perform a 10 min, 25 °C HOLD prior to the 12 min, 42 °C reverse transcription for primer annealing. For more information, refer to “Reagent Optimization Guidelines” on page 13 and “Cycle Optimization Guidelines” on page 17.

b. The cycle number can be reduced when amplifying more than 10,000 copies of target RNA. For more information refer to “Cycle Optimization Guidelines” on page 17.

Protocol for Two-Step RNA PCR

Overview **Note** The following procedure is optimized for two-step RNA PCR using the pAW109 control reagents to generate the 308-bp amplification product.

To design RNA PCR experiments for your samples and primers, refer to “Reagent Optimization Guidelines” on page 13 and “Cycle Optimization Guidelines” on page 17.

Preparation of Reagents The kit is stored at –15 to –25 °C. Prior to use, thaw all reagents except the enzymes. When the reagents are thawed, keep them on ice. Keep the enzymes (AmpliTaq Gold DNA Polymerase, MultiScribe Reverse Transcriptase, and RNase Inhibitor) in a freezer until immediately prior to use. Mix kit components by vortexing (or gentle mixing for enzymes), and using a microcentrifuge, briefly spin down the tube contents. When finished with the kit, return it to the –15 to –25 °C freezer.

Reaction Mix Preparation Preparing a reaction mix of RT and PCR components for multiple samples is recommended in order to increase the precision of the results. The use of a reaction mix reduces the number of reagent transfers and minimizes volume loss due to pipetting.

General steps for Reaction Mix preparation

Step	Action
1	Prepare reaction mix by combining all the nonenzymatic components listed in the appropriate table.
2	Mix the components by pipetting up and down.
3	Vortex briefly.
4	Add the enzymatic components listed for the appropriate Reaction Mix.
5	Mix the components by inverting the microcentrifuge tube or by vortexing gently.

Two-Step Reaction Mix Preparation The two-step RT-PCR reaction requires two reaction mixes:

- ◆ RT Reaction Mix
- ◆ PCR Reaction Mix

RT Reaction Mix Preparation

The ingredients of a 20 μL , RT Reaction Mix are listed in the table below. To make the RT Reaction Mix, follow the instructions described in “Reaction Mix Preparation” on page 22.

The reaction volume for the RT step can be varied from 10–100 μL . A 10- μL reaction volume is sufficient to reverse transcribe up to 200 ng of total RNA. Increase the reaction volume when >200 ng of total RNA is reverse transcribed. For quantitation, a larger RT reaction volume can improve accuracy. When changing the reaction volume, make sure the final concentration is consistent with that described in the table below.

IMPORTANT MultiScribe Reverse Transcriptase and RNase Inhibitor are sensitive to air oxidation. Keep these reagents at -20°C until just prior to use.

Two-Step RT Reaction Mix

Component	Volume/Tube μL	Final Concentration
pAW109 RNA ^a	1.0	10^6 copies
RNase-free water	7.7	–
5X RT-PCR Buffer	4.0	1X
25 mM Magnesium Chloride	2.0	2.5 mM
10 mM dNTP Blend	2.0	1 mM (250 μM of each dNTP)
RNase Inhibitor	0.5	10 Units/20 μL
100 mM DTT	2.0	10 mM
Random Hexamer ^b	0.5	1.25 μM
MultiScribe Reverse Transcriptase (50 Units/ μL)	0.3	15 Units/20 μL
Total	20.0	

a. The sensitivity of the Control Reagent Kit reagents allows amplification with as little as 1,000 copies of pAW109 RNA (refer to “Performance Characteristics” on page 29). When testing the kit at low copy number, make dilutions in the presence of 0.1 $\mu\text{g}/\mu\text{L}$ of carrier (such as *E.coli* ribosomal or transfer) RNA.

b. Random hexamers, oligo d(T)₁₆, or sequence-specific DM 152 downstream primer can be used for primers of cDNA synthesis. When using an oligo d(T)₁₆ primer, use 0.5 μL of primer in a 20 μL reaction. This ensures that the final concentration of oligo d(T)₁₆ in the Reaction Mix is 1.25 μM . When using DM 152 primer use 0.5 μL in a 20 μL reaction (0.375 μM final concentration).

**RT Thermal
Cycling
Parameters for
Two-Step RNA
PCR**

The thermal cycler parameters described here are optimized for the GeneAmp Gold RNA PCR Reagent Kit pAW109 control reagents to generate a 308-bp amplification product.

The parameters described here are only for the GeneAmp PCR Systems 2400, 9600, and 9700. See thermal cycler manuals for details on operation.

Note All GeneAmp PCR Systems display the Calculated Sample Temperature.

Cycling parameters for the RT step of the two-step process

Step	Hybridization ^a	Reverse Transcription ^b
	HOLD	HOLD
Time	10 min	12 min
Temperature	25 °C	42 °C

a. If using random hexamers or oligo d(T)₁₆ primers for first-strand cDNA synthesis, incubate the primers at 25 °C for ten minutes. If using a sequence-specific reverse primer, this step is not necessary.

b. For information on optimizing reverse transcriptase reactions refer to “Cycle Optimization Guidelines” on page 17.

PCR Reaction Mix Preparation There are two possible reaction mixes for the second cDNA amplification step, reaction mix one and reaction mix two. Use reaction mix one if you do not need to perform multiple PCRs or to archive cDNA. Use reaction mix two if you have multiple targets for PCR or if you need to archive a portion of the cDNA.

Reaction Mix One

Use all 20 μL of the first step (RT) and add the following for a 50 μL reaction:

Two-Step PCR Reaction Mix

Component	Volume/Tube (μL)	Final Concentration
RT reaction	20.0	–
RNase-free water	19.0	–
5X RT-PCR Buffer	6.0	1X
25 mM Magnesium Chloride	1.5	1.75 mM
10 mM dNTP Blend	2.0	0.8 mM (200 μM of each dNTP)
15 μM Gene Specific Upstream Primer DM151	0.5	0.15 μM
15 μM Gene Specific Downstream Primer DM152 ^a	0.5	0.15 μM
AmpliTaq Gold DNA Polymerase (5.0 Units/ μL)	0.5	2.5 Units/50 μL
Total	50.0	

a. Omit primer DM152 if it was included in the RT step as the reverse transcription primer and add 19.5 μL of RNase-free water instead of 19.0 μL .

Reaction Mix Two

Use an aliquot (3 μL) of the first step (RT) and add the following for a 50 μL reaction:

Two-Step PCR Reaction Mix (two)

Component	Volume/Tube (μL)	Final Concentration
RT reaction	3.0	–
RNase-free water	29.2	–
5X RT-PCR Buffer	9.4	1X
25 mM Magnesium Chloride	3.2	1.75 mM
10 mM dNTP Blend	3.7	0.8 mM (200 μM of each dNTP)
15 μM Gene Specific Upstream Primer DM151	0.5	0.15 μM
15 μM Gene Specific Downstream Primer DM152	0.5	0.15 μM
AmpliTaq Gold DNA Polymerase (5.0 Units/ μL)	0.5	2.5 Units/50 μL
Total	50.0	

**PCR Thermal
Cycling
Parameters for
Two-Step RNA
PCR**

The thermal cycler parameters described here are optimized for the GeneAmp Gold RNA PCR Reagent Kit pAW109 control reagents to generate a 308-bp amplification product.

Some thermal cyclers may require longer (~60 sec) denaturation times. Refer to the manufacturers operating manuals.

The parameters described here are only for the GeneAmp PCR Systems 2400, 9600, and 9700. See thermal cycler manuals for details on operation.

Note All GeneAmp PCR Systems display the Calculated Sample Temperature.

Cycling parameters for the PCR step of the two-step process

Step	AmpliAq Gold Activation	PCR CYCLE (43 cycles) ^a		PCR (Final Step)
	HOLD	Denature	Anneal/ Extend	HOLD
				Anneal/Extend
Time	10 min	20 sec	1 min	7 min
Temperature	95 °C	94 °C	62 °C	72 °C

a. The cycle number can be reduced when amplifying more than 5,000 copies of target RNA. For more information refer to "Cycle Optimization Guidelines" on page 17.

Appendix A. Additional Primer and PCR Product Information

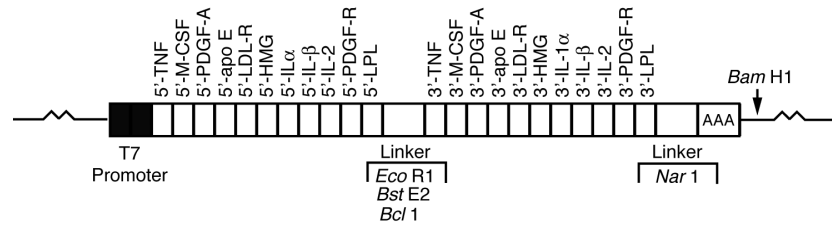
Additional Primers and PCR Products from pAW109 The DNA sequences for the other GeneAmplicer® primers in the control pAW109 sequences are shown below. Refer to “pAW109 Transcribed RNA” on page 29 to see a map of the pAW109 polylinker.

GeneAmplicer primers for amplicons in pAW109 RNA

Gene Name	Primer Information		PCR Product Size (bp) From	
	Name	Sequence	pAW109	mRNA
TNF	AW112	5'-CAGAGGGAAGAGTTCCCCAG-3'	301	325
	AW113	5'-CCTTGGTCTGGTAGGAGACG-3'		
M-CSF	AW111	5'-GAACAGTTGAAAGATCCAGTG-3'	302	192
	GM20	5'-TCGGACGCAGGCCTTGTCATG-3'		
PDGF-A	AW116	5'-CCTGCCCATTCGGAGGAAGAG-3'	301	225
	AW117	5'-TTGGCCACCTTGACGCTGCG-3'		
PDGF-B	AW118	5'-GAAGGAGCCTGGGTTCCCTG-3'	300	226
	AW119	5'-TTTCTCACCTGGACAGGTCG-3'		
LDL-R	AW125	5'-CAATGTCTCACCAAGCTCTG-3'	301	258
	AW126	5'-TCTGTCTCGAGGGGTAGCTG-3'		
HMG	AW102	5'-TACCATGTCAGGGGTACGTC-3'	303	247
	AW104	5'-CAAGCCTAGAGACATAATCATC-3'		
IL-1 α	DM151	5'-GTCTCTGAATCAGAAATCCTTCTATC-3'	308	421
	DM152	5'-CATGTCAAATTTCACTGCTTCATCC-3'		
IL-1 β	DM155	5'-AAACAGATGAAGTGCTCCTTCCAGG-3'	306	391
	DM156	5'-TGGAGAACACCACTTGTTGCTCCA-3'		
IL-2	DM135	5'-GAATGGAATTAATAATTACAAGAATCCC-3'	305	229
	DM136	5'-TGTTTCAGATCCCTTTAGTTCCAG-3'		
PDGF-R	AW158	5'-TGACCACCCAGCCATCCTTC-3'	300	228
	AW159	5'-GAGGAGGTGTTGACTTCATTC-3'		
LPL	AW155	5'-GAGATTTCTCTGTATGGCACC-3'	301	276
	AW157	5'-CTGCAAATGAGACACTTCTC-3'		

**pAW109
Transcribed RNA**

A schematic map of transcribed pAW109 RNA is shown below.



A schematic diagram of the general arrangement of the RNA transcribed from plasmid pAW109.

**Performance
Characteristics**

Each lot of the GeneAmp Gold RNA PCR Reagent Kit has been shown, under the conditions stated for One-Step RNA PCR (page 19), to yield a visible 308-bp band on an agarose gel stained with ethidium bromide starting with 1000 copies of pAW109 RNA.

Appendix B. References

- Chou, Q., Russell, M., Birch, D., Raymond, J., and Bloch, W. 1992. Prevention of pre-PCR mis-priming and primer dimerization improves low-copy-number amplifications. *Nucleic Acids Res.* 20:1717–1723.
- Faloona, F., Weiss, S., Ferre, F., and Mullis, K. 1990. Direct detection of HIV sequences in blood high-gain polymerase chain reaction [abstract]. In: *6th International Conference on AIDS*, University of California, San Francisco: San Francisco (CA). Abstract 1019.
- Innis, M.A., Myambo, K. B., Gelfand, D.H., and Brow, M.A.D. 1988. DNA sequencing with *Thermus aquaticus* DNA polymerase and direct sequencing of polymerase chain reaction-amplified DNA. *Proc. Natl. Acad. Sci. USA* 85:9436–9440.
- Kwok, S. and Higuchi, R. 1989. Avoiding false positives with PCR. *Nature* 339:237–238.
- Lawyer, F.C., Stoffel, S., Saiki, R.K., *et al.* 1989. Isolation, characterization, and expression in *Escherichia coli* of the DNA polymerase gene from *Thermus aquaticus*. *J. Biol. Chem.* 264:6427–6437.
- Mullis, K.B., and Faloona, F.A. 1987. Specific synthesis of DNA *in vitro* via a polymerase-catalyzed chain reaction. *Methods in Enzymol.* 155:335–350.
- Richardson, C.C. 1966. DNA polymerase from *Escherichia coli*. In: *Procedures in Nucleic Acid Research*. Cantoni, G.L. and Davies, D.R., eds. New York: Harper & Row. 263–276.
- Rychlik, W., Spencer, W.J., and Rhoads, R.E. 1990. Optimization of the annealing temperature for DNA amplification *in vitro* [published erratum appears in *Nucleic Acids Res* 1991 Feb 11;19(3):698]. *Nucleic Acids Res.* 18:6409–6412.
- Saiki, R.K., Scharf, S., Faloona, F., *et al.* 1985. Enzymatic amplification of b-globin genomic sequences and restriction site analysis for diagnosis of sickle cell anemia. *Science* 230:1350–1354.
- Saiki, R.K., Gelfand, D.H., Stoffel, S., *et al.* 1988. Primer-directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* 239:487–491.

Wang, A.M., Doyle, M.V., and Mark, D.F. 1989. Quantitation of mRNA by the polymerase chain reaction [published erratum appears in Proc. Natl. Acad. Sci. USA 1990 Apr;87(7):2865]. *Proc. Natl. Acad. Sci. USA* 86:9717–9721.

Appendix C. Technical Support

Contacting Technical Support You can contact Applied Biosystems for technical support by telephone or fax, by e-mail, or through the Internet. You can order Applied Biosystems user documents, MSDSs, certificates of analysis, and other related documents 24 hours a day. In addition, you can download documents in PDF format from the Applied Biosystems Web site (please see the section “To Obtain Documents on Demand” following the telephone information below).

To Contact Technical Support by E-Mail Contact technical support by e-mail for help in the following product areas:

Product Area	E-mail address
Genetic Analysis (DNA Sequencing)	galab@appliedbiosystems.com
Sequence Detection Systems and PCR	pcrlab@appliedbiosystems.com
Protein Sequencing, Peptide and DNA Synthesis	corelab@appliedbiosystems.com
Biochromatography, PerSeptive DNA, PNA and Peptide Synthesis systems, CytoFluor®, FMat™, Voyager™, and Mariner™ Mass Spectrometers	tsupport@appliedbiosystems.com
Applied Biosystems/MDS Sciex	api3-support@sciex.com
Chemiluminescence (Tropix)	tropix@appliedbiosystems.com

Hours for Telephone Technical Support In the United States and Canada, technical support is available at the following times:

Product	Hours
Chemiluminescence	8:30 a.m. to 5:30 p.m. Eastern Time
Framingham support	8:00 a.m. to 6:00 p.m. Eastern Time
All Other Products	5:30 a.m. to 5:00 p.m. Pacific Time

**To Contact
Technical Support
by Telephone or
Fax**

In North America

To contact Applied Biosystems Technical Support, use the telephone or fax numbers given below. (To open a service call for other support needs, or in case of an emergency, dial **1-800-831-6844** and press **1**.)

Product or Product Area	Telephone Dial...	Fax Dial...
ABI PRISM® 3700 DNA Analyzer	1-800-831-6844 , then press 8	1-650-638-5981
DNA Synthesis	1-800-831-6844 , then press 21	1-650-638-5981
Fluorescent DNA Sequencing	1-800-831-6844 , then press 22	1-650-638-5981
Fluorescent Fragment Analysis (includes GeneScan® applications)	1-800-831-6844 , then press 23	1-650-638-5981
Integrated Thermal Cyclers (ABI PRISM®877 and Catalyst 800 instruments)	1-800-831-6844 , then press 24	1-650-638-5981
ABI PRISM® 3100 Genetic Analyzer	1-800-831-6844 , then press 26	1-650-638-5981
Bioinformatics (includes BioLIMS™, BioMerge™, and SQL GT™ applications)	1-800-831-6844 , then press 25	1-505-982-7690
Peptide Synthesis (433 and 43X Systems)	1-800-831-6844 , then press 31	1-650-638-5981
Protein Sequencing (Procise® Protein Sequencing Systems)	1-800-831-6844 , then press 32	1-650-638-5981
PCR and Sequence Detection	1-800-762-4001 , then press 1 for PCR, 2 for the 7700 or 5700, 6 for the 6700 or dial 1-800-831-6844 , then press 5	1-240-453-4613

Product or Product Area	Telephone Dial...	Fax Dial...
Voyager™ MALDI-TOF Biospectrometry and Mariner™ ESI-TOF Mass Spectrometry Workstations	1-800-899-5858 , then press 13	1-508-383-7855
Biochromatography (BioCAD® Workstations and Poros® Perfusion Chromatography Products)	1-800-899-5858 , then press 14	1-508-383-7855
Expedite™ Nucleic acid Synthesis Systems	1-800-899-5858 , then press 15	1-508-383-7855
Peptide Synthesis (Pioneer™ and 9050 Plus Peptide Synthesizers)	1-800-899-5858 , then press 15	1-508-383-7855
PNA Custom and Synthesis	1-800-899-5858 , then press 15	1-508-383-7855
FMAT™ 8100 HTS System and Cytofluor® 4000 Fluorescence Plate Reader	1-800-899-5858 , then press 16	1-508-383-7855
Chemiluminescence (Tropix)	1-800-542-2369 (U.S. only), or 1-781-271-0045	1-781-275-8581
Applied Biosystems/MDS Sciex	1-800-952-4716	1-650-638-6223

Outside North America

Region	Telephone Dial...	Fax Dial...
Africa and the Middle East		
Africa (English Speaking) and West Asia (Fairlands, South Africa)	27 11 478 0411	27 11 478 0349
South Africa (Johannesburg)	27 11 478 0411	27 11 478 0349
Middle Eastern Countries and North Africa (Monza, Italia)	39 (0)39 8389 481	39 (0)39 8389 493

Region	Telephone Dial...	Fax Dial...
Eastern Asia, China, Oceania		
Australia (Scoresby, Victoria)	61 3 9730 8600	61 3 9730 8799
China (Beijing)	86 10 64106608	86 10 64106617
Hong Kong	852 2756 6928	852 2756 6968
Korea (Seoul)	82 2 593 6470/6471	82 2 593 6472
Malaysia (Petaling Jaya)	60 3 758 8268	60 3 754 9043
Singapore	65 896 2168	65 896 2147
Taiwan (Taipei Hsien)	886 2 2358 2838	886 2 2358 2839
Thailand (Bangkok)	66 2 719 6405	66 2 319 9788
Europe		
Austria (Wien)	43 (0)1 867 35 75 0	43 (0)1 867 35 75 11
Belgium	32 (0)2 712 5555	32 (0)2 712 5516
Czech Republic and Slovakia (Praha)	420 2 61 222 164	420 2 61 222 168
Denmark (Naerum)	45 45 58 60 00	45 45 58 60 01
Finland (Espoo)	358 (0)9 251 24 250	358 (0)9 251 24 243
France (Paris)	33 (0)1 69 59 85 85	33 (0)1 69 59 85 00
Germany (Weiterstadt)	49 (0) 6150 101 0	49 (0) 6150 101 101
Hungary (Budapest)	36 (0)1 270 8398	36 (0)1 270 8288
Italy (Milano)	39 (0)39 83891	39 (0)39 838 9492
Norway (Oslo)	47 23 12 06 05	47 23 12 05 75
Poland, Lithuania, Latvia, and Estonia (Warszawa)	48 (22) 866 40 10	48 (22) 866 40 20
Portugal (Lisboa)	351 (0)22 605 33 14	351 (0)22 605 33 15
Russia (Moskva)	7 095 935 8888	7 095 564 8787
South East Europe (Zagreb, Croatia)	385 1 34 91 927	385 1 34 91 840
Spain (Tres Cantos)	34 (0)91 806 1210	34 (0)91 806 1206
Sweden (Stockholm)	46 (0)8 619 4400	46 (0)8 619 4401
Switzerland (Rotkreuz)	41 (0)41 799 7777	41 (0)41 790 0676
The Netherlands (Nieuwerkerk a/d IJssel)	31 (0)180 331400	31 (0)180 331409

Region	Telephone Dial...	Fax Dial...
United Kingdom (Warrington, Cheshire)	44 (0)1925 825650	44 (0)1925 282502
All other countries not listed (Warrington, UK)	44 (0)1925 282481	44 (0)1925 282509
Japan		
Japan (Hacchobori, Chuo-Ku, Tokyo)	81 3 5566 6230	81 3 5566 6507
Latin America		
Del.A. Obregon, Mexico	305-670-4350	305-670-4349

**To Reach
Technical Support
Through the
Internet**

We strongly encourage you to visit our Web site for answers to frequently asked questions and for more information about our products. You can also order technical documents or an index of available documents and have them faxed or e-mailed to you through our site. The Applied Biosystems Web site address is

<http://www.appliedbiosystems.com/techsupp>

To submit technical questions from North America or Europe:

Step	Action
1	Access the Applied Biosystems Technical Support Web site.
2	Under the Troubleshooting heading, click Support Request Forms , then select the relevant support region for the product area of interest.
3	Enter the requested information and your question in the displayed form, then click Ask Us RIGHT NOW (blue button with yellow text).
4	Enter the required information in the next form (if you have not already done so), then click Ask Us RIGHT NOW . You will receive an e-mail reply to your question from one of our technical experts within 24 to 48 hours.

To Obtain Documents on Demand

Free, 24-hour access to Applied Biosystems technical documents, including MSDSs, is available by fax or e-mail or by download from our Web site.

To order documents...	Then...
by index number	a. Access the Applied Biosystems Technical Support Web site at http://www.appliedbiosystems.com/techsupp b. Click the Index link for the document type you want, then find the document you want and record the index number. c. Use the index number when requesting documents following the procedures below.
by phone for fax delivery	a. From the U.S. or Canada, call 1-800-487-6809 , or from outside the U.S. and Canada, call 1-858-712-0317 . b. Follow the voice instructions to order the documents you want. Note There is a limit of five documents per request.
through the Internet for fax or e-mail delivery	a. Access the Applied Biosystems Technical Support Web site at http://www.appliedbiosystems.com/techsupp b. Under Resource Libraries , click the type of document you want. c. Enter or select the requested information in the displayed form, then click Search . d. In the displayed search results, select a check box for the method of delivery for each document that matches your criteria, then click Deliver Selected Documents Now (or click the PDF icon for the document to download it immediately). e. Fill in the information form (if you have not previously done so), then click Deliver Selected Documents Now to submit your order. Note There is a limit of five documents per request for fax delivery but no limit on the number of documents you can order for e-mail delivery.

Headquarters

850 Lincoln Centre Drive
Foster City, CA 94404 USA
Phone: +1 650.638.5800
Toll Free: +1 800.345.5224
Fax: +1 650.638.5884

Worldwide Sales Offices

Applied Biosystems vast distribution and service network, composed of highly trained support and applications personnel, reaches into 150 countries on six continents. For international office locations, please call our local office or refer to our web site at www.appliedbiosystems.com.

www.appliedbiosystems.com



Applied Biosystems is committed to providing the world's leading technology and information for life scientists.

Printed in the USA, 07/2010
Part Number 4310253D