

Custom TaqMan[®] SNP Genotyping Assays

Protocol

© Copyright 2006, 2010 Applied Biosystems. All rights reserved.

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

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

NOTICE TO PURCHASER: LIMITED LICENSE

A license to perform the patented 5' nuclease process for research is obtained by the use of a Licensed 5' Nuclease Kit (containing Licensed Probe), or the combination of an Authorized 5' Nuclease Core Kit plus Licensed Probe, or by license rights that may be purchased from Applied Biosystems.

The Custom TaqMan® SNP Genotyping Assay contains Licensed Probe. Use of this product is covered by U.S. patent claims and corresponding patent claims outside the U.S. 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. Separate purchase of an Authorized 5' Nuclease Core Kit would convey rights under the applicable claims of US patents, and corresponding patent claims outside the United States, which claim 5' nuclease methods. No right under any other patent claims 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 from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Dr., Foster City, California 94404, USA.

TRADEMARKS:

AB (Design), ABI PRISM, Applied Biosystems, GeneAmp, and VIC are registered trademarks and Applera, FAM, MicroAmp, ROX, and TAMRA are trademarks of Applied Biosystems or its subsidiaries in the U.S. and/or certain other countries.

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

Microsoft is a registered trademark of Microsoft Corporation.

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

Part Number 4334431 Rev. F
07/2010

Contents

Preface	v
Safety	v
How to Obtain Support	ix
Product Overview	1
Custom TaqMan Genomic Assays Service	1
Custom TaqMan SNP Genotyping Assays	3
SNP Genotyping Assay Storage Recommendations	5
Assay Workflow	7
Sample Preparation	8
PCR Amplification	10
Allelic Discrimination Plate Read and Analysis	17
Troubleshooting	18
Appendix A: Assay Information in the AIF	25
Appendix B: Ordering Custom TaqMan SNP Genotyping Assays	31
Appendix C: User-Supplied Materials and Equipment .	37
Appendix D: Chemistry Overview	43
Appendix E: PCR Laboratory Practices	47
Bibliography	49

Preface

This preface covers:

Safety.	v
How to Obtain Support	xi


Safety


Safety Alert Words


Four safety alert words appear in Applied Biosystems user documentation at points in the document where you need to be aware of relevant hazards. Each alert word—**IMPORTANT**, **CAUTION**, **WARNING**, **DANGER**—implies a particular level of observation or action, as defined below.

Definitions


IMPORTANT! – Indicates information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.

 **CAUTION** – Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.

 **WARNING** – Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.

 **DANGER** – Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury. This signal word is to be limited to the most extreme situations.

Chemical Hazard Warning

 **WARNING CHEMICAL HAZARD.** Some of the chemicals used with Applied Biosystems instruments and protocols are potentially hazardous and can cause injury, illness, or death.

Chemical Safety Guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. (See “About MSDSs” on page vi.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer’s cleanup procedures as recommended in the MSDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

About MSDSs

Chemical manufacturers supply current Material Safety Data Sheets (MSDSs) with shipments of hazardous chemicals to *new* customers. They also provide MSDSs with the first shipment of a hazardous chemical to a customer after an MSDS has been updated. MSDSs provide the safety information you need to store, handle, transport, and dispose of the chemicals safely.

Each time you receive a new MSDS packaged with a hazardous chemical, be sure to replace the appropriate MSDS in your files.

Obtaining MSDSs

The MSDS for any chemical supplied by Applied Biosystems is available to you free 24 hours a day. To obtain MSDSs:

1. Go to <https://docs.appliedbiosystems.com/msdssearch.html>
2. In the Search field of the MSDS Search page:
 - a. Type in the chemical name, part number, or other information that you expect to appear in the MSDS of interest.
 - b. Select the language of your choice.
 - c. Click **Search**.
3. To view, download, or print the document of interest:

- a. Right-click the document title.
- b. Select:
 - **Open** – To view the document
 - **Save Target As** – To download a PDF version of the document to a destination that you choose
 - **Print Target** – To print the document
4. To have a copy of an MSDS sent by fax or e-mail, in the Search Results page:
 - a. Select **Fax** or **Email** below the document title.
 - b. Click **RETRIEVE DOCUMENTS** at the end of the document list.
 - c. Enter the required information.
 - d. Click **View/Deliver Selected Documents Now**.

Note: For the MSDSs of chemicals not distributed by Applied Biosystems, contact the chemical manufacturer.

Chemical Waste Hazards



CAUTION HAZARDOUS WASTE. Refer to Material Safety Data Sheets and local regulations for handling and disposal.



WARNING CHEMICAL WASTE HAZARD. Wastes produced by Applied Biosystems instruments are potentially hazardous and can cause injury, illness, or death.



WARNING CHEMICAL STORAGE HAZARD. Never collect or store waste in a glass container because of the risk of breaking or shattering. Reagent and waste bottles can crack and leak. Each waste bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.

Chemical Waste Safety Guidelines

To minimize the hazards of chemical waste:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the manufacturers of the chemicals in the waste container before you store, handle, or dispose of chemical waste.

- Provide primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Handle chemical wastes in a fume hood.
- After emptying the waste container, seal it with the cap provided.
- Dispose of the contents of the waste tray and waste bottle in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations.

Waste Disposal

If potentially hazardous waste is generated when you operate the instrument, you must:

- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure the health and safety of all personnel in your laboratory.
- Ensure that the instrument waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.

IMPORTANT! Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Biological Hazard Safety



WARNING

BIOHAZARD.

Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to: protective eyewear, face shield, clothing/lab coat, and gloves. All

work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

- U.S. Department of Health and Human Services guidelines published in *Biosafety in Microbiological and Biomedical Laboratories* (stock no. 017-040-00547-4; <http://bmbi.od.nih.gov>)
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030; http://www.access.gpo.gov/nara/cfr/waisidx_01/29cfr1910a_01.html).
- Your company's/institution's Biosafety Program protocols for working with/handling potentially infectious materials.

Additional information about biohazard guidelines is available at:

<http://www.cdc.gov>

Send Us Your Comments

Applied Biosystems welcomes your comments and suggestions for improving its user documents. You can e-mail your comments to:

techpubs@appliedbiosystems.com

IMPORTANT! The e-mail address above is only for submitting comments and suggestions relating to documentation. To order documents, download PDF files, or for help with a technical question, go to <http://www.appliedbiosystems.com>, then click the link for **Support**. (See “How to Obtain Support” below).

How to Obtain Support

For the latest services and support information for all locations, go to <http://www.appliedbiosystems.com>, then click the link for **Support**.

At the Support page, you can:

- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support

- Order Applied Biosystems user documents, MSDSs, certificates of analysis, and other related documents
- Download PDF documents
- Obtain information about customer training
- Download software updates and patches

In addition, the Support page provides access to worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities.

Product Overview

The Custom TaqMan[®] SNP Genotyping Assays are part of the Custom TaqMan[®] Genomic Assays service. This section provides a brief overview of the assays service (below) and the assays ([page 3](#)).

Custom TaqMan Genomic Assays Service

Product Description	<p>The Custom TaqMan[®] Genomic Assays service is an assay development service that designs, synthesizes, formulates, and delivers analytically quality-controlled primer and probe sets based on sequence information you submit. All information you submit is kept secure and confidential.</p> <p>Use the Custom TaqMan[®] Genomic Assays service to order:</p> <ul style="list-style-type: none">• Custom TaqMan[®] SNP Genotyping Assays – For performing genotyping studies with single-nucleotide polymorphisms (SNPs), detecting multinucleotide polymorphisms (MNPs), and detecting insertions/deletions (in/dels).• Custom TaqMan[®] Gene Expression Assays – For quantitative gene expression analysis and DNA sequence detection. <p>Note: This protocol describes the Custom TaqMan SNP Genotyping Assays. For information about the Custom TaqMan Gene Expression Assays, see protocol PN 4334429.</p>
Service Advantages	<p>The following are advantages of the Custom TaqMan Genomic Assays service:</p> <ul style="list-style-type: none">• Use of the service allows you to save the time, money, and effort involved in primer design, synthesis, and formulation of the gene-specific assays.• Quality-control tested assays reduce the need to test assays in your laboratory, saving you time and labor.• Orders are processed in a secure and confidential manner.• Assays are delivered ready-to-use under universal conditions.

Assay Quality

Assays are shipped only if they pass all analytical quality control checks below:

- Every primer and probe is individually tested by mass spectroscopy to verify the accuracy of the resulting synthesized oligonucleotide.
- Assays must meet probe and primer yield specifications, as performed by UV/Vis spectrophotometry.
- Human SNP Genotyping Assays are tested using 20 human genomic DNA samples. To pass this functional test, amplification must occur and at least one allelic discrimination cluster (heterozygous or homozygous, compared to No Template Controls) must be generated. Non-human SNP Genotyping Assays are not functionally tested.

Submission Guidelines

Refer to the *Custom TaqMan® Genomic Assays Protocol: Submission Guidelines* (PN 4367671) for guidelines and procedures on:

- Selecting, evaluating, and annotating a target sequence to create a submission file
- Ordering and reordering the custom assays

Custom TaqMan SNP Genotyping Assays

Product Capabilities

Custom TaqMan SNP Genotyping Assays allow you to:

Action	Example
Perform genotyping studies with any possible single-nucleotide polymorphism (SNP) in any organism	AGTTCATCCATGGTCA --> ACTTCATACATGGTCA Annotated as: AGTTCAT[C/A]CATGGTCA
Detect insertions/deletions (in/dels) of up to six bases for genotyping studies	AGTTCATCCATGGTCA --> AGTTCATGGTCA Annotated as: AGTTCAT[CCAT/*]GGTCA
Detect multinucleotide polymorphisms (MNPs) of up to six bases for genotyping studies	AGTTCATCCGGTCA --> AGTTCATATGGTCA Annotated as: AGTTCAT[CC/AT]GGTCA

The assays use the TaqMan[®] assay-based (5' nuclease) chemistry for amplifying and detecting specific polymorphisms in purified genomic DNA (gDNA) or complementary DNA (cDNA) samples.

IMPORTANT! To ensure optimal assay performance, Applied Biosystems recommends that you quality-check your target sequence through bioinformatic analysis (for example, perform a BLAST[®] search of the target sequence against the genome for specificity and masking of repetitive sequences). For more information, refer to the:

- *Bioinformatic Evaluation of a Sequence for Custom TaqMan[®] SNP Genotyping Assays* tutorial (available from <http://www.appliedbiosystems.com>)
- *Custom TaqMan[®] Genomic Assays Protocol: Submission Guidelines* (PN 4367671)
- Sequence Evaluation Information section in File Builder Software v3.0

Product Properties

All Custom TaqMan SNP Genotyping Assays:

- Are designed and optimized to work with TaqMan[®] Universal PCR Master Mix (with or without AmpErase[®] UNG) using the same thermal cycling conditions.

-
- Require only three components:
 - 1 to 20 ng of purified gDNA sample per well
 - Note:** You can use cDNA samples; however, Applied Biosystems does not currently provide recommendations for using cDNA with the Custom TaqMan SNP Genotyping Assays.
 - 40X SNP Genotyping Assay or 80X SNP Genotyping Assay (specific for each polymorphism)
 - TaqMan[®] Universal PCR Master Mix (with or without AmpErase[®] UNG)
 - Require only one PCR amplification step and an endpoint reading to obtain results.

SNP Genotyping Assay Contents

The 40X or 80X SNP Genotyping Assay contains:

- Sequence-specific forward and reverse primers to amplify the polymorphic sequence of interest.
- Two TaqMan[®] MGB probes:
 - One probe labeled with VIC[®] dye detects the Allele 1 sequence
 - One probe labeled with FAM[™] dye detects the Allele 2 sequence

Shipment Contents

Each Custom TaqMan SNP Genotyping Assay shipment consists of:

- One tube for each assay ordered, containing the Custom TaqMan SNP Genotyping Assay. Each tube is identified with a label on the side and a 2-D bar code on the bottom.
- Printed copy of the Data Sheet, which includes information about each assay ordered.
- Printed copy of the product insert, which includes an overview of the assays and PCR amplification protocols.
- CD-ROM containing:
 - A pdf file of the *Custom TaqMan[®] SNP Genotyping Assays Protocol* (PN 4334431)
 - A pdf file of the *Custom TaqMan[®] Genomic Assays Protocol: Submission Guidelines* (PN 4367671)
 - A pdf file of the product insert
 - The assay information file (AIF)

- A text file titled “Understanding Your Assay Information File” (PN 4343831)

About the Assay Information File

The assay information file (AIF) contains information about your Custom TaqMan SNP Genotyping Assay, including your order number and the primer and probe sequences.

For more information on the AIF, see [“Assay Information in the AIF” on page 25](#)

SNP Genotyping Assay Storage Recommendations

Storage and Stability

- Store the SNP Genotyping Assays at -15 to -25 °C in the dark.
IMPORTANT! Protect all Custom TaqMan SNP Genotyping Assays from direct exposure to light. Excessive exposure to light may affect the fluorescent probes.
- Do not perform more than 10 freeze-thaw cycles. If you expect to freeze-thaw the SNP Genotyping Assays more than three times, consider sub-aliquoting the SNP Genotyping Assays to minimize the number of freeze-thaw cycles.
- Applied Biosystems recommends diluting the SNP Genotyping Assays to a 20X working stock; see [“Diluting SNP Genotyping Assays” on page 6](#) below.
- It may be possible to store an assay for longer than a year. For long term storage, dilute assays to a 20X stock for best stability.

Diluting SNP Genotyping Assays

Applied Biosystems recommends that you dilute the SNP Genotyping Assays to a 20× working stock, then aliquot for routine use. By having aliquots, you can minimize freeze-thaw cycles and protect the SNP Genotyping Assays from exposure to light.



WARNING

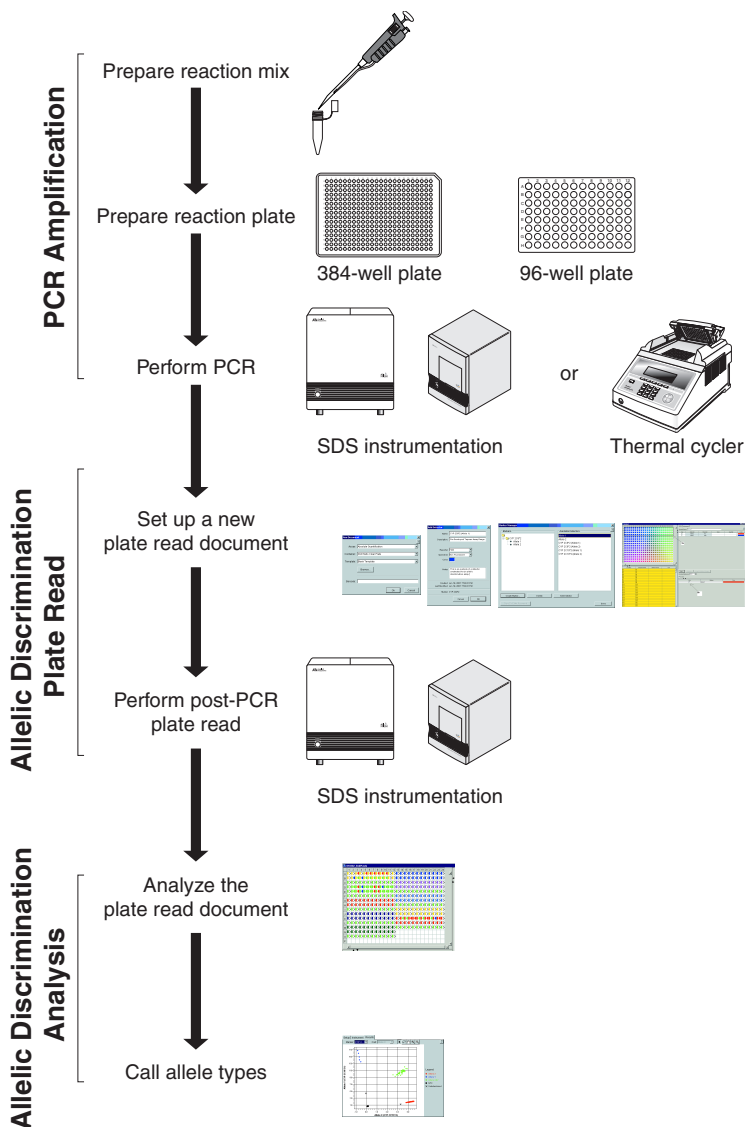
CHEMICAL HAZARD. SNP Genotyping Assay (<2% formamide). Exposure causes eye, skin, and respiratory tract irritation. It is a possible developmental and birth defect hazard. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

To dilute SNP Genotyping Assays:

1.	Dilute the 40× or 80× SNP Genotyping Assay to a 20× working stock with 1× TE buffer. Note: The 1× TE buffer should be 10 mM Tris-HCl, 1 mM EDTA, pH 8.0, and made using DNase-free, sterile-filtered water.
2.	Vortex, then centrifuge the mixture.
3.	Store multiple aliquots of the SNP Genotyping Assays at -15 to -25 °C.

Assay Workflow

The figure below shows the workflow for Custom TaqMan SNP Genotyping Assays. See [“Ordering Custom TaqMan SNP Genotyping Assays”](#) on page 31 and [“User-Supplied Materials and Equipment”](#) on page 37 for the materials needed to perform the assay.



Sample Preparation

Recommended DNA Template

The recommended DNA template for Custom TaqMan SNP Genotyping Assays is 1 to 20 ng of purified gDNA sample per well, with all wells in the same study containing the same amount of gDNA.

Note: Although you can use cDNA template, Applied Biosystems does not currently provide recommendations for using cDNA with the Custom TaqMan SNP Genotyping Assays.

IMPORTANT! If your assay is designed for gDNA, use gDNA template. If your assay is designed for cDNA, use cDNA template.

Quantifying gDNA

Applied Biosystems recommends quantifying the amount of gDNA in samples before using Custom TaqMan SNP Genotyping Assays. Use a reliable method, such as those listed below.

- UV/Vis spectrophotometry (A_{260}/A_{280} measurement) (Haque *et al.* 2003)
- Real-time quantification by RNase P
If you use the RNase P method, generate a standard curve using the DNA template standards in the TaqMan[®] DNA Template Reagents Kit (PN 401970) and the RNase P gene primers and probe provided in the TaqMan[®] RNase P Detection Reagents Kit (PN 4316831).

Note: Refer to the appropriate instrument user guide for details on performing and analyzing real-time runs using RNase P reagents.

Preparing the DNA Reaction Plate

Methods for Adding DNA

The Custom TaqMan SNP Genotyping Assays allow you to use either wet or dried-down DNA. If your experiment requires multiple plates that use the same DNA, or if you plan to use the same DNA in several experiments, it is convenient to dry down the DNA in the plates, which are then ready for use at any time. Both methods are described below.

Note: For both methods, use a calibrated, positive-displacement pipettor to minimize contamination and error.

Preparing a Plate with Wet DNA:

1.	<p>Dilute each DNA sample with DNase-free water to deliver a final DNA mass in the range of 1 to 20 ng per well.</p> <p>IMPORTANT! All wells in the same assay must contain the same amount of sample or control.</p> <table border="1" data-bbox="521 413 1193 645"> <thead> <tr> <th data-bbox="521 413 809 534">If you prepare a...</th> <th data-bbox="809 413 1193 534">Then the volume of DNA sample and DNase-free water per reaction should be...</th> </tr> </thead> <tbody> <tr> <td data-bbox="521 534 809 590">384-well reaction plate</td> <td data-bbox="809 534 1193 590">2.25 μL</td> </tr> <tr> <td data-bbox="521 590 809 645">96-well reaction plate</td> <td data-bbox="809 590 1193 645">11.25 μL</td> </tr> </tbody> </table> <p>Note: Multiple assays can be run on one reaction plate, but they must be analyzed separately.</p>	If you prepare a...	Then the volume of DNA sample and DNase-free water per reaction should be...	384-well reaction plate	2.25 μ L	96-well reaction plate	11.25 μ L
If you prepare a...	Then the volume of DNA sample and DNase-free water per reaction should be...						
384-well reaction plate	2.25 μ L						
96-well reaction plate	11.25 μ L						
2.	<p>Into each well of the 96-well or 384-well optical reaction plate, pipette one control or sample aliquot of the volume appropriate for the plate type (indicated in step 1).</p>						

Preparing a Plate with Dried-Down DNA:

1.	<p>Into each well of a 96-well or 384-well optical reaction plate, pipette one control or sample aliquot (1 to 20 ng of DNA).</p> <p>IMPORTANT! All wells belonging to the same assay must contain the same amount of sample or control.</p>
2.	<p>Dry-down the samples completely by evaporation at room temperature in a dark, amplicon-free location. Cover the plate with a lint-free tissue while drying.</p> <p>IMPORTANT! If you use gDNA, do not accelerate drying by heating the plate. Heating the plate may cause problems with gDNA recovery.</p> <p>Note: Multiple assays can be run on one reaction plate, but must be analyzed separately.</p>

PCR Amplification

Overview During the first step of a Custom TaqMan® SNP Genotyping Assay experiment, AmpliTaq Gold® DNA polymerase from the TaqMan Universal PCR Master Mix, No AmpErase® UNG, amplifies target DNA using sequence-specific primers. TaqMan MGB probes from the SNP Genotyping Assay provide a fluorescence readout of the amplification of each allele.

Note: Applied Biosystems recommends using TaqMan Universal PCR Master Mix, No AmpErase UNG, which is the master mix referred to in this protocol. However, you can use TaqMan Universal PCR Master Mix, which contains AmpErase UNG. For more information on UNG, see “[PCR Laboratory Practices](#)” on [page 47](#).

General Process PCR amplification requires that you:

- Prepare the reaction mix ([page 12](#)).
- Add the reaction mix to the prepared DNA reaction plate ([page 14](#)).
- Perform PCR ([page 16](#)).

See “[PCR Laboratory Practices](#)” on [page 47](#) for general instructions on avoiding false positive amplifications.

Assay Setup Guidelines

To ensure optimal analysis and troubleshooting of the Custom TaqMan SNP Genotyping Assays, prepare an optical reaction plate containing the following for each assay:

- DNA samples with unknown genotype at the SNP of interest.
- (Strongly recommended) Two no template controls (NTCs); use DNase-free water. Applied Biosystems strongly recommends using at least two NTCs per assay to:
 - Orient the VIC-dye and/or FAM-dye clusters to an origin.
 - Enable the detection of DNA contamination on a given set of plates.
- (Optional) DNA controls with known genotype at the SNP of interest.

Reagent and Sample Preparation Guidelines

- Keep all Custom TaqMan SNP Genotyping Assays protected from light, in the freezer, until you are ready to use them. Excessive exposure to light may affect the fluorescent probes.
- Minimize freeze-thaw cycles.
- To dilute the 40X or 80X SNP Genotyping Assay to a 20X working stock solution, see [“Diluting SNP Genotyping Assays” on page 6](#).
- Prior to use:
 - Thoroughly mix the TaqMan Universal PCR Master Mix, No AmpErase UNG, by swirling the bottle.
 - Resuspend the SNP Genotyping Assay by vortexing, then centrifuge the tube briefly.
 - (For wet DNA only) After thawing frozen DNA samples, resuspend the samples by vortexing, then centrifuge the tubes briefly.
- Prepare the reaction mix for each assay before transferring it to the optical reaction plate for thermal cycling.
- After adding the reagents to the DNA samples, mix thoroughly to avoid stratification of the reagents and/or air bubbles in the well. Stratification can lead to “stringy” clusters (see [page 20](#) in the [“Troubleshooting”](#) section).

Preparing the Reaction Mix

The reaction mix is made from 40X or 80X SNP Genotyping Assay, TaqMan Universal PCR Master Mix, No AmpErase UNG, and DNase-free water. The recommended final reaction volume per well is 5 µL for a 384-well plate and 25 µL for a 96-well plate.

Note: For instructions on diluting the 40X or 80X SNP Genotyping Assay to a 20X working stock solution, see [“Diluting SNP Genotyping Assays” on page 6.](#)



WARNING CHEMICAL HAZARD. SNP Genotyping Assay (<2% formamide). Exposure causes eye, skin, and respiratory tract irritation. It is a possible developmental and birth defect hazard. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.



CAUTION CHEMICAL HAZARD. TaqMan® Universal PCR Master Mix (2X), No AmpErase® UNG may cause eye and skin irritation. Exposure may cause discomfort if swallowed or inhaled. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

To prepare the reaction mix :

1. Calculate the number of reactions to be performed for each assay.

Applied Biosystems recommends including:

- At least two NTCs on each plate.
- If available, at least one known DNA control on each plate.

To prepare the reaction mix (*continued*):

2. Calculate the total volume of each component needed for each assay, using the table below. Be sure to choose the appropriate DNA delivery method for your experiment.

Component	Wet DNA Method		Dry-Down DNA Method	
	Volume ($\mu\text{L}/\text{Well}$)			
	384-well plate	96-well plate	384-well plate	96-well plate
TaqMan Universal PCR Master Mix (2X), No AmpErase UNG	2.50	12.50	2.50	12.50
20X working stock of SNP Genotyping Assay	0.25	1.25	0.25	1.25
DNase-free water	(none)	(none)	2.25	11.25
Total Volume per Well	2.75	13.75	5.00	25.00

Note: In your calculations, include some extra reactions to compensate for the volume loss that occurs during pipetting.

3. Gently swirl the bottle of TaqMan Universal PCR Master Mix (2X), No AmpErase UNG (abbreviated as “UMM” in subsequent steps). Ensure the UMM is well mixed before use.
4. Vortex and centrifuge the 20X SNP Genotyping Assay briefly.
5. Pipette the required total volumes of UMM and 20X SNP Genotyping Assay into a sterile microcentrifuge tube.
6. Cap the tube and invert several times to mix.

To prepare the reaction mix (continued):

- | | |
|----|---|
| 7. | Centrifuge the tube briefly to spin down the contents and to eliminate any air bubbles from the reaction mix. |
|----|---|

Adding the Reaction Mix

To add the reaction mix to the prepared DNA reaction plate :

- | 1. | Into each well of your DNA reaction plate (as prepared per the procedures on page 8), pipette the reaction mix as indicated below. | | | | | | | | | | | |
|------------|--|---------------------|--|--|-------------------------|---------------------|----------|------|---|---------|-------|----|
| | <table border="1" style="width: 100%; text-align: center;"> <thead> <tr> <th rowspan="2">Plate Type</th> <th colspan="2">Volume of Reaction Mix ($\mu\text{L}/\text{Well}$)</th> </tr> <tr> <th>Wet DNA Delivery Method</th> <th>Dry-Down DNA Method</th> </tr> </thead> <tbody> <tr> <td>384-well</td> <td>2.75</td> <td>5</td> </tr> <tr> <td>96-well</td> <td>13.75</td> <td>25</td> </tr> </tbody> </table> <p>IMPORTANT! Be sure that no cross-contamination occurs from well to well during pipetting.</p> | Plate Type | Volume of Reaction Mix ($\mu\text{L}/\text{Well}$) | | Wet DNA Delivery Method | Dry-Down DNA Method | 384-well | 2.75 | 5 | 96-well | 13.75 | 25 |
| Plate Type | Volume of Reaction Mix ($\mu\text{L}/\text{Well}$) | | | | | | | | | | | |
| | Wet DNA Delivery Method | Dry-Down DNA Method | | | | | | | | | | |
| 384-well | 2.75 | 5 | | | | | | | | | | |
| 96-well | 13.75 | 25 | | | | | | | | | | |
| 2. | Inspect all the wells for uniformity of volume, and note which wells do not appear to contain the proper volume. (See page 19 in the “Troubleshooting” section.) | | | | | | | | | | | |
| 3. | Seal the plate with the appropriate cover. (See page 38 for a list of covers compatible with your instrument.) | | | | | | | | | | | |
| 4. | Vortex the plate to mix the wells. | | | | | | | | | | | |
| 5. | Centrifuge the plate briefly to spin down the contents and eliminate any air bubbles. | | | | | | | | | | | |

Selecting a Thermal Cycler

Applied Biosystems instruments that can be used for PCR amplification are shown in the table below. See [page 38](#) for a list of reaction plates and covers that can be used on each instrument.

Instrument Type	Instrument Name
Thermal Cycler	Applied Biosystems 9800 Fast Thermal Cycler, using the 9700/9600 emulation mode Note: Custom TaqMan SNP Genotyping Assays can be performed on a 9800 Fast Thermal Cycler using <i>standard</i> reagents and <i>standard</i> cycling protocols. Custom TaqMan SNP Genotyping Assays are not supported using Fast reagents or Fast protocols.
	GeneAmp® PCR System 9700
Real-Time PCR System These instruments allow real-time analysis of PCR, which is helpful for troubleshooting. If using a Real-Time PCR System for PCR amplification, perform the endpoint plate read separately.	Applied Biosystems 7900HT Fast Real-Time PCR System [‡] Note: Custom TaqMan SNP Genotyping Assays can be performed on a 7900HT Fast System using <i>standard</i> reagents and <i>standard</i> cycling protocols. Custom TaqMan SNP Genotyping Assays are not supported using Fast reagents or Fast protocols.
	Applied Biosystems 7500 Fast Real-Time PCR System Note: Custom TaqMan SNP Genotyping Assays can be performed on a 7500 Fast System using <i>standard</i> reagents and <i>standard</i> cycling protocols. Custom TaqMan SNP Genotyping Assays are not supported using Fast reagents or Fast protocols.
	Applied Biosystems 7500 Real-Time PCR System
	Applied Biosystems 7300 Real-Time PCR System
	ABI PRISM® 7000 Sequence Detection System

[‡] *Applied Biosystems 7900HT Fast Real-Time PCR System* refers to all 7900HT instruments, regardless of the sample block being used (Fast or standard).

IMPORTANT! Because of differences in ramp rates and thermal accuracy, you may need to adjust the settings if you use thermal cyclers other than those indicated above.

IMPORTANT! Use of thermal cyclers from manufacturers other than Applied Biosystems is not supported by Applied Biosystems.

Performing PCR To perform PCR :

- Specify the thermal cycling conditions.

IMPORTANT! These conditions are optimized for use only with Custom TaqMan SNP Genotyping Assays on the instruments specified on [page 15](#).

Standard Protocol		
AmpliTaq Gold Enzyme Activation	PCR (40 Cycles)	
HOLD	Denature	Anneal/Extend
10 min at 95 °C	15 sec at 92 °C	1 min at 60 °C
Alternate Protocol [‡]		
AmpliTaq Gold Enzyme Activation	PCR (50 Cycles)	
HOLD	Denature	Anneal/Extend
10 min at 95 °C	15 sec at 92 °C	90 sec at 60 °C

Note: Refer to the appropriate instrument user guide for help with programming your thermal cycler.

- Specify the reaction volume according to the table below.

Plate Type	Reaction Volume (µL/Well)
384-well	5
96-well	25

- Load the reaction plate into the thermal cycler, then start the run.

Allelic Discrimination Plate Read and Analysis

Overview After PCR amplification, perform an endpoint plate read using an Applied Biosystems Real-Time PCR System. The Sequence Detection System (SDS) Software uses the fluorescence measurements made during the plate read to plot fluorescence (Rn) values based on the signals from each well. The plotted fluorescence signals indicate which alleles are in each sample.

General Process Performing a plate read and analyzing the data from Custom TaqMan SNP Genotyping Assays requires that you:

- Create and set up an allelic discrimination plate read document in the SDS software.
- Perform an allelic discrimination plate read on an Applied Biosystems Real-Time PCR System.
- Analyze the plate read document.
- Make manual allele calls or review automatic allele calls.
- Convert allele calls to genotypes.

For More Information For information about analyzing your data, refer to the following documents:

- 7900HT Fast System
 - *Applied Biosystems 7900HT Fast Real-Time PCR System Allelic Discrimination Getting Started Guide* (PN 4364015)
 - *Sequence Detection Systems Software Online Help for the Applied Biosystems 7900HT Fast Real-Time PCR System*
- 7300/7500/7500 Fast Systems
 - *Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Allelic Discrimination Getting Started Guide* (PN 4347822)
 - *Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Online Help*
- 7000 System – *ABI PRISM® 7000 Sequence Detection System User Guide* (PN 4330228)

Troubleshooting

Troubleshooting Table

Observation	Possible Cause	Recommended Action
NTCs generated fluorescence signals that cluster with DNA samples rather than close to the origin.	DNA contamination of the NTC wells and also other wells may have occurred.	Test your reagents for the presence of contaminating nucleic acid.
NTCs generated high fluorescence signal but did not cluster with DNA samples.	Some assays have high NTC fluorescence.	Measure Rn-NTC values for each cluster. If clusters are well-separated from NTCs, make allele calls as usual.
Distinct FAM- or VIC-dye homozygote clusters or heterozygote clusters were not observed.	Reporter dyes were not appropriately assigned.	1. Verify that reporter dyes are assigned to the correct allele.
	Quencher dye was improperly selected.	2. Reanalyze the plate read.

Observation	Possible Cause	Recommended Action
<p>A sample did not cluster with one specific allele type.</p>	<p>The sample may contain:</p> <ul style="list-style-type: none"> • More or less DNA than other samples • A rare allelic variation or sequence duplication • Mixtures of multiple alleles from DNA contamination 	<ol style="list-style-type: none"> 1. Recheck the DNA concentrations of the samples. 2. Retest the samples to verify true DNA outlier versus one-time PCR artifact. 3. Test the sample using a different SNP genotyping assay to see if the sample itself is contaminated.
	<p>Inaccurate reagent delivery or evaporation occurred.</p>	<p>Check all wells for uniformity in volume, noting which wells do not appear to contain the proper volume. Redo any reactions that did not contain the proper volume.</p>
	<p>Air bubbles are present in the wells during the PCR.</p>	<p>Remove air bubbles by swinging, tapping, or briefly centrifuging the reaction plate.</p> <p>For future assays, use a pipetting technique that does not form air bubbles. See “Avoiding Pipetting Errors” on page 23.</p>
	<p>A second SNP exists under the probe.</p>	<p>Perform bioinformatics. For more information, refer to the:</p> <ul style="list-style-type: none"> • <i>Custom TaqMan® Genomic Assays Protocol: Submission Guidelines</i> (PN 4367671) • <i>Bioinformatic Evaluation of a Sequence for Custom TaqMan® SNP Genotyping Assays</i> tutorial (Applied Biosystems Web site)
	<p>A second SNP exists under the primer.</p>	
<p>Unknown DNA samples did not generate fluorescence signals.</p>	<p>The sample may:</p> <ul style="list-style-type: none"> • Not contain DNA • Contain PCR inhibitors 	<ol style="list-style-type: none"> 1. Recheck the DNA concentrations of the samples using one of the methods listed in “Quantifying gDNA” on page 8. 2. Retest the sample to verify the result. 3. Test the sample using a different SNP genotyping assay.

Observation	Possible Cause	Recommended Action
<p>Clusters appear “stringy,” that is, in long stringy clusters on the three vectors from the NTCs. The Rn of the points varies widely, and the cluster may string all the way back to the NTCs.</p>	<p>Inefficient mixing of reagents.</p>	<ul style="list-style-type: none"> • Prepare the reaction mix first (page 12), then add the reaction mix to the prepared DNA reaction plate (page 14). • Mix the contents of each well after adding the reaction mix to the DNA samples (especially to wet DNA) by briefly vortexing the reaction plate and then centrifuging the plate prior to thermal cycling.
	<p>DNA samples are not of equal concentration due to:</p> <ul style="list-style-type: none"> • Poor or inaccurate DNA quantitation • Poor DNA quality 	<ul style="list-style-type: none"> • All the DNA samples on the plate should be of equivalent concentration, within the recommended range of 1 to 20 ng. If 1 ng of DNA is selected, then all samples on the plate should be at 1 ng. Do not mix 1-ng samples with 20-ng samples on the same plate. In addition, low DNA quantity (<1 ng/well) can cause poor amplification results. • Make sure DNA samples do not contain varying levels of DNA inhibitors. • Look at the ROX™ dye signal to ensure uniformity.
	<p>Poor thermal cycler performance.</p>	<p>Ensure that all wells of the thermal cycler are performing equally, then calibrate if necessary.</p>
	<p>Incorrect reagents were used (for example, master mix, assay, or water).</p>	<p>Repeat the experiment, making sure to use the correct reagents. (The reagents are listed on page 13).</p>

Observation	Possible Cause	Recommended Action
<p>All samples cluster with the NTCs (PCR amplification did not occur)</p>	<p>One or more of the reaction components was not added.</p>	<p>Make sure the DNA, Custom TaqMan SNP Genotyping Assay, and UMM were added to the reaction plates.</p>
	<p>AmpliAq Gold DNA polymerase was not activated efficiently.</p>	<p>Make sure AmpliTaq Gold DNA polymerase was activated by implementing the initial 10-minute cycle at 95 °C.</p>
	<p>The annealing temperature on the thermal cycler was too high or too low for the primers and/or probes.</p>	<p>Make sure the thermal cycler is set to the correct annealing and extension temperatures and times. Ensure that thermal cyclers are calibrated and maintained regularly.</p>
	<p>DNA was impure or of insufficient quantity.</p>	<p>Make sure that 1 to 20 ng of high-quality DNA sample was added to each well.</p>
	<p>A PCR inhibitor is present in the reaction. Impure DNA samples, as well as other inhibitors such as high concentrations of EDTA, can inhibit the PCR process.</p>	<p>Make sure you are following correct laboratory practices when preparing your samples. See "PCR Laboratory Practices" on page 47.</p>
<p>Some samples cluster with the NTCs.</p>	<p>The samples that did not amplify may have a known null allele (for example, GSTM1 or GSTT1 genes are known to be absent in a significant percentage of the population).</p>	<ul style="list-style-type: none"> • Check the literature for the possibility of a null genotype. • Try amplifying the gene with another set of PCR primers. If another primer set within the gene also does not amplify, it suggests the presence of a null genotype.
	<p>The samples may have evaporated.</p>	<p>Repeat the experiment, making sure to:</p>
	<p>The samples may be missing DNA template.</p>	<ul style="list-style-type: none"> • Use the correct reagents. (The reagents are listed on page 13).
	<p>The samples may be missing a reagent (for example, master mix, SNP Genotyping Assay, or water).</p>	<ul style="list-style-type: none"> • Use the recommended volumes: 5 µL for 384-well reaction plates and 25 µL for 96-well reaction plates.
<p>DNA was impure or of insufficient quantity.</p>	<p>Make sure that 1 to 20 ng of high-quality DNA sample was added to each well.</p>	

Observation	Possible Cause	Recommended Action
<p>Samples appear in a single cluster</p>	<p>The primer and probe sequences target non-specific sequences.</p>	<p>Perform bioinformatics. For more information, refer to the:</p> <ul style="list-style-type: none"> • <i>Custom TaqMan® Genomic Assays Protocol: Submission Guidelines</i> (PN 4367671) • <i>Bioinformatic Evaluation of a Sequence for Custom TaqMan® SNP Genotyping Assays</i> tutorial (http://www.appliedbiosystems.com)
	<p>The minor allele frequency (MAF) is low.</p>	<p>Inspect the MAF for the SNP. You may need a larger sample size to see the minor allele.</p>
	<p>The SNP is a pseudo-SNP, non-polymorphic SNP, or non-informative SNP for the population.</p>	<ul style="list-style-type: none"> • Verify that it is a SNP. • Check for a population-specific MAF.
<p>Scattered points creating “cloudy” or diffuse clusters.</p>	<p>Plate contains sample DNA of varying concentrations.</p>	<ul style="list-style-type: none"> • Requantitate sample DNA plate. • Ensure accurate DNA delivery to each well of the plate, so that each well contains the same amount of DNA (for example, all samples are at 1 ng).
	<p>Incorrect reagents were used (for example, master mix, assay, or water).</p>	<p>Repeat the experiment, making sure to use the correct reagents. (The reagents are listed on page 13).</p>
	<p>Varied amounts of reagent were dispensed to the plate.</p>	<p>Inspect the level of ROX™ dye across the plate to verify this cause.</p> <p>For future assays, use a pipetting technique that ensures uniform amounts of reagent are dispensed to the plate. See “Avoiding Pipetting Errors” on page 23.</p>

Avoiding Pipetting Errors

- Improve pipetting precision, as follows:
 - Calibrate and service the pipettors regularly.
 - Pipette larger volumes.
 - Reduce the number of pipetting steps whenever possible.
 - Increase the consistency of the pipetting method
 - Consult the manufacturer about the correct method of dispensing liquid volumes accurately from the pipettor. For example, some pipettors are designed to deliver the designated volume at the first plunger stop, so “blowing out” the residue may cause error.
- Use master reaction mixes. Applied Biosystems highly recommends using a master reaction mix, as follows:
 - a. Mix all common components (including the same template) to a set of reactions together.

Note: When you make each master reaction mix, add 5 to 10% additional volume to compensate for pipetting losses.
 - b. Dispense the mix to the replicate wells of the plate.

Appendix A: Assay Information in the AIF

This appendix describes the assay information that you receive with your shipment of a Custom TaqMan[®] SNP Genotyping Assay.

About the Assay Information File (AIF)

With each Custom TaqMan SNP Genotyping Assay order, you receive a CD-ROM on which is an assay information file (AIF). The AIF:

- Is identical to other AIFs for TaqMan[®]-based assays.
- Includes the number from the bar code on the box in which the assays were shipped.
- Is provided in tab-delimited format.
- Includes 55 data fields; this information includes 55 columns and 2 rows (see [Table on page 26](#)).
- Is most easily read when opened in Microsoft[®] Excel software or a similar spreadsheet program.

You can use the AIF to:

- Identify which assay is included in each assay tube.
- Associate the 2-D bar code on each assay tube with the Vial ID.
- Determine assay IDs.
- Determine the primer and probe sequences.

For LIMS Users

The AIF on your CD-ROM is named so that your LIMS system automatically uploads it. The AIF file name has the format: AbD_SNP_xx.yy.txt, where:

- AbD_SNP refers to the Custom TaqMan SNP Genotyping Assays.
- xx is the order number (up to 10 characters).
- yy is the plate ID number (up to 10 characters).

Determining Tube Contents

To determine the contents of each assay tube, match the Assay ID on the tube label with values in the Assay ID and Well Loc columns in the AIF.

AIF Columns

The table below describes the columns in the AIF.

Note: Because the information in the AIF varies by product line, all the fields are not filled for the Custom TaqMan SNP Genotyping Assays. In the table below, the Example column indicates “Blank” for those cases.

Contents of the assay information file:

Column	Example	Description
Customer Name	Blank	The name of your organization or institution
Order Number	123456789	Your sales order number
Ship Date	15-JAN-06	The date that the product is packaged for shipment
Delivery Number	See packing slip	A unique number for shipment (see packing slip for details)
Part Number	4331349	Product number used for ordering the assay
Product Type	Custom TaqMan® SNP Genotyping Assay Service	The type of product, which is indicated by the part number (part numbers are listed on page 31)
Assay ID	MyAssay	A unique identifier for the assay
Lot Number	Blank	A unique identifier for the manufacturing batch
Plate Type	96-position tube rack v1	The type of container in which the assay is shipped
Plate ID	1234567	Barcode ID of the box in which the assay is shipped
Vial Type	10-digit barcoded tube	The type of tube containing the assay
Vial ID	0004696076	A unique 2-D barcode of the assay tube
Well Location	B02	The well location of the assay in the associated barcoded plate
Assay Mix Conc.	40X	The concentration of the SNP Genotyping Assay [primers and probe(s)]

Contents of the assay information file: (continued)

Column	Example	Description
Forward Primer Name	MyAssayF	The name for the forward primer assigned by the design software; F designates the forward primer
Forward Primer Seq.	GGATCCTGAGCAATCA	The nucleotide sequence of the forward primer
Forward Primer Conc.	36	The concentration (μM) of the forward primer
Reverse Primer Name	MyAssayR	The name for the reverse primer assigned by the design software; R designates the reverse primer
Reverse Primer Seq.	TTACCGATCTTAATAA	The nucleotide sequence of the reverse primer
Reverse Primer Conc.	36	The concentration (μM) of the reverse primer
Reporter 1 Name	MyAssayV1	The name for the reporter 1 oligo (probe) assigned by the design software; V1 indicates the dye label and the design strand
Reporter 1 Dye	VIC [®] dye	The dye label for reporter 1
Reporter 1 Sequence	TTTCGCTGGCGTGAAG	The nucleotide sequence of reporter 1
Reporter 1 Conc.	8	The concentration (μM) of reporter 1
Reporter 1 Quencher	NFQ	A quencher used for reporter 1 of the assay
Reporter 2 Name	MyAssayM1	The name for the reporter 2 oligo (probe) assigned by the design software; M1 indicates the dye label and the design strand
Reporter 2 Dye	FAM [™] dye	The dye label for reporter 2
Reporter 2 Sequence	TCGCTGGCATGAAG	The nucleotide sequence of reporter 2
Reporter 2 Conc.	8	The concentration (μM) of reporter 2
Reporter 2 Quencher	NFQ	A quencher used for reporter 2 of the assay
Context Sequence	Blank	The nucleotide sequence surrounding the probe

Contents of the assay information file: *(continued)*

Column	Example	Description
Design Strand	Forward	<ul style="list-style-type: none"> • Forward – The probe binds to the forward strand • Reverse – The probe binds to the reverse strand
Category	Blank	Chromosome location of the SNP
Category ID	Blank	Chromosome location of the SNP
Group	Blank	Microsatellite markers associated with the SNP
Group ID	Blank	Microsatellite markers associated with the SNP
Gene Symbol	Blank	LocusLink symbol for the associated gene
Gene Name	Blank	LocusLink gene name
Chromosome	Blank	Chromosome on which gene/SNP is located
Species	Blank	Organism for which the assay was designed
Target Exons	Blank	The exon or exons (as referenced by the public accession number) that are spanned by the probe
NCBI Gene Reference	Blank	NCBI transcript ID that is detected by the assay
NCBI SNP Reference	Blank	Reference ID from the NCBI-dbSNP database
Medline Reference	Blank	PubMed references for the gene
Celera ID	Blank	Unique transcript or SNP ID in the Celera Discovery System (CDS)
Cytogenetic Band	Blank	Chromosomal band location of the gene (if not available, then chromosome number is provided)

Contents of the assay information file: *(continued)*

Column	Example	Description
SNP Type	Blank	Type of SNP (based on Celera Assembly); Acceptor Splice Site, Donor Splice Site, Intergenic/Unknown, Intron, Mis-sense Mutation, Nonsense Mutation, Putative UTR 5', Silent Mutation, UTR 3', UTR 5'
Minor Allele Freq - Caucasian	Blank	As calculated by SNP genotyping at Applied Biosystems - Caucasian
Minor Allele Freq - African-American	Blank	As calculated by SNP genotyping at Applied Biosystems - African-American
Minor Allele Freq - Japanese	Blank	As calculated by SNP genotyping at Applied Biosystems - Japanese
Minor Allele Freq - Chinese	Blank	As calculated by SNP genotyping at Applied Biosystems - Chinese
Celera Build Assembly Number	Blank	Version of the Celera Assembly from which the coordinate position is obtained
Location on Celera Assembly	Blank	Nucleotide location on the Celera human genome assembly (as referenced)
NCBI Assembly Number	Blank	Version of the NCBI assembly from which the coordinate position is obtained
Location on NCBI Assembly	Blank	Nucleotide location on the NCBI human genome assembly (as referenced)

Appendix B: Ordering Custom TaqMan SNP Genotyping Assays

Assay Part Numbers The Custom TaqMan® SNP Genotyping Assays part numbers indicate the type of assay and the number of reactions to order.

SNP human assay part numbers

Scale	Conc.	Number of Reactions		Part Number
		25- μ L Reaction 96-Well	5- μ L Reaction 384-Well	
Small	40X	300	1,500	4331349
Medium	40X	1,000	5,000	4332072
Large	80X	2,400	12,000	4332073

SNP non-human assay part numbers

Scale	Conc.	Number of Reactions		Part Number
		25- μ L Reaction 96-Well	5- μ L Reaction 384-Well	
Small	40X	300	1,500	4332077
Medium	40X	1,000	5,000	4332075
Large	80X	2,400	12,000	4332076

Considerations for Choosing Human SNP Genotyping Assays

Applied Biosystems performs a functional test on all human assays, in which genomic DNAs (gDNAs) from 20 unrelated individuals (from 4 populations and both sexes) are amplified under universal conditions to test for assay amplification and clustering. Assays that fail this test are not shipped. Consequently, assays to human targets that are expected to fail the functional test should be ordered as non-human assays, including the following:

- Human cDNA sequences.
The test fails because intronic sequences prevent primer or probe binding or separate assay component binding sites and prevent efficient amplification.
- Human Y-chromosome-specific sequences.
The test fails because >90% of the samples in the test must amplify to pass and the female samples in the functional test will not amplify.

Ordering Assays

To order Custom TaqMan SNP Genotyping Assays, create a submission file, then send it to Applied Biosystems.

Brief procedures are provided below for ordering new assays and for reordering previously received assays. For detailed procedures, please refer to the *Custom TaqMan® Genomic Assays Protocol: Submission Guidelines* (PN 4367671).

Ordering New Assays

Follow this procedure to order any Custom TaqMan SNP Genotyping Assay for the first time. The next time you want to order the *same* assay, see “[Reordering Assays](#)” on page 34.

To order a new assay:

1.	<p>From your computer, launch the File Builder software:</p> <ol style="list-style-type: none"> a. Go to: http://www.appliedbiosystems.com/filebuilder. b. Click the appropriate software version, then save the application file to your computer. c. Follow the onscreen instructions to install the software.
2.	<p>In File Builder, create a submission file with the following information:</p> <ul style="list-style-type: none"> • Name • Email address • Phone number • DNA sequence name • DNA sequence (~300 base pairs) • Target site coordinate containing target site, equal sign, and target site name

To order a new assay: *(continued)*

3. Send the submission file to Applied Biosystems using one of the following methods:

- Using the File Builder software, upload the file to the Applied Biosystems store.
- Send the file on a CD-ROM. For shipping addresses, contact your local Applied Biosystems sales office.
- Email the file to your regional Custom TaqMan Genomic Assays sales office as an attachment. In the Subject line, type **Custom TaqMan Genomic Assays Order Information**. Use the following email addresses:

Region	Email Address
Africa, CIS, Europe, Middle East, and West Asia	assays.europe@eur.appliedbiosystems.com
Australia and New Zealand	abozorders@appliedbiosystems.com
Japan	JPOrderAdminCS@appliedbiosystems.com
North America	genomics@appliedbiosystems.com

Note: For other regional offices, go <http://www.appliedbiosystems.com>.

Reordering Assays

Follow this procedure to order a Custom TaqMan SNP Genotyping Assay that you have previously ordered. To order a *new* assay, see [“Ordering New Assays” on page 32](#).

To reorder an assay:

1. Provide the following information:
 - Part number of the assay that you want to order
 - Original Assay ID (on the assay tube and on the Data Sheet)
 - Original sales order number(s) (on the plate, packing slip, AIF on the CD-ROM, and invoice)

Note: To ensure that you receive the identical sequences for the primers and probe, Applied Biosystems uses the assay ID and sale order number to retrieve your assay design information.

IMPORTANT! Do not submit the target DNA sequence unless you want to redesign your assay. New versions of Applied Biosystems design software may create assays with different primer and probe sequences, which may lead to a difference in assay performance.

To reorder an assay: *(continued)*

2. Email the information to your regional Custom TaqMan Genomic Assays sales office. In the Subject line, type **Custom TaqMan Genomic Assays Reorder Information**. Use the following email addresses:

Region	Email Address
Africa, CIS, Europe, Middle East, and West Asia	assays.europe@eur.appliedbiosystems.com
Australia and New Zealand	abozorders@appliedbiosystems.com
Japan	JPOrderAdminCS@appliedbiosystems.com
North America	genomics@appliedbiosystems.com

Note: For other regional offices, go to <http://www.appliedbiosystems.com>.

Note: You can reorder multiple assays in a single email.

Appendix C: User-Supplied Materials and Equipment

The following tables list materials and equipment required for using the Custom TaqMan[®] SNP Genotyping Assays.

These items are not supplied with the Custom TaqMan SNP Genotyping Assays. Unless otherwise noted, the listed items are available from major laboratory suppliers (MLSs).

Instruments

Instrument	Source
Thermal Cyclers	
Applied Biosystems 9800 Fast Thermal Cycler	Contact your local Applied Biosystems sales office.
GeneAmp [®] PCR System 9700	
Real-Time PCR Systems	
Applied Biosystems 7900HT Fast Real-Time PCR System [‡]	Contact your local Applied Biosystems sales office.
Applied Biosystems 7500 Fast Real-Time PCR System	
Applied Biosystems 7500 Real-Time PCR System	
Applied Biosystems 7300 Real-Time PCR System	
ABI PRISM [®] 7000 Sequence Detection System	

[‡] *Applied Biosystems 7900HT Fast Real-Time PCR System* refers to all 7900HT instruments, regardless of the sample block (Fast or standard).

Reaction Plates and Covers

The table below lists the reaction plates and covers you can use to perform Custom Taqman SNP Genotyping Assays. The instruments on which each reaction plate and cover can be used are shown at right.

Reaction Plate and Cover	Applied Biosystems Part Number	Compatible Instrument
96-Well Plates		
MicroAmp™ Optical 96-Well Reaction Plate with Barcode, 20 plates	4306737	<ul style="list-style-type: none"> • 9700 instrument • 7900HT Fast instrument, standard blocks • 7500 instrument • 7300 instrument • 7000 instrument
MicroAmp™ Optical 96-Well Reaction Plate with Barcode, 500 plates	4326659	
MicroAmp™ Optical 96-Well Reaction Plate with Barcode and Optical Caps, 20 plates/2400 caps	403012	
MicroAmp™ Optical 96-Well Reaction Plate with Barcode and Optical Adhesive Films, 100 plates/100 covers	4314320	
MicroAmp™ Optical 96-Well Reaction Plate, 10 plates	N8010560	
MicroAmp™ Optical 96-Well Reaction Plate, 500 plates	4316813	
MicroAmp™ Fast Optical 96-Well Thermal Cycling Plate with Barcode, 20 plates	4346906	<ul style="list-style-type: none"> • 9800 instrument • 7900HT Fast instrument • 7500 Fast instrument
MicroAmp™ Fast Optical 96-Well Thermal Cycling Plate with Barcode, 200 plates	4366932	
384-Well Plates		
MicroAmp™ Optical 384-Well Reaction Plate with Barcode, 50 plates	4309849	<ul style="list-style-type: none"> • 9700 instrument • 7900HT Fast instrument, standard blocks
MicroAmp™ Optical 384-Well Reaction Plate with Barcode, 500 plates	4326270	
MicroAmp™ Optical 384-Well Reaction Plate, No Barcode, 1000 plates	4343370	
MicroAmp™ Optical 384-Well Reaction Plate with Barcode, 1000 plates	4343814	

Reaction Plate and Cover	Applied Biosystems Part Number	Compatible Instrument
Covers		
MicroAmp™ Optical Adhesive Film Kit, 20 covers	4313663	<ul style="list-style-type: none"> • 9800 instrument • 9700 instrument
MicroAmp™ Optical Adhesive Film, 100 covers	4311971	<ul style="list-style-type: none"> • 7900HT Fast instrument
MicroAmp™ Optical Adhesive Films, 25 films	4360954	<ul style="list-style-type: none"> • 7500/7500 Fast instrument • 7300 instrument • 7000 instrument
MicroAmp™ Optical Caps, 8 caps/strip, 300 strips	4323032	<ul style="list-style-type: none"> • 9800 instrument • 9700 instrument • 7900HT Fast instrument, standard blocks • 7500 instrument • 7300 instrument
MicroAmp™ Clear Adhesive Films, 100 films	4306311	<ul style="list-style-type: none"> • 9800 instrument • 9700 instrument
MicroAmp™ Optical 8-Tube Strip (0.2 mL), 125 strips	4316567	<ul style="list-style-type: none"> • 9700 instrument • 7500 instrument • 7300 instrument

Reagents

Reagent	Source
Dilution of 40X or 80X SNP Genotyping Assay	
TE buffer (10 mM Tris-HCl, 0.1 mM EDTA, pH 8.0, made using DNase-free, sterile-filtered water)	MLS (Major Laboratory Supplier)
Sample Preparation: gDNA Quantification	
TaqMan [®] DNA Template Reagents Kit	Applied Biosystems (PN 401970)
TaqMan [®] RNase P Detection Reagents Kit	Applied Biosystems (PN 4316831)
PCR Amplification	
TaqMan [®] Universal PCR Master Mix, No AmpErase [®] UNG, 200 reactions	Applied Biosystems (PN 4324018)
TaqMan [®] Universal PCR Master Mix, 200 reactions	Applied Biosystems (PN 4304437)
TaqMan [®] Universal PCR Master Mix, No AmpErase [®] UNG, 2000 reactions	Applied Biosystems (PN 4326614)
TaqMan [®] Universal PCR Master Mix, 2000 reactions	Applied Biosystems (PN 4326708)
10-Pack, TaqMan [®] Universal PCR Master Mix, No AmpErase [®] UNG	Applied Biosystems (PN 4324020)
10-Pack, TaqMan [®] Universal PCR Master Mix	Applied Biosystems (PN 4305719)
Various Procedures	
DNase-free, sterile-filtered water	MLS

**Other Equipment
and
Consumables**

Item	Source
Centrifuge with plate adapter	MLS
Microcentrifuge	MLS
Vortexer	MLS
2-D barcode reader	Consumables suppliers
MicroAmp™ Cap Installing Tool	Applied Biosystems (PN 4330015)
MicroAmp™ Multi Removal Tool	Applied Biosystems (PN 4313950)
Polypropylene tubes	MLS
Pipettors: <ul style="list-style-type: none">• Positive-displacement• Air-displacement• Multichannel• Robotic pipetting station	MLS
Pipette tips, aerosol-resistant	MLS
Disposable gloves	MLS

Documents

Document	Applied Biosystems Part Number
<i>ABI PRISM® 7000 Sequence Detection System User Guide</i>	4330228
<i>Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Allelic Discrimination Getting Started Guide</i>	4347822
<i>Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Online Help</i>	NA
<i>Applied Biosystems 7900HT Fast Real-Time PCR System Allelic Discrimination Getting Started Guide</i>	4364015
<p><i>Bioinformatic Evaluation of a Sequence for Custom TaqMan® SNP Genotyping Assays tutorial</i></p> <p>To obtain this document:</p> <ol style="list-style-type: none"> Go to: http://www.appliedbiosystems.com/support/apptech/. Click Real-Time PCR/TaqMan® Genomic Assays. Under the TaqMan® Genomic Assays subheading, click the document link to open the pdf file. 	NA
<i>Custom TaqMan® Genomic Assays Protocol: Submission Guidelines</i>	4367671
<i>Real-Time PCR Systems Chemistry Guide</i>	4348358
<i>Sequence Detection Systems Software Online Help for the Applied Biosystems 7900HT Fast Real-Time PCR System</i>	NA
<i>TaqMan® Universal PCR Master Mix Protocol</i>	4304449

Software

Software	Source
File Builder Software v3.0	Contact your local Applied Biosystems sales office.
SDS Software v1.2.3 or later for the 7000 System	
SDS Software v1.3.1 or later for the 7300/7500/7500 Fast Systems	
SDS Software v2.2.2 or later for the 7900HT Fast System	
Microsoft® Excel software or equivalent spreadsheet software	Software suppliers

Appendix D: Chemistry Overview

- Assay Components** Each Custom TaqMan[®] SNP Genotyping Assay consists of a single tube containing:
- Two primers for amplifying the polymorphic sequence of interest
 - Two TaqMan[®] MGB probes for distinguishing between the two alleles

- About the Probes** Each TaqMan MGB probe contains:
- A reporter dye at the 5' end of each probe
 - VIC[®] dye is linked to the 5' end of the Allele 1 probe.
 - FAM[™] dye is linked to the 5' end of the Allele 2 probe.
 - A minor groove binder (MGB) at the 3' end of each probe. This modification increases the melting temperature (T_m) for a given probe length (Afonina *et al.*, 1997; Kutuyavin *et al.*, 1997), which allows the design of shorter probes. Shorter probes result in greater differences in T_m values between matched and mismatched probes, producing robust allelic discrimination. Even single nucleotide mismatches between a probe and the target sequence reduce the efficiency of probe hybridization, which in turn reduces the amount of reporter dye cleaved from a quenched probe. Furthermore, AmpliTaq Gold[®] DNA polymerase is more likely to displace a mismatched probe without cleaving it. Each of these factors minimizes the production of nonspecific fluorescence signals.
 - A nonfluorescent quencher (NFQ) at the 3' end of each probe. Because the quencher does not fluoresce, Applied Biosystems Real-Time PCR Systems can measure reporter dye contributions with greater sensitivity than with TAMRA[™] dye quencher-based probes.

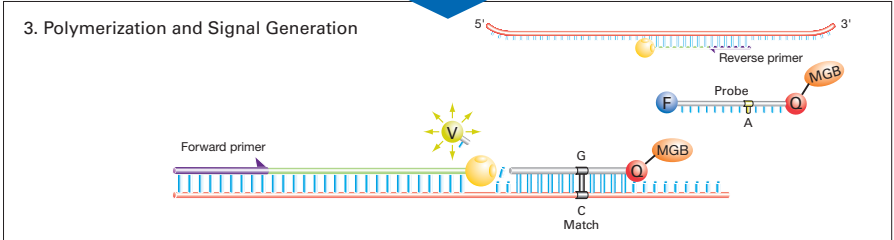
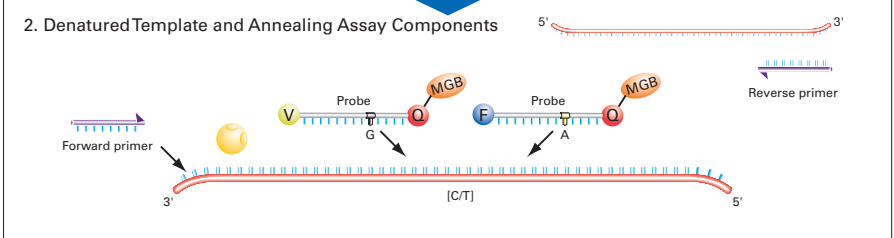
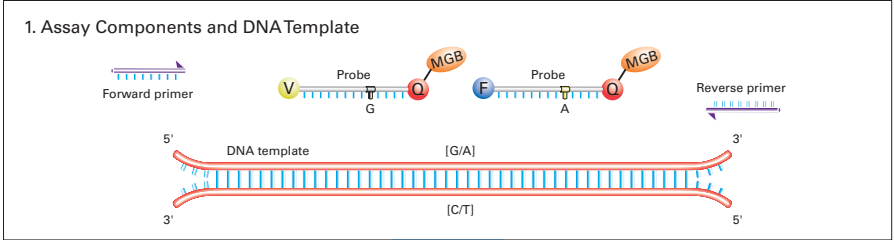
- 5' Nuclease Assay** During PCR, the following steps occur:

- Each TaqMan MGB probe anneals specifically to its complementary sequence between the forward and reverse primer sites.
- When the oligonucleotide probe is intact, the proximity of the reporter dye to the quencher dye results in quenching of the reporter fluorescence primarily by Förster-type energy transfer (FRET; Förster, 1948; Lakowicz, 1983).
- AmpliTaq Gold[®] DNA polymerase extends the primers bound to the template DNA.
- AmpliTaq Gold DNA polymerase cleaves only probes that are hybridized to the target.
- Cleavage separates the reporter dye from the quencher dye, which results in increased fluorescence by the reporter.
- The increase in fluorescence signal occurs when probes that have hybridized to the complementary sequence are cleaved. Thus, the fluorescence signal generated by PCR amplification indicates which alleles are present in the sample.

The table below shows the correlation between fluorescence signals and sequences in a sample.

A substantial increase in...	Indicates...
VIC-dye fluorescence only	Homozygosity for Allele 1
FAM-dye fluorescence only	Homozygosity for Allele 2
Both VIC- and FAM-dye fluorescence	Allele 1-Allele 2 heterozygosity

The figure on [page 45](#) illustrates the 5' nuclease assay process.



LEGEND	
	VIC® dye
	FAM™ dye
	Quencher
	Minor Groove Binder
	AmpliTaq Gold® DNA Polymerase
	Probe
	Primer
	Template
	Extended Primer

5' Nuclease assay process

Appendix E: PCR Laboratory Practices

Introduction PCR assays require special laboratory practices to avoid false positive amplifications (Kwok and Higuchi, 1989). The high throughput and repetition of these assays can lead to amplification of a single DNA molecule (Saiki *et al.*, 1985; Mullis and Faloona, 1987).

General PCR Practices

- 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, dedicated equipment, and supplies for:
 - Sample preparation
 - PCR setup
 - PCR amplification
 - Analysis of PCR products
- Never bring amplified PCR products into the PCR setup area.
- Open and close all sample tubes and reaction plates carefully. Try not to splash or spray PCR samples.
- Do not open sealed reaction plates.
- Keep reactions and components sealed as much as possible.
- Use positive-displacement pipettes or aerosol-resistant pipette tips.
- Clean lab benches and equipment periodically with freshly diluted 10% bleach solution.
- If you must open a sealed reaction plate, use TaqMan® Universal PCR Master Mix containing AmpErase® UNG (uracil-N-glycosylase) to minimize the risk of carryover contamination.

About AmpErase UNG

AmpErase uracil-N-glycosylase (UNG) is a 26-kDa recombinant enzyme encoded by the *Escherichia coli* uracil-N-glycosylase gene. This gene has been inserted into an *E. coli* host to direct expression of the native form of the enzyme (Kwok and Higuchi, 1989).

UNG acts on single- and double-stranded dU-containing DNA. It acts by hydrolyzing uracil-glycosidic bonds at dU-containing sites. The enzyme causes the release of uracil, thereby creating an alkali-sensitive apyrimidic site in the DNA. The enzyme has no activity on RNA or dT-containing DNA (Longo *et al.* 1990).

For 5' nuclease activities, AmpErase UNG treatment can prevent the reamplification of carryover PCR products. When dUTP replaces dTTP in PCR amplification, AmpErase UNG treatment can remove up to 200,000 copies of amplicon per 50- μ L reaction.

Bibliography

Afonina, I., Zivarts, M., Kutyavin, I., *et al.* 1997. Efficient priming of PCR with short oligonucleotides conjugated to a minor groove binder. *Nucleic Acids Res.* 25:2657–2660.

Förster, V. T. 1948. Zwischenmolekulare Energiewanderung und Fluoreszenz. *Annals of Physics (Leipzig)* 2:55–75.

Haque, K.A., Pfeiffer, R.M., Beerman M.B., *et al.* 2003. Performance of high-throughput DNA quantification methods. *BMC Biotechnology* 3:20.

Kutyavin, I.V., Lukhtanov, E.A., Gamper, H.B., and Meyer, R.B. 1997. Oligonucleotides with conjugated dihydropyrroloindole tripeptides: base composition and backbone effects on hybridization. *Nucleic Acids Res.* 25:3718–3723.

Kwok, S. and Higuchi, R. 1989. Avoiding false positives with PCR. *Nature* 339:237–238.

Lakowicz, J.R. 1983. Energy Transfer. In: *Principles of Fluorescence Spectroscopy*, New York: Plenum Press 303–339.

Longo, M.C., Berninger, M.S., and Hartley, J.L. 1990. Use of uracil DNA glycosylase to control carry-over contamination in polymerase chain reactions. *Gene* 93:125–128.

Mullis, K.B. and Faloona, F.A. 1987. Specific synthesis of DNA *in vitro* via a polymerase-catalyzed chain reaction. *Methods Enzymol.* 155:335–350.

Saiki, R.K., Scharf, S., Faloona, F., *et al.* 1985. Enzymatic amplification of β -globin genomic sequences and restriction site analysis for diagnosis of sickle cell anemia. *Science* 230:1350–1354.

Worldwide Sales and Support

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

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

Headquarters

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

07/2010