

USER GUIDE

applied
biosystems®
by *life* technologies™

TaqMan® *Salmonella enterica* Detection Kit

Catalog Number 4366104

Publication Number 4381939

Revision E



ABI 29/01 – 09/07

[End of validity: refer to certificate available at
www.afnor-validation.com]

ALTERNATIVE ANALYTICAL METHODS FOR AGRIBUSINESS

Certified by AFNOR Certification

For testing of Food and Environmental samples only.

life
technologies™

The information in this guide is subject to change without notice.

Disclaimer

LIFE TECHNOLOGIES CORPORATION 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 LIFE TECHNOLOGIES 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.

Limited Use Label License: environmental testing, quality control/quality assurance testing, food and agricultural testing

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product (a) to perform internal research for the sole benefit of the purchaser; and (b) for environmental testing, quality control/quality assurance testing, food and agricultural testing, including reporting results of purchaser's activities in environmental testing, quality control/quality assurance testing, food and agricultural testing for a fee or other commercial consideration. No other right is hereby granted expressly, by implication, or by estoppel. This product is for environmental testing, quality control/ quality assurance testing, food and agricultural testing and research purposes only.

The purchase of this product does not grant the purchaser any additional rights, including (without limitation) the right to transfer or resell the product in any form, the right to use the product as a therapeutic agent or diagnostics test component, or to use the product to perform other tests on a contract or fee per test basis for or on behalf of third parties. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008. Licensed to Life Technologies, Inc. for use in research only.

Human diagnostic uses require a separate license from Roche.

TRADEMARKS

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

AOAC is a registered trademark and Performance Tested Methods is a service mark of AOAC International.

AmpliQaq Gold and TaqMan are registered trademarks of Roche Molecular Systems, Inc., used under permission and license.

© 2012 Life Technologies Corporation. All rights reserved.

Contents

TaqMan® <i>Salmonella enterica</i> Detection Kit	5
Overview	5
About this protocol	5
Product description	5
Description of target microorganisms	5
Kit sensitivity	5
Kit specificity	6
Intended user	6
AOAC Performance Tested Methods sm Certification	6
ISO 16140 Validation	6
Expiration date	7
Operational conditions	7
Definitions of terms	8
Chemistry overview	9
Reaction components	9
Polymerase Chain Reaction (PCR)	9
Fluorescence detection	9
Internal Positive Control (IPC)	9
Materials and equipment	10
Kit contents	10
Storage	10
Shelf life	10
Equipment and materials not included	10
Kit workflow	12
Perform pre-enrichment	12
Overview	12
Pre-enrichment process	12
Prepare samples	13
Overview	13
Sample preparation process	13
Prepare PCR	14
Overview	14
Preparing PCR (without RapidFinder™ software)	14
Run the reactions	16
Overview	16
Running a plate without RapidFinder™ Software	16
View results	16

Overview	16
General process without RapidFinder™ Express Software	16
Resources for viewing results	17
■ APPENDIX A Specificity	19
Inclusivity	19
Exclusivity	21
■ APPENDIX B Troubleshooting	23
■ APPENDIX C References	25
■ APPENDIX D Good PCR Practices	27
Prevent contamination and nonspecific amplification	27
Overview	27
PCR good laboratory practices	27
Plate layout suggestions	28
■ APPENDIX E Safety	29
Chemical safety	30
Specific chemical handling	31
Biological hazard safety	31
Documentation and Support	33
Obtaining SDSs	33
Obtaining Certificates of Analysis	33
Obtaining support	33
Food safety support	33
Limited product warranty	34



TaqMan[®] *Salmonella enterica* Detection Kit

IMPORTANT! Before using this product, read and understand the information in [Appendix E, “Safety” on page 29](#) in this document.

Overview

About this protocol

This protocol provides:

- Background material about the TaqMan[®] *Salmonella enterica* Detection Kit
- A list of materials and equipment that can be used with the TaqMan[®] *Salmonella enterica* Detection Kit
- Guidelines for sample enrichment and preparation
- Instructions for preparing PCR using the TaqMan[®] *Salmonella enterica* Detection Kit on an Applied Biosystems Sequence Detection System (SDS) or Real-Time PCR System
- General troubleshooting guidelines

Product description

TaqMan[®] Pathogen Detection Kits provide a simple, reliable, and rapid procedure for detecting the presence of a specific bacterial pathogen. The assay uses the polymerase chain reaction (PCR) to amplify a target unique to that microorganism, and TaqMan[®] probes to detect the presence of the specific organism.

Description of target microorganisms

The TaqMan[®] *Salmonella enterica* Detection Kit detects all subspecies of *Salmonella enterica* including *enterica*, *salamae*, *arizonae*, *diarizonae*, *houtenae*, and *indica* only. *Salmonella enterica* causes disease in humans and other warm-blooded animals and therefore incorporates the most important clinical serovars. Hence, this method permits detection of *Salmonella enterica* (Groups I to IV) and is not suitable to detect *Salmonella bongori* (Group V). The genus *Salmonella*, consisting of the two species *Salmonella enterica* and *Salmonella bongori*, is a member of the *Enterobacteriaceae* family. More than 2,400 *Salmonella* serotypes have been reported, all of which are potentially pathogenic. *Salmonella enterica* with its six subspecies is of clinical relevance for humans. *Salmonella* is a frequently reported cause of foodborne illness, occurring in both food poisoning-triggered epidemics and isolated cases. *Salmonella* bacteria are the causative agent for Salmonellosis. Outbreaks have been associated with raw meats and poultry, eggs, milk and dairy products, seafood, coconut sauces, salad dressings, cocoa and chocolate, spices, frozen products, and vegetables such as hot peppers.

Kit sensitivity

TaqMan[®] Pathogen Detection Kits work on enriched samples, and detect as little as 1 CFU in 25 grams of food. The sensitivity of the assay in real culture samples depends on the quality of the sample preparation method that is used.

Kit specificity

The TaqMan® *Salmonella enterica* Detection Kit has demonstrated 100% inclusivity for more than 100 strains of *Salmonella enterica*. It has also demonstrated 100% exclusivity for 30 other non-*Salmonella* strains. For more information, refer to [Appendix A, “Specificity” on page 19](#).

Intended user

The intended users of the TaqMan® *Salmonella enterica* Detection Kit are investigators who need to detect the presence of *Salmonella enterica* in food. The TaqMan® *Salmonella enterica* Detection Kit is for use in food and environmental testing only. Not for any animal or human therapeutic or diagnostic use.

AOAC Performance Tested Methodssm Certification



The TaqMan® *Salmonella enterica* Detection Kit earned the Performance Tested Methodssm Certification from the AOAC Research Institute. The validation was conducted using ISO 6579 as the reference method. The validated workflow includes:

- The PrepMan® Ultra Sample Preparation Reagent Kit
- The TaqMan® *Salmonella enterica* Detection Kit
- The Applied Biosystems® 7300/7500 Real-Time PCR Systems
- RapidFinder™ Software

The workflow was certified for use with the following matrices:

Reference method	Matrix
ISO 6579	Food matrices: raw ground beef, chicken wings, cheddar cheese, and dry pet food

ISO 16140 Validation



ABI 29/01 – 09/07
[End of validity: refer to certificate at www.afnor-validation.com]

ALTERNATIVE ANALYTICAL METHODS FOR AGRIBUSINESS
Certified by AFNOR Certification

The TaqMan® *Salmonella enterica* Detection Kit has been certified “NF Validation”. The certification uses the ISO 16140 standard for the validation of alternative methods (Alternative Analytical Methods for Agribusiness. Certified by NF Validation; www.afnor-validation.com). This kit was compared and found equivalent to the ISO 6579 reference method. The validated workflow includes:

- The TaqMan® *Salmonella enterica* Detection Kit
- The PrepMan® Ultra Sample Preparation Reagent Kit
- The Applied Biosystems® 7000/7300/7500/7500 Fast/7700/StepOnePlus® Real-Time PCR Systems
- RapidFinder™ Software, Sequence Detection System (SDS) Software, and StepOne® Software

Reference method	Matrix
ISO 6579	All food and feed categories The tested product categories are meat products (processed and unprocessed), dairy products, egg products, seafood, vegetables, and feeding stuffs.



In the context of NF Validation, all samples identified as positive by the TaqMan® *Salmonella enterica* method must be confirmed, from the BPW enrichment broth, by performing one of the following tests:

- By performing an enrichment step in RVS (0.1 mL BPW in 10 mL of RVS broth): incubate at $41.5 \pm 1^\circ\text{C}$ for 24 ± 3 hours and streak onto XLD agar or another selective agar plate. Perform a Latex test on the observed characteristic colonies. In the case of a negative Latex test, perform a biochemical gallery on purified colonies characteristic of *Salmonella*. If the confirmatory result remains negative, proceed to a selective enrichment in MKTTn broth, as described in the CEN ISO standardized methods.
- According to classical tests described in methods standardized by CEN ISO from colonies (including purification step).
- Any other method certified by “NF Validation” that is based on a different principle than the TaqMan® *Salmonella enterica* method. It is necessary that the complete protocol of the validated method be performed entirely, which means that all steps that precede the confirmation step must be common to both methods.

In the event of discordant results (positive with the alternative method, non-confirmed by one of the means described above), the laboratory must follow the necessary steps to guarantee the validity of the obtained result.

General recommendations:

- Comply with Good Laboratory Practices (GLP; refer to EN ISO 7218 standard).
- We recommend that ISO 6579 and ISO 6887 standards be followed for the preparation of “master suspensions”.
- In the context of NF Validation, samples of more than 25 grams have not been tested.

Expiration date

For more information about the expiration date of the “NF Validation” certification, please refer to the certificate, ABI 29/01 – 09/07, available on the website at www.afnor-validation.com.

Operational conditions

Altitude

The Applied Biosystems® 7000/7300/7500/7500 Fast/7700/StepOnePlus® Real-Time PCR Systems are for indoor use only and for altitudes not exceeding 2,000 m (6,500 ft) above sea level.

Temperature and humidity requirements

Condition	Acceptable range
Temperature	15 to 30°C (50 to 90°F) Maximum change of less than 15°C (59°F) per 24 hours
Humidity	20 to 80% relative humidity, noncondensing

Definitions of terms

This protocol uses the following terms:

- **Amplification** – The process of making copies of, and thereby increasing the amount of, a specific DNA sequence.
- **Environmental Master Mix (EMM)** – A common reagent for all pathogen detection assays. The EMM is used to prepare the premix solution. It contains the polymerase enzyme that initiates PCR in the presence of the necessary primers and DNA sample.
- **Internal Positive Control (IPC)** – A control present in all reaction wells (contained in the Target Assay Mix). The IPC should always yield a positive result. If it does not, there may be a problem with amplification.
- **Negative control** – Monitors for contamination (unexpected amplification in the absence of a target) and reagent integrity (IPC signal should be present). One negative control is required for each target assay.
- **Polymerase Chain Reaction (PCR)** – Technology used to amplify, or increase the amount of a DNA sequence.
- **Positive control** – Monitors for the expected amplification of a target. Target signal not detected in a positive control well indicates a pipetting error or a problem with amplification. A positive control is optional for each target assay and is not recommended because it can cause cross-contamination.
- **Premix solution** – A solution you prepare that contains Environmental Master Mix (EMM) and Target Assay Mix (TAM).
- **Primer** – A segment of DNA that is complementary to the target DNA sequence or Internal Positive Control DNA sequence. Primer is needed to initiate amplification.
- **Probe** – A segment of DNA that is complementary to the target DNA sequence or Internal Positive Control DNA sequence. The probe is labeled with a reporter dye. When the probe binds to the target or Internal Positive Control, a reaction detected by the Sequence Detection System (SDS) or Real-Time PCR System indicates the presence of the target or Internal Positive Control.
- **Target** – The pathogen being tested.
- **Target Assay Mix (TAM)** – Target-specific reaction for the pathogen detection assay. TAM is used to prepare the premix solution. TAM contains specific primers and probes for the target and the IPC.
- **Unknown sample** – A DNA sample from a food substance that you are testing for the presence of one or more food pathogens.



Chemistry overview

Reaction components

Reaction components include:

- Target Assay Mix
- Environmental Master Mix
- DNA isolated from enrichment culture of raw materials or food samples, which you supply

Polymerase Chain Reaction (PCR)

PCR is a method used to amplify, or increase the amount of, a specific DNA sequence. Typically, the target DNA sequence is amplified using a reaction containing DNA polymerase, nucleotides, and primers complementary to that DNA sequence. When the reaction is heated, the DNA sequence denatures, separating into separate strands. As the solution cools, the primers anneal, or bind, to the target sequences in the separated DNA strand. The DNA polymerase then creates a new strand by extending the primers with nucleotides, creating a copy of the DNA sequence. When repeated, this cycle of denaturing, annealing, and extending exponentially increases the number of target DNA sequences. Ideally, no amplification occurs if the target DNA sequence is not present.

Fluorescence detection

The TaqMan® probe contains a fluorescent dye on one end and a quencher (that suppresses fluorescence) on the other. If the target sequence is present during the PCR, then the probe is degraded, resulting in an increase in fluorescence. The instrument detects accumulation of PCR products by monitoring the increase in fluorescence. The increase in fluorescence signal is detected only if the target sequence is complementary to the probe and is amplified during PCR. Because of these requirements, separate nonspecific amplification is not detected.




Internal Positive Control (IPC)

Applied Biosystems includes an IPC in the TaqMan® Pathogen Detection Kits. An IPC in each reaction avoids false negatives due to the presence of substances that inhibit PCR. The IPC also demonstrates whether or not PCR reagents are working and amplifying properly. This IPC eliminates the need for a positive control, reducing the risk of cross-contamination in unknown samples.

Materials and equipment

Kit contents

The TaqMan® *Salmonella enterica* Detection Kit (Part no. 4366104) contains reagents for 100 reactions. Components are shown in the table below.

Cap Color	Component name	Volume (µL)
	<i>Salmonella enterica</i> Target Assay Mix, 1 tube	300
	Negative Control, 1 tube	1000
	2X Environmental Master Mix, 2 tubes	750

Storage

- On receipt, store the 10X *Salmonella enterica* Target Assay Mix at –20°C or below and protect from light. Excessive exposure to light may affect the fluorescent probes.
- Store the Negative Control and 2X Environmental Master Mix at 5 ± 3°C.
- Minimize freeze-thaw cycles.

Shelf life

See expiration date on the outer box label.

Equipment and materials not included

The following table includes materials that are required for using (but not included in) the TaqMan® *Salmonella enterica* Detection Kit. Unless otherwise indicated, many of the listed items are available from major laboratory suppliers (MLS).

Instruments, equipment, consumables, and reagents	
Item	Source
Instruments	
Applied Biosystems® 7000 Real-Time PCR System	Contact your local Life Technologies sales office.
Applied Biosystems® 7300 Real-Time PCR System	
Applied Biosystems® 7500 Real-Time PCR System	
Applied Biosystems® 7500 Fast Real-Time PCR System	
Applied Biosystems® 7700 Real-Time PCR System	
StepOne® Real-Time PCR System	
Equipment	
Benchtop microcentrifuge	MLS
Block heater	MLS
Incubator shaker	MLS
Consumables	
Aerosol-resistant pipette tips	MLS
Disposable gloves	MLS

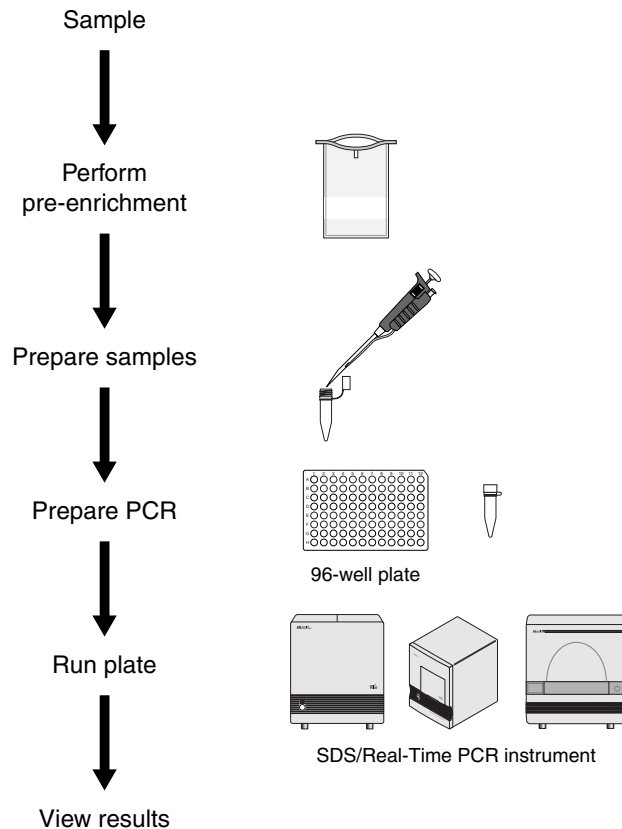


Instruments, equipment, consumables, and reagents	
Item	Source
Pipettors: <ul style="list-style-type: none"> • Positive-displacement • Air-displacement • Multichannel 	MLS
Applied Biosystems® Optical Adhesive Covers, 25 covers	Life Technologies Cat. no. 4360954
MicroAmp® Optical Caps, (8 caps/strip), 300 strips	Life Technologies Cat. no. 4323032
MicroAmp® Optical Adhesive Covers and MicroAmp® 96-Well Optical Reaction Plate with Barcode (code 128), 100 plates with covers	Life Technologies Cat. no. 4314320
MicroAmp® Optical Adhesive Cover Starter Kit	Life Technologies Cat. no. 4313663
MicroAmp® 96-Well Optical Reaction Plate with Barcode (code 128), 20 plates	Life Technologies Cat. no. 4306737
Sterile microcentrifuge tubes with attached screw-cap lid	MLS
Stomacher	MLS
Stomacher bags	MLS
Reagents	
PrepMan® Ultra Sample Preparation Reagent	Life Technologies Cat. no. 4318930
Buffered Peptone Water (BPW)	MLS
RNase-free, sterile-filtered water	MLS
<i>Salmonella enterica</i> subsp. <i>enterica</i> serovar <i>Typhi</i> (Source strain Ty2)‡	ATCC Cat. no. 700931D

‡ Provided for use as a positive control. We do not recommend the use of positive controls with TaqMan® Pathogen Detection Kits. Refer to “[Internal Positive Control \(IPC\)](#)” on page 9 for more information.

Kit workflow

The TaqMan® *Salmonella enterica* Detection Kit workflow is shown below.



Perform pre-enrichment

Overview

Pre-enrichment is the first step in using the TaqMan® *Salmonella enterica* Detection Kit. Pre-enrichment consists of growing the specific pathogen from a selected sample. The recommended enrichment procedure is based on ISO method 6579. For more information, refer to [Appendix C, “References” on page 25](#).

Pre-enrichment process

To perform pre-enrichment:

1. (Optional) Prepare an enrichment control, or an incubated pre-enrichment broth without any sample, as a negative control.
2. Warm the Buffered Peptone Water (BPW) broth to $37 \pm 1^\circ\text{C}$.
3. Perform pre-enrichment:
For ground beef, cheddar cheese, and dry pet food:
 - a. Combine the sample with BPW broth in a 1:9 ratio; up to 25 grams of sample in 225 mL of BPW broth.



- b. Close the stomacher bag, then place in a stomacher for at least one minute and homogenize at normal speed.
- c. Incubate the stomacher bag at $37 \pm 1^\circ\text{C}$ for 18 ± 2 hours.

For chicken wings:

- a. Add whole chicken wings to a filtered stomacher bag followed by 225 mL of BPW broth.
- b. Close the stomacher bag, then hand-massage it for 30 ± 15 seconds.
- c. Incubate the stomacher bag at $37 \pm 1^\circ\text{C}$ for 18 ± 2 hours.

Note: We recommend the use of filtered stomacher bags to prevent pipetting inaccuracy due to particulants.

Prepare samples

Overview

Sample preparation is the second step in using the TaqMan® *Salmonella enterica* Detection Kit. Sample preparation consists of preparing the pre-enriched sample for inclusion in your PCR.

Sample preparation process



WARNING! CHEMICAL HAZARD. PrepMan® Ultra Sample Preparation Reagent contains a material that can cause eye, skin, and respiratory tract irritation; and adverse effects on the kidneys, blood, and central nervous system. Read the SDS and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

To prepare the pre-enriched sample:

1. Transfer 1 mL of pre-enriched sample from the sample stomacher bag to a sterile microcentrifuge tube, then cap the tube. Use caution when handling enriched samples. Treat samples as biohazard waste.
Note: Sterile microcentrifuge tubes with attached screw-cap lids are recommended to avoid:
 - Lids opening during sample heating
 - Sample cross-contamination (resulting from the removed lids mistakenly placed on the wrong sample tubes)
2. Centrifuge the tube at maximum speed in a microcentrifuge for 4 ± 1 minutes to spin down the contents.
3. Remove the supernatant from the tube, leaving the pellet behind. Dispose of the supernatant according to biohazardous material disposal procedures.
Note: Do not disturb the pellet during supernatant extraction.
4. Add 100 μL of PrepMan® Ultra Sample Preparation Reagent, resuspend the pellet, then vortex the tube for 10 ± 5 seconds to mix the contents.
5. Heat the tube in a heat block at $100 \pm 5^\circ\text{C}$ for 12 ± 2 minutes, then vortex the tube for 10 ± 5 seconds to mix the contents.

Note: Incubate at $95 \pm 5^\circ\text{C}$ to reduce the frequency of the Eppendorf lids from popping open during this step.

6. Centrifuge the tube at maximum speed in a microcentrifuge for 4 ±1 minute to spin down the contents.
7. Transfer 50 µL of the supernatant into a new microcentrifuge tube. The sample can be stored at 5 ±3°C for up to a month.
Note: Do not disturb the pellet during supernatant extraction.
8. Dilute the sample DNA with RNase-free water by combining 10 µL of sample DNA with 90 µL of RNase-free water.
9. Vortex the new tube for 10 ±5 seconds to mix the contents. The sample DNA is now ready for PCR.


Prepare PCR

Overview

Preparing PCR is the third step in using the TaqMan® *Salmonella enterica* Detection Kit. Preparing PCR consists of creating the plate document and the reaction mix. The procedure for preparing PCR changes depending on whether or not you use RapidFinder™ Software.

Note: All TaqMan® Pathogen Detection Kits run with a single standard PCR protocol, allowing them to be combined on the same plate. This feature allows screening for multiple pathogens in the same PCR run.

IMPORTANT! If you use RapidFinder™ Software, refer to the “Create or Edit Plate and Pipette a Plate” instructions in the *RapidFinder™ Express Software Online Help*.

 **CAUTION!** CHEMICAL HAZARD. PrepMan® Ultra Sample Preparation Reagent contains a material that can cause eye, skin, and respiratory tract irritation; and adverse effects on the kidneys, blood, and central nervous system. Read the SDS and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Preparing PCR (without RapidFinder™ software)

To prepare the reaction mix:

1. Create and set up a plate document with thermal cycling conditions specified in the following table. Refer to the appropriate instrument user guide for more details.

Step	AmpliTaq Gold® enzyme activation	PCR	
		Cycle (45 cycles)	
	HOLD	Denature	Anneal/extend
Temp.	95°C	95°C	60°C
Time	10 min	15 sec	1 min

2. Thaw all reagents completely.
3. Label a tube with the target name.



4. Create the Premix Solution according to the following table.

Component	Vol. (µL) for one 30-µL reaction	Vol. (µL) for four 30-µL reactions‡
2X Environmental Master Mix (EMM)	15.0	66.0
10X Target Assay Mix (TAM)	3.0	13.2
Total Volume	18.0	79.2

‡ This example is illustrative for 4 reactions. Volumes shown include 10% overage to compensate for pipetting errors.

Note: Use a new tip when pipetting EMM and TAM.

Note: At least one negative control is required per target organism tested. Inclusion of a positive control is optional.

5. Mix the solution by gently pipetting up and down, then cap the tube.
6. Repeat steps 3 through 5 for each target assay.
7. Transfer 18 µL of Premix Solution into each well to be used, gently pipetting at the bottom of the well.

Note: Use a new tip for each Target Premix Solution.

Note: Refer to “Plate layout suggestions” on page 28 for more information.

IMPORTANT! Steps 8–10: Mix very gently with the pipette tip at the bottom of the tube to minimize aerosol formation and cross-contamination. Use a new tip for each well.

8. Transfer 12 µL of unknown sample (from [step 9 on page 14](#)) into each sample well, gently pipetting up and down to mix the solution.

IMPORTANT! Steps 8–10: Mix very gently with the pipette tip at the bottom of the tube to minimize aerosol formation and cross-contamination.

9. Transfer 12 µL of negative control into each negative-control well, gently pipetting up and down to mix the solution.
10. (*Optional*) Close all unknown sample and negative control tubes. Transfer 12 µL of positive control into each positive-control well, gently pipetting up and down to mix the solution.
- Note:** Use a new tip for each well, even when pipetting the same positive control.
- Note:** We recommend diluting the positive control with RNase-free water to 1 pg/mL, resulting in approximately 2,000 copies.
11. Close the tubes or apply an optical cover to the plate.
12. Make sure that the reagents are in the bottom of the wells. If available, use a centrifuge with a plate adapter to briefly centrifuge the plate.

Run the reactions

Overview

Running a plate is the fourth step in using the TaqMan® *Salmonella enterica* Detection Kit. Running a plate involves using an Applied Biosystems Sequence Detection System (SDS) or Real-Time PCR System to analyze your sample.

IMPORTANT! If you use RapidFinder™ Software, refer to the “Running a Plate” instructions in the *RapidFinder™ Express Software Online Help* (Pub. no. 4401842).

Running a plate without RapidFinder™ Software

To run the plate:

1. Open the plate document that corresponds to the reaction plate.
2. Load the reaction plate into the SDS or Real-Time PCR System.
3. Start the run.

IMPORTANT! To avoid false positives due to amplified material in your work area, do not open tubes after amplification.

View results

Overview

Viewing results varies depending on the instrument you use. Refer to the appropriate instrument user guide of your real-time PCR instrument or the *RapidFinder™ Software Online Help* (Pub. no. 4401842) for instructions on how to analyze data and view your results.

IMPORTANT! If you use RapidFinder™ Express Software, refer to the “Viewing Results” instructions in the *RapidFinder™ Express Software Online Help* (Pub. no. 4401842).

General process without RapidFinder™ Express Software

The general process for viewing results from TaqMan® Pathogen Detection Kits involves:

- Viewing the amplification plots for the entire plate.
- Setting the baseline and threshold values.
- Checking each sample for a FAM™ dye (target-specific) signal and a VIC® dye signal (IPC). The following table provides a basic guide for interpreting the results:

FAM™ dye signal (target)	VIC® dye signal (IPC)	Result
+	+, -	Positive
-	+	Negative
-	-	See Appendix B, “Troubleshooting” on page 23.



Resources for viewing results

For more information about analyzing your data, refer to:

- The appropriate instrument user guide
- *RapidFinder™ Express Software Online Help* (Pub. no. 4401842)

We do not recommend using the same method to both screen samples and confirm the results. The confirmation methods that can be used (for confirmation of positive results) in the context of NF Validation are described on [pages 6 and 7](#).





Specificity

Inclusivity

Genus serotypes	Source[‡]
<i>Salmonella</i> <i>Abaetetuba</i>	Abaetetuba Food Company
<i>Salmonella</i> <i>Abony</i>	NCTC 6017
<i>Salmonella</i> <i>Adelaide</i>	Silliker
<i>Salmonella</i> <i>Agona</i>	Silliker
<i>Salmonella</i> <i>Ajiobo</i>	Ajiobo Food Company
<i>Salmonella</i> <i>Alachua</i>	Silliker
<i>Salmonella</i> <i>Albany</i>	CDC 151-86
<i>Salmonella</i> <i>Anatum</i>	Silliker
<i>Salmonella</i> <i>Anatum</i>	ATCC 9270
<i>Salmonella</i> <i>Arizonae</i>	CDC 346-86
<i>Salmonella</i> <i>Bareilly</i>	Silliker
<i>Salmonella</i> <i>Bere</i>	CDC 315-86
<i>Salmonella</i> <i>Berta</i>	CDC 212-86
<i>Salmonella</i> <i>Betioky</i>	CDC 946
<i>Salmonella</i> <i>Binza</i>	Silliker
<i>Salmonella</i> <i>Bockenheim</i>	CDC 1097
<i>Salmonella</i> <i>Bovismorbificans</i>	CDC 889-83
<i>Salmonella</i> <i>Brandenburg</i>	Brandenburg Food Company
<i>Salmonella</i> <i>Bredeney</i>	Bredeney Feed Company
<i>Salmonella</i> <i>California</i>	ATCC 23201
<i>Salmonella</i> <i>Cerro</i>	ATCC 10723
<i>Salmonella</i> <i>Choleraesuis</i>	ATCC 10708
<i>Salmonella</i> <i>Choleraesuis</i> <i>Kunzendorf</i>	ATCC 12011
<i>Salmonella</i> <i>Crossness</i>	CDC 1795
<i>Salmonella</i> <i>Cubana</i>	Silliker
<i>Salmonella</i> <i>Derby</i>	Derby Feed Company
<i>Salmonella</i> <i>Drypool</i>	Silliker
<i>Salmonella</i> <i>Eastbourne</i>	Silliker
<i>Salmonella</i> <i>Enteritidis</i>	ATCC 4931
<i>Salmonella</i> <i>Ferlac</i>	ATCC 43976
<i>Salmonella</i> <i>Flint</i>	CDC 16-86
<i>Salmonella</i> <i>Gallinarum</i>	CDC 2950-79
<i>Salmonella</i> <i>Gaminara</i>	ATCC 8324
<i>Salmonella</i> <i>Glostrup</i>	ATCC BAA 556
<i>Salmonella</i> <i>Hadar</i>	ATCC 51956
<i>Salmonella</i> <i>Havana</i>	Silliker
<i>Salmonella</i> <i>Heidelberg</i>	ATCC 8326

Genus serotypes	Source†
<i>Salmonella Hilversum</i>	ATCC 15784
<i>Salmonella Harmelen</i>	ATCC 15783
<i>Salmonella Houton</i>	ATCC 29834
<i>Salmonella Illinois</i>	ATCC 11646
<i>Salmonella Indiana</i>	CDC 343-86
<i>Salmonella Infantis</i>	CDC 139-84
<i>Salmonella Inverness</i>	CDC 133-87
<i>Salmonella Iaviana</i>	Silliker
<i>Salmonella Jerusalem</i>	Silliker
<i>Salmonella Johannesburg</i>	CDC 500-86
<i>Salmonella Kentucky</i>	Kentucky Food Company
<i>Salmonella Kottbus</i>	CDC 292-85
<i>Salmonella Krefeld</i>	CDC 939-84
<i>Salmonella Livingstone</i>	CDC 110-85
<i>Salmonella Locarno</i>	CDC 886
<i>Salmonella London</i>	CDC 382-84
<i>Salmonella Madelia</i>	CDC 288-85
<i>Salmonella Malawi</i>	CDC 327-80
<i>Salmonella Manila</i>	Manila Feed Company
<i>Salmonella Maregrosso</i>	CDC 235-77
<i>Salmonella Mbandaka</i>	Silliker
<i>Salmonella Miami</i>	CDC 114-84
<i>Salmonella Michigan</i>	CDC 47-85
<i>Salmonella Minnesota</i>	CDC 245-86
<i>Salmonella Montevideo</i>	Silliker
<i>Salmonella Muenchen</i>	CDC 427-85
<i>Salmonella Newbrunswick</i>	Silliker
<i>Salmonella Newport</i>	CDC 520-86
<i>Salmonella Nienstedten</i>	Silliker
<i>Salmonella Ohio</i>	CDC 299-86
<i>Salmonella Oranienburg</i>	CDC 447-86
<i>Salmonella Ouakam</i>	CDC 32-87
<i>Salmonella Panama</i>	CDC 551-86
<i>Salmonella Paratyphi A</i>	CDC 1072-82
<i>Salmonella Paratyphi B</i>	ATCC 8759
<i>Salmonella Paratyphi C</i>	ATCC 13428
<i>Salmonella Phoenix</i>	CDC 696-84
<i>Salmonella Poona</i>	CDC 251-86
<i>Salmonella Potsdam</i>	ATCC 25957
<i>Salmonella Putten</i>	ATCC 15787
<i>Salmonella Rissen</i>	Silliker
<i>Salmonella Roodepoort</i>	Roodepoort Pet Food Company
<i>Salmonella Rubislaw</i>	Silliker
<i>Salmonella Saintpaul</i>	Silliker
<i>Salmonella Salamae</i>	ATCC 15786
<i>Salmonella Sandiego</i>	ATCC 23199

Genus serotypes	Source [‡]
<i>Salmonella Saphra</i>	CDC 479-85
<i>Salmonella Schwarzengrund</i>	Schwarzengrund Feed Company
<i>Salmonella Senftenberg</i>	Silliker
<i>Salmonella Simsbury</i>	ATCC 12004
<i>Salmonella Sundsvall</i>	CDC 668-84
<i>Salmonella Tallahasee</i>	ATCC 12002
<i>Salmonella Tennessee</i>	Tennessee Bakery
<i>Salmonella Thompson</i>	Thompson Food Company
<i>Salmonella Tranora</i>	CDC 857
<i>Salmonella Treforest</i>	CDC 657
<i>Salmonella Typhi</i>	CDC 2049-87
<i>Salmonella Typhimurium</i>	ATCC 6994
<i>Salmonella Urbana</i>	Silliker
<i>Salmonella Vellore</i>	ATCC 15611
<i>Salmonella Wassenaar</i>	CDC 121-87
<i>Salmonella Weltevreden</i>	CDC 110-83
<i>Salmonella Weslasco</i>	CDC 185-87
<i>Salmonella Worthington</i>	CDC 247-85

[‡] NCTC: National Collection of Type Cultures, USA; ATCC: American Type Culture Collection, USA; Silliker: Silliker Food Science Center Culture Collection, USA; CDC: Centers for Disease Control, USA.

Exclusivity

Genus serotypes	Source [‡]
<i>Acetobacter aceti</i>	ATCC 23746
<i>Citrobacter amalonaticus</i>	Testing Kit Company
<i>Citrobacter farmieri</i>	ATCC 51112
<i>Citrobacter freundii</i>	ATCC 6879
<i>Citrobacter braakii</i>	ATCC 29063
<i>Citrobacter youngae</i>	ATCC 29935
<i>Enterobacter aerogenes</i>	ATCC 35028
<i>Enterobacter agglomerans</i>	ATCC 29917
<i>Enterobacter cloacae</i>	ATCC 35030
<i>Enterobacter cloacae</i>	ATCC 7256
<i>Enterobacter sakazakii</i>	ATCC 51329
<i>Enterococcus faecalis</i>	ATCC 49452
<i>Enterococcus faecalis</i>	ATCC 33186
<i>Escherichia blattae</i>	ATCC 29907
<i>Escherichia coli</i>	ATCC 11229
<i>Escherichia coli O157:H7</i>	ATCC 35150
<i>Hafnia alvei</i>	ATCC 25927
<i>Klebsiella oxytoca</i>	ATCC 13182
<i>Klebsiella pneumoniae</i>	ATCC 9591
<i>Morganella morganii</i> subsp. <i>Morganii</i>	ATCC 29853

Genus serotypes	Source [‡]
<i>Providencia stuartii</i>	Testing Kit Company
<i>Proteus mirabilis</i>	ATCC 25933
<i>Proteus vulgaris</i>	ATCC 8427
<i>Serratia liquefaciens</i>	ATCC 27592
<i>Serratia marcescens</i>	ATCC 14041
<i>Shigella boydii</i>	ATCC 9207
<i>Shigella dysenteriae</i>	ATCC 11456b
<i>Shigella flexneri</i>	ATCC 9199
<i>Shigella sonnei</i>	ATCC 25931
<i>Staphylococcus aureus</i>	ATCC 13565
<i>Yersinia enterocolitica</i>	ATCC 29913
<i>Yersinia frederksensii</i>	Silliker
<i>Yersinia pseudotuberculosis</i>	Testing Kit Company

[‡] ATCC: American Type Culture Collection, USA; Silliker: Silliker Food Science Center Culture Collection, USA; CDC: Centers for Disease Control, USA.

B

Troubleshooting

Observation	Possible cause	Recommended action
No IPC or target-specific signal is detected in unknown wells.	Inhibition of PCR occurred.	Repeat sample preparation, then repeat assay. If PCR is still inhibited, dilute the sample (for example, 1:5 or 1:10) to dilute inhibitors. Alternatively, use a Bacterial Genomic DNA Purification Kit (Major Laboratory Supplier) to remove inhibitors.
	Environmental Master Mix not stored properly	Repeat the assay using properly stored assay components.
	Target-specific 10X Assay Mix not stored properly	Avoid freezing and thawing assay components. Protect the Assay Mix from light.
	Pipetting error (no premix solution added)	Repeat the assay. Make sure to pipette premix solution into all wells.
No IPC signal is detected, but target-specific signal is detected in positive control wells.	A high copy number of target DNA exists in samples, resulting in preferential amplification of the target-specific DNA.	No action is required.
No IPC signal is detected, but target-specific signal is detected in negative control wells.	Carryover contamination and one of the following: <ul style="list-style-type: none"> A high copy number of target DNA exists in samples, resulting in preferential amplification of the target-specific DNA. A problem occurred with IPC amplification. 	Examine unknowns to determine if an IPC signal is present. If an IPC signal is present in unknown wells, IPC amplification is not a problem. Repeat the assay using fresh aliquots of all reagents and clean pipetting equipment.
No IPC signal is detected, but target-specific signal is detected in unknown wells.	A high copy number of target DNA exists in samples, resulting in preferential amplification of the target-specific DNA.	No action required.
No IPC or target-specific signal is detected in positive control wells.	<ul style="list-style-type: none"> Environmental Master Mix was not stored properly. Target specific 10X Assay Mix not stored properly. 	Repeat the assay using properly stored assay components. Avoid freezing and thawing assay components. Protect the Assay Mix from light.

Observation	Possible cause	Recommended action
No target-specific signal is detected in positive control wells.	A pipetting error occurred (no positive control added).	Repeat the assay. Make sure to pipet the positive control into all positive control wells.
Target-specific signal is detected in negative control well.	Carryover contamination occurred.	<p>Repeat the assay using fresh aliquots of all reagents and clean pipetting equipment.</p> <p>If the negative control still shows contamination, repeat the assay using a new kit.</p> <p>If the negative control continues to show contamination, contact Life Technologies Technical Support.</p>
Replicate results for this sample are inconsistent.	All replicate wells for a sample do not have the same result. Replicates may have positive, negative, and/or warning results.	<p>If more than two replicates yield the same result (for example, you ran 3 replicates and 2 replicates are negative, but 1 replicate is positive), it is probable that the result with the larger number of replicates is accurate. However, your laboratory protocol may require that you repeat the assay using fresh samples and reagents.</p> <p>If you ran only two replicates and results are not consistent, repeat the assay using fresh samples and reagents.</p>

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

Microbiology of food and animal feeding stuffs — Horizontal method for the detection of *Salmonella* spp. Reference number ISO 6579:2002 (E).

Microbiology of food and animal feeding stuffs — General requirements and guidance for microbiological examinations. Reference number ISO 7218:2007.

Microbiology of food and animal feeding stuffs — Preparation of test samples, initial suspension and decimal dilutions for microbiological examination. Reference number ISO 6887.

Microbiology of food and animal feeding stuffs — Protocol for the validation of alternative methods. Reference number ISO 16140:2003.

2010. Association of Analytical Chemists (AOAC). <http://www.aoac.org>. Accessed November 2010.

2010. Association Francaise pour la Normalisation (AFNOR). <http://www.afnor-validation.com>. Accessed November 2010.



Prevent contamination and nonspecific amplification

Overview

PCR assays require special laboratory practices to avoid false positive amplifications. The high throughput and repetition of these assays can lead to amplification of a single DNA molecule.

PCR good laboratory practices

When preparing samples for PCR amplification:

- Maintain separate work areas, dedicated equipment, and supplies for:
 - Sample preparation
 - PCR setup
 - PCR amplification
 - Analysis of PCR products

Note: Rooms can be simulated using a clean bench or PCR bench available from major laboratory suppliers.

- Wear clean gloves and a clean lab coat (not previously worn while handling amplified PCR products or used during sample preparation).
- Change gloves whenever you suspect that they are contaminated and before leaving the work area.
- 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.
- Keep reactions and components capped as much as possible.
- Use positive-displacement pipettes or aerosol-resistant pipette tips.
- Clean lab benches and equipment after use with 10% bleach solution, followed by a 70% ethanol rinse to remove residual bleach.

IMPORTANT! To avoid false positives due to cross-contamination:

- Prepare and close all negative-control and unknown sample tubes before pipetting the positive control.
 - Do not open tubes after amplification.
 - Use different sets of pipettors when pipetting negative-control, unknown, and positive-control samples.
-

Plate layout suggestions

- Separate different targets by a row if enough space is available.
- Put at least one well between unknown samples and controls if possible.
- Separate negative and positive controls by one well if possible.
- Place replicates of one sample for the same target next to each other.
- Start with the unknown samples and put controls at the end of the row or column.
- Put positive controls in one of the outer rows or columns if possible.
- Consider that caps come in strips of 8 or 12.



WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
 - Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, etc). To obtain SDSs, see [“Obtaining SDSs” on page 33](#).
-

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below, and consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see “[Obtaining SDSs](#)” on page 33.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of 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.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the 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.



WARNING! CHEMICAL HAZARD. All chemicals in the instrument, including liquid in the lines, are potentially hazardous. Always determine what chemicals have been used in the instrument before changing reagents or instrument components. Wear appropriate eyewear, protective clothing, and gloves when using the instrument.



WARNING! CHEMICAL HAZARD. CHEMICAL STORAGE HAZARD. Never collect or store waste in a glass container because of the risk of breaking or shattering. Four-liter reagent and waste bottles can crack and leak. Each 4-liter 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.

Specific chemical handling



WARNING! CHEMICAL HAZARD. MicroSEQ® *Salmonella enterica* Detection Kit. Harmful if swallowed or inhaled. Causes eye, skin, and respiratory tract irritation. Do not taste or swallow. Avoid breathing vapor. Read the SDS and follow the handling instructions. Use with adequate ventilation. Avoid contact with eyes, mouth, and skin. Wear appropriate protective eyewear, clothing, and gloves.

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:

In the U.S.:

- U.S. Department of Health and Human Services guidelines published in Biosafety in Microbiological and Biomedical Laboratories found at: www.cdc.gov/biosafety
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030), found at: 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: www.cdc.gov

In the EU:

Check local guidelines and legislation on biohazard and biosafety precaution and refer to the best practices published in the World Health Organization (WHO) Laboratory Biosafety Manual, third edition, found at: www.who.int/csr/resources/publications/biosafety/WHO_CDS_CSR_LYO_2004_11/en/



Documentation and Support

Obtaining SDSs

Safety Data Sheets (SDSs) are available from www.lifetechnologies.com/support.

Note: For the SDSs of chemicals not distributed by Life Technologies, contact the chemical manufacturer.

Obtaining Certificates of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

Obtaining support

For the latest services and support information for all locations, go to:

www.lifetechnologies.com/support

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

Food safety support

Website: www.lifetechnologies.com/foodsafety

Support email: foodsafety@lifetech.com

Phone number (Inside North America): 1-800-500-6885

Phone number (Outside North America): Go to www.lifetechnologies.com/contactus.html and select the appropriate country from the drop-down menu.

Limited product warranty

Life Technologies and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

December 2012

