

Understanding Your Shipment

For use with single tube TaqMan® Assays

Pub. No. MAN0017153 Rev. A.0

Purpose

This document is provided to help you better understand the items included with your TaqMan® Assays shipment.

Items included in your shipment

Items included with every shipment:

- Dispatch Note (Packing List)—Contains your order information, a list of the assays that were ordered, and the quantity of the assays shipped.
- Box containing a storage rack with assay tubes.
Note: Assays from different TaqMan® product lines are grouped onto separate racks.
- Data Sheet—Provides a summary of the assays in your shipment, the assay details, and the storage conditions.

For further information about the assays, go to www.thermofisher.com/taqmanfiles. One or more of the following documents will be available for download, depending on the product ordered.

- Assay Information File
- User Instruction Documents (Protocols, User Guides, and Quick Reference Cards)
- Certificates of Analysis
- Safety Data Sheets
- Data Sheets

Obtain your documents online

1. Go to thermofisher.com/taqmanfiles.
2. Select **TaqMan® Assays**

Download TaqMan™ Assay / Plates & Cards Files

Select Type:

- TaqMan™ Assays ⓘ
- Fixed TaqMan™ Array Cards (TAC / TLDA) ⓘ
- Custom/Flexible TaqMan™ Cards (TAC / TLDA) ⓘ
- Fixed TaqMan™ Array Plates (TAP) ⓘ
- Custom/Flexible TaqMan™ Array Plates (TAP) ⓘ
- OpenArray™ Plates ⓘ

3. Enter your **Sales Order #** and **Rack/Plate ID**, then click **Submit**.

Download TaqMan™ Assay / Plates & Cards Files

Enter Sales Order # ⓘ

Enter Rack/Plate ID ⓘ [Clear All](#)

Select Document Type:

- Download the current Rack/Plate ID ⓘ
- Show all the Rack/Plate IDs ⓘ

[I do not have this information](#)

4. Select the type of files to download for your order.


TaqMan™ Assays

[Your Rack ID \(Assay type\)](#) [Reference Documents for \(your assay type\)](#)

Note: Files are available for download in PC and MAC formats.

Dispatch Note

The **Dispatch Note** contains information specific to your shipment. This document is in an envelope that is labeled **Packing List Enclosed**. The envelope is attached to the outside of your shipping container.

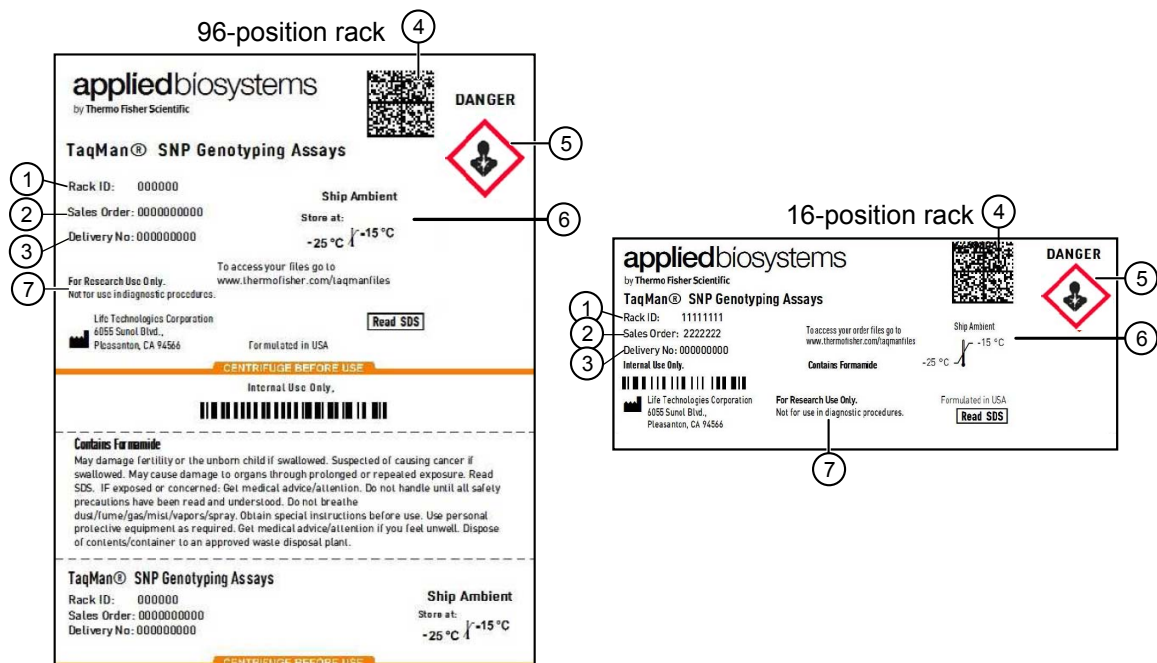
DISPATCH NOTE		Life Technologies Corporation			
Dispatch Note No : 8306917650 Manufacturing Sales Order No: 6415125		5781 Van Allen Way 92008 Carlsbad United States T: 18009556288 F: 18003312286			
					
① Delivery No : 862581857 Page : 1 of 2	DISPATCH TO : 215458				
ACCOUNT TO : 1 Accounts payable		ORDERING ENTITY AND ADDRESS Ordering entity and address			
CONTRACT NO: RELEASE NO:		ORDER PLACED BY: J. Doe PHONE NO : XXXXXXXXX			
FREIGHT TERMS : F02-FOB FACTORY - FRT PPD & ADD					
② Order Date Received 04/25/2017	Purchase Order Number 61491639	Dispatched On 04/28/2017	③ Carrier Federal Express SO		
④ Line No.	Catalogue No.	Description & Size	Lot Number	Quantity Shipped	⑧
0001	4448892	Hs00944495_g1 TAQMAN GENE EX A		1.000	
⑤ 0002	4451936	SSAYS MTO, FG, TAQMAN® GENE EX ASSAYS MTO,	2936626	1.000	
		⑥	⑦		

- ① Delivery number—A unique number that is assigned by Thermo Fisher Scientific to each customer shipment. This number matches the Delivery number that is listed on your box of assays.
- ② Purchase order number—Customer-provided Purchase Order (P.O.) Number.
- ③ Dispatched on—The date the product is packaged for shipment (mm/dd/yyyy).
- ④ Catalog No. 001 (top)—The number that is used for reordering the product, sometimes called the **Part Number**. This catalog number matches the catalog number on your assay tube.
- ⑤ Catalog No. 002 (bottom)—An internal number that is used for tracking purposes. This catalog number does not match the catalog number on your assay tube.
- ⑥ Description & Size—The name and size of the item ordered.
- ⑦ Lot Number—Manufacturing tracking number that is assigned by Thermo Fisher Scientific. This lot number does not match the lot number on your assay tube.
- ⑧ Quantity Shipped—The number of items in this shipment.

Note: Each item that is ordered displays two line items (duplicated under the **Quantity Shipped** heading). Only 1 item is included with your order.

TaqMan® Assays storage racks

Your shipment contains either 96- or 16-position racks. Racks are labeled with order-specific and product-specific information.



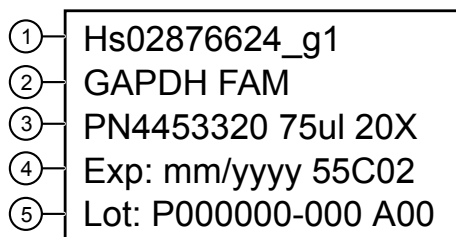
- ① Rack ID—A unique number that is assigned to each rack, either 96- or 16-position.
- ② Sales Order Number—A unique number that is assigned to each customer Sales Order.
- ③ Delivery Number—A unique number that is assigned to each customer shipment.
- ④ Barcode—Used by Thermo Fisher Scientific to track your order and shipment.
- ⑤ Warning—Information on potential hazards.
- ⑥ Storage Conditions—Instructions for product storage.

Note: Products are shipped at room temperature.

- ⑦ Regulatory Statement—For Research Use Only. Not for use in diagnostic procedures.

Tube labels


All tube labels contain five lines of text.



- ① Assay ID
- ② Product-specific information [for example, gene symbol, RS number, assay name, dye label, primer limited (PL)]
- ③ Part Number, volume, and assay concentration
- ④ Expiration Date and Internal Processing Code
- ⑤ Lot Number and Manufacturing Order Group Well Location

Data Sheets

The Data Sheet provides a summary of the assays in a specific rack, and supporting details for those assays.



① — **TaqMan™ Gene Expression Assays Data Sheet**

Sales Order: 36356650S8 **Order Date:** dd-mm-yyyy

Rack ID: P01966855 **Customer P.O.:** 10033082

Delivery No: 862633330 **Ship-To Country:** US

Customer: Jane Doe **Contact:**

Contact Email: jane.doe@email.com

Contact Phone: 000-000-0000

Number of Assays in this Rack: 1 **Number of Assays in this Order**
Made to Order: 1
Inventoried: 1

② — **Assays in this Shipment**

Tube Position	Assay ID*	Gene Symbol*	Part Number
A01	Hs02786624_g1	GAPDH	4453320

* "+" sign appears for entries greater than 20 characters. Full entry recorded in Assay Information File (AIF)

Figure 1 Data Sheet—Page 1

- ① A summary of your order, including the order date, customer P.O. and contact information, Sales Order and Rack ID Numbers, number of assays in the rack, and the number of assays ordered.
- ② Assays in this Shipment—Identification of the assays in your shipment by the Assay ID, target name, tube position in the rack, and Part Number.

Other Details

① Assay Details

Part Number	Product	Reporter / Quencher	Volume (μL)	Formulation	No. of Reactions (20μL Reaction Size)
4331182	TaqMan™ Gene Expression Assays, Inventored, SM	FAM/MGB-NFQ	250	20x Mix (TE)	250
4448892	TaqMan™ Gene Expression Assays, XS	FAM/MGB-NFQ	75	20x Mix (TE)	75
4351372	TaqMan™ Gene Expression Assays, SM	FAM/MGB-NFQ	360	20x Mix (TE)	360
4351370	TaqMan™ Gene Expression Assays, MED	FAM/MGB-NFQ	750	20x Mix (TE)	750
4351368	TaqMan™ Gene Expression Assays, LG	FAM/MGB-NFQ	967	60x Mix (TE)	2,900
4448489	TaqMan™ Gene Expression Assays, SM	VIC/MGB-NFQ	360	20x Mix (TE)	360
4448490	TaqMan™ Gene Expression Assays, MED	VIC/MGB-NFQ	750	20x Mix (TE)	750
4448491	TaqMan™ Gene Expression Assays, LG	VIC/MGB-NFQ	967	60x Mix (TE)	2,900
4448484	TaqMan™ Gene Expression Assays, Primer-Limited, SM	VIC/MGB-NFQ	360	20x Mix (TE)	360
4448485	TaqMan™ Gene Expression Assays, Primer-Limited, MED	VIC/MGB-NFQ	750	20x Mix (TE)	750
4448486	TaqMan™ Gene Expression Assays, Primer-Limited, LG	VIC/MGB-NFQ	967	60x Mix (TE)	2,900

② Storage Conditions

-25°C to -15°C

③ Supporting Documentation

Supporting documentation and files associated with this product are available for download at thermofisher.com/taqmanfiles:

- User Guide: "How to" guide for performing the application
- Assay Information File (AIF): Technical details for all assays in the shipment
- Certificate of Analysis: Validation results for your assays
- Data Sheet: Electronic version of this document
- Safety Data Sheet (SDS): Safety guidelines
- Quick Reference Card (QRC): Summary of the User Guide
- Understanding Your Shipment: Directory for understanding the contents of your shipment

④ Additional Quality Information

A standard certificate of analysis is included in the supporting documentation for this product. For additional quality information, visit thermofisher.com/support.

⑤ Technical Support

Visit www.thermofisher.com/contactus.

Limited product warranty

Life Technologies Corporation 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.thermofisher.com/us/en/home/global/terms-and-conditions. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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

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

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Limited Use Label License No. 537: Minor Groove Binder (MGB) in Real-Time PCR for Research Use

Figure 2 Data Sheet—Pages 2 and 3

- ① Other Details—The formulation, product description, reporter/quencher, fill volume, concentration, and number of reactions.
- ② Storage Conditions—Instructions for product storage.
- ③ Supporting Documentation—Additional assay-related documentation that can be downloaded from thermofisher.com/taqmanfiles.
- ④ Additional Quality Information—Instructions for obtaining additional quality-related information.
- ⑤ Technical Support—Instructions for obtaining Thermo Fisher Scientific Technical Support.

Assay Information File

The Assay Information File is provided in TXT format, and/or in XML and HTML formats, depending on the product line and order date.

- HTML format files are provided as a reference for users; open them in a web browser.
- XML and TXT format files are provided primarily for electronic data importation and manipulation.

Table 1 Definitions for Assay Information File fields (alphabetical)

Field	Description
AB Minor Allele Frequency—African-American	Minor allele frequency in a collection of 45 African American samples, as calculated by Applied Biosystems™
AB Minor Allele Frequency—Caucasian	Minor allele frequency in a collection of 45 Caucasian samples, as calculated by Applied Biosystems™
AB Minor Allele Frequency—Chinese	Minor allele frequency in a collection of 45 Chinese samples, as calculated by Applied Biosystems™
AB Minor Allele Frequency—Japanese	Minor allele frequency in a collection of 45 Japanese samples, as calculated by Applied Biosystems™
Amino Acid Change	(TaqMan® SNP Genotyping Assays only) The amino acid change associated with a coding SNP
Amplicon Size	Size of the PCR amplification product
Annotation Source	Public annotation source used for a specific product family
Assay ID	Alphanumeric string that identifies the assay
Assay Mix Concentration	Concentration of the assay mix (including all primers and probe components)
Assay Name	For custom products, the customer-designated name of the sequence submitted; for predesigned assays, the Life Technologies–designated name assigned at the time of assay design
Category	(TaqMan® Copy Number Assays only) Chromosome(s) targeted by the assay, based on Celera* Panther Protein Classification, Level 1
Category ID	(TaqMan® Copy Number Assays only) Unique 10-digit ID for the category name
Celera ^[1] Assembly Build Number	(TaqMan® SNP Genotyping Assays only) The version of the Celera assembly from which the coordinate position is taken
Celera ID (for Predesigned assays)	For Predesigned TaqMan ^{fi} Gene Expression and SNP Genotyping Assays, this refers to the Celera ID—a unique transcript or SNP ID, respectively, in the Celera Discovery System (CDS)
Celera SNP ID	(TaqMan® SNP Genotyping Assays only) The hCV SNP number in the Celera Discovery System (CDS)
Chromosome	Chromosome on which the assay target (e.g., gene, SNP, intergenic region) is found
Contact Name	Customer contact name provided with order submission
Component Type	For kit products, this differentiates TM (TaqMan® Assay) or RT (Reverse Transcription primer) component
Context Sequence	Nucleotide sequence surrounding the TaqMan® probe. For TaqMan® SNP Genotyping Assays, the sequence surrounding the target SNP. The first nucleotide in the brackets corresponds to the VIC™ allele and the second to the FAM™ allele. For TaqMan® MicroRNA Assay, the mature microRNA sequence itself.
Customer Name	Name of the customer's organization or institution
Cytogenetic Band	Cytoband/chromosomal band location of assay target sequence
Delivery Number	Unique number assigned to each customer shipment
Design Strand	Strand used to design the TaqMan® probe
DGV ^[2] Database Version	(TaqMan® Copy Number Assays only) Reference genome assembly version used by the Database of Genomic Variants (DGV)
DGV** Locus Location	(TaqMan® Copy Number Assays only) Position on the reference genome of the DGV Locus
DGV** Variation ID	(TaqMan® Copy Number Assays only) Target DGV Variation ID
DGV** Variation Type	(TaqMan® Copy Number Assays) Type of genomic variation; for example, Copy Number
Distance to Stem-Loop (TaqMan® MicroRNA Assays only)	Distance from TaqMan® MicroRNA Assay context sequence to the base of the stem-loop sequence
Entrez Gene ID	The Entrez Gene database numeric identifier for the interrogated gene (www.ncbi.nlm.nih.gov/gene)

Field	Description
Expiration Date	Expiration date of the assay
Forward Primer Concentration	Concentration (in μM) of the forward PCR primer
Forward Primer Name	Name of the forward PCR primer, given by the design software; consists of the Assay ID or Assay Name, with an F suffix
Forward Primer Sequence	Nucleotide sequence of the forward PCR primer
Gene Name	Gene name in the Entrez Gene database
Gene Symbol	Entrez Gene symbol for the interrogated gene
Group	(Copy number assays only) Based on Celera* Panther Protein Classification, Level 2
Group ID (TaqMan® Copy Number Assays only)	Unique 10-digit ID for the group name
HapMap Minor Allele Freq—CEU (TaqMan® SNP Genotyping Assays only)	Minor allele frequency in CEPH (Utah residents with ancestry from northern and western Europe), as calculated by the HapMap Project
HapMap Minor Allele Freq—CHB (TaqMan® SNP Genotyping Assays only)	Minor allele frequency in Han Chinese in Beijing, China, as calculated by the HapMap Project
HapMap Minor Allele Freq—JPT (TaqMan® SNP Genotyping Assays only)	Minor allele frequency in Japanese in Tokyo, Japan, as calculated by the HapMap Project
HapMap Minor Allele Freq—YRI (TaqMan® SNP Genotyping Assays only)	Minor allele frequency in Yorubans in Ibadan, Nigeria, as calculated by the HapMap Project
Location on Celera* Assembly (SNP genotyping assays only)	The nucleotide location on the Celera* human genome assembly (as referenced)
Location on NCBI Assembly	Nucleotide location on the NCBI genome assembly of the assay target. It is near the base location in the middle of the context sequence
Lot Number	Unique identifier assigned to each manufacturing lot
Location on Transcript or Gene	The assay's nucleotide location on the target transcript or gene. For TaqMan® Gene Expression Assays, it is the target location on the first RNA accession that is listed. For TaqMan® Copy Number Assays, it is the exon and/or intron gene location
Mapping Date	Date that Thermo Fisher Scientific completed assay mapping
Mature MicroRNA Accession	(TaqMan® MicroRNA Assays only) miRBase ^[3] microRNA name
Medline Reference	PubMed references for the gene
Minor Allele Frequency—African-American	(TaqMan® SNP Genotyping Assays only) Minor allele frequency in African-Americans, as calculated by SNP genotyping at Thermo Fisher Scientific
Minor Allele Frequency—Caucasian	(TaqMan® SNP Genotyping Assays only) Minor allele frequency in Caucasians, as calculated by SNP genotyping at Thermo Fisher Scientific
Minor Allele Frequency—Chinese	(TaqMan® SNP Genotyping Assays only) Minor allele frequency in Chinese people, as calculated by SNP genotyping at Thermo Fisher Scientific
Minor Allele Frequency—Japanese	(TaqMan® SNP Genotyping Assays only) Minor allele frequency in Japanese people, as calculated by SNP genotyping at Thermo Fisher Scientific
NCBI Assembly Build Number	The version of the NCBI reference genome assembly used for assay target location
NCBI Gene Reference	NCBI transcript ID that is detected by the assay
NCBI Reference SNP ID	(TaqMan® SNP Genotyping Assays only) The reference SNP number (RS Number) in the NCBI-dbSNP database
NCBI SNP Reference	(TaqMan® SNP Genotyping Assays only) Reference from the NCBI-dbSNP database
Order Number	A customer sales order number assigned by Thermo Fisher Scientific, unique to an order
Part Number ^[4]	Product number used in conjunction with the Assay ID to order the assay
Product Type	Thermo Fisher Scientific product category or family (corresponds with the part number)
Rack or Plate ID	Unique number on the label of each rack or plate; also referred to as Rack ID

Field	Description
Rack or Plate Type	Type of container in which the assay is shipped (e.g., 96- or 16-position rack)
Reporter 1 Concentration	Concentration (in μM) of reporter 1 TaqMan [®] probe
Reporter 1 Dye	Dye label for reporter 1 TaqMan [®] probe
Reporter 1 Name	Name for the reporter 1 TaqMan [®] probe, determined by the design software. It consists of the Assay ID or Assay Name, with a suffix that contains a letter to indicate the dye label (M for FAM [™] dye and V for VIC [™] dye). Some assays also include a number to indicate which DNA strand was used to design the probe: 1 indicates that the forward strand was used to design the probe; 2 that it was the reverse strand; and 3 that the design strand is not identified.
Reporter 1 Quencher	Quencher used for reporter 1 TaqMan [®] probe, for example, nonfluorescent quencher (NFQ)
Reporter 1 Sequence	Nucleotide sequence of reporter 1 TaqMan [®] probe
Reporter 2 Concentration	(TaqMan [®] SNP Genotyping Assays only) Concentration (in μM) of reporter 2 TaqMan [®] probe
Reporter 2 Dye	(TaqMan [®] SNP Genotyping Assays only) Dye label for reporter 2 TaqMan [®] probe
Reporter 2 Name	Name for the reporter 2 TaqMan [®] probe, determined by the design software. It consists of the Assay ID or Assay Name, with a suffix that contains a letter to indicate the dye label (M for FAM [™] dye). Some assays also include a number to indicate which DNA strand was used to design the probe: 1 indicates that the forward strand was used to design the probe; 2 that it was the reverse strand; and 3 that the design strand is not identified.
Reporter 2 Quencher	(TaqMan [®] SNP Genotyping Assays only) Quencher used for reporter 2 TaqMan [®] probe
Reporter 2 Sequence	(TaqMan [®] SNP Genotyping Assays only) Nucleotide sequence of reporter 2 TaqMan [®] probe
Reverse Primer Concentration	Concentration (μM) of the reverse PCR primer
Reverse Primer Name	Name for the reverse primer, given by the design software; consists of the Assay ID or Assay Name, with an R suffix
Reverse Primer Sequence	Nucleotide sequence of the reverse PCR primer
RNA Database ^[5] Gene Symbol	The RNA database gene symbol for the interrogated gene product
RNA Database ID	The RNA database identifier for the interrogated gene
RT Primer Sequence	Sequence of the RT primer for the interrogated transcript
Sanger dB Link	(TaqMan [®] MicroRNA Assays only) Link to microRNA page on the miRBase ^[3] database website (based on accession #)
Sales Order Number	A customer sales order number assigned by Thermo Fisher Scientific, unique to an order
Ship Date	Date that the product is shipped
Shipment ID	Unique number assigned to each customer shipment
(Shipping) Rack or Plate ID	Unique number on the label of each rack or plate; also referred to as Rack ID
(Shipping) Rack or Plate Type	Type of container in which the assay is shipped (e.g., 96- or 16-position rack)
SNP Type	(TaqMan [®] SNP Genotyping Assays only) Type of SNP, e.g., intron, acceptor splice site, donor splice site, intergenic/unknown, etc.
Species	Organism for which the assay was designed
Stem-Loop Chromosome Location	(TaqMan [®] MicroRNA Assays only) Genomic position of stem-loop sequence
Stem-Loop ID	(TaqMan [®] MicroRNA Assays only) miRBase ^[3] stem-loop alphanumeric identifier
Target Exons	Exons (as referenced by the public accession number) that are spanned by the probe
Target Type	Target sequence identifier (for example, Gene Symbol, Reference SNP (RS) Number, etc.)
Target Value	Publicly available gene or RNA name, or customer-provided identifier
Transcript Accession	The NCBI alphanumeric sequence identifier for the interrogated transcript
Tube/Well Position	Well location of the assay in the rack or plate
Vial (or Tube) ID	10-digit number that uniquely identifies each tube

Field	Description
Vial (or Tube) Type	Type of vial or tube containing the assay
Well Location	Well location of the assay in the rack or plate
XML Template Version	Version of the XML template used for the Assay Information File

^[1] Historical reference to Celera Discovery System

^[2] Database of Genomic Variants (DGV), <http://projects.tcag.ca/variation/>. Iafrate AJ, Feuk L, Rivera MN et al. (2004) Detection of large-scale variation in the human genome. *Nat Genet* 36(9):949–51.

^[3] miRBase microRNA database, <http://www.mirbase.org>. Griffiths-Jones S, Saini HK, van Dongen S, Enright AJ (2008) *Nucl Acids Res* 36(Database Issue):D154–D158.

^[4] Part numbers for Predesigned, Custom, and Custom Plus Assays reflect assay scale and whether it is a human or nonhuman target.

^[5] Mammalian noncoding RNA database (RNADB), <http://research.imb.uq.edu.au/rnadb/>. Pang KC, Stephen S, Engstrom PG et al. (2005) RNADB - a comprehensive mammalian noncoding RNA database. *Nucl Acids Res* 33 (Database Issue): D125–D130.


User Instruction Documents

A specific set of user instruction documents is provided with each TaqMan[®] Assay product. These documents can include protocols, quick reference cards, and user guides for products and software. For example, user documents that are provided with TaqMan[®] Gene Expression Assays and Custom TaqMan[®] Gene Expression Assays include the following:

- User Protocol—A detailed guide that provides step-by-step instructions for using the product.
- Quick Reference Card—A shorter summary of the key steps in the User Protocol. This guide is designed for use by experienced users.
- Design and Ordering Guide—A guide that provides step-by-step instructions for designing and ordering Custom TaqMan[®] Gene Expression Assays and Custom TaqMan[®] SNP Genotyping Assays.

Certificates of Analysis

The Certificate of Analysis contains assay quality information.



applied biosystems
by Thermo Fisher Scientific

Certificate of Analysis

TaqMan® Gene Expression Assays

Sales Order: 36356650S8
Delivery Number: 862633330
Ship Date: 07-JUN-2017
Page: 1 of 1

①
②
③

④ Part Number: 4453320 ⑤ Part Description: TaqMan® Gene Expression Assays, Inventoried, XS

⑥ Assay ID: Hs02786624_g1

⑦ Gene Symbol: GAPDH

⑧ Lot Number: P170507-001A02

⑨ Expiration Date: 06-MAY-2022

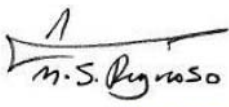
Vial Type	Quality Control Test	Result
TaqMan (TM)	Fill Volume	PASS
TaqMan (TM)	Mass Spectrometry	PASS
TaqMan (TM)	No Template Control	PASS

Applied Biosystems' oligo manufacturing process requires a fill volume check on each TM assay lot to verify that the volume is consistent with the required volume printed on the product label. Applied Biosystems' oligo manufacturing process also requires a mass spectrometry test for each component of the TM assay to verify that the identity of each oligo meets set specifications. A no template control (NTC) test is conducted on the assay level with a pass = Ct > 35 (18S rRNA assays are an exception). A "*" sign appears for Assay IDs and Gene Symbols greater than 20 characters. Full entry recorded in Assay Information File (AIF).

⑪ For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

ISO 9001
ISO 13485
REGISTERED

Pleasanton, CA



Quality Assurance

To obtain a Certificate of Analysis on-line go to www.thermofisher.com/coa or email us at cofarequest@lifetech.com

Doc p/n: 4440784 rev.D

- ① Sales Order—A unique number that is assigned by Thermo Fisher Scientific to each customer Sales Order.
- ② Delivery Number—A unique number that is assigned by Thermo Fisher Scientific to each customer shipment.
- ③ Ship Date—The date the product is packaged for shipment (mm/dd/yyyy).
- ④ Part Number—The number that is used for (re)ordering the product; sometimes called the Product Number.
- ⑤ Part Description—The unique name for part and scale; based on Part Number.
- ⑥ Assay ID—A unique identifier that is assigned by Thermo Fisher Scientific to each assay.
- ⑦ Assay-Related Target—Field heading and content varies by TaqMan® Assay type (for example, Gene Symbol for gene expression assays, RS Number for SNP assays, Gene Identifier for non-coding RNA, or Assay Name for custom, small RNA, or pri-miRNA assays).
- ⑧ Lot Number—A unique identifier that is assigned to each manufacturing batch; allows Thermo Fisher Scientific to identify the manufacturing site and date of manufacture.
- ⑨ Expiration Date—The date after which the material should no longer be used by the customer.
- ⑩ Quality Control—List of quality control tests, confirmation that they were completed, and results [pass/fail/previously passed (TaqMan® SNP Genotyping Assays only)].
- ⑪ Regulatory Statement—For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

Safety Data Sheets

The material Safety Data Sheet (SDS) provides important safety information about your shipment. It includes information on the composition and properties of hazardous materials in the shipment; handling, storage, and disposal of these materials; and regulatory information by geographic location.


ThermoFisher SCIENTIFIC		Safety Data Sheet	
SECTION 1: Identification of the substance/mixture and of the company/undertaking			
<u>Identification of the substance or mixture</u>			
Product code	4331349, 4332072		
Product name	TaqMan Assays		
Synonyms	4332075, 4332077, 4332079, 4399989, 4400293, 4400296, 4426963, 4427014, 4440880, 4440888, 4441118, 4442488, 4459411, 4351368, 4448481, 4448486, 4351379, 4351384, 4351376, 4351382, 4448492, 4448510, 4448513, 4448516, 4362691		
<u>Company/undertaking identification</u>			
Life Technologies Corporation 5781 Van Allen Way PO Box 6482 Carlsbad, CA 92008 +1 760 603 7200	Life Technologies 5250 Mainway Drive Burlington, ONT CANADA L7L 6A4 800/263-6236		
24 hour Emergency Response for Hazardous Materials Within the USA + Canada: 1-800-424-9300 and +1 [or Dangerous Goods] Incident. Spill, Leak, Fire, 703-527-3887 Exposure, or Accident. Call CHEMTREC Outside the USA + Canada: +1 703-741-5970			
Country specific Emergency Number (if available): CHEMTREC Brazil (Rio De Janeiro) +(55)-2139581449 (português)			
Research Use Only. Not for use in diagnostic procedures.			
SECTION 2: Hazards identification			
<u>GHS - Classification</u>			
Signal Word DANGER			
			
<u>Health hazards</u>			
Specific target organ systemic toxicity (repeated exposure)	Category 2		
Carcinogenicity	Category 2		
Reproductive Toxicity	Category 1B		
Revision date	22-Aug-2016	Page	1 / 7
Product code	4331349, 4332072	Product name	TaqMan Assays
www.thermofisher.com			

Figure 3 Sample Safety Data Sheet

Limited product warranty

Life Technologies Corporation 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.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Manufacturer: Multiple Life Technologies Corporation manufacturing sites are responsible for manufacturing the products associated with the workflow covered in this guide.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision History: Pub. No. MAN0017153

Revision	Date	Description
A.0	13 July 2017	New document. Supersedes publication number C016152 1210 Rev E.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.