

QuantiGene™ Singleplex Assay Kit

Catalog Number QS0012, QS0013, QS0014, QS0016, and QS0384

Pub. No. MAN0019612 Rev. C



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

The QuantiGene Singleplex Assay uses branched DNA technology to quantify RNA directly from a variety of sample types without RNA purification or target amplification. The QuantiGene Singleplex Assay is comprised of 3 modules:

- QuantiGene Singleplex Assay Kit
- QuantiGene Singleplex target-specific probe sets
- QuantiGene Sample Processing Kit (not required for purified RNA samples)

Contents and storage

Refer to box label for expiration date.

Item	Description	Cat. No. QS0012	Cat. No. QS0013	Cat. No. QS0014 , QS0016	Cat. No. QS0384	Storage
Blocking Reagent	Aqueous buffered solution containing a preservative	400 µL	1.8 mL	10 mL	2 x 10 mL	2–8°C
Pre-Amplifier Solution	DNA in an aqueous buffered solution	23 mL	120 mL	550 mL	3 x 400 mL	
Amplifier Solution	DNA in an aqueous buffered solution	23 mL	120 mL	550 mL	3 x 400 mL	
Capture Plate	96-well polystyrene plate coated with capture probes	2 each ^[1]	10 each ^[1]	50 each ^[2]	50 each ^[1]	
Label Probe Solution	Oligonucleotide-alkaline phosphatase conjugate in aqueous buffered solution	23 mL	120 mL	550 mL	3 x 400 mL	
Substrate	Chemiluminescent substrate ^[3]	24 mL	120 mL	530 mL	4 x 300 mL	15–30°C
Lysis Mixture	Aqueous buffered solution containing a preservative	18 mL	93 mL	410 mL	510 mL	
Wash Buffer Component 1	Aqueous solution	3 mL	15 mL	75 mL	2 x 75 mL	
Wash Buffer Component 2	Aqueous buffered solution	5 mL	25 mL	125 mL	2 x 125 mL	
Plate Seals Day 1	Adhesive backed foil plate seal	2 ea	10 ea	50 ea	260 ea	
Plate Seals Day 2	Adhesive backed foil plate seal	10 ea	50 ea	210 ea	–	

^[1] Captures plates in QS0012, QS0013 and QS0384 are packaged individually.

^[2] Capture plates in QS0014 are packaged individually. QS0016 has 5 packs of 10 plates.

^[3] Lumigen™ APS-5

Safety warnings and precautions

All chemicals should be considered potentially hazardous. This product and its components should be handled by those trained in laboratory techniques and be used according to the principles of good laboratory practice.

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates

This QuantiGene™ Singleplex Assay Kit contains all of the general reagents required to run 192, 960, or 4800 QuantiGene Singleplex assays, depending on the kit. Sample processing kits and target-specific probe sets are sold separately.

Product use

QuantiGene™ Singleplex Assay Kit is designed for use with QuantiGene Singleplex Probe Sets and QuantiGene Sample Processing Kits for quantitation of target-specific RNA or DNA from cultured cell or whole blood lysates, fresh, frozen, or FFPE tissue homogenates, or purified RNA. For more information, refer to the *QuantiGene Singleplex Assay User Manual*. For research use only. Not for use in diagnosis of disease in human or animals.

- Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions



Revision history: Pub. No. MAN0019612 C

Revision	Date	Description
C	30 August 2024	Created Rev C, added new catalog number QS0384.
B.0	16 February 2021	Created Rev B.0, updated quantity volumes in Contents and storage table.
A.0	30 September 2020	New document for <i>QuantiGene™ Singleplex Assay Kit Product Information Sheet</i> .

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

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

©2020-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.