

QuantiGene Singleplex Assay Kits

Cat. Nos. QS0008, QS0009, QS0010, QS 0011, QS0018, and QS0019

About QuantiGene Singleplex Assay

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 reagent system is comprised of 3 modules:

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

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

Intended 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, please refer to the *QuantiGene Singleplex Assay User Manual*. For research use only. Not for use in diagnosis of disease in human or animals.

Contents and Storage

Component	Description	Quantity QS0008 QS0018	Quantity QS0009 QS0019	Quantity QS0010 QS0011	Storage
Blocking Reagent	Aqueous buffered solution containing a preservative	400 µL	1.8 mL	10 mL	-20 °C
Pre-Amplifier	DNA in an aqueous buffered solution	24 µL	120 µL	550 µL	-20 °C
Amplifier	DNA in an aqueous buffered solution	24 µL	120 µL	550 µL	-20 °C
Capture Plate	96-well polystyrene plate coated with capture probes	2 ea*	10 ea*	50 ea†	2-8 °C
Label Probe	Oligonucleotide-alkaline phosphatase conjugate in aqueous buffered solution	24 µL	120 µL	550 µL	2-8 °C
Substrate	Chemiluminescent substrate‡	24 mL	120 mL	530 mL	2-8 °C
Amplifier/Label Probe Diluent	Aqueous buffered solution with a protein-containing preservative	72 mL	360 mL	1,800 mL	15-30 °C
Lysis Mixture	Aqueous buffered solution containing a preservative	18 mL	93 mL	410 mL	15-30 °C
Wash Buffer Component 1	Aqueous solution	3 mL	15 mL	75 mL	15-30 °C
Wash Buffer Component 2	Aqueous buffered solution	5 mL	25 mL	125 mL	15-30 °C
Plate Seals	Adhesive backed foil plate seal	12 ea	60 ea	260 ea	15-30 °C

* Captures plates in QS0008, QS0018, QS0009, & QS0019 are packaged individually.

† Capture plates in QS0010 are packaged individually. QS0011 has 1 pack of 50 plates.

‡ Lumigen® APS-5

Safety Warnings and Precautions

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

www.affymetrix.com/contact Please visit our website for international distributor contact information.
For research use only. Not for use in diagnostic procedures.

P/N 13069, 13070, 13071 Rev. C 021716

© 2016 Affymetrix, Inc. All rights reserved. Affymetrix® is a registered trademark of Affymetrix, Inc. QuantiGene® is a registered trademark exclusively licensed to Affymetrix, Inc. All other trademarks are the property of their respective owners.