

QuantiGene ViewRNA ISH Tissue Positive Control Kit (2-Plex)

Cat. No. QVT0516

Product Overview

The QuantiGene ViewRNA ISH Tissue Positive Control Kit is designed for use with QuantiGene ViewRNA ISH Tissue 2-Plex Assays. This kit includes 5 ± 1 μm thick rat kidney FFPE sections attached to positively charged slides as well as TYPE 1 rat Spp1 and TYPE 6 rat Ubc Probe Sets. Refer to the appropriate QuantiGene ViewRNA ISH Tissue Assay User Manual, based on equipment used, for instructions on performing *in situ* hybridization assays.

Contents and Storage

The QuantiGene ViewRNA ISH Tissue Positive Control Kit contains the following components. The Rat Kidney Tissue Slides have a shelf life of 3 months and the QuantiGene ViewRNA Probe Sets have a shelf life of 1 year from date of receipt when stored as recommended.

Table 1 QuantiGene ViewRNA ISH Tissue Positive Control Kit Contents and Storage

Component	Description	Quantity	Storage
Rat Kidney FFPE Tissue Slide	5 μm -thick rat kidney FFPE section attached to a positively, charged slide.	4	-20 °C
QuantiGene ViewRNA TYPE 1 Rat Spp1 Probe Set (VC1-10477)	120 μL of a TYPE 1 rat Spp1 Probe Set. See individual Product Insert for Probe Set specificity.	1	-20 °C
QuantiGene ViewRNA TYPE 6 Rat Ubc Probe Set (VC6-10447)	120 μL of a TYPE 6 rat Ubc Probe Set. See individual Product Insert for Probe Set specificity.	1	-20 °C

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.

Recommendations for Use

Allow slides to thaw for 30 minutes at room temperature. Perform one positive (rat Spp1 and Ubc probe sets) and one negative (omit probe sets) control assay using a total of 2 rat kidney tissue sections. For these control FFPE slides, use a 10 minute pretreatment boiling time and a 20 minute Protease incubation time as described in the QuantiGene ViewRNA ISH Tissue Assay User Manual.

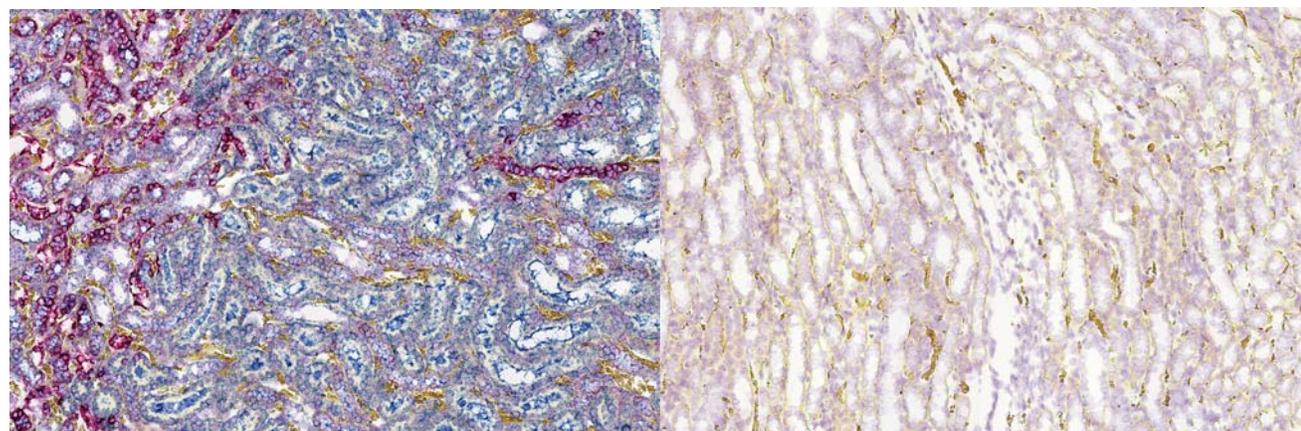
Expected Results

Figure 1 shows the expected results from the control assay using Spp1 and Ubc probe sets and rat kidney FFPE tissue sections. Images are the 20X bright-field image of the rat kidney staining with the Spp1 (red) and Ubc (blue) probes (positive control) and with no probe (negative control). Red color is Fast Red chromogenic precipitate of the labeled Spp1 mRNA, aqua blue color is Fast Blue chromogenic precipitate of the labeled Ubc mRNA, purplish-blue color is the hematoxylin staining of the nuclei, and yellow color is the red blood cells in the sample.

The positive control slide should exhibit aqua blue signal for Ubc and a saturating red signal in a subpopulation of cells for Spp1. The nucleus, stained with hematoxylin, should be uniformly stained purplish-blue and exhibit good morphology.

The negative control slide should exhibit a clean background of less than 1 dot per 10 cells.

Figure 1 Expected results from a positive control assay using Spp1 and Ubc probe sets. Left image is the positive control with Spp1 (red) and Ubc (blue) probes. Right image is the negative control which has gone through the assay procedure without target probes.



US Headquarters

Affymetrix, Inc.
3420 Central Expressway
Santa Clara, CA 95051
Tel: 1-888-362-2447
Direct: 1-408-731-5000
Fax: 1-408-731-5380
Email: info@affymetrix.com
Email: orders_fremont@affymetrix.com
Email: pqbhelp@affymetrix.com

European Headquarters

Panomics Srl
Via Sardegna 1
20060 Vignate-Milano (Italy)
Tel: +39-02-95-360-250
Fax: +39-02-95-360-992
Email: info_europe@affymetrix.com
Email: order_europe@affymetrix.com
Email: techsupport_europe@affymetrix.com

Asia Pacific Headquarters

Affymetrix Singapore Pte Ltd.
No.7 Gul Circle #2M-01/08
Keppel Logistics Building
Singapore 629563
Tel: +65-6395-7200
Fax: +65-6395-7300
Email: info_asia@affymetrix.com
Email: order_asia@affymetrix.com
Email: techsupport_asia@affymetrix.com

www.affymetrix.com/panomics

For research use only. Not for use in diagnostic procedures.

P/N 17405 Rev.1 121016

© 2012 Affymetrix, Inc. All rights reserved. Affymetrix® and QuantiGene® are registered trademarks of Affymetrix, Inc. All other trademarks are the property of their respective owners.