



Agilent Gene Expression Wash Buffer Kit

Preparation Guide

Add Triton X-102 detergent to the Gene Expression Wash Buffer 1 and 2 to reduce the chance of wash-related background artifacts. This preparation step is *required* for miRNA profiling applications and *optional* for gene expression applications.

To add Triton X-102 to the wash buffers

Add Triton X-102 to both wash buffer 1 and 2 when the cubitainer is first opened.

- 1 Open the cardboard box with the cubitainer of wash buffer and carefully remove the outer and inner caps of the cubitainer.
- 2 Use a pipette to put 2 mL of the provided 10% Triton X-102 into the wash buffer in the cubitainer.
- 3 Replace the original inner and outer caps and mix the buffer carefully but thoroughly by inverting the cubitainer 5 to 6 times.
- 4 Carefully remove the outer and inner caps and install the spigot provided with the wash buffer.
- 5 Prominently label the wash buffer box to indicate that Triton X-102 has been added and indicate the date the Triton X-102 was added.

To add Triton X-102 to small amounts of wash buffer

- 1 Add the desired amount of wash buffer to a clean graduated cylinder.
- 2 Add 10% Triton X-102 to final concentration of 0.005%. (For example, for 500 mL, add 250 μ L 10% Triton X-102.)
- 3 Transfer the buffer to a clean sterile bottle and mix thoroughly.



NOTE

Follow good laboratory practices. Wear appropriate personal protective equipment when using these products. Maintain a clean and uncluttered work area. Read protocols prior to initiating procedures. Calibrate pipettes at least yearly to avoid over-drawing.

Certain wash related background artifacts have been observed when using Agilent's gene expression wash buffer. These background artifacts can cause degraded assay performance. The addition of Triton X-102 to the wash buffer significantly reduces the likelihood of such artifacts with no loss of data quality or change to buffer stringency.

Artifacts are more pronounced in miRNA profiling experiments and less pronounced when used in gene expression profiling experiments. Agilent recommends for the addition of Triton X-102 to gene expression buffers:

- miRNA profiling application – mandatory
- gene expression application – optional

See www.agilent.com/chem/washbufferupdate for details.

Limited Use License

This product is distributed and sold to the End-User for life science research and commercial applications, but *not for diagnostic use*. End user does not have a right to resell or transfer the nucleotide analog components of this product either alone or as components of another product. Any use of this product other than for life science research and commercial applications is strictly prohibited.

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

© Agilent Technologies, Inc. 2007, 2015

Revision A1



5190-0448



Agilent Technologies