



Agilent Low-Input QuickAmp Labeling Kit and Agilent Low-Input QuickAmp WT Labeling Kit

Product Information

The Agilent Low-Input QuickAmp (LIQA) and Agilent Low-Input QuickAmp WT (LIQA WT) Labeling Kits include reagents to amplify and label RNA samples for hybridization to gene expression (for LIQA) or exon (for LIQA WT) oligo microarrays. This product insert applies to:

- Low-Input QuickAmp Labeling Kit (p/n 5190-2305, 5190-2306, 5190-2307 and 5190-2308). Refer to the *Agilent One Color Microarray-Based Gene Expression Analysis Protocol* or the *Agilent Two Color Microarray-Based Gene Expression Analysis Protocol* versions 6.0 or later.
- Low-Input QuickAmp WT Labeling Kit (p/n 5190-2942, 5190-2943, 5190-2944, and 5190-3386). Refer to the *Agilent One Color Microarray-Based Exon Analysis Protocol* or the *Agilent Two Color Microarray-Based Exon Analysis Protocol* versions 1.0 or later.

Protocols are available at www.agilent.com/chem/dnamanuals-protocols and included on the Microarray Kit Data Disk.

Storage

Store kit containing enzyme and WT Primer Mix (when included) in a non-frostfree -20°C freezer. Keep all components on ice while in use, minimizing the amount of time on ice. Store Cyanine CTPs at 4°C after initial use to avoid multiple freeze-thaw cycles. For long-term storage, store Cyanine CTPs at -20°C.

Components

All kits contain the components below. The 5190-2306 and 5190-2944 two-color kits include 2 vials of each component. Some kits also contain the dyes listed in [Table 1](#).



Agilent Technologies

T7 Primer	24 µL
5x First Strand Buffer	100 µL
0.1M DTT	70 µL
10 mM dNTP Mix	20 µL
AffinityScript RNase Block Mix	36 µL
5x Transcription Buffer	160 µL
NTP Mix	35 µL
T7 RNA Polymerase Blend	10 µL
Nuclease-Free Water	250 µL

Table 1 Additional components for selected kits

Component	5190-2305	5190-2306	5190-2307	5190-2942	5190-2943	5190-2944	5190-3386
Cyanine 3-CTP	8 µL	8 µL	N/A	N/A	8 µL	8 µL	N/A
Cyanine 5-CTP	N/A	8 µL	8 µL	N/A	N/A	8 µL	8 µL
WT Primer Mix	N/A	N/A	N/A	30 µL	30 µL	2 × 30 µL	30 µL

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.

WARNING

Avoid contact with skin or eyes. In case of contact with skin or eyes, wash immediately with water.

Limited Use License for End Users - For Research Use Only

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.

© Agilent Technologies, Inc. 2009-2011
 Printed in USA
 Revision C, August 2011



5190-2331-p