

## Turbo Labeling™ Kit - Biotin

### Quick Start Guide

KIT0608

This Quick Start Guide is intended to provide a quick reference once the complete User Guide has been reviewed and understood. This is not a stand-alone document. For optimal kit performance, please read the entire User Guide to review licensing, labeling restrictions and the labeling procedure before using the kit.

The user guide is downloadable at [www.appliedbiosystems.com](http://www.appliedbiosystems.com) or a printed copy is available free of charge by contacting technical support at 1-800-831-6844 option 5.

#### STEP Biotin labeling procedure

- Starting with 1-20 µg of aRNA in 13 µl of nuclease-free water, add 5 µl Turbo Biotin reagent and 2 µl Turbo 10x labeling buffer:

Component	Volume (µl)
aRNA (1-20 µg) + water	13
Turbo Biotin Reagent	5
10x Labeling Buffer	2
<b>Total</b>	<b>20</b>

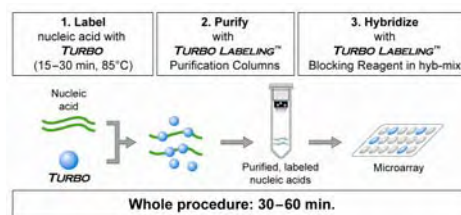
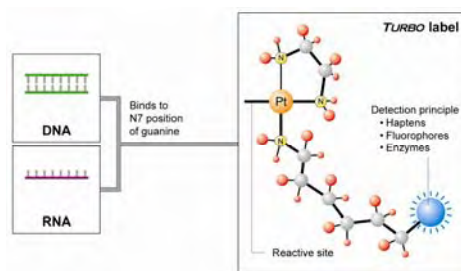
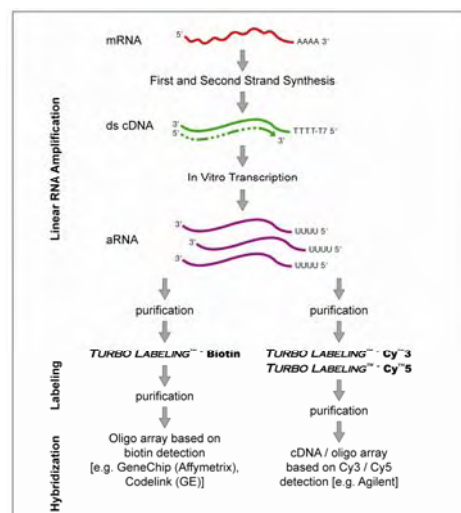
- Mix by flicking or pipetting gently and incubate on a thermocycler:
  - 85 °C for 30 minutes
  - 4 °C for 1-30 minutes
- Briefly centrifuge to collect tube contents

#### STEP Purification procedure

- Briefly vortex purification column, loosen cap 1/4 turn and snap off bottom closure
- Place tube in provided 2 ml collection tube
- Centrifuge column at 16,000 x g for 1 minute
- Discard flow-through and retain collection tube
- Wash column by pipetting 300 µl of nuclease-free water onto the column and centrifuge at 16,000 x g for 1 minute
- Discard flow-through and collection tube
- Transfer column to a new 1.5 ml micro-centrifuge collection tube
- Pipette biotin-labeled aRNA onto column bed
- Centrifuge at 16,000 x g for 1 minute
- Discard column and retain flow-through which is purified labeled aRNA (20 µl)

### Array Hybridization

Hybridization recommendations are detailed in the user guide.



## Turbo Labeling™ Kit - Cy<sup>®</sup>3/Cy<sup>®</sup>5

### Quick Start Guide

Cy3 KIT0609

Cy5 KIT0610

This Quick Start Guide is intended to provide a quick reference once the complete User Guide has been reviewed and understood. This is not a stand-alone document. For optimal kit performance, please read the entire User Guide to review licensing, labeling restrictions and the labeling procedure before using the kit.

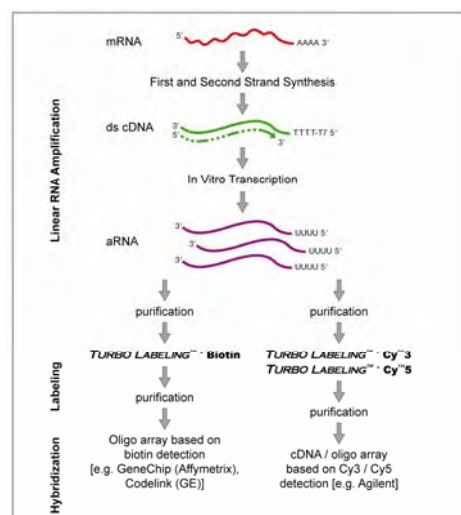
The user guide is downloadable at [www.appliedbiosystems.com](http://www.appliedbiosystems.com) or a printed copy is available free of charge by contacting technical support at 1-800-831-6844 option 5.

#### STEP Cy3/Cy5 labeling procedure

- Starting with 1-15 µg of aRNA in 40 µl of nuclease-free water, add 5 µl Turbo Cy3 or Cy5 reagent and 5 µl Turbo 10x labeling buffer:

Component	Volume (µl)
aRNA (1-15 µg) + water	40
Turbo Cy3 or Cy5 Reagent	5
10x Labeling Buffer	5
<b>Total</b>	<b>50</b>

- Mix by flicking or pipetting gently and incubate in thermocycler:
  - 85° C for 15 minutes
  - 4° C for 1-30 minutes
- Briefly centrifuge to collect tube contents



#### STEP Purification procedure

- Briefly vortex purification column, loosen cap 1/4 turn and snap off bottom closure
- Place tube in provided 2 ml collection tube
- Centrifuge column at 16,000 x g for 1 minute
- Discard flow-through and retain collection tube
- Wash column by pipetting 300 µl of nuclease-free water onto the column and centrifuge at 16,000 x g for 1 minute
- Discard flow-through and collection tube
- Transfer column to a new 1.5 ml micro-centrifuge collection tube
- Pipette Cy3- or Cy5-labeled aRNA onto column bed
- Centrifuge at 16,000 x g for 1 minute
- Discard column and retain flow-through which is purified labeled aRNA (50 µl)

#### Array Hybridization

Hybridization recommendations are detailed in the user guide.

