

# User Bulletin

## TaqMan<sup>®</sup> Reverse Transcription Reagents Kit

---

June 2008

**SUBJECT:** How to Find the *TaqMan<sup>®</sup> Reverse Transcription Reagents Kit Protocol*

<b>In This User Bulletin</b>	Use this user bulletin to:	
	Find a PDF copy of the protocol by using the document part number or the product description . . . . .	2
	Find a PDF copy of the protocol by using the product part number	2
	Order a print copy of the protocol . . . . .	3
<b>Overview</b>	You can use the Applied Biosystems web site to find the protocol for the TaqMan <sup>®</sup> Reverse Transcription Reagents Kit.	

## Find a PDF copy of the protocol by using the *document* part number or the product description

1. Go to [www.appliedbiosystems.com](http://www.appliedbiosystems.com).
2. Click **Support** in the top menu bar. The Support web page is displayed.
3. In the Product & Technical Literature box, enter the part number **402876** or the description **Gold RT-PCR** in the Keyword Search field.
4. Ensure that the drop-down menu says “Search all documents,” then click **Search**. The Product & Technical Literature page is displayed.
5. Scroll down to the Protocols section, then click **TaqMan® Gold RT-PCR: Protocol: Rev D**. The PDF copy of the protocol opens.

Refer to chapters 2 and 3 in the protocol for instructions on how to use the TaqMan Reverse Transcription Reagents Kit.

## Find a PDF copy of the protocol by using the *product* part number

1. Go to [www.appliedbiosystems.com](http://www.appliedbiosystems.com).
2. Enter the appropriate part number (for example, **N8080234**) in the Search box that is just below the top blue menu bar. You can enter other part numbers of products that require the TaqMan® Reverse Transcription Reagents Kit, including part numbers 402869, 4304133, 4304134, N8080232, or N8080233.
3. Ensure that “All Categories” is selected in the drop-down list to the right of the Search box, then click the orange arrow. The search results are displayed.
4. Click **Ordering Information** to the right of the appropriate part number (for example, N8080234). The TaqMan® Reverse Transcription Reagents web page is displayed.
5. Select the **Literature/Resources** tab.

6. Click **Protocol: TaqMan<sup>®</sup> Gold RT-PCR Kit: Protocol**. The PDF copy of the protocol opens.

Refer to chapters 2 and 3 in the protocol for instructions on how to use the TaqMan Reverse Transcription Reagents Kit.

## Order a print copy of the protocol

1. Go to [www.appliedbiosystems.com](http://www.appliedbiosystems.com).
2. Click **Log In** just below the top blue menu bar.
3. Enter your username and password to enter the Applied Biosystems Store, then click **Login**. The Products web page is displayed.
4. Enter the part number **402876** in the Search box just below the top blue menu bar.
5. Ensure that “All Categories” is selected in the drop-down list to the right of the Search box, then click the orange arrow. The search results are displayed.
6. Click **TaqMan<sup>®</sup> Gold RT-PCR Reagents Protocol**. The TaqMan<sup>®</sup> Gold RT-PCR Reagents Protocol web page is displayed.
7. Click **Add to Basket**.
8. In the gray bar across the top of the page, click **Shopping Basket**. The Shopping Basket: Applied Biosystems Store web page is displayed.
9. Confirm your order. then click **Continue**. The Shipping and Billing page is displayed.
10. Enter the required shipping and billing information, then click **Continue**. The Summary page is displayed.
11. Click **Place Order**. The page confirming your order is displayed.

Refer to chapters 2 and 3 in the protocol for the instructions on how to use the TaqMan<sup>®</sup> Reverse Transcription Reagents Kit.

© Copyright 2008, Applied Biosystems. All rights reserved.

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

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights (such as 5' Nuclease Process patent rights) are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

### **TRADEMARKS:**

Applera, Applied Biosystems, and AB (Design) are registered trademarks of Applera Corporation or its subsidiaries in the U.S. and/or certain other countries.

TaqMan is a registered trademarks of Roche Molecular Systems, Inc.

All other trademarks are the sole property of their respective owners.

### **Worldwide Sales and Support**

Applied Biosystems vast distribution and service network, composed of highly trained support and applications personnel, reaches 150 countries on six continents. For sales office locations and technical support, please call our local office or refer to our Web site at [www.appliedbiosystems.com](http://www.appliedbiosystems.com).

Applera is committed to providing the world's leading technology and information for life scientists. Applera Corporation consists of the Applied Biosystems and Celera Genomics businesses.

### **Headquarters**

850 Lincoln Centre Drive  
Foster City, CA 94404 USA  
Phone: +1 650.638.5800  
Toll Free (In North America): +1 800.345.5224  
Fax: +1 650.638.5884

06/2008

