

# TaqMan® AIV-Matrix Reagents

Do not store in a frost-free freezer.

Catalog # (P/N): 4405543

Materials Provided: Reagents for 100, 25 µL RT-PCRs are supplied. Store the reagents in a non-frost-free freezer.

Component	Amount	Storage
2X RT-PCR Buffer	1.375 mL	-20°C
25X RT-PCR Enzyme Mix	110 µL	-20°C
25X AIV-Matrix Primer Probe Mix	110 µL	-20°C
Detection Enhancer	190 µL	-20°C
Nuclease-free Water	1.75 mL	-20°C, 4°C, or room temp

## TaqMan® Probe Reporter/Quencher:

Target	Reporter	Quencher
AIV-Matrix	FAM™ dye	BHQ <sup>®</sup> -1 dye
Xeno™ RNA Control	CF560 dye*	BHQ-1 dye

\*Cal Fluor<sup>®</sup> Orange 560 (CF560) dye has an absorbance maximum of 538 nm and an emission maximum of 559 nm. When using an Applied Biosystems Real-Time PCR System, it is best to calibrate with Cal Fluor Orange 560 dye. If that is not possible, use the VIC<sup>®</sup> dye detector for Cal Fluor Orange 560 dye.

## Safety Information:

Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

**WARNING! CHEMICAL HAZARD. 2X RT-PCR Buffer** causes eye, skin, and respiratory tract irritation. May be harmful if swallowed, inhaled, or absorbed through skin. Avoid breathing vapor. Use with adequate ventilation. Avoid contact with eyes and skin.

**WARNING! CHEMICAL HAZARD. 25X RT-PCR Enzyme Mix** may cause eye, skin, and respiratory tract irritation. May be harmful if swallowed. Avoid breathing vapor. Use with adequate ventilation. Avoid contact with eyes and skin.

**WARNING! CHEMICAL HAZARD. Detection Enhancer** may cause cancer! Causes eye, skin, respiratory tract, and gastrointestinal irritation. Avoid breathing vapor. Use with adequate ventilation. Avoid contact with eyes and skin.

**WARNING!** This product contains a chemical known to the State of California to cause cancer: Acetamide.

## USER INFORMATION

### General Information:

TaqMan<sup>®</sup> AIV-Matrix Reagents enable one-step, real-time, RT-PCR targeting the Avian Influenza Virus matrix gene RNA and Xeno RNA Control. The RT-PCR tolerates a wide range of input RNA amounts, but we typically do not exceed 100 ng of input RNA.

### RNA isolation

- Add Xeno RNA Control to the lysis solution used for RNA isolation, to serve as a positive control both for recovery of RNA and for the RT-PCR. Add 5,000–10,000 copies of Xeno RNA Control per isolation (0.5–1.0 µL at 10,000 copies/µL).
- When isolating RNA from cell-free sample sources such as serum, add carrier RNA to the lysis solution to maximize RNA recovery. Include carrier RNA even if you have added Xeno RNA Control to the lysis solution.
- It is important to use RNA that is free of RT-PCR inhibitors in the procedure.
- To isolate RNA for use in this procedure, we recommend using the Ambion<sup>®</sup> MagMAX™ RNA Isolation Kit appropriate for your sample type. Search for "MagMAX" at [www.appliedbiosystems.com](http://www.appliedbiosystems.com).

### Applications:

#### Reaction setup

1. Program the real-time PCR instrument according to the manufacturer's recommendation.

ROX™ passive reference dye is included in the RT-PCR Buffer.

2. Assemble RT-PCR master mix(es), for 25 µL reactions, on ice. Include 5–10% extra to ensure that you will have enough master mix for all the samples.

Component	Amount
2X RT-PCR Buffer	12.5 µL
25X AIV-Matrix Primer Probe Mix	1 µL
25X RT-PCR Enzyme Mix	1 µL
Detection Enhancer	1.67 µL
Total volume master mix per reaction	16.17 µL

3. Distribute RT-PCR master mix to the wells of a PCR plate or to PCR tubes.
4. Add sample or Nuclease-free Water for a final reaction volume of 25 µL.
5. Run the thermal cycler program.

---

## QUALITY CONTROL

---

<b>Functional Testing:</b>	Each lot of supplied reagent is tested for detection of RNA target using the instructions in this protocol. Relevant kit components are tested in the following nuclease assays.
<b>RNase Activity:</b>	A sample is incubated with labeled RNA, followed by PAGE analysis.
<b>Nonspecific Endonuclease Activity:</b>	A sample is incubated with supercoiled plasmid DNA and analyzed by agarose gel electrophoresis.
<b>Exonuclease Activity:</b>	A sample is incubated with labeled double-stranded DNA, followed by PAGE analysis.

---

## OTHER INFORMATION

---

**Material Safety Data Sheets:** Material Safety Data Sheets (MSDSs) for any chemical product supplied by Applied Biosystems or Ambion are available 24 hours a day. At [www.appliedbiosystems.com](http://www.appliedbiosystems.com), select Support, then MSDS. Search by chemical name, product name, product part number, or MSDS part number. Right-click to print or download the MSDS of interest. At [www.ambion.com](http://www.ambion.com), go to the web catalog page for the product of interest. Select MSDS, then right-click to print or download. Or, e-mail ([MSDS\\_Inquiry\\_CCRM@appliedbiosystems.com](mailto:MSDS_Inquiry_CCRM@appliedbiosystems.com)), telephone (650-554-2756; USA), or fax (650-554-2252; USA) your request, specifying the catalog or part number(s) and the name of the product(s). We will e-mail the associated MSDSs unless you request fax or postal delivery. Requests for postal delivery require 1-2 weeks for processing.

**Warranty and Liability:** *For research use only. Not for use in diagnostic procedures.*

Applied Biosystems is committed to delivering superior product quality and performance, supported by industry-leading global service and technical support teams. Warranty information for the accompanying consumable product is available at [www.ambion.com/info/warranty](http://www.ambion.com/info/warranty) in "Limited Warranty for Consumables," which is subject to the exclusions, conditions, exceptions, and limitations set forth under the caption "EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS" in the full warranty statement. Please contact Applied Biosystems if you have any questions about our warranties or would like information about post-warranty support.

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

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, Patents, and Licensing:**

Trademarks of Life Technologies Corporation and its affiliated companies: AB Logo™, Ambion®, Applied Biosystems®, FAM™, MagMAX™, ROX™, VIC®, and Xeno™. TaqMan is a registered trademark of Roche Molecular Systems, Inc. Black Hole Quencher®, 'BHQ' and 'CAL Fluor' are registered trademarks of Biosearch Technologies, Inc., Novato, CA. All other trademarks are the sole property of their respective owners.

**Limited Use Label License: Veterinary Diagnostics**

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform veterinary diagnostic services including reporting the results of purchaser's activities for a fee and for the purchaser's own internal research. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. For information on obtaining additional rights, please contact [outlicensing@lifetech.com](mailto:outlicensing@lifetech.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008. No right under any other patent claim (such as claims to methods, apparatus or reagents) is conveyed expressly, by implication, or by estoppel. Human diagnostic uses require a separate license from Roche.

© 2009, 2011 Life Technologies Corporation. All rights reserved. 4412867C