

VetMAX™ Rotavirus A/B/C Controls

Catalog Number A57451

Pub. No. MAN0029472 Rev. A.0



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

Applied Biosystems™ VetMAX™ Rotavirus A/B/C Controls (Cat. No. [A57451](#)) are designed for use in multiplex real-time reverse transcription quantitative PCR (RT-qPCR) workflows for identification of porcine rotavirus species A, B, and C (RVA, RVB, and RVC) RNA and Xeno™ RNA. VetMAX™ Rotavirus A/B/C Controls contain single- and double-stranded RVA, RVB, RVC, and Xeno™ RNA controls in a ready-to-use, premixed format.

Xeno™ RNA Control is included as an internal positive control. Add Xeno™ RNA Control to each sample during RNA isolation to monitor RNA recovery and identify PCR inhibitors in molecular detection workflows.

Contents and storage

Item	Concentration	Amount ^[1]	Storage
VetMAX™ Rotavirus A/B/C Controls	<ul style="list-style-type: none"> RVA: 1,000 copies/μL RVB: 1,000 copies/μL RVC: 1,000 copies/μL Xeno™ RNA: 1,000 copies/μL 	80 μL	–30°C to –10°C
Xeno™ RNA	10,000 copies/μL	250 μL	

^[1] Sufficient for 10 positive-control reactions. We recommend 2 positive control reactions per run.

Guidelines for use

- Follow good laboratory practices to prevent false positives and contamination of test samples with PCR products (see “Good laboratory practices for PCR and RT-PCR”).
- Include two positive-control reactions in each RT-qPCR run.

Table 1 Control and experimental RT-qPCR (20-μL reactions)

Component	Sample	Volume per reaction
Test sample	RNA	8 μL
No-template control (NTC)	Nuclease-free water	8 μL
Positive control	VetMAX™ Rotavirus A/B/C Controls	8 μL

Good laboratory practices for PCR and RT-PCR

- Wear clean gloves and a clean lab coat.
 - Do not wear the same gloves and lab coat that you have previously used when handling amplified products or preparing samples.
- Change gloves if you suspect that they are contaminated.
- Maintain separate areas and dedicated equipment and supplies for:
 - Sample preparation and reaction setup.
 - Amplification and analysis of products.
- Do not bring amplified products into the reaction setup area.
- Open and close all sample tubes carefully. Avoid splashing or spraying samples.
- Keep reactions and components capped as much as possible.
- Use a positive-displacement pipettor or aerosol-resistant barrier pipette tips.
- Clean lab benches and equipment periodically with 10% bleach solution or DNA decontamination solution.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms

and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 2130 Woodward Street | Austin, Texas 78744 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN0029472 A.0

Revision	Date	Description
A.0	16 May 2023	New document for VetMAX™ Rotavirus A/B/C Controls.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.