



Research use only. Not for use in diagnostic procedures.

NeoMDx™ cCMV real-time PCR assay (CMV-96 and CMV-384)

NeoMDx™ Elution Solution

Product number: CMV-ELU-384 Lot Number: 3330144

Expiration date: 31-May-2026

Store ambient.

For research use only. Not for use in diagnostic procedures.

The above listed component lot of NeoMDx™ cCMV real-time PCR assay (CMV-96 and CMV-384) has been tested and approved by Quality.

Approved by: 

DocuSigned by:

Sarah Parker

B6C893CB69B9479...

  
Date: 20 September 2024 | 11:50 AM EDT

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

CMV-ELU-R Rev 01

