



Research use only. Not for use in diagnostic procedures.

# HostDetect™ CHO DNA Quantification Kit

Product number: DXMDX-RGT-1003

Lot Number: 3342495

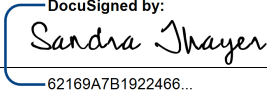
Kit Expiration date: 15-OCT-2025

PARAMETER	SPECIFICATION	RESULT
CHO Sensitivity Panel	Sensitivity panel passes QC.	Pass
Internal Control Sensitivity Panel	Sensitivity panel passes QC.	Pass
Kit IC	Kit IC passes QC.	Pass
Negative Control	Kit Negative control passes QC.	Pass

Store at -20°C or below.

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

The above listed lot of the CHO DNA Quantification kit has been tested and approved by Quality.

Approved by:  62169A7B1922466...

Date: 12 November 2024 | 9:08 AM EST

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.

DXMDX-RGT-1003 Rev 01

