



Research use only. Not for use in diagnostic procedures.

## Candida auris Real-time PCR Reagents

**Product number:** DXMDX-RGT-1001 **Lot Number:** 3186331

**Kit Expiration date:** 04-FEB-2024

**Kit Contents:**

Component	Lot No.	Expiration Date
Reagent A	CC20220203	04-FEB-2024
CAU Reagent B	3186315	15-JUN-2024
Reagent C	3186317	10-JUL-2024
Negative Control	3186257	19-JUL-2024
Internal Control	3186322	16-JUN-2024
CAU Positive Control	3186323	15-JUN-2024

Store at -20°C or below.

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

The above listed component lot of Candida auris Real-time PCR Kit has been tested and approved by Quality.

Approved by: DocuSigned by:  
Sandra Thayer  
93C1E01C92BB4AE...

Date: 20 September 2023 | 11:00 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.

DXMDX-RGT-1001-R Rev 01

