



Product Manual

AMPIPROBE[®] Candida 1 Assay Kit

Catalog #: ENZ-GEN202-0100

100 Assays



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

TRADEMARKS

ENZO, AMPIPROBE, and AMPIGENE are trademarks of Enzo Biochem, Inc. Qiagen and Rotor-Gene are registered trademarks of Qiagen GmbH.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

TABLE OF CONTENTS



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

Description.....	2
Introduction	2
AMPIPROBE® Technology.....	3
Principle of AMPIPROBE® Candida 1 Assay.....	3
AMPIPROBE® Technology Schematic	4
Safety Warnings & Precautions.....	5
Materials Supplied.....	6
Storage	6
Other Materials Required But Not Supplied.....	6
Procedural Notes	7
Sample Compatibility	7
Reagent Preparation	7
PCR Setup.....	7
PCR Run Profile.....	9
Additional Instrument Settings.....	9
Data Analysis	10
Interpretation of Results	10
Performance Characteristics	11
Frequently Asked Questions	12
References.....	13
Contact Information.....	14



DESCRIPTION

The AMPIPROBE[®] Candida 1 Assay is a real-time polymerase chain reaction (qPCR) assay for the qualitative detection of *Candida albicans*, *Candida glabrata* and *Candida krusei* genomic DNA. The kit uses the AMPIPROBE[®] assay platform which takes advantage of paired fluorophore- and quencher-labeled primers specific for each of the target species as well as an internal control. The kit contains all reagents necessary for PCR-based detection of *C. albicans*, *C. glabrata* and *C. krusei* DNA. A positive PCR control consisting of a mixture of Candida target templates and a negative PCR control that results negative for *Candida spp*, but positive for the internal control (human β -globin). Please read the complete kit insert before performing this assay.

INTRODUCTION

Vulvovaginal candidiasis (VVC) is second, after bacterial vaginosis, among the many causes of vaginitis, and is diagnosed in up to 40% of women with vaginal complaints in the primary care setting¹. Approximately 70-75% of women will have at least one symptomatic *Candida* vulvovaginal infection during their lives^{1,2}. Accurate detection and differentiation of *Candida* species that cause VVC leads to a significant improvement in clinical management of women with vaginal complaints, especially in women with recurrent VVC (RVVC).

The current assay provides rapid and accurate results for the detection of *C. albicans*, *C. glabrata*, and *C. krusei* DNA in a user-supplied sample of interest.

AMPIPROBE® TECHNOLOGY

Enzo's AMPIPROBE® technology incorporates probe detection technology in primer design. It employs a combination of fluorescent reporter-labeled primers and quencher-labeled primers to amplify DNA, akin to traditional PCR. When free in solution, fluorescent primers generate a signal. However, as the primers are incorporated into amplified DNA, the quencher and the fluorophore are brought within close proximity and exhibit Förster resonance energy transfer (FRET). This causes a logarithmic decay of signal with respect to the number of PCR amplification cycles. The cycle number at which fluorescence signal drops below a defined threshold, is indicative of the amount of target nucleic acid present in the sample.

Enzo's AMPIPROBE® Assay kits provide the following benefits:

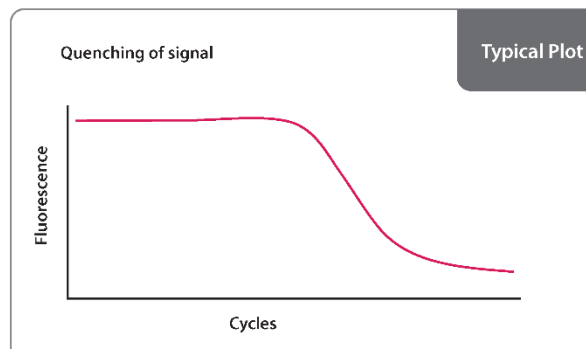
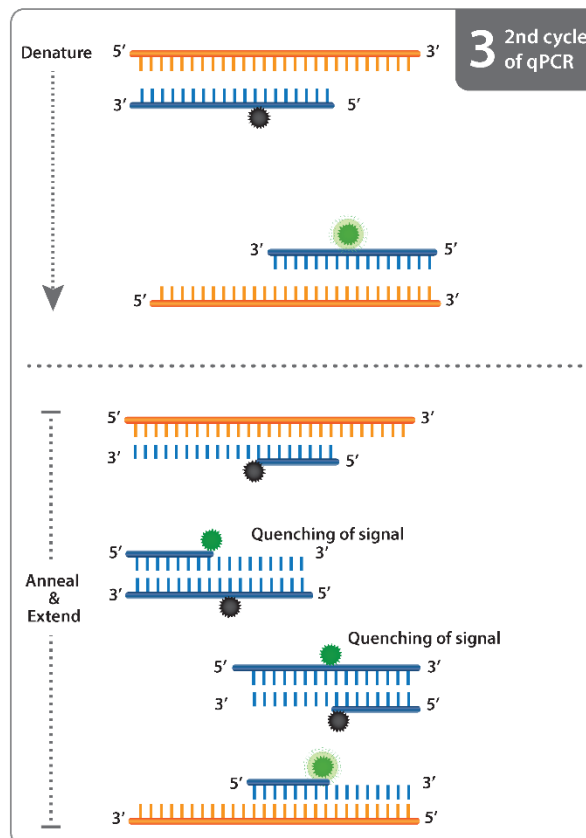
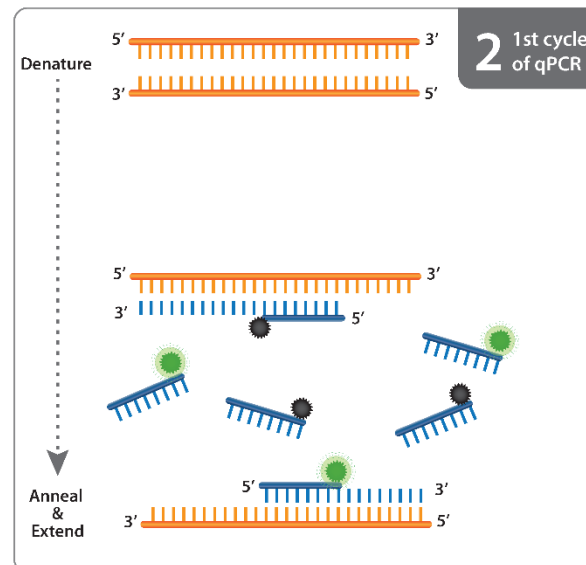
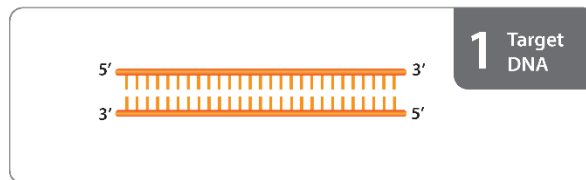
- High sensitivity
- Low sample input
- Adaptable to multiplexing
- Compatible with open qPCR instruments

PRINCIPLE OF AMPIPROBE® CANDIDA 1 ASSAY

The AMPIPROBE® Candida 1 Assay is a qualitative real-time PCR assay for the detection of *C. albicans*, *C. glabrata* and *C. krusei* genomic DNA. The nucleotide sequences of the PCR primers have been optimized to target sequences within species-specific regions of the *Candida* genome. The *Candida* primers have been designed using AMPIPROBE® technology which incorporates reporter and quencher dyes into the primers. Target detection is based on fluorescence decay when successive rounds of amplification bring fluorophore and quencher-labeled primer pairs in close proximity resulting in FRET. Loss of fluorescence below a defined threshold in a particular channel indicates sample positivity for the corresponding *Candida* species.

The AMPIPROBE® Candida 1 Primer Mix includes primers for a ubiquitously conserved human housekeeping gene (human β -globin). Successful amplification of the internal control serves as an indicator of sample adequacy, extraction efficiency and successful amplification in each individual sample.

AMPIPROBE® TECHNOLOGY SCHEMATIC



SAFETY WARNINGS & PRECAUTIONS



Avoid freeze /
thaw cycles



Handle with
care

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

- Kit components should be stored at or below -20°C . Care should be taken to limit the number of freeze-thaw cycles.
- The AMPIPROBE[®] Candida 1 Primer Mix contains fluorescently labeled primers. To avoid photobleaching, protect from prolonged exposure to light.
- Some components of this kit may contain hazardous substances. Reagents can be harmful if ingested or absorbed through the skin and may cause irritation to the eyes. Reagents should be treated as possible mutagens and should be handled with care and disposed of properly.
- Practice aseptic technique when handling reagents to avoid introduction of contaminants that might interfere with assay interpretation.
- The use of screw-cap tubes and barrier pipette tips is strongly encouraged to prevent samples and reagents from becoming aerosolized which might lead to contamination.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas. Any biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

MATERIALS SUPPLIED

Reagent	Quantity
AMPIGENE[®] HS Taq DNA Polymerase	100 µL
AMPIGENE[®] dNTP Mix	50 µL
AMPIPROBE[®] 5X Assay Buffer	500 µL
AMPIPROBE[®] Candida 1 Primer Mix	500 µL
Candida 1 Positive PCR Control	100 µL
Negative PCR Control	100 µL
Nuclease-free Water	1 mL

STORAGE

All components of this kit are stable at -20°C until the kit's expiration date.

OTHER MATERIALS REQUIRED BUT NOT SUPPLIED

1. Vortex mixer
2. Mini centrifuge
3. Calibrated pipettes capable of delivering volumes between 1 and 1000 µL
4. DNase/RNase free barrier pipette tips
5. DNase/RNase free 1.5 mL Screw Cap Micro tube (for preparation of master mix)
6. qPCR instrument and compatible accessories

PROCEDURAL NOTES

Do not mix components from different kit lots or use reagents beyond the kit expiration date.

SAMPLE COMPATIBILITY

Sample types and DNA extraction methods should be evaluated for compatibility with the AMPIPROBE[®] Candida 1 Assay Kit as part of the user's validation process.

REAGENT PREPARATION

1. Pre-cool Loading Block at 4°C for at least one hour prior to setup.
2. Thaw the following kit reagents at room temperature (15-25°C): Nuclease-free Water, AMPIGENE[®] dNTP Mix, AMPIPROBE 5X Assay Buffer, AMPIPROBE[®] Candida 1 Primer Mix, Candida 1 Positive PCR Control, and Negative PCR Control.

Once thawed, vortex the reagents to mix. Spin briefly to collect the contents at the bottom of tubes. Keep the reagents on ice until ready to use. Protect the AMPIPROBE[®] Candida 1 Primer Mix from prolonged exposure to light.

Note: Some precipitation may occur in the AMPIPROBE[®] 5X Assay Buffer. Vortex and warm (37°C) vial to dissolve precipitate prior to use.

3. Remove the AMPIGENE[®] HS Taq DNA Polymerase from storage and briefly spin to collect contents at the bottom of tube. **Do not vortex.** Keep tube on ice until ready to use.

PCR SETUP

The procedure described below uses the Qiagen Rotor-Gene Q instrument, but other real-time PCR machines capable of detecting fluorescence quenching may also be used. Refer to the manufacturer's instrument manual for details regarding operation of the PCR machine.

For each assay run, include the following controls: (a) Candida 1 Positive PCR Control, (b) Negative PCR Control, and (c) No Template Control (not included in kit). Depending upon the user's requirements, additional controls may be included.

1. Immediately prior to use, prepare sufficient **PCR Master Mix** for the DNA samples and controls in a labeled, DNase/RNase free screw-cap tube, according to the volumes given in Table 1. Add the reagents to the tube in the order listed. Include sufficient volume for 2 extra reactions to allow for any pipetting variance.

The **PCR Master Mix** contains all of the components needed for the reaction except the DNA template (sample).

2. Thoroughly mix the **PCR Master Mix** by either inverting the tube or by pipetting up and down until swirls are no longer visible (10-20 times). **Do not vortex**. Centrifuge briefly to bring contents to the bottom of the tube. Avoid prolonged exposure to light.
3. Dispense 15 μL of **PCR Master Mix** into each of the 0.1 mL PCR tubes, taking care to deliver the solution to the bottom of the tube.
4. Add 10 μL sample or control to the PCR tube and mix by pipetting up and down at least 3 times.
5. Cap the tubes and visually inspect to confirm a tight seal and correct volume.
6. Place the tubes in the thermal cycler and run the method. See **PCR RUN PROFILE** section for the method details.

Table 1. Preparation of PCR Master Mix

Reagent	Volume per reaction	Total volume required ^{a,b}
Nuclease-free Water	3.5 μL	3.5 μL x (N+2)
AMPIPROBE™ 5X Assay Buffer	5 μL	5 μL x (N+2)
AMPIPROBE™ dNTP Mix	0.5 μL	0.5 μL x (N+2)
AMPIPROBE™ Candida 1 Primer Mix	5.0 μL	5.0 μL x (N+2)
AMPIGENE™ HS Taq DNA Polymerase	1 μL	1 μL x (N+2)
TOTAL	15 μL	15 μL x (N+2)

^a Total volume required for **N + 2** reactions, where **N** is the number of samples plus controls. When preparing the Master Mix, prepare enough for 2 extra reactions to allow for any pipetting variance.

^b If the number of samples plus controls exceeds 24, more than 2 extra reaction volumes may be needed.

PCR RUN PROFILE

Create a temperature profile on your PCR instrument as indicated in Table 2. The method described has been verified using a Qiagen Rotor-Gene Q instrument. Use of other instruments may require minor modifications.

Table 2. PCR Cycling Parameters

Step	Parameter	Temp [°C]	Cycles	Hold [mm:ss]
Hold	Enzyme Activation	95°	1	3:00
Cycling	Amplification	95°	40	00:10
		68°		00:30 (fluorescent signal detection)

Fluorescence is detected in the **FAM**/Green, **HEX**/Yellow, **ROX**/Orange and **Cy5**/Red channels at the end of the 2nd segment of the cycling step (68°C).

ADDITIONAL INSTRUMENT SETTINGS

The settings described in Table 3 are recommended for use with the Qiagen Rotor-Gene Q instrument. Perform gain optimization on a tube corresponding to No Template Control and set the instrument to perform optimization before first acquisition.

Table 3. Gain Optimization Settings

Channel	Min Reading	Max Reading	Min Gain	Max Gain
FAM /Green	70FI	80FI	-10	10
HEX /Yellow	70FI	80FI	-10	10
ROX /Orange	70FI	80FI	-10	10
Cy5 /Red	70FI	80FI	-10	10

DATA ANALYSIS

Fluorescent signal intensity is detected in four channels:

- The signal for *C. albicans* amplification is detected in the **FAM**/Green channel.
- The signal for *C. glabrata* amplification is detected in the **HEX**/Yellow channel.
- The signal for *C. krusei* amplification is detected in the **ROX**/Orange channel.
- The signal for Internal Control amplification is detected in the **Cy5**/Red channel.

Results are interpreted using the instrument software by determining the cycle number, **Ct**, corresponding to the point where fluorescence signal drops below a defined threshold.

Using the Rotor-Gene Q software, the data for each of the four channels must be analyzed separately. For all four channels convert the data to display on a linear scale, select “Dynamic Tube” and ensure that “Slope Correct” is OFF. The settings recommended in Table 4 can be used to generate **Ct** values.

Table 4. Recommended Data Analysis Parameters

Channel	Ignore First	Threshold
FAM /Green	5	-0.09
HEX /Yellow	10	-0.03
ROX /Orange	10	-0.09
Cy5 /Red	10	-0.03

INTERPRETATION OF RESULTS

- The generation of a **Ct** value using the above analysis parameters is indicative of a positive result.
- The absence of a **Ct** value when the above analysis parameters are applied is indicative of a negative result.
- The Candida 1 Positive PCR control must produce a positive result in all channels (see Table 5). If this condition is not met, the entire run is invalid and must be repeated.
- The Negative PCR control must produce negative results in the green, yellow and orange channels AND a positive result in the red channel (see Table 5). If these conditions are not met, the entire run is invalid and must be repeated.

- If included in the assay run, any No Template Control must test negative in the green, yellow and orange channels. If amplification is observed in the red channel the **Ct** must be >35 (see Table 5).
- For any individual samples of human origin there must be a positive result in the red channel with **Ct** < 35. If this condition is not met, results for the individual sample are considered invalid and the sample should be rerun.

Table 5. Expected Results for Assay Controls

Channel	Candida 1 Positive PCR Control	Negative PCR Control	No Template Control
FAM/Green	POS	NEG	NEG
HEX/Yellow	POS	NEG	NEG
ROX/Orange	POS	NEG	NEG
Cy5/Red	POS	POS	NEG or Ct >35

PERFORMANCE CHARACTERISTICS

Overall assay detection sensitivity will vary based on the methods used for nucleic acid extraction and should be verified independently.

Analytical sensitivity of the PCR reaction, based on results obtained using plasmid controls, is estimated to be in the range of 1 to 50 target copies per reaction.

FREQUENTLY ASKED QUESTIONS

Can this assay be run on an ABI 7500?

Most thermal cyclers can be used, provided the software is equipped to detect fluorescence decay. However, the PCR method profile and subsequent data analysis parameters may require instrument-specific modification. The software accompanying the ABI 7500 instrument is not capable of generating Ct values for a fluorescence quenching assay and although the assay can be run, data must be analyzed independently.

What is the composition of the Candida 1 Positive PCR Control and the Negative PCR Control?

The PCR controls supplied with the AMPIPROBE™ Candida 1 Assay Kit are composed of linearized plasmids containing the specific sequences targeted by the primers in this assay. The amount of target in each mixture is formulated to produce a positive result falling within the dynamic range of the assay when it is carried out according to the procedure described in this user manual.

Can the AMPIPROBE Candida 1 Assay be run together with the AMPIPROBE Candida 2 Assay?

Due to differences in primer fluorescence, Candida 1 and Candida 2 assays should not be run simultaneously. The gain should be set on each assay individually.



Product Manual

REFERENCES

1. Sobel, J. D. 2007. Vulvovaginal candidosis. *Lancet* **369**:1961-1971.
2. Anderson, M. R., K. Klink, and A. Cohrssen. 2004. Evaluation of vaginal complaints. *JAMA* **291**:1368-1379.



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com