



AMPIPROBE[®] Bacterial Vaginosis Extraction Controls

Catalog Number: ENZ-GEN206

For the latest product information, including support documentation, visit us online:

www.enzolifesciences.com

Technical Support (US): 800-942-0430
Technical Support (EU): +41 61 926 8989



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 use only. Not for use in diagnostic procedures. 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, Rotor-Gene and QIASymphony SP are registered trademarks of Qiagen GmbH. APTIMA is registered trademark of Hologic, Inc.

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

INTENDED USE

The AMPIPROBE® Bacterial Vaginosis Extraction Controls are designed to be used during the genomic DNA extraction of *Atopobium vaginae*, *Gardnerella vaginalis*, *Lactobacillus spp.*, *Megasphaera spp.*, and BVAB2.



Avoid freeze /
thaw cycles



Handle with
care

SAFETY WARNINGS & PRECAUTIONS

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.
- 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
Negative Extraction Control	300 µL
BV Positive Extraction Control	300 µL
No BV Target Control	300 µL

NOTE

All Performance specifications must be validated by end-user and will vary based on chosen extraction methods. Eluent from extraction can be used as extraction control for AMPIPROBE[®] Bacterial Vaginosis Assay kit (ENZ-GEN 205-0100).

Plasmids in Extraction Controls are present at higher concentrations compared to assay controls included in the qPCR kit. Extra care should be taken to avoid contamination.

STORAGE

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

PROCEDURAL NOTES

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

REAGENT PREPARATION

1. Thaw the kit reagents at room temperature (15-25°C).
2. Vortex the reagents to mix.
3. Spin briefly to collect the contents at the bottom of tubes.

SETUP OF DNA EXTRACTION CONTROL

The procedure described below is optimized for the QIASymphony SP instrument (Qiagen) starting from an APTIMA[®] Multitest Swab Specimen collection tube.

Once thawed at room temperature, the AMPIPROBE[®] Extraction Control tubes may be prepared. Label three 2.0 mL sample tubes with the appropriate control type (PC, NC, No BV Target), and then add 1.0 mL of blank APTIMA[®] media to each tube.

Next, add 25 µL of each AMPIPROBE[®] Extraction Control into their associated, labelled 2.0 mL tube. The extraction controls are now prepared and ready for use. Continue extraction procedure same as samples.

PCR SETUP

Please refer to the PCR SETUP protocol present in Product Manual **AMPIPROBE® Bacterial Vaginosis Assay Kit** (Catalog #: ENZ-GEN205-0100, 100 Assays) and treat BV Positive Extraction Control, No BV Target Control and Negative Extraction controls as Samples.

Here below a brief summary of the PCR Setup protocol:

1. 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.
2. Dispense 15 μ L 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.
3. Add 10 μ L of control to the PCR tube and mix by pipetting up and down at least 3 times.
4. Cap the tubes and visually inspect to confirm a tight seal and correct volume.
5. Place the tubes in the thermal cycler and run the method. See PCR RUN PROFILE section for the method details in the AMPIPROBE® Bacterial Vaginosis Assay kit.

Note: The reagents required for the PCR Master Mix provided in the AMPIPROBE® Bacterial Vaginosis Assay kit will be sufficient to run also the Bacterial Vaginosis Extraction Controls.



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