# AMPIPROBE<sup>®</sup> Bacterial Vaginosis Extraction Controls

## The AMPIPROBE<sup>®</sup> Bacterial Vaginosis Extraction Controls are designed to ensure

The AMPLIPADE Candida Extraction Controls are designed to be used during the genomic DNA extraction of Atopobium vaginae, Gardnerella and detection of Atopobium vaginae, Gardnerella vaginalis, Lactobacillus spp., Megasphaera spp., and BVAB2.

And the formance specifications, must be validated by Endouser and will vary thanks of the property of the property of the extraction can be used as extraction control for AMPIPROBE® Bacterial Vaginosis Assay kit (Prod. No. ENZ-GEN205).

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.

Bacterial vaginosis (BV) is the most common vaginal infection in the United States and affects approximately 29% of women. BV complications include an increased risk to develop other sexually transmitted infections and an increased risk of infection after pelvic surgery. In addition, BV during pregnancy has been linked with adverse pregnancy outcomes including preterm labor and delivery, low birth weight, premature rupture of membranes, postpartum metritis and intra-amniotic infection.

Clinical signs and symptoms (Amsel criteria) are often used to diagnose BV but the gold standard method for diagnosis is enumeration of bacterial morphotypes on a Gram-stained smear.

Although the exact etiology of BV is still undetermined, the consensus is that it occurs when *Lactobacillus spp..*, which normally comprising 90-95% of total bacteria in healthy vaginal flora, is replaced by anaerobes, mainly but not restricted to *Gardnerella vaginalis*. More recently, molecular detection of other microbes, that are mostly uncultureable, have been associated with BV including *Atopobium vaginae*, BVAB2, and *Megasphaera spp.*.

#### **Ordering Information**

Order Online »

**ENZ-GEN206-0500** 500μl

Manuals, SDS & CofA

**View Online** »

- Compatible with manual and automated DNA extraction methods
- Compatible with most open qPCR platforms

#### **Handling & Storage**

**Use/Stability** All components are stable at -20°C until the kit's expiration date.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Dry Ice

### Regulatory Status RUO - Research Use Only

#### **Product Details**

**Application** qPCR

**Application Notes**The AMPIPROBE<sup>®</sup> Bacterial Vaginosis Extraction

Controls is compatible with manual and automated DNA extraction methods as well as any properly calibrated qPCR thermal cycler capable to detect fluorescence decay. It has been validated for use on the QIAGEN

QIAsymphony SP and Rotor-Gene Q.

Contents Negative Extraction Control

**BV Positive Extraction Control** 

No BV Target Control

**Species Reactivity** Atopobium vaginae, BVAB-2, Gardnerella vaginalis,

Lactobacillus spp., Megasphaera-1

Last modified: May 29, 2024

