

AMPIPROBE®

Bacterial Vaginosis Extraction Controls

The AMPIPROBE® Bacterial Vaginosis Extraction Controls are designed to ensure

an efficient genomic DNA extraction process during the genomic DNA extraction of *Atopobium vaginae*, *Gardnerella vaginalis*, *Lactobacillus spp.*, *Megasphaera spp.*, and BVAB2.

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 (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.*

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

Ordering Information

[Order Online »](#)

ENZ-GEN206-0500	500µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

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[®] 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 QIA Symphony 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)