AMPIPROBE[®] Ureaplasma spp. / M. genitalium / M. hominis Assay Kit

Special Speci

Ureaplasma and Mycoplasma, which belong to the same Mycoplasmataceae family and Mollicutes class, are the smallest self-replicating organisms and are characterized by their lack of a cell wall. These characteristics and limited biosynthetic capabilities contribute to the parasitic nature of Ureaplasma and Mycoplasma.

The class of pathogens is present in healthy individuals but has been associated with many adverse conditions affecting the reproductive tract. *Mycoplasma genitalium* is associated with urethritis, cervical inflammation, and pelvic inflammatory disease. *M. hominis* is often present concurrently with *Ureaplasma* species and is associated with a variety of conditions ranging from pelvic inflammatory diseases, chorioamnionitis, postpartum endometritis bacterial vaginosis, arthritis, osteoarthritis, wound infections, and several conditions in neonates.

Ureaplasma and *Mycoplasma*, especially in combination with other conditions such as bacterial vaginosis or cervical incompetence, have been associated with adverse pregnancy outcomes, such as chorioamnionitis, spontaneous preterm labor and preterm premature rupture of membranes. The AMPIPROBE® UMM Assay provides rapid and accurate results for the qualitative detection of *Ureaplasma spp.* (*Ureaplasma parvum* and *Ureaplasma urealyticum*), *Mycoplasma genitalium*, and *Mycoplasma hominis* DNA in a user-supplied sample of interest.

- Highly sensitive and specific assay
- Low-cost alternative to other methods of Ureaplasma spp. / M. genitalium / M. hominis detection
- Compatible with most open qPCR platforms
- Smaller sample input that allows remaining extracted samples to be used in other tests

Ordering Information

Order Online »

ENZ-GEN209-0100

100 tests

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[®] Ureaplasma spp. / M. genitalium / M.

hominis Assay is compatible with any properly calibrated qPCR thermal cycler capable to detect

fluorescence decay. It has been validated for use on the

QIAGEN Rotor-Gene Q

Contents AMPIGENE[®] HS Taq DNA Polymerase

AMPIGENE® dNTP Mix

AMPIPROBE® 5X Assay Buffer
AMPIPROBE® UMM Primer Mix
UMM Positive PCR Control
Negative PCR Control
Nuclease-free Water

Sensitivity In a validation study approved by the New York State

Department of Health, the AMPIPROBE® Ureaplasma spp. / M. genitalium / M. hominis Assay was determined to have the following percentage of Sensitivity:Ureaplasma spp. – 97.80%M. genitalium – 87.30M. hominis – 79.70%

Species Reactivity Mycoplasma genitalium, Mycoplasma hominis,

Ureaplasma spp.

Specificity

In a validation study approved by the New York State Department of Health, the AMPIPROBE® Ureaplasma spp. / M. genitalium / M. hominis Assay was determined to have the following percentage of Specificity:

Ureaplasma spp. - 95.00% M. genitalium - 100.00 M. hominis - 100.00%

Last modified: May 29, 2024



info-