

AMPIPROBE®

Ureaplasma spp. / M. genitalium / M. hominis Assay Kit

The AMPIPROBE® *Ureaplasma* spp. / *M. genitalium* / *M. hominis* (UMM) Assay is a real-time polymerase chain reaction (qPCR) assay for the qualitative detection of *Ureaplasma* spp. (*Ureaplasma parvum* and *Ureaplasma urealyticum*), *Mycoplasma genitalium*, and *Mycoplasma hominis* 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 *Ureaplasma* spp., *Mycoplasma genitalium*, and *Mycoplasma hominis* DNA. A positive PCR control consisting of a mixture of the target templates and a negative PCR control that results negative for UMM, but positive for the internal control (human β -globin). Please read the complete kit insert before performing this assay.

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 – 87.30% M. 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%



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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