

AMPIPROBE®

UP/Mg/Mh Extraction Controls

The AMPIPROBE® UP/Mg/Mh Extraction Controls are designed to ensure an efficient genomic DNA extraction process and detection of *Ureaplasma* spp. (*Ureaplasma parvum* and *Ureaplasma urealyticum*), *Mycoplasma genitalium*, and *Mycoplasma hominis*.

The AMPIPROBE® UP/Mg/Mh Extraction Controls are designed to be used during the genomic DNA extraction of *Ureaplasma* spp. (*Ureaplasma parvum* and *Ureaplasma urealyticum*), *Mycoplasma genitalium*, and *Mycoplasma hominis*.

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® UP/Mg/Mh Assay kit (Prod. No. ENZ-GEN209).

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.

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.

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

Ordering Information

[Order Online »](#)

ENZ-GEN210-0500

500µl

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® UP/Mg/Mh 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 QIAasympyony SP and Rotor-Gene Q.

Contents Negative Extraction Control
UP/Mg/MH Positive Extraction Control

Species Reactivity Mycoplasma genitalium, Mycoplasma hominis, Ureaplasma spp.



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)