

Bacteria, Fungi, and Mycoplasma Detection Kits

DNA extraction and real-time PCR based quality control

DNA Extraction

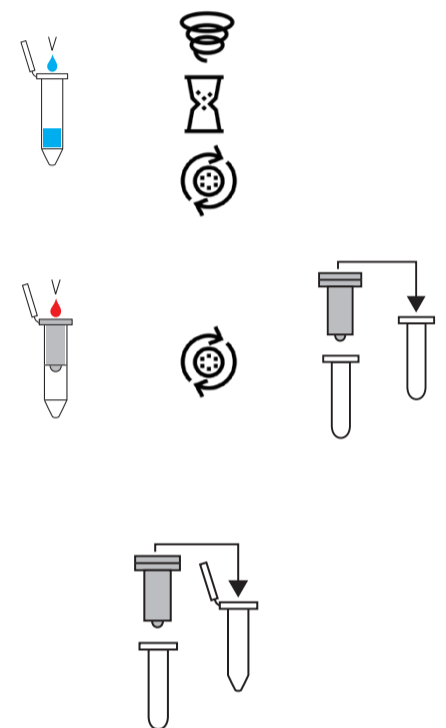
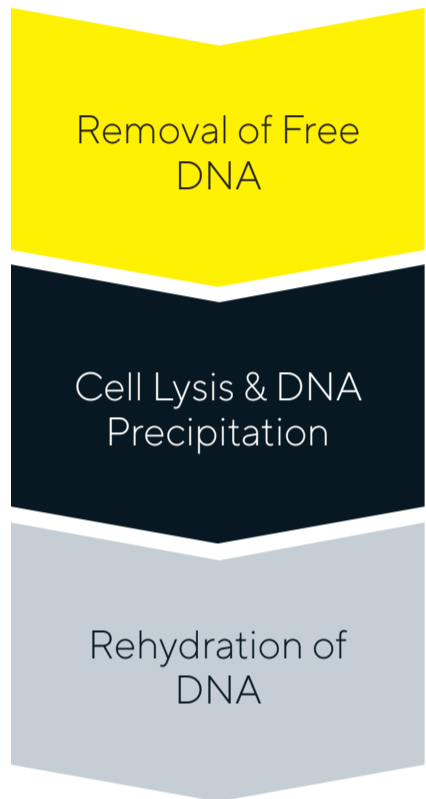
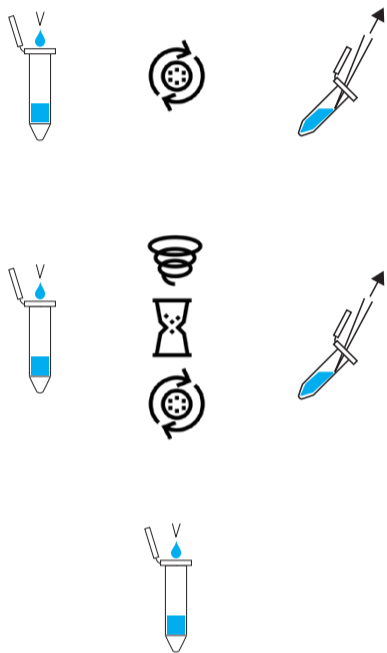
Microsart® ATMP Extraction
for efficient extraction of total (gram-positive & negative) bacterial and fungal DNA.

A DNA isolation with only a few handling steps to reduce contamination and DNA loss.



Microsart® AMP Extraction
for efficient extraction of Mycoplasma.

A column-based extraction designed to capture all DNA of Mycoplasma.



DNA Ready for PCR

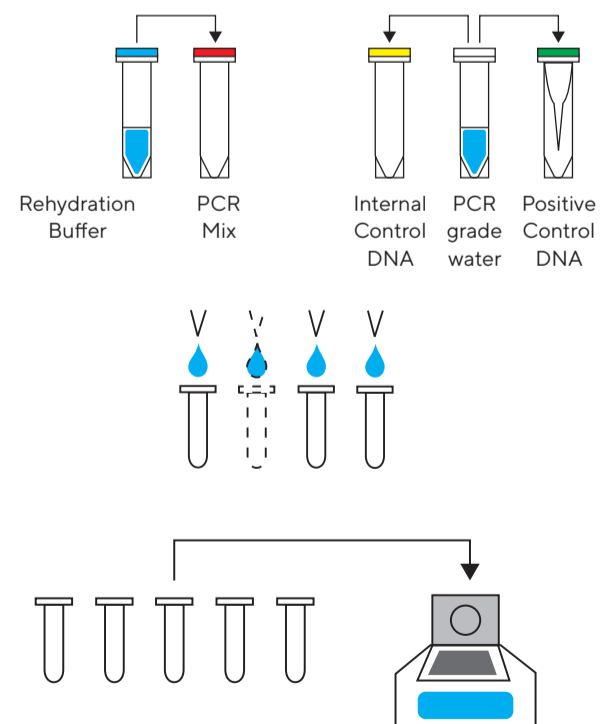
PCR Setup



Microsart® ATMP Bacteria & Microsart® ATMP Fungi
for comprehensive amplification of DNA derived from bacterial and fungal contamination.



Microsart® ATMP Mycoplasma
for specific amplification of DNA derived from Mycoplasma contamination.



Rapid Sterile Release/
Contamination Detection