

# Suspect Mycoplasma Contamination?



## Find out with the easy-to-use EZ-PCR™ Mycoplasma Detection Kit

BI's EZ -PCR Mycoplasma detection kit is a rapid, highly specific PCR-based platform, designed for the detection of a broad range of Mycoplasma species directly from the culture medium.

The kit contains an optimized complete PCR mix, which includes Mycoplasma-specific primers (to amplify the conserved 16S ribosomal RNA coding region within the Mycoplasma genome), dNTP's and Taq polymerase.

The kit also includes Internal Control (IC) DNA template (results in a 357bp PCR product, to verify a successful PCR run) as well as positive control DNA template (results in a 270bp PCR product on an agarose gel).

Samples can be prepared in less than 10 minutes, with accurate and rapid results, no special equipment is required.

### New and improved version with internal control to eliminate the possibility of PCR inhibition by the test samples

- Easy to use - direct addition of cell culture supernatant to PCR reaction - no DNA isolation/ purification steps required
- Reliable results - internal and positive control
- Detects over 100 Mycoplasma species\*
- Rapid results - less than 5 hours from sampling to final results
- Highly sensitive - detection of 10 CFU/ml



\* The primer set allows detection of various mycoplasma species common in cell culture: *M. fermentans*, *M. hyorhinis*, *M. arginini*, *M. orale*, *M. salivarium*, *M. hominis*, *M. pulmonis*, *M. arthritis*, *M. bovis*, *M. pneumoniae*, *M. pirum* and *M. capricolum*, as well as *Acholeplasma* and *Spiroplasma* species, with high sensitivity and specificity.

### Order Information

Product Name	Cat. No.	Size
EZ-PCR Mycoplasma Detection Kit	20-700-20	20 assays



**BI**  
Biological Industries  
Culture of Excellence

Biological Industries Israel Beit Haemek Ltd.  
Kibbutz Beit Haemek 2511500, Israel  
T.+972.4.9960595 F.+972.4.9968896  
Email: info@bioind.com  
[www.bioind.com](http://www.bioind.com)



**cGMP**  
Manufacturing  
Facility