

MYCOPLASMA DETECTION KIT

MycoStrip™

Mycoplasma

NEW

A new way to detect mycoplasma in your cell culture

- ❖ Fast and easy-to-use
- ❖ No special equipment needed
- ❖ Highly sensitive and reliable
- ❖ No false positives

Mycoplasma contamination is a major problem in cell culture that can greatly impact the reproducibility and validity of experimental results. InvivoGen's MycoStrip™ detects mycoplasma contamination using isothermal PCR and the results are rapidly visualized using a lateral flow detection cassette within 5 minutes.

MycoStrip™ Detection Kit



WWW.INVIVOGEN.COM/MYCOSTRIP

