

# VisionArray® MYCO

## PreCise Master Mix

For the amplification of  
mycobacterialspecific sequences

The VisionArray MYCO PreCise Master Mix is a ready to use PCR Master Mix for amplifying ITS and, in case of *M. tuberculosis* complex, IS6110 region of mycobacterial genomes including but not limited to those detected by the VisionArray MYCO Chip. The VisionArray MYCO PreCise Master Mix contains the VisionArray MYCO Primer Mix 1.0, dNTP/dUTP Solution, the VisionArray PreCise Taq DNA Polymerase, the PCR-Buffer, MgCl<sub>2</sub>, and the VisionArray Uracil-DNA Glycosylase. During PCR, specific sections of the ITS and IS6110 region of the mycobacterial genomes will be biotinylated. The hybridization and detection of the amplified sequences is subsequently performed using the VisionArray Detection Kit and the VisionArray MYCO Chip.

### Ordering Information

[Order Online »](#)

ZTV-ES-0008-50	50 tests
----------------	----------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Technical Info / Product Notes      For the Manual please [click here](#).



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)