

VisionArray® MYCO

Chip 2.0

For the specific detection of genus Mycobacterium and several clinically relevant mycobacterial species that have been produced with the help of the Vision

Array MYCO PreCise Master Mix 2.0. The VisionArray MYCO Chip 2.0 is intended to be used with a VisionArray Analysis Package for the qualitative detection and identification of PCR amplicons of the genera Mycobacterium, Mycobacteriodes, Mycolicibacillus, Mycolicibacter, and Mycolicibacterium as well as several additional clinically relevant mycobacterial species that have been produced with the help of the VisionArray MYCO PreCise Master Mix 2.0.

Ordering Information

[Order Online »](#)

ZTV-VA-0005-10	10ea
----------------	------

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 For the Manual please [click here](#).
Notes



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)