

VisionArray® FUNGI

PreCise Master Mix

1.0

For the amplification of fungi specific sequences

The VisionArray® FUNGI PreCise Master Mix 1.0 is intended to be used to amplify and biotinylate specific sections of the 16S-23S ITS region of fungi genomes by polymerase chain reaction (PCR). The VisionArray® FUNGI PreCise Master Mix 1.0 is designed to amplify fungi types including but not limited to those detected by the corresponding VisionArray® FUNGI Chips and genomic sequences of the human HLA-DQA1 gene as a PCR positive control.

Ordering Information

[Order Online »](#)

ZTV-ES-0009-50	0.75ml
----------------	--------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Quantity 50 tests

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com