

VisionArray® FUNGI

Chip 1.0

For the specific detection of 30 fungi types that have been produced with the help of the VisionArray FUNGI PreCise Master Mix

The VisionArray® FUNGI Chip 1.0 is intended to be used for the qualitative detection and genotyping of PCR-amplificates of 30 clinically relevant fungi genotypes that have been produced with the help of the VisionArray® FUNGI PreCise Master Mix 1.0 (Prod. No. ES-0009-50) from formalin-fixed, paraffin-embedded specimens. The chip is intended to be used in combination with a VisionArray® Software.

Ordering Information

[Order Online »](#)

ZTV-VA-0006-10	10ea
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

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

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