

# T. vaginalis native antigen

Cell lysate ideal for use as a broad positive control

## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-630-130-1000 | 1ml |
|------------------|-----|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Trichomonas vaginalis native antigen

**Appearance** Cloudy liquid.

**Formulation** Liquid. In sterile distilled water.

**Purity** ≥99%

**Purity Detail** Lysed suspension of whole protozoa.

**Source** From *Trichomonas vaginalis*, strain C-1:NIH. Inactivated by freezing/thawing in water.

**Last modified: May 29, 2024**



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)