

# Microcystins (Adda specific) ELISA kit

Patented ELISA for rapid detection of microcystins and nodularins.

Polyclonal antibody prepared against the Adda moiety binds to microcystins and nodularins, allowing the congener-independent determination of these toxins and many of its congeners, and does not cross-react with other non-related toxins or compounds. No pre-sample preparation required. Total time for measurement is less than 2.5 hours. Enables simultaneous measurement of multiple samples at reasonable costs. *May require a license for import, please [contact us](#) for more information.*

Microcystins are extremely toxic compounds produced by cyanobacteria (blue-green algae), belonging to species of *Microcystis*, *Oscillatoria*, *Anabaena* and *Nostoc*. The contamination of drinking water or water of recreational areas can cause severe health problems to exposed humans and animals. Microcystins possess a cyclic heptapeptide structure of the general composition cyclo(-D-Ala-L-X-D-*erythro*- $\beta$ -methylisoAsp-L-Y-Adda-D-iso-Glu-N-methyldehydroAla), where Adda is the unusual C20 aa 3-amino-9-methoxy-2,6,8-trimethyl-10-phenyldeca-4,6-dienoic acid and X and Y are variable L-aa.

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

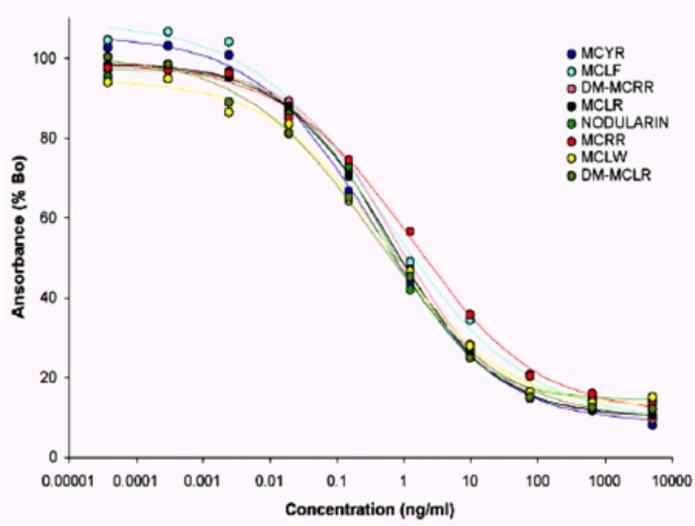
ALX-850-319-KI01

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Specific – antibody to Adda moiety specifically binds microcystin & nodularin toxins
- Sensitive – measure as little as 0.1ng/ml toxin
- Convenient – no pre-sample preparation required
- Rapid – analyze 41 duplicate samples in <2.5 hours



Cross-reactivity pattern against microcystins and nodularin congeners

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application ELISA

Application Notes For the quantitative and sensitive congener-independent detection of microcystins and nodularin in water samples.

Assay Time 2 hour 30 minutes

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Sensitivity 0.1ng/ml (range 0.15 to 5ng/ml)

Species Reactivity Species independent

Technical Info / Product Notes Patented technology. *U.S. Patent No. 6,967,240.*  
*Worldwide Patent PCT WO 01/18059 A2.*

Wavelength 450 nm

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)