

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*- β -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 »](#)

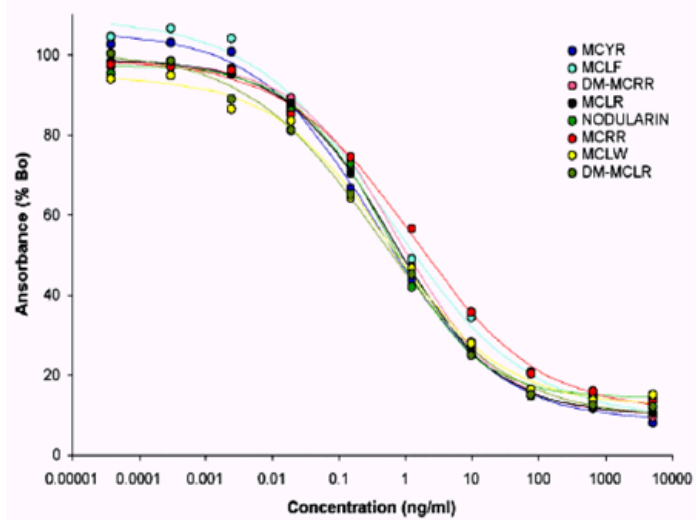
ALX-850-319-KI01

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

Long Term Storage +4°C

Regulatory Status RUO - Research Use Only



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

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.