

Hepatotox Set 1

May require a license for import, please [contact us](#) for more information.

Ordering Information

[Order Online »](#)

ALX-850-325-KI01	1Set
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Contents Set contains separate vials of: Microcystin-LA (25µg); Microcystin-LF (25µg); Microcystin-LR (100µg); Microcystin-LW (25µg); Microcystin-LY (25µg); Microcystin-RR (50µg); Microcystin-YR (25µg); Nodularin (100µg).

Purity ≥95% (HPLC)

Source Isolated from *Microcystis aeruginosa* and *Nodularia spumigena*.

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)