

Kinetoplast DNA (catenated)

KDNA is an aggregate of interlocked DNA minicircles (mostly 2.5kb) from the kinetoplast of the insect trypanosome *Crithidia fasciculata*, that form extremely large networks of high molecular weight. As a result, these networks fail to enter an agarose gel.

Ordering Information

[Order Online »](#)

TOP-TG2013-3	250µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name KDNA (catenated)



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)