

DNA, (purified) (DNAse I activated)

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-840-040-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

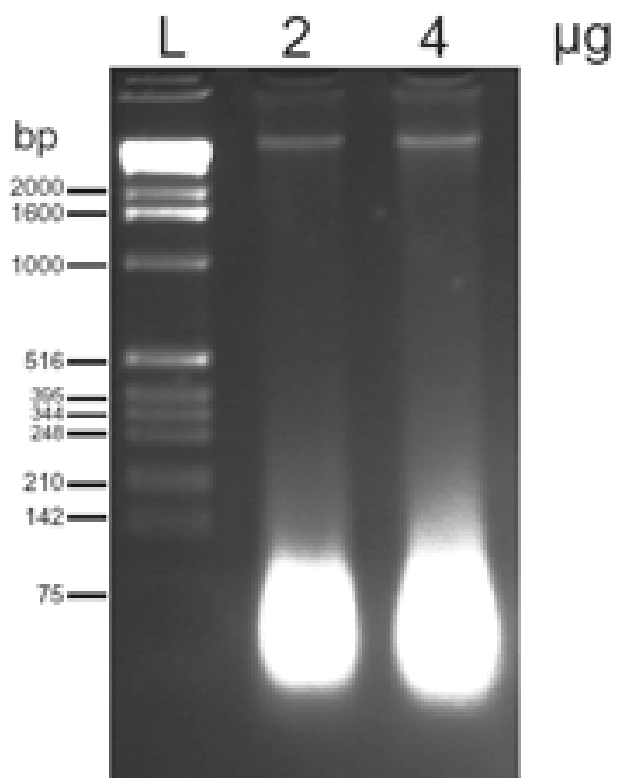


Figure: Agarose gel of DNA (purified) (DNase I activated) (Prod. No. [ALX-840-040](#)) (Lot L20796).",DNA

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes	For use with PARP-1 (Prod. No. ALX-201-063) or PARP-2 (Prod. No. ALX-201-064) <i>in vitro</i> activity tests. For details please see protocol.
Concentration	0.2mg/ml
Formulation	Liquid. In TRIS-EDTA.
Purity	≥99% (SDS-PAGE)
Source	Purified from calf thymus and activated by digestion with DNase I.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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