

AMPIGENE[®] DNA

Ladder 250-10,000 bp

Ready to use DNA standards for quantifying PCR yield

AMPIGENE[®] DNA Ladder 250-10,000 bp contains DNA fragments ranging from 250bp to 10,000bp that can be used as molecular weight standards for agarose gel electrophoresis. The bands are evenly spaced and are intensified at 1,000bp and 3,000bp to serve as easily identifiable reference bands. The approximate mass of DNA in each of the bands is provided, assuming 5µl load, for determining the approximate mass of DNA in samples with comparable intensity and size.

Ordering Information

[Order Online »](#)

ENZ-GEN103-0100	100lanes
-----------------	----------

Manuals, SDS & CofA

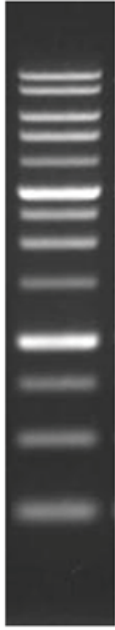
[View Online »](#)

- Ready to use format can be loaded straight onto a gel
- Evenly spaced bands with intensified reference bands
- Room temperature stable

DNA Mass
(ng/5 μ l)

28
28
28
28
18
92
34
34
20

92
23
30
45



Base Pairs

10,000
8,000
6,000
5,000
4,000
3,000
2,500
2,000
1,500

1,000
750
500
250

1% TAE agarose gel

Handling & Storage

Short Term Storage Ambient

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

For size determination and DNA quantification using agarose electrophoresis

Contents

Ready-to-use AMPIGENE[®] DNA Ladder 250-10,000 bp
Sample Loading Buffer

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