

AMPIGENE[®] DNA

Ladder 50-1,500 bp

Ready to use DNA standards for quantifying
PCR yield

AMPIGENE[®] DNA Ladder 50-1,500 bp contains DNA fragments ranging from 50bp to 1,500bp that can be used as molecular weight standards for agarose gel electrophoresis. The bands are evenly spaced and are intensified at 200bp, 500bp and 1,200bp 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-GEN102-0100	100lanes
-----------------	----------

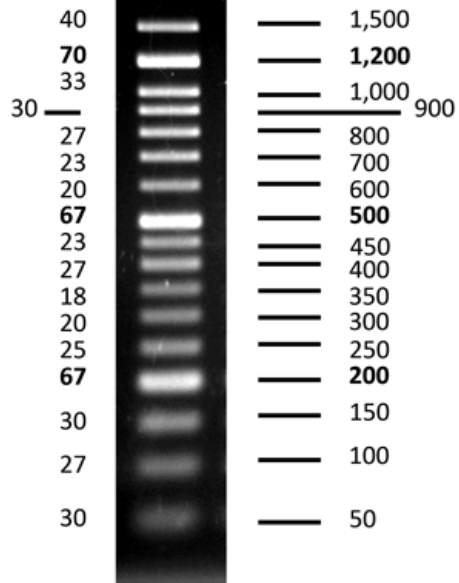
Manuals, SDS & CofA

[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/5μl)

Base Pairs



2% TAE agarose gel

Short Term Storage	Ambient
--------------------	---------

Long Term Storage -20°C

Shipping Ambient Temperature

Product Details

Application Notes

For size determination and DNA quantification using agarose electrophoresis

Contents

Ready-to-use AMPIGENE[®] DNA Ladder 50-1,500 bp
Sample Loading Buffer

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