

AMPIGENE® DNA

Ladder 100-1,500 bp

Ready to use DNA standards for quantifying
PCR yield

AMPIGENE® DNA Ladder 100-1,500 bp contains DNA fragments ranging from 100bp to 1,500bp that can be used as molecular weight standards for agarose gel electrophoresis. The bands are evenly spaced and are intensified at 500bp and 1,500bp 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-GEN101-0100	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)

Base Pairs

72.5		1,500
50		1,000
40		900
40		800
27.5		700
30		600
85		500
40		400
35		300
40		200
40		100



1.7% TAE agarose gel

Short Term Storage	Ambient
Long Term Storage	-20°C
Shipping	Ambient Temperature

Product Details

For size determination and DNA quantification using agarose electrophoresis

Ready-to-use AMPIGENE® DNA Ladder 100-1,500 bp
Sample Loading Buffer



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