

Product Information

Ready-to-Use DNA Ladder

Cat. No.	Product	Unit Size
D001	1kb DNA Ladder	250 -L
D002	1kb plus DNA Ladder	250 -L
D003	100bp plus DNA Ladder	250 -L

Storage upon receipt:

Store at 4°C for 6 months, or -20 °C for 24 months.

Product Description

The 1 kb DNA Ladder is suitable for sizing linear double-stranded DNA fragments from 1 kb to 10 kb. The ladder contains a total of eight bands: 1000, 2000, 3000, 4000, 5000, 6000, 8000, 10000bp. The 4 kb band has a concentration of 100 ng/5 μ L with increased intensity to provide internal orientation. The other bands have a concentration of 50 ng/5 μ L. The 1 kb DNA Ladder is provided in 1X loading buffer.

The 1 kb plus DNA Ladder is suitable for sizing linear double-stranded DNA fragments from 300 bp to 10 kb. The ladder contains a total of twelve bands: 300, 500, 800, 1000, 1500, 2000, 3000, 4000, 5000, 6000, 8000, 10000bp. The 2 kb band has a concentration of 100 ng/5 μ L with increased intensity to provide internal orientation. The other bands have a concentration of 50 ng/5 mL. The 1 kb plus DNA Ladder is provided in 1X loading buffer.

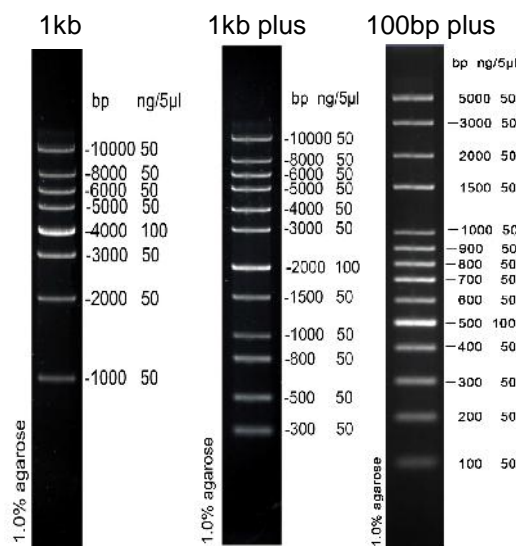
The 100bp plus DNA Ladder is suitable for sizing linear double-stranded DNA fragments from 100 bp to 5000 bp. The ladder contains a total of fourteen bands: 100, 200, 300, 400, 500, 600, 700, 800, 900, 1000, 1500, 2000, 3000, 5000bp.

The 500 bp band has a concentration of 100 ng/5 μ L with increased intensity to provide internal orientation. The other bands have a concentration of 50 ng/5 μ L. The 100 bp plus DNA Ladder is provided in 1X loading buffer.

The ladders are generated from PCR and restriction enzyme digestion of double stranded DNA. The DNA is purified by phenol extraction, and equilibrated to 10 mM Tris-HCl (pH 8.0) and 1 mM EDTA.

Product Protocols

The Ready-to-Use DNA Ladders are supplied in a ready-to-load format. For agarose gel electrophoresis, load ~1 μ L of DNA ladder per mm lane.



Related Products

Cat. No	Product Name	Unit Size
N100	GreenView DNA Gel Stain, 10,000X in H ₂ O	500 μ L
N101	GreenView Plus DNA Gel Stain, 10,000X in DMSO	500 μ L
N102	RedView DNA Gel Stain, 10,000X in DMSO	500 μ L
N103	GreenView™ Ultra DNA Gel Stain, 10,000X in DMSO	500 μ L