

Product Information

1 KB DNA Ladder

Catalog Number: 31021

Unit Size: 150 applications

Kit Contents

Component A: 1 KB DNA Ladder (30 µg/300 µl)
Component B: 6X Loading buffer with two tracking dyes (1.5 ml)

Storage

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

Storage Buffer

10 mM Tris-HCl (pH 8.0), 1 mM EDTA

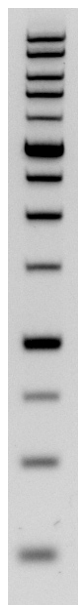
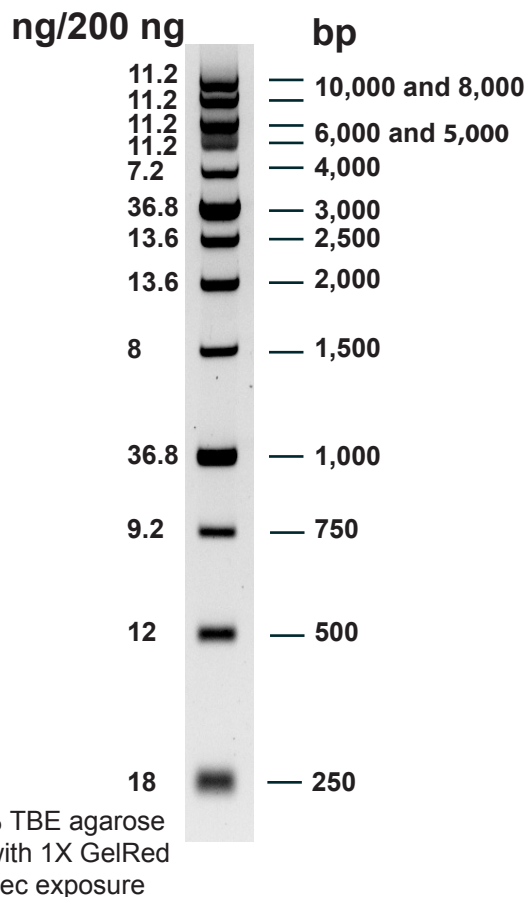
Product Description

The 1 KB DNA Ladder is suitable for sizing linear double-stranded DNA fragments from 250 bp to 10 kb. The ladder consists of 13 bands that are generated from PCR followed by restriction enzyme digestion, phenol extraction, and equilibration to 10 mM Tris-HCl (pH 8.0) and 1 mM EDTA. The 1 kb and 3 kb bands have increased intensity to provide internal orientation.

Recommended Procedure

For agarose gel electrophoresis, it is recommended the 1 KB DNA ladder be used in combination with the included 6X loading buffer (component B), at a final concentration of 1X. Load 0.2 µg of DNA ladder per mm lane.

The 1 KB DNA Ladder has not been tested with polyacrylamide gels and was not designed for quantification of unknown DNA electrophoresed on the same gel.



**200 ng 1 KB DNA ladder
1% TAE agarose gel with
1X GelGreen
0.5 second exposure**

Related Products

Catalog number	Product Description
41001	GelRed™ Nucleic Acid Gel Stain, 3X in water
41002	GelRed™ Nucleic Acid Gel Stain, 10,000X in DMSO
41003	GelRed™ Nucleic Acid Gel Stain, 10,000X in water
41004	GelGreen™ Nucleic Acid Gel Stain, 10,000X in DMSO
41005	GelGreen™ Nucleic Acid Gel Stain, 10,000X in water
41006	TBE, 5X
31000	EvaGreen® dye, 20X in water
31003	Fast EvaGreen® qPCR Master Mix (200 rxn)
31014	Fast Plus EvaGreen® qPCR Master Mix - Low Rox (200 rxn)
31015	Fast Plus EvaGreen® qPCR Master Mix - High Rox (200 rxn)
31020	Fast Plus EvaGreen® qPCR Master Mix (200 rxn)
31022	Ready-to-Use 1 KB DNA Ladder

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.