

ReadiUse™ GeneRuler 1 kb DNA Ladder

Catalog number: 60055, 60056 Unit size: 100 ul, 2X250 ul

Component	Storage	Amount (Cat No. 60055)	Amount (Cat No. 60056)
ReadiUse™ GeneRuler 1 kb DNA Ladder	Freeze (< -15 °C), Minimize light exposure	100 μL	2X250 μL

OVERVIEW

ReadiUse™ GeneRuler 1 kb DNA ladder is a ready-to-use DNA ladder, which is pre-mixed with loading dye for direct gel loading. It is composed of 14 individual DNA fragments: 25k, 10k, 8k, 6k, 5k, 4k, 3k, 2.5k, 2k, 1.5k, 1k, 750, 500, and 250 bp derived from a mixture of PCR products and specifically digested plasmid DNA. This product contains two enhanced bands (3 kb and 1 kb) for easy reference. In addition, two tracking dyes, Xylene cyanol FF and Bromophenol blue which mimic the migration of 4,000 bp and 500 bp dsDNA during electrophoresis are added for real time monitoring.

SAMPLE EXPERIMENTAL PROTOCOL

ReadiUseTM GeneRuler 1 kb DNA Ladder is pre-formulated and ready to use. 5 μ L per well is recommended to use for each gel.

EXAMPLE DATA ANALYSIS AND FIGURES

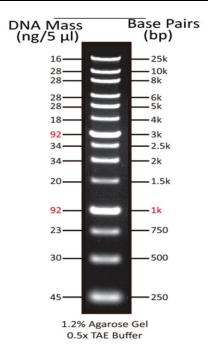


Figure 1. ReadiUse™ GeneRuler 1 kb DNA Ladder (5 uL/well) was run on 1.2% agarose gel with 0.5X TAE Buffer.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.