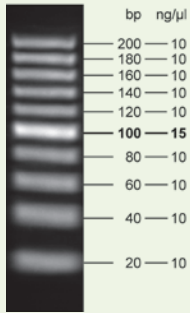
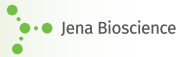


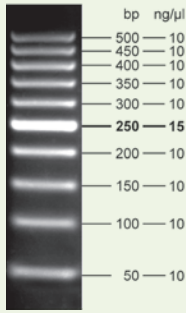
20 bp DNA Ladder



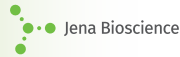
3.6% Agarose



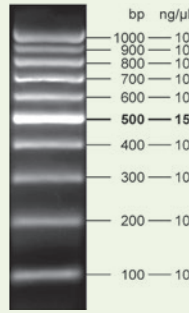
50 bp DNA Ladder



3.0% Agarose



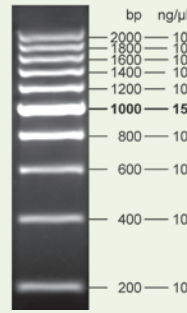
100 bp DNA Ladder



2.4% Agarose



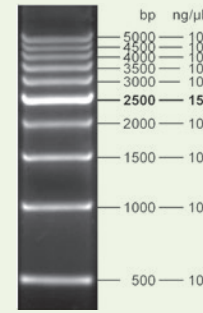
200 bp DNA Ladder



1.8% Agarose



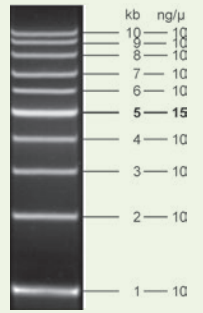
500 bp DNA Ladder



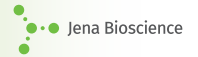
1.2% Agarose



1 kb DNA Ladder



0.6% Agarose



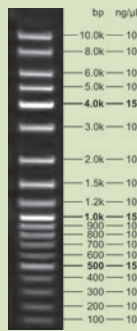
Log Scale DNA Ladders

are well suited for standard gel electrophoresis applications. The ladders can be combined if small and large fragments need to be analyzed on the same gel.

Linear Scale DNA Ladders

show virtually uniform spacing over a large fragment size range. →

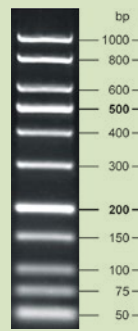
Long Range DNA Ladder



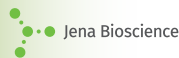
1.0% Agarose



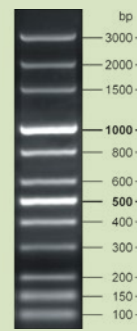
Low Range DNA Ladder



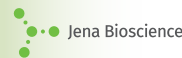
3.0% Agarose



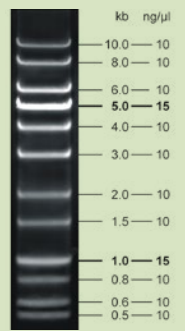
Mid Range DNA Ladder



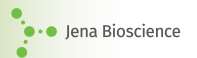
1.8% Agarose



High Range DNA Ladder



0.6% Agarose



Jena Bioscience DNA Ladders for easy-to-track sample sizing and quantification after PCR or restriction digest. All DNA Ladders are provided ready-to-use containing tracking dyes (orange/blue). In addition, fluorescent DNA Ladders are supplied with EveGreen™ fluorescent intercalator dye.

Log Scale DNA Ladders, ready-to-load

Product	Length	Tracking dye	Concentration	S-Pack [μl lanes]	L-Pack [μl lanes]	Unstained DNA Ladder	Fluorescent DNA Ladders
20 bp DNA Ladder	20 – 200 bp	orange	105 ng/μl	500 100	5x 500 500	M-212	M-232
50 bp DNA Ladder	50 – 500 bp	orange	105 ng/μl	500 100	5x 500 500	M-213	M-233
100 bp DNA Ladder	100 – 1000 bp	orange / blue	105 ng/μl	500 100	5x 500 500	M-214	M-234
200 bp DNA Ladder	200 – 2000 bp	blue / blue	105 ng/μl	500 100	5x 500 500	M-215	M-235
500 bp DNA Ladder	500 – 5000 bp	blue / blue	105 ng/μl	500 100	5x 500 500	M-216	M-236
1 kb DNA Ladder	1 – 10 kb	blue / blue	105 ng/μl	500 100	5x 500 500	M-217	M-237

Linear Scale DNA Ladders, ready-to-load

Product	Length	Tracking dye	Concentration	S-Pack [μl lanes]	L-Pack [μl lanes]	Unstained DNA Ladder	Fluorescent DNA Ladders
Long Range DNA Ladder	100 bp – 10 kb	orange / blue	205 ng/μl	500 100	5x 500 500	M-205	M-225
Low Range DNA Ladder	50 bp – 1 kb	orange / blue	120 ng/μl	500 100	5x 500 500	M-202	M-222
Mid Range DNA Ladder	100 bp – 3 kb	orange / blue	130 ng/μl	500 100	5x 500 500	M-203	M-223
High Range DNA Ladder	500 bp – 10 kb	blue / blue	130 ng/μl	500 100	5x 500 500	M-204	M-224

