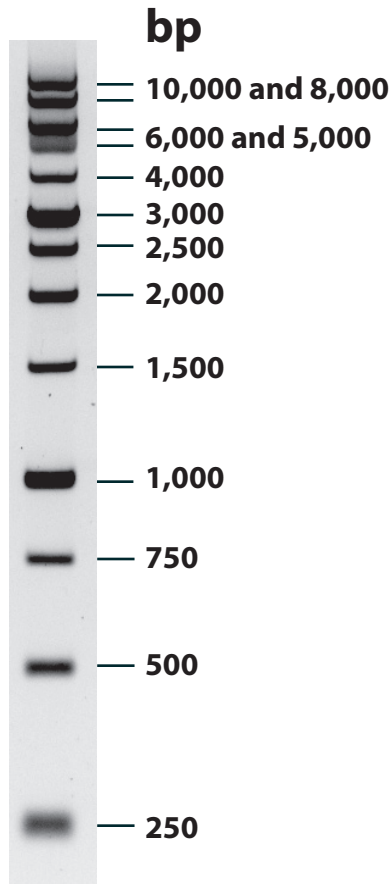


1 KB DNA Ladder: Compatible with GelRed™ and GelGreen™ Dyes

200 ng



0.8% TBE
agarose gel with
1X GelRed

Biotium now introduces DNA ladders for use in agarose gel electrophoresis.

- **Compatible with GelRed™ and GelGreen™ Products!**
- **Range:** 250 bp to 10 kb
- **Number of bands:** 13
- **Concentration:** 100 ng/μl
- **Recommended use:** 200 ng per well
- **Available in 2 formats:** ladder alone and a ready-to-load format

Cat.#	Product Name	Size	Price
31021	1 KB DNA Ladder	30 μg	\$55
31022	Ready-to-Use 1 KB DNA Ladder	150 applications	\$65