

---

## EpiQuik 1 Kb DNA Ladder

(Catalog No. R12006)

---

### Description

EpiQuik 1 Kb DNA Ladder consists of 13 fragments: 300 bp, 500 bp, 700 bp, 1000 bp, 1500 bp, 2000 bp, 2500 bp, 3000 bp, 4000 bp, 5000 bp, 6000 bp, 8000 bp and 10000 bp. EpiQuik 1 kb DNA Ladder is a ready-to-use molecular weight marker suitable for DNA fragment size determination on gel electrophoresis. EpiQuik 1 kb DNA Ladder is formulated to run accurately and to provide crisp band patterns. It contains two dyes: bromophenol blue and xylene cyanole which serve as visual aid to monitor the progress of migration during agarose gel electrophoresis.

### Applications

For precise sizing of double-stranded DNA fragments from 300 to 10000 bp on agarose gels.

### Storage Conditions

-20°C up to one year or 4°C for one month.

### Safety Warnings and Precautions

This product is designed for research purposes and in vitro use only. According to common laboratory safety practice, it is recommended to wear protective clothing, gloves and safety glasses. Please refer to [www.epigentek.com](http://www.epigentek.com) for Material Safety Data Sheet of the product. Some applications this product is used in may require a license which is not provided by the purchase of this product. Users should obtain the license if required.

### Ordering Information

Products	Size	Cat. No.
EpiQuik 1 Kb DNA Ladder	500 µl	R12006-1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*