
EpiQuik Mid-Range DNA Ladders Express

(Catalog No. R12004)

Description

EpiQuik Mid-Range DNA Ladder consists of ten fragments: 300 bp, 500 bp, 500 bp, 1000 bp, 1500 bp, 2000 bp, 2500 bp, 3000 bp, 4000 bp and 5000 bp. Supplied in sufficient quantity for 100 loadings using 4-5 μ l per loading. Not designed for quantifying of DNA concentration in a sample. Ladders available premixed as ready-to-use format premixed with loading-dye-solution (EXPRESS) or plain ladder to be mixed with loading dye of choice.

Applications

For precise sizing of double-stranded DNA fragments from 300 to 5000 bp on agarose gels.

Storage Conditions

Stable at room temperature for 12 months.

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 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 Mid-Range DNA Ladders Express	500 μ l	R12004-1

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