

EpiQuik 100 bp DNA Ladder Express

(Catalog No. R12002)

Description

The EpiQuik 100 bp DNA Ladder Express can be used for precise visual sizing of double-stranded DNA fragments from 100 to 1000 bp on agarose gels. Note: Not designed for quantifying DNA concentration in a sample.

Composition

Consists of multiple repeats of 100 bp fragments from 100 bp up to 1000 bp. Ready-to-use: the ladder is premixed with Loading Dye Solution for direct loading onto agarose gels.

Applications

For precise sizing of double-stranded DNA fragments from 100 to 1000 bp on agarose gels.

Storage Conditions

Stable at room temperature for 12 months.

Recommendations

For best results, please load 10-20 μ l of the EpiQuik 100 bp DNA Ladder per well.

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

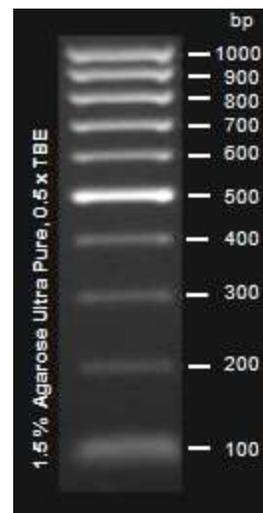
EpiQuik 100 bp DNA Ladder Express

Size

500 μ l

Cat. No.

R12002-1



▲ Agarose gel depicting the EpiQuik 100 bp DNA Ladder Express.

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