



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

DNA Gel Extraction Kit

A spin column-based DNA extraction kit for purifying DNA from agarose gels.



Product attributes

Product Description

Biotium's DNA Gel Extraction Kit is a silica-gel spin column based DNA extraction kit designed to purify DNA fragments from agarose gels.

- Purify DNA fragments from 40 bp to 40 kb
- Maximum binding capacity of each column is 10 ug DNA
- Purify DNA from agarose gels made with TAE or TBE buffer

The DNA Gel Extraction Kit removes GelRed® and GelGreen® Nucleic Acid Gel Stains, along with other commonly used nucleic acid gel stains. The purified DNA is compatible with common downstream applications such as transformation, sequencing, PCR, and restriction digests.

This datasheet was generated on June 15, 2026 at 02:40:24 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dna-gel-extraction-kit/>