## **DNA Gel Extraction Kit**

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



## **Product attributes**

Call us: 800-304-5357 Email: btinfo@biotium.com

## **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 November 11, 2025 at 11:47:36 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dna-gel-extraction-kit/