

ReadiPrep™ Spin Column for PCR Product Cleanup *Compatible with Qiagen QIAQuick PCR Cleanup Kit*

Catalog number: 60510
Unit size: 10 Columns

Component	Storage	Amount (Cat No. 60510)
ReadiPrep™ Spin Column for PCR Product Cleanup	Room temperature (10-25 °C)	10 Columns

OVERVIEW

The ReadiPrep™ Spin Column for PCR Product Cleanup is a quick and reliable tool designed to purify DNA from PCR reactions and other enzymatic processes. This column removes primers, unused nucleotides, salts, and other contaminants that might be present in the sample to yield high-quality DNA suitable for a variety of downstream applications.

These columns have a pre-packed silica-based membrane to selectively bind DNA for the removal of impurities. The protocol is straightforward and requires minimal hands-on time, making it ideal for quick and reproducible workflows.

It is compatible with a various sample type and is optimized for use in downstream applications such as DNA sequencing, cloning, ligation, and quantitative PCR (qPCR).

AT A GLANCE

1. Add 5 volumes of binding buffer into PCR sample.
2. Apply the mix to ReadiPrep™ Spin Column and centrifuge the column.
3. Wash the column with wash buffer.
4. Elute the sample from the column with elution buffer.

Note: Binding, wash and elution buffers are not provided in #60510.

SAMPLE EXPERIMENTAL PROTOCOL

Refer the protocol of the ReadiPrep™ PCR Purification Kit (cat #60511)
<https://www.aatbio.com/products/readiprep-pcr-purification-kit?unit=60511>

EXAMPLE DATA ANALYSIS AND FIGURES

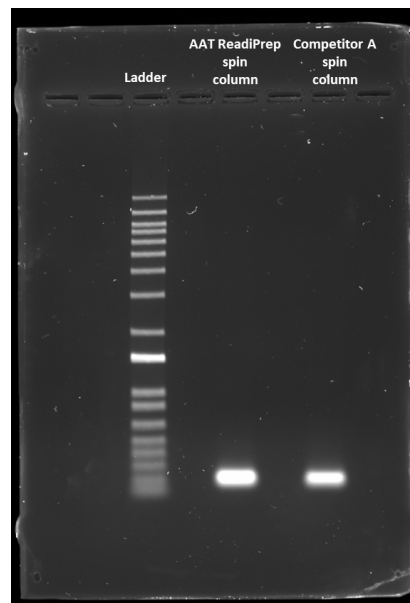


Figure 1. The 18S rRNA PCR product was purified using the ReadiPrep™ Spin Column for PCR Cleanup (Cat# 60510) or a comparable column from Competitor A. The purified products were electrophoresed on a 1% agarose gel in 1X TAE buffer and subsequently stained with Gelite Safe DNA Gel Stain (Cat# 17700).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.