

ReadiPrep™ PCR Purification Kit

 Catalog number: 60511
 Unit size: 25 Preps

Component	Storage	Amount (Cat No. 60511)
Component A: ReadiPrep™ Binding Buffer	Refrigerated (2-8 °C)	1 bottle (15 mL)
Component B: ReadiPrep™ Spin Column	Room temperature (10-25 °C)	25 columns
Component C: ReadiPrep™ Wash Buffer	Refrigerated (2-8 °C)	1 bottle (10 mL)
Component D: ReadiPrep™ Elution Buffer	Refrigerated (2-8 °C)	1 bottle (5 mL)

OVERVIEW

The ReadiPrep™ PCR Purification Kit is a fast and reliable DNA cleanup system designed for the purification of amplified DNA from PCR reactions. Using a silica membrane-based spin column, the kit efficiently removes primers, unused dNTPs, salts, enzymes, and other contaminants to deliver high-purity DNA ready for downstream applications. It is equivalent to QIAquick PCR purification kit.

This PCR cleanup kit is optimized for a wide range of molecular biology workflows including DNA sequencing, cloning, ligation, and quantitative PCR (qPCR). Its user-friendly protocol minimizes hands-on time and ensures consistent, reproducible results—ideal for both routine and high-throughput labs. The ReadiPrep™ PCR Purification Kit offers the speed, simplicity, and performance required for efficient and high-quality DNA recovery.

AT A GLANCE

1. Add 5 volumes of ReadiPrep™ Binding Buffer into PCR sample.
2. Apply the mix to ReadiPrep™ Spin Column and centrifuge the column.
3. Wash the column with ReadiPrep™ Wash Buffer working solution.
4. Elute the sample from the column with ReadiPrep™ Elution Buffer.

PREPARATION OF WORKING SOLUTION

1. Prepare ReadiPrep™ Wash Buffer working solution by adding 200 μ L of ReadiPrep™ Wash Buffer (Component C) into 800 μ L of Ethanol (96-100%, not provided).

Note: 1 mL of working solution is enough for 1 prep.

SAMPLE EXPERIMENTAL PROTOCOL

1. Add 5 volumes of ReadiPrep™ Binding Buffer (Component A) to 1 volume of the PCR sample, and then mix well. E.g. Add 500 μ L of ReadiPrep™ Binding Buffer to 100 μ L of PCR sample.
Note: For samples lower than 100 μ L volume, add molecular grade water to make up the volume to 100 μ L.
2. Apply the sample to ReadiPrep™ Spin Column (Component B) and centrifuge the column at 13,000 rpm for 60 seconds.
Note: For better binding, samples in columns can be incubated for 1 minute before centrifugation.
3. Discard the flow-through and place the column back into the collection tube.
4. Add 600 μ L of ReadiPrep™ Wash Buffer working solution and centrifuge the column at 13,000 rpm for 60 seconds.
Note: For additional salt removal, wash buffer in columns can be incubated for 1 minute before performing centrifuge.

Note: To prevent salt carryover for sensitive procedures, additional wash steps can be performed with 350 μ L of ReadiPrep™ Wash Buffer working solution and repeat step-4.

5. Remove the wash buffer and perform an additional centrifuge at 13,000 rpm for 2 minutes to dry the columns completely.
6. Add 30 μ L of ReadiPrep™ Elution Buffer (Component D) to the center of the column and centrifuge the column at 13,000 rpm for 60 seconds. To increase yield, perform another elution step with 30 μ L of ReadiPrep™ Elution Buffer and centrifuge.

Note: For better yield, elution buffer in columns can be incubated for 1 minute before performing centrifuge.

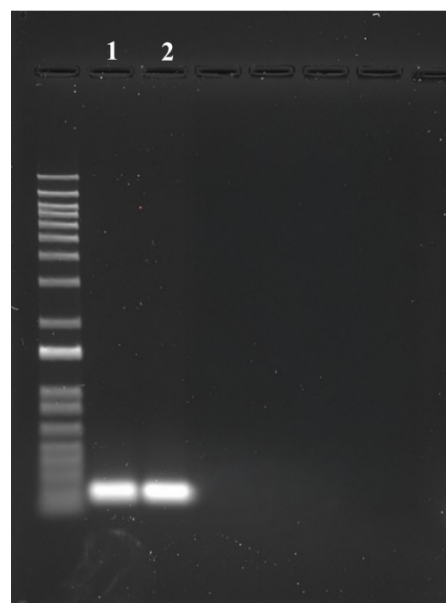
EXAMPLE DATA ANALYSIS AND FIGURES


Figure 1. Comparison of 18S rRNA PCR product purified with Competitor A's purification kit (Labeled as 1) and ReadiPrep™ PCR clean up kit (cat# 60511, labeled as 2). Purified products were run on 1% agarose gel with 1X TAE buffer followed by Gelite safe DNA gel staining (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.