
QuickChiP™ DNA Purification Kit (100 purification)

Catalog No.: NBP2-29498

Kit Contents:

Membrane Binding Buffer -----	50 ml
Wash Buffer -----	40 ml
Nuclease Free water -----	20 ml
Binding Columns -----	100
Collection Tubes -----	100

Storage: Kit components should be stored at room temperature.

Applications: DNA purification from PCR samples.

Description: This protocol is designed to purify single or double stranded DNA fragments from PCR and other enzymatic reactions. Fragments ranging from 100 bp to 10 kb can be purified from primers, nucleotides and salt using Novus QuickChiP™ DNA purification kit.

Preparation of Wash Buffer: Add 68 ml of 95% ethanol before using Wash Buffer (**Ethanol is not supplied in the kit**).

Procedure:

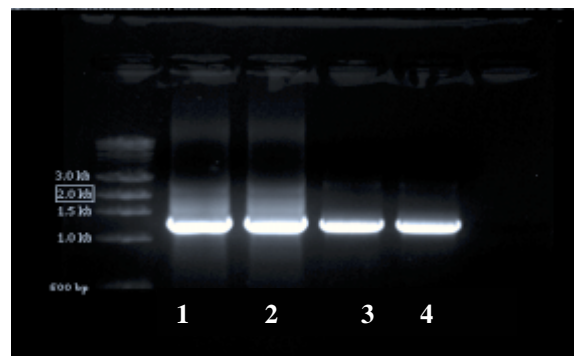
1. Amplify the target of choice using standard amplification condition in the total volume up to 200 µl.
2. After the PCR reaction is completed, transfer them to 1.5 ml centrifuge tube and add 300 µl of Membrane Binding Solution to the PCR reaction.
3. Place a spin column in a 2 ml collection tube (provided in the kit).
4. Transfer prepared PCR product into the spin column and incubate for 1 minute at room temperature.
5. Centrifuge at 16,000 x g (14,000rpm) in a microcentrifuge for 1 minute at room temperature. Remove the spin column from the collection tube and discard the flow-through. Reinsert the spin column into the collection tube.
7. Add 700 µl of Column Wash Solution with 95% Ethanol. Centrifuge at 16,000 x g (14,000 rpm) in a microcentrifuge for 30 seconds at room temperature. Remove the

spin column from the collection tube and discard the flow-through and place the spin column back into the same tube.

8. Centrifuge again at 16,000 x g (14,000 rpm) in microcentrifuge for 30 seconds at room temperature to dry out the spin column.
9. Place the spin column into a clean 1.5 ml microcentrifuge tube. Apply 50 µl of Elution buffer directly to the center of the column without touching the membrane with pipette tip. Incubate at room temperature for 1 minute. Centrifuge at 16,000 x g (14,000 rpm) in a microcentrifuge for 1 minute at room temperature. The collected sample contains PCR products free of primers, free nucleotides, and salt. The eluted PCR products can be analyzed by agarose gel electrophoresis.

Note: If the eluted DNA need to be more concentrated, the DNA can be eluted in as little as 20 µl of nuclease-free water.

Figure 1. Agarose gel electrophoresis of PCR products.



Lane #1 and 2 are unpurified PCR samples.
Lane #3 and 4 are purified PCR samples using DNA purification kit.