PCR & Gel Clean-up Kit

Quick, easy kit for rapid concentration and clean-up of PCR products

PCR & Gel clean-up columns use silica membrane technology in a mini spin column format to deliver easy and efficient DNA purification (removal of unused, single stranded primers) from PCR products, agarose gels, and labeling reactions.

Sample material: <400 µl PCR reaction mixture, <400mg agarose gel

Fragment size: 50bp - ~20kbp

Recovery: 70 - 95%

Elution volume: 15 - 50 μl

Preparation time: 10 min for 6 preps

Binding capacity: 25 µg

Handling: Centrifugation, vacuum Typical purity (A260/A280): 1.8 – 1.9

The PCR & Gel Clean-up Kits can be used downstream of DNA labeling with Enzo's CYTAG[®] CGH labeling Kit (Prod. No. ENZ-42671) and are included with the CYTAG[®] TotalCGH labeling kit (Prod. No. ENZ-42674) to clean up excess nucleotides and enzymes.

A rapid and simple mini kit for purifying DNA from PCR reactions, agarose and polyacrylamide gels, and labeling reactions. Our silica membrane PCR & Clean-up mini spin columns reliably remove primers, enzymes, mineral oil, PCR additives (e.g., salts, betaine, DMSO), detergents (e.g., Tween 20, Triton X-100), dyes, SDS, unbound labels, and free dNTPs, delivering high quality and pure DNA for use in CGH arrays, PCR, ligation, sequencing, etc.

Ordering Information

Order Online »

ENZ-GEN100-0050	50 tests
ENZ-GEN100-0200	200 tests

Manuals, SDS & CofA

View Online »

Key Features

High performance: Recovers up to 95% of DNA fragments down to 50 bp with binding capacity up to 25 μg

Advanced: Binding buffer with pH indicator ensuring optimal binding conditions (pH < 7.0)

Concentrated recovery: elution volumes as small 15 µl

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application NotesFor the purification of PCR products, extraction of DNA

from agarose gels, and removal of nucleotides, primers,

enzymes, mineral oil, detergents, and dyes.

Contents Binding buffer (w/ pH indicator)

Wash Buffer 5x Elution Buffer

PCR & gel clean-up columns (yellow ring)

Collection tubes (2 mL)

Technical Info / Product Notes Application Notes:

Compatibility of CYTAG® CGH Labeling Kit from Enzo Life

Sciences with SurePrint® G3 Human CGH 1x1M

Microarrays from Agilent

Last modified: May 29, 2024

