

# 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<sup>®</sup> CGH labeling Kit (Prod. No. ENZ-42671) and are included with the CYTAG<sup>®</sup> 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-GEN100-0020	20 tests
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 Notes**      For the purification of PCR products, extraction of DNA from agarose gels, and removal of nucleotides, primers, enzymes, mineral oil, detergents, and dyes.

**Assay Time**      <10 min

**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](#)



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)