

## Nucleotide/Dye Removal Kit

### Fast removal of unincorporated dye-nucleotides or dye-terminators

#### DNA Cleanup

	Cat.-No.	Amount
	PP-216XS	10 preparations
	PP-216S	50 preparations
	PP-216L	250 preparations

For *in vitro* use only  
 Quality guaranteed for 12 months  
 Store at room temperature

#### Kit contents

Gel Filtration Spin Columns  
 2 ml Collection Tubes

#### To be provided by you

1.5-2 ml Microcentrifuge Tubes

#### Description

Nucleotide/Dye Removal Kit is designed for fast and easy separation of unincorporated dye-labeled, marker-labeled or unlabeled nucleotides from DNA. The kit contains ready-to-use spin columns preloaded with a gel filtration resin. The simple cleanup procedure is based on gel filtration technology and guarantees highest recovery rates.

The kit has been optimized for cleanup of enzymatic DNA labeling reactions like PCR probe labeling, DNA sequencing, Nick Translation or DNA end-labeling. It allows the removal of labeled dNTPs or ddNTPs in a single step within a few minutes. The purified DNA is applicable in down-stream applications like FISH (fluorescence in situ hybridization) or Southern - / Northern blotting.

#### Specification

Recommended sample volume: 10-20  $\mu$ l  
 Recommended centrifugal force: 1,000 g  
 DNA fragment cut-off: 10-20 bp

#### Preparation procedure

##### 1 Column set-up

- Gently vortex the spin column to resuspend the resin.
- Loosen the cap of the column a quarter turn to avoid vacuum inside the spin column.
- Snap off the bottom closure of the spin column and place the spin column in one of the provided 2 ml collection tubes.
- Centrifuge for 1 min at 1,000 g.
- Discard the flow-through, remove the cap and transfer the spin column to a new 1.5-2 ml microcentrifuge tube.

##### 2 DNA cleanup

- Carefully load the DNA sample to the column.
- Pipet the sample (10-20  $\mu$ l) to the center of the gel surface without touching the gel-bed with the pipet tip.
- The sample must be absorbed from the gel. Avoid pipeting of the sample directly on the tube inside or into the space between gel and tube.
- Centrifuge for 1 min at 1,000 g.
- Remove the spin column from the microcentrifuge tube.
- The eluate contains the purified DNA and can directly be applied in down-stream applications, dried in a vacuum centrifuge or stored at -20°C.