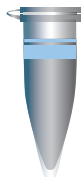


Sample Preparation



1. Add 1-1.5 mL overnight **bacteria culture** to 1.5 mL microcentrifuge tube.
2. Centrifuge at $10\,000 \times g$ for 30 sec
3. Discard the supernatant.
- » Repeat these steps to collect more cells.

Cell Lysis



4. Resuspend completely the pellet of bacterial cells in 250 μL of **Resuspension Buffer**.
5. Add 250 μL **Lysis Buffer** and gently invert the tube 4-6 times to mix.
- » Do not allow this step for more than 5 min
- » Please close the bottle of the Lysis buffer immediately after usage to avoid acidification
- » Do not vortex to avoid shearing of genomic DNA.

Neutralization



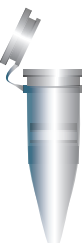
6. Add 350 μL of **Neutralization Buffer** and mix by gently inverting the tube 4-6 times. The solution should become cloudy and no local precipitate should be visible.
7. Centrifuge 10 min at $>14\,000 \times g$ until a compact white pellet form.

Binding



8. Place a SmartPure column into a collection tube and apply the **supernatant** to the SmartPure column
9. Centrifuge for 30-60 sec at $6\,000 \times g$.
10. Discard the flow-through.

Washing



11. Add 650 μL of **Wash Buffer** to the Smart Pure column.
12. Centrifuge for 30-60 sec at $12\,000 \times g$.
13. Discard the flow-through.
14. Repeat washing step once.
15. Centrifuge for an additional 1 min at $12\,000 \times g$ to remove residual liquid

Elution



16. Transfer the SmartPure column to a sterile 1.5 mL micro centrifuge tube.
17. Add 50 μL of **Elution Buffer**, ddH_2O or TE Buffer to the SmartPure column and let it stand for 1 min at room temperature.

Purified DNA

18. Centrifuge for 1 min at $12\,000 \times g$ to recover the purified plasmid DNA in the microcentrifuge tube.
- » The purified plasmid DNA can be used directly or stored at -20°C for long term storage.

The SmartPure Plasmid Kit ensures an easy, fast and effective purification of up to 40 µg Plasmid DNA from 1-5 mL overnight culture of *E. coli*.

First use

- Add all the RNase solution to the SmartPure Resuspension Buffer, mix and store at 2–8°C.
- Add required ethanol to the SmartPure Wash Buffer (see label on the bottle) and mix well.

→ If precipitates appear in the SmartPure Lysis Buffer and the Neutralization Buffer, please incubate the solutions at 37°C for several minutes.

Kit components (100 preps)

| Component | Amount |
|---------------------------------|-----------|
| SmartPure Resuspension Buffer | 25 mL |
| SmartPure Lysis Buffer | 25 mL |
| SmartPure Neutralization Buffer | 35 mL |
| SmartPure Wash Buffer | 2 x 15 mL |
| SmartPure Elution Buffer | 20 mL |
| RNase solution | 1 |
| SmartPure Column | 100 |

Material not supplied

- Sterile 1.5 mL microcentrifuge tubes
- 10/100/1000 µL tips
- Microcentrifuge (14 000 x g)
- Vortex mixer
- Absolute ethanol

Shipping & Storage conditions

The SmartPure Plasmid is shipped at room temperature. The Kit is stable for 18 months if stored in a dry environment at room temperature (15°C-25°C).

The RNase solution must be stored at 2-8°C.

After addition of the RNase solution, the SmartPure Resuspension buffer must be stored at 2-8°C

Troubleshooting

No plasmid recovery

Please check whether the ethanol had been added to the Wash Buffer bottle.

Low plasmid yield

- 1] Please make sure that the cell culture is fresh and cell density is optimized
- 2] Please ensure that the bacteria pellet is completely resuspended in the Resuspension buffer and adequately mixed.
- 3] Incubating Elution Buffer at 30–60°C may increase elution yields.
- 4] Please note that low copy number plasmids and large insert give lower yield.

Electrophoresis problem

- 1] Presence of Genomic DNA: the sample was mixed too vigorously in the Lysis Buffer. Invert the tubes gently 5-6 times when Lysis Buffer and Neutralization Buffer are added.
- 2] Presence of RNA: Add RNase A to the Resuspension buffer (100 µg/mL).

Absorbance problem

Please use the Elution Buffer to adjust zero value and dilute the sample.

→ **If the OD260/OD230 ratio is low:** wash the SmartPure column one more time.

→ **If the OD260-320/OD280-320 ratio is low:** there is a protein contamination. Please add Neutralization Buffer and centrifuge the mixture with a sufficient rotating speed to compact the precipitate. Be careful to pipette the supernatant only.

→ **If the OD260-320/OD280-320 ratio is high:** add RNase A to the Resuspension Buffer (100 µg/mL).

EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM

• Tel.: +32 4 372 74 00

• Fax: +32 4 372 75 00

• Toll-free: + 800 666 00 123 • info@eurogentec.com

• www.eurogentec.com

NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA

• Tel.: +1-510-791-9560

• Toll-free: +1 800-452-5530

• Fax: +1 510-791-9572 • service@anaspec.com

• www.anaspec.com