

# Thermo Scientific GeneJET Plasmid Miniprep Kit

K0502, K0503

**Note.** All steps should be carried out at room temperature. All centrifugations should be carried out in a microcentrifuge at  $\geq 12\ 000 \times g$  (10 000-14 000 rpm, depending on the rotor type).

**Harvest bacteria.** Harvest the bacterial culture by centrifugation at 8000 rpm (6800  $\times g$ ) in a microcentrifuge for 2 minutes at room temperature. Decant the supernatant and remove all remaining medium.

## 1 Ressuspend Cells, Lyse and Neutralize



Add to the pelleted cells:

250  $\mu\text{L}$  of **Resuspension Solution** and vortex.

250  $\mu\text{L}$  of **Lysis Solution** and invert the tube 4-6 times.

350  $\mu\text{L}$  of **Neutralization Solution** and invert the tube 4-6 times.

Centrifuge 5 minutes.



## 2 Bind DNA



Transfer the supernatant to the **Thermo Scientific GeneJET Spin Column**.

Centrifuge 1 minute.



## 3 Wash the column



Add 500  $\mu\text{L}$  of **Wash Solution** and centrifuge for 30-60 s. }  $\times 2$  times

Discard the flow-through.

Centrifuge empty column for 1 minute.



## 4 Elute purified DNA



Transfer the column into a new tube.

Add 50  $\mu\text{L}$  of **Elution Buffer** to the column and incubate 2 minutes.

Centrifuge 2 minutes.

Collect the flow-through.

[thermoscientific.com/onebio](http://thermoscientific.com/onebio)

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

**Technical Support**

**North America:**

TS.molbio@thermofisher.com

**Europe and Asia:**

TS.molbio.eu@thermofisher.com

**Thermo**  
SCIENTIFIC

Part of Thermo Fisher Scientific