



## QIAGEN Supplementary Protocol:

### Purification of total DNA from insects using the DNeasy® Blood & Tissue Kit

This protocol is designed for purification of DNA from up to 50 mg of insects, such as drosophila.

#### Introduction

In this protocol, insects are ground using liquid nitrogen and a mortar and pestle or homogenized using the TissueRuptor, an equivalent electric homogenizer, or a disposable microtube pestle. Samples are then processed according to the standard DNeasy protocols.

**IMPORTANT:** Please read the *DNeasy Blood & Tissue Handbook*, paying careful attention to the "Safety Information" and "Important Notes" sections, before beginning this procedure. DNeasy Blood & Tissue Kits are intended for research use. No claim or representation is intended to provide information for the diagnosis, prevention, or treatment of a disease.

#### Equipment and reagents required

When working with chemicals, always wear a suitable lab coat, disposable gloves, and protective goggles. For more information, consult the appropriate material safety data sheets (MSDSs), available from the product supplier.

- DNeasy Blood & Tissue Kit (cat. no. 69504 or 69506)
- Pipets and pipet tips
- Vortexer
- Microcentrifuge tubes (1.5 ml)
- Microcentrifuge with rotor for 1.5 ml and 2 ml tubes
- Thermomixer, shaking water bath, or rocking platform for heating at 56°C
- Ethanol (96–100%)\*

#### Using a mortar and pestle

- Mortar and pestle
- Liquid nitrogen

\* Do not use denatured alcohol, which contains other substances such as methanol or methylethylketone.

### Using an electric homogenizer or disposable microtube pestle

- TissueRuptor with disposable probes or an equivalent electric homogenizer; or a disposable microtube pestle
- PBS, pH 7.2 (50 mM potassium phosphate, 150 mM NaCl)

### Important points before starting

- If using the DNeasy Blood & Tissue Kit for the first time, read “Important Notes” in the *DNeasy Blood & Tissue Handbook*.
- All centrifugation steps are carried out at room temperature (15–25°C) in a microcentrifuge.
- Vortexing should be performed by pulse-vortexing for 5–10 s.
- Two separate protocols are given. Samples can be ground using a mortar and pestle (see below) or homogenized using the TissueRuptor, an equivalent electric homogenizer, or a microtube pestle (page 4). Do not use a mixture of the two protocols.
- PBS is required for use in the protocol using an electric homogenizer or a microtube pestle. Buffer ATL is not required in this protocol.

### Things to do before starting

- Buffer ATL and Buffer AL may form precipitates upon storage. If necessary, warm to 56°C until the precipitates have fully dissolved.
- Buffer AW1 and Buffer AW2 are supplied as concentrates. Before using for the first time, add the appropriate amount of ethanol (96–100%) as indicated on the bottle to obtain a working solution.
- Preheat a thermomixer, shaking water bath, or rocking platform to 56°C for use in step 3.
- If using frozen insects, equilibrate the sample to room temperature. Avoid repeated thawing and freezing of samples since this will lead to reduced DNA size.

### Procedure — using a mortar and pestle

- 1. Grind up to 50 mg insects in liquid nitrogen using a mortar and pestle. Place the powder in a 1.5 ml microcentrifuge tube.**

To prevent cross-contamination, thoroughly clean the mortar and pestle between samples.

- 2. Add 180 µl Buffer ATL.**

- 3. Add 20 µl proteinase K. Mix thoroughly by vortexing, and incubate at 56°C until the insects are completely lysed. Vortex occasionally during incubation to disperse the sample, or place in a thermomixer, shaking water bath, or on a rocking platform.**

Lysis time varies depending on the type of insect processed. Lysis is usually complete in 1–3 h. If it is more convenient, samples can be lysed overnight; this will not affect them adversely.

After incubation the lysate may appear viscous, but should not be gelatinous as it may clog the DNeasy Mini spin column. If the lysate appears very gelatinous, see the “Troubleshooting Guide”, in the *DNeasy Blood & Tissue Handbook*, for recommendations.

- 4. Vortex for 15 s. Add 200  $\mu$ l Buffer AL to the sample, and mix thoroughly by vortexing. Then add 200  $\mu$ l ethanol (96–100%), and mix again thoroughly by vortexing.**

It is essential that the sample, Buffer AL, and ethanol are mixed immediately and thoroughly by vortexing or pipetting to yield a homogeneous solution. Buffer AL and ethanol can be premixed and added together in one step to save time when processing multiple samples.

A white precipitate may form on addition of Buffer AL and ethanol. This precipitate does not interfere with the DNeasy procedure. Some insect types may form a gelatinous lysate after addition of Buffer AL and ethanol. In this case, vigorously shaking or vortexing the preparation is recommended.

- 5. Pipet the mixture from step 4 (including any precipitate) into the DNeasy Mini spin column placed in a 2 ml collection tube (provided). Centrifuge at  $\geq 6000 \times g$  (8000 rpm) for 1 min. Discard flow-through and collection tube.\***
- 6. Place the DNeasy Mini spin column in a new 2 ml collection tube (provided), add 500  $\mu$ l Buffer AW1, and centrifuge for 1 min at  $\geq 6000 \times g$  (8000 rpm). Discard flow-through and collection tube.\***
- 7. Place the DNeasy Mini spin column in a new 2 ml collection tube (provided), add 500  $\mu$ l Buffer AW2, and centrifuge for 3 min at 20,000  $\times g$  (14,000 rpm) to dry the DNeasy membrane. Discard flow-through and collection tube.**

It is important to dry the membrane of the DNeasy Mini spin column, since residual ethanol may interfere with subsequent reactions. This centrifugation step ensures that no residual ethanol will be carried over during the following elution.

Following the centrifugation step, remove the DNeasy Mini spin column carefully so that the column does not come into contact with the flow-through, since this will result in carryover of ethanol. If carryover of ethanol occurs, empty the collection tube, then reuse it in another centrifugation for 1 min at 20,000  $\times g$  (14,000 rpm).

- 8. Place the DNeasy Mini spin column in a clean 1.5 ml or 2 ml microcentrifuge tube (not provided), and pipet 200  $\mu$ l Buffer AE directly onto the DNeasy membrane. Incubate at room temperature for 1 min, and then centrifuge for 1 min at  $\geq 6000 \times g$  (8000 rpm) to elute.**

Elution with 100  $\mu$ l (instead of 200  $\mu$ l) increases the final DNA concentration in the eluate, but also decreases the overall DNA yield (see *DNeasy Blood & Tissue Handbook*).

- 9. Recommended: For maximum DNA yield, repeat elution once as described in step 8.**

This step leads to increased overall DNA yield.

A new microcentrifuge tube can be used for the second elution step to prevent dilution of the first eluate. Alternatively, to combine the eluates, the microcentrifuge tube from step 8 can be reused for the second elution step.

**Note:** Do not elute more than 200  $\mu$ l into a 1.5 ml microcentrifuge tube because the DNeasy Mini spin column will come into contact with the eluate.

\* Flow-through contains Buffer AL or Buffer AW1 and is therefore not compatible with bleach. See *DNeasy Blood & Tissue Handbook* for safety information.

## Procedure — using an electric homogenizer or disposable microtube pestle

1. Place up to 50 mg insects in a 1.5 ml microcentrifuge tube.
2. Add 180  $\mu$ l PBS and homogenize the sample using the TissueRuptor, an equivalent electric homogenizer, or a disposable microtube pestle.

The TissueRuptor uses disposable probes to help prevent cross-contamination. We recommend using disposable probes or disposable microtube pestles only once. With other electric homogenizers, the probe should be thoroughly cleaned between samples.

3. Add 20  $\mu$ l proteinase K and 200  $\mu$ l Buffer AL (without added ethanol). Mix thoroughly by vortexing, and incubate at 56°C for 10 min.

Ensure that ethanol has not been added to Buffer AL (see “Buffer AL” in the *DNeasy Blood & Tissue Handbook*). Buffer AL can be purchased separately.

It is essential that the sample and Buffer AL are mixed immediately and thoroughly by vortexing or pipetting to yield a homogeneous solution.

4. Add 200  $\mu$ l ethanol (96–100%) to the sample, and mix thoroughly by vortexing.

It is important that the sample and the ethanol are mixed thoroughly to yield a homogeneous solution.

5. Pipet the mixture from step 4 (including any precipitate) into the DNeasy Mini spin column placed in a 2 ml collection tube (provided). Centrifuge at  $\geq 6000 \times g$  (8000 rpm) for 1 min. Discard flow-through and collection tube.\*

6. Place the DNeasy Mini spin column in a new 2 ml collection tube (provided), add 500  $\mu$ l Buffer AW1, and centrifuge for 1 min at  $\geq 6000 \times g$  (8000 rpm). Discard flow-through and collection tube.\*

7. Place the DNeasy Mini spin column in a new 2 ml collection tube (provided), add 500  $\mu$ l Buffer AW2, and centrifuge for 3 min at 20,000  $\times g$  (14,000 rpm) to dry the DNeasy membrane. Discard flow-through and collection tube.

It is important to dry the membrane of the DNeasy Mini spin column, since residual ethanol may interfere with subsequent reactions. This centrifugation step ensures that no residual ethanol will be carried over during the following elution.

Following the centrifugation step, remove the DNeasy Mini spin column carefully so that the column does not come into contact with the flow-through, since this will result in carryover of ethanol. If carryover of ethanol occurs, empty the collection tube, then reuse it in another centrifugation for 1 min at 20,000  $\times g$  (14,000 rpm).

8. Place the DNeasy Mini spin column in a clean 1.5 ml or 2 ml microcentrifuge tube (not provided), and pipet 200  $\mu$ l Buffer AE directly onto the DNeasy membrane. Incubate at room temperature for 1 min, and then centrifuge for 1 min at  $\geq 6000 \times g$  (8000 rpm) to elute.

Elution with 100  $\mu$ l (instead of 200  $\mu$ l) increases the final DNA concentration in the eluate, but also decreases the overall DNA yield (see *DNeasy Blood & Tissue Handbook*).

\* Flow-through contains Buffer AL or Buffer AW1 and is therefore not compatible with bleach. See *DNeasy Blood & Tissue Handbook* for safety information.

**9. Recommended: For maximum DNA yield, repeat elution once as described in step 8.**

This step leads to increased overall DNA yield.

A new microcentrifuge tube can be used for the second elution step to prevent dilution of the first eluate. Alternatively, to combine the eluates, the microcentrifuge tube from step 8 can be reused for the second elution step.

**Note:** Do not elute more than 200  $\mu$ l into a 1.5 ml microcentrifuge tube because the DNeasy Mini spin column will come into contact with the eluate.

## Troubleshooting

For general troubleshooting, please consult the Troubleshooting Guide in the *DNeasy Blood & Tissue Handbook*.

QIAGEN kit handbooks can be requested from QIAGEN Technical Services or your local QIAGEN distributor.

Selected kit handbooks can be downloaded from [www.qiagen.com/literature/handbooks/default.aspx](http://www.qiagen.com/literature/handbooks/default.aspx).

Material safety data sheets (MSDS) for any QIAGEN product can be downloaded from [www.qiagen.com/ts/msds.asp](http://www.qiagen.com/ts/msds.asp).

Trademarks: QIAGEN®, DNeasy® (QIAGEN Group); DU® (Beckman Instruments, Inc.).

© 2006 QIAGEN, all rights reserved.