

Important Note

The kit contains alkaline reagents.

Please handle it with care: wear gloves and goggles...

Sample Preparation



1. Transfer the sample in a microtube.

Sample	Amount
Mouse tail	5 mm (~10 mg)
Blood	1-3 μ L
Seed	1-2
Leaf	Piece of 2-3 mm

Lysis



2. Add 100 μ L of Reagent A.

3. Incubate for 10 min at 98°C.

Recovery



4. Add 14 μ L of Reagent B.

→ The recovered genomic DNA can be used directly in (q)PCR or stored at -20°C for long term storage.

PCR Setup

For a subsequent PCR reaction please use:

1 μ L of extracted solution from blood, leaves or seeds

1-5 μ L of extracted solution from mouse tail

The SmartExtract DNA Kit is optimized to perform in a single tube an easy, fast and effective extraction of (q)PCR-ready genomic DNA from almost any crude sample type (e.g. blood, human & animal tissues, plant seeds & leaves...)

Important Note

The kit contains alkaline reagents.

Please handle it with care: wear gloves and goggles...

Kit components (100 preps)

Component	Amount
Reagent A (Blue Cap)	5 x 2 mL
Reagent B (Yellow Cap)	2 x 700 μ L

Material not supplied

- Sterile microtubes
- 10/100/1000 μ L tips
- Microcentrifuge
- Thermocycler

Shipping & Storage conditions

The SmartExtract DNA Kit is shipped at room temperature. The Kit is stable for 24 months if stored protected from light in a dry environment at room temperature (15°C-25°C).

Troubleshooting

No or low yield

Please adjust the sample amount.

In case of poor PCR results, dilution of the extracted sample can improve subsequent amplifications.

EUROPE

KANEKA EUROGENTEC S.A.
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