

FitAmp[™] Gel DNA Isolation Kit

Base Catalog # P-1007

PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The FitAmpTM Gel DNA Isolation Kit allows isolating DNA fragment from low melting point agarose gel. DNA size can be from 200 bp to 1 kb; DNA quantity from 1 ng to 1 μ g, optimal at between 10 ng and 500 ng. Recovery rate of DNA is greater than 70%.

Cleaned DNA fragment is ready for direct use in the following applications: ligation/transformation, sequencing, restriction digestion, microarray analysis, labeling, microinjection, and in vitro transcription.

110 Bi County Blvd. Ste. 122, Farmingdale, NY 11735 Tel: 1-877-374-4368 ■ Fax: 1-718-484-3956 ■ E-mail: info@epigentek.com ■ Web: www.epigentek.com © Epigentek Group Inc. All rights reserved. Products are for research use only.

Page 1 Printed 2014-05-29 P-1007



KIT CONTENTS

Components	50 samples P-1007-1	100 samples P-1007-2
K1 (Lysis & Binding Buffer) K2 (Wash Buffer) K3 (DNA Elution Solution)	21 ml 2.5 ml 1 ml	2 x 21 ml 2 x 2.5 ml 2 ml
F-Spin Column	50	100
F-Spin Collection Tube	50	100
User Guide	1	1

SHIPPING & STORAGE

Upon receipt, store **K1** at 2-8°C away from light. Store **all other components** at room temperature (15-25°C). The kit is stable for up to 1 year from the date of shipment when stored properly.

GENERAL PRODUCT INFORMATION

Usage Limitation: The *FitAmp*[™] kit is for research use only and is not intended for diagnostic or therapeutic application.

Quality Control: Epigentek guarantees the performance of all products in the manner described in our product instructions.

Product Updates: Epigentek reserves the right to change or modify any product to enhance its performance and design.

Intellectual Property: *FitAmp*[™] is a trademark of Epigentek Group, Inc.

A BRIEF OVERVIEW

The *FitAmp*[™] Gel DNA Isolation Kit uses a unique procedure and composition to efficiently isolate DNA from low melting point agarose gel. The kit has the following features:

- Fast procedure (< 15 minutes).
- Recovery of ultra-pure DNA from agarose gel at high concentration in minimal volume.
- Use of proprietary binding solution without chaotropic or other irritant reagents.



PRINCIPLE & PROCEDURE

The FitAmpTM Gel DNA Isolation Kit applies our proprietary lysis and binding buffer to gel-DNA sample. After transferring the sample to the specifically designed Fast-Spin Column and wash with washing buffer, the DNA is easily recovered in 8-20 μ l with elution solution. Isolated DNA is ready for down-stream application.



Schematic Procedure for Using the *FitAmp™* Gel DNA Isolation Kit

PROTOCOL

Note: Always cap spin columns before placing them in the microcentrifuge.

Before starting, add 20 ml of 100% ethanol to each bottle of K2 to make the final wash buffer.

- 1. Cut the agarose gel slice that contains DNA fragment and weigh the gel slice. In general, it should not be over 150 mg.
- Transfer the gel slice to a 1.5 ml microcentrifuge tube. Add 3 volumes of K1 to each volume of gel slice and break it down into several pieces using a pipette tip. Incubate at 60-65°C for 10-12 minutes (during this period, briefly vortex 2-3 times or until the gel slice is dissolved at regular intervals).
- 3. Transfer the dissolved gel solution to the column. Centrifuge at 12,000 rpm for 45-60 seconds. Discard the flowthrough. Replace the column to the collection tube (Note: maximum volume of the column is 600μ L)
- 4. Add 200 μ l of **K2** (final wash buffer) to the column and centrifuge at 12,000 rpm for 30 seconds. Discard the flowthrough and replace the column to the collection tube. Add 200 μ l of **K2** (final wash buffer) to the column and centrifuge at 12,000 rpm for 45 seconds.

110 Bi County Blvd. Ste. <mark>1</mark>22, Farmingdale, NY 11735 Tel: 1-877-374-4368 ■ Fax: 1-718-484-3956 ■ E-mail: info@epigentek.com ■ Web: www.epigentek.com © Epigentek Group Inc. All rights reserved. Products are for research use only.



5. Place the column in a new 1.5 ml vial. Add 8-18 μ l of **K3** directly to the column filter, centrifuge at 12,000 rpm for 30 seconds to elute cleaned DNA.

DNA is now ready for use or storage at -20° C.

RELATED PRODUCTS

- P-1003 FitAmp[™] General Tissue Section DNA Isolation Kit
- P-1004 FitAmp[™] Plasma/Serum DNA Isolation Kit

110 Bi County Blvd. Ste. 122, Farmingdale, NY 11735 Tel: 1-877-374-4368 ■ Fax: 1-718-484-3956 ■ E-mail: info@epigentek.com ■ Web: www.epigentek.com © Epigentek Group Inc. All rights reserved. Products are for research use only.