



Plasmid Screening *Toothpick*[™]

INTRODUCTION

Plasmid Screening *Toothpick*[™] allows rapid analysis of plasmid DNA by restriction enzyme(s) from a bacterial colony itself, without growing an overnight culture. Simply pick a bacterial colony and screen with Plasmid Screening *Toothpick*[™] to see if you have the right construct. The solution is enough for 300 Preps.

KIT COMPONENTS Cat# 786-026

Plasmid Screening <i>Toothpick</i> [™]	4.5ml
-------------------------------------------------	-------

STORAGE CONDITIONS

It is shipped at ambient temp. Upon arrival, store at 4⁰C and is stable for one year, if stored and used properly.

PROTOCOL

1. Pick a freshly plated bacterial colony (1-2mm in diameter) with a sterile toothpick or pipette tip and suspend it in 5µl fresh growth media (e.g. LB medium). Mix it thoroughly with gentle hand.

Note: *If the colony size is smaller, incubate the tube at room temperature for 60-90 minutes.*

OPTIONAL - *Remove 1µl culture from the tube and suspend in a 10µl growth medium containing an appropriate selection of antibiotic. Incubate it for scaling up the culture. Alternatively, freeze the culture for later use.*

2. Add 15µl of the Plasmid Screening *Toothpick*[™] solution to the bacterial suspension and mix by pipetting up and down 4-5 times. Incubate at room temperature for 1-2 minutes and mix again.

3. Heat the sample at 100⁰C in a boiling water bath for 45-50 seconds.

4. Cool the sample to room temperature.

5. Transfer 10µl sample to a tube containing 2µl of restriction digestion mix (i.e. 1µl of restriction enzyme; 1-10 units and 1µl of 10X restriction buffer).

6. Incubate the sample for 30-60 minutes at the appropriate temperature for the restriction enzyme used.

7. After incubation, add 2µl of loading buffer to the sample. Mix and load into the wells of an agarose gel. Once electrophoresis is complete, stain the gel in 0.2mg/ml ethidium bromide solution for 10 minutes and visualize the DNA under UV. **Note:** *Ethidium bromide may be added in the agarose gel, before run.*

RELATED PRODUCTS

1. *GET*[™] Plasmid Maxiprep (Cat# 786-363 to 786-364):

G-Biosciences has developed pinkRESIN[™], a superior binding matrix for purification of plasmid DNA. It has an enhanced binding affinity for plasmids and resuspend with ease, thus eliminates the loss of plasmid.

2. *Dot-Blot Enhancer*[™] System (Cat# 786-163): *The device is useful for Nucleic acid and Protein dot-blot procedures. It concentrates the sample in a small area at the point of application and enhances the signal strength.*

Note: For other related products, visit our web site at www.GBiosciences.com or contact us.

Rev 11.07.08-SA/MM/IA

