



379PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Plasmid Screening Toothpick™

(Cat. # 786-026)



think proteins! think G-Biosciences www.GBiosciences.com

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.

ITEM(S) SUPPLIED (Cat. # 786-026)

Description	Size
Plasmid Screening Toothpick™	4.5ml

STORAGE CONDITIONS

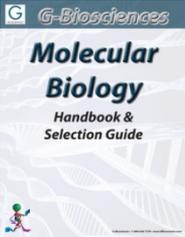
It is shipped at ambient temp. Upon arrival, store at 4°C and is stable for one year, if stored and used properly.

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 100°C 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 with LabSafe™ Nucleic Acid Stain (Cat. # 786-409) or 0.2mg/ml ethidium bromide solution for 10 minutes and visualize the DNA under UV.
NOTE: *LabSafe™ Nucleic Acid Stain and ethidium bromide may be added in the agarose gel, before electrophoresis.*

RELATED PRODUCTS

Download our Molecular Biology Handbook.



<http://info.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 10/23/2012 CMH



www.GBiosciences.com