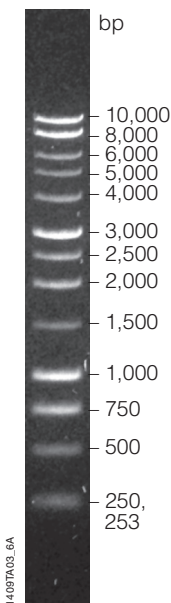


Quality and Stability of DNA Ladders Shipped at Ambient Temperature

Promega Corporation
2800 Woods Hollow Rd., Madison, WI 53711



BenchTop 1kb DNA Ladder (Cat.# G7541)

By removing gel or dry ice for shipping DNA markers, we will eliminate the need for 17,000 kilograms of dry ice a year.

Introduction

We are committed to preserving and improving the natural environment by focusing on a long-term view of what is needed to thrive for the next 100 years and beyond. One of the initial steps in that process is the reducing the amount of dry ice or cold gel used for shipping reagents.

Due to demonstrated stability of DNA, we decided to evaluate the need for any added refrigerant during shipping of several DNA markers. By removing gel or dry ice for these shipments, we will eliminate the need for 17,000 kilograms of dry ice a year.

In this paper, we describe how each DNA ladder (Table 1) was evaluated by an outside testing lab (Gaynes Lab Incorporated) and Promega under simulated shipping conditions. To replicate temperatures experienced during shipping samples, we followed recommendations outlined in ISTA 7D (Figure 1). All testing was conducted in standard Promega corrugated fiberboard boxes. Functional gel analysis was accomplished using inventory material stored at recommended conditions. Fill volume evaporation tests were also carried out and compared to inventory controls.

Materials and Methods

Samples

Table 1. DNA Markers Evaluated at Gaynes Lab Inc.

| Cat.# | Product |
|-------|-----------------------------------------|
| G2101 | 100bp DNA Ladder |
| G5711 | 1kb DNA Ladder |
| G7511 | BenchTop Φ X174 DNA/HaeIII Markers |
| G7521 | BenchTop pGEM [®] DNA Markers |
| G7531 | BenchTop PCR Markers |
| G7541 | BenchTop 1kb DNA Ladder |
| G8291 | BenchTop 100bp DNA Ladder |

ISTA 7D Testing Method

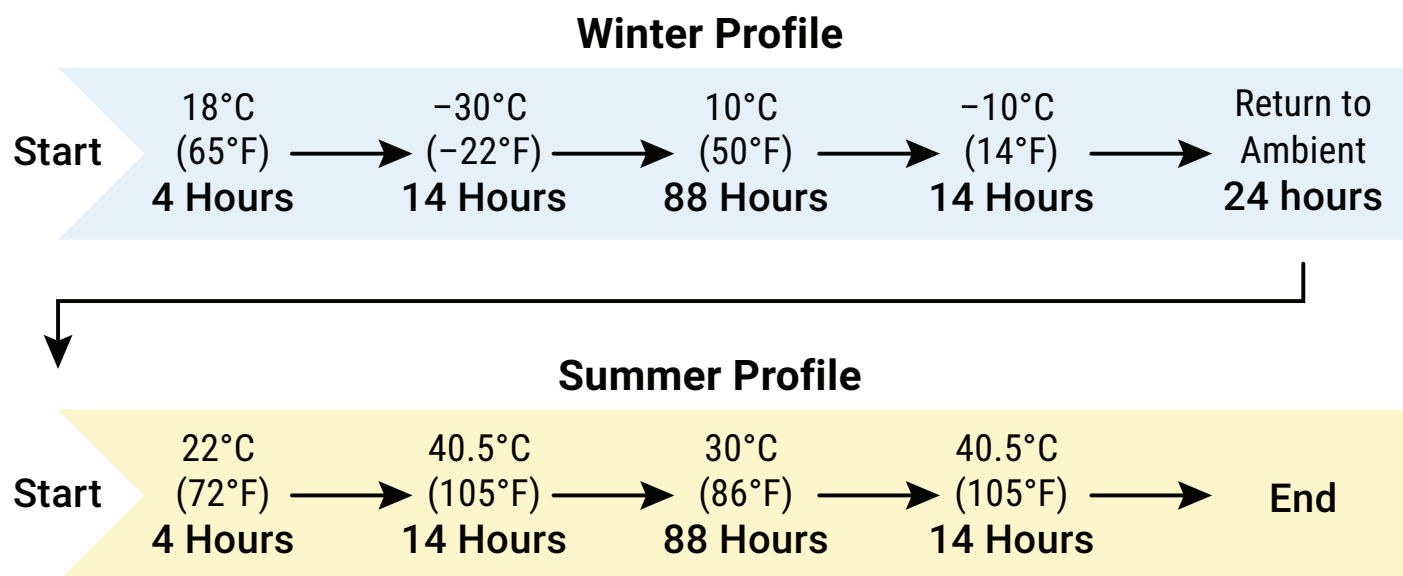


Figure 1. Temperature and time cycles used at the Gaynes Lab for ambient testing of DNA markers.

Additional Internal Testing

In addition to the markers noted in Table 1, testing was conducted on several additional DNA markers by Promega. This testing consisted of samples being exposed to 37°C for 120 hours. Fill volume evaporation and gel analysis were conducted and compared to controls.

Results and Discussion

In both studies, evaporation analysis results showed no significant loss of material. Sample gel results were consistent with inventory samples stored at recommended storage conditions (Figure 2).

For additional information regarding individual DNA markers, please contact Promega Technical Services.

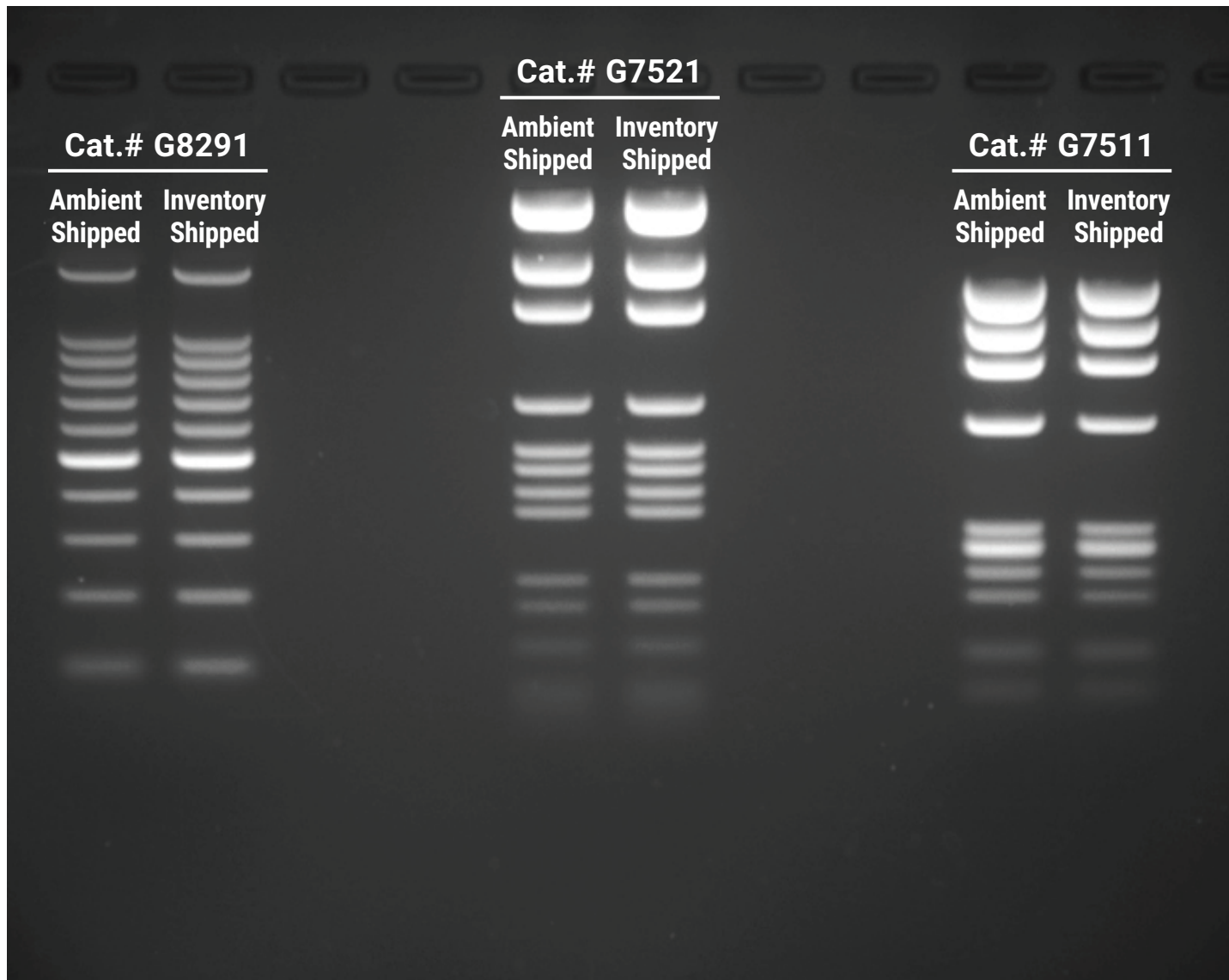


Figure 2. Comparison of samples loaded on a 2% agarose gel at 80 Volts for 3 hours. Lanes from left to right: BenchTop 100bp DNA Ladder (Cat.# G8291), BenchTop pGEM[®] DNA Markers (Cat.# G7521) and BenchTop Φ X174 DNA/HaeIII Markers (Cat.# G7511).

pGEM is a registered trademark of Promega Corporation.
 Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

