

## Product Information

### GeneSnare™ Arbitrary Primer, Set 1

Catalog Number **GS 0040**  
Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

The GeneSnare™ Arbitrary Primer, Set 1 contains the first set of Annealing Control Primer (ACP)™ arbitrary primers, A1 to A24, from a total of 72 arbitrary primers that Sigma offers for the GeneSnare differential expression system. The GeneSnare system is a simple and efficient mRNA differential display method using the proprietary ACP technology. Sigma offers the GeneSnare Differential Expression Kit (Product Code GS 0100) as a complete system for performing differential display. GeneSnare Arbitrary Primer, Set 1 is a critical component of the kit. All arbitrary primers provided in this set are validated in *Arabidopsis*, corn, soybean, CHO, and HeLa cell cultures, and should be suitable for any organism.

### Components

The arbitrary primers are provided in a lyophilized format in a 96 well plate. 24 primers are arrayed in the first three columns of the plate (see Figure 1).

**Figure 1.** Primer array

	1	2	3	4	5	6	7	8	9	10	11	12
A	A1	A9	A17									
B	A2	A10	A18									
C	A3	A11	A19									
D	A4	A12	A20									
E	A5	A13	A21									
F	A6	A14	A22									
G	A7	A15	A23									
H	A8	A16	A24									

This product provides only the arbitrary primers needed for a PCR reaction. To perform a complete differential display, ACP anchor primer set (Product Code GS 0020) and RT-PCR kit (Product Code GS 0030) are required. To increase genome coverage for gene discovery, more than one set of arbitrary primers may be used for differential display.

### Precautions and Disclaimer

This product is for R&D use only. Not for drug, household or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Storage/Stability

The primer set should be stored at  $-20\text{ }^{\circ}\text{C}$ . Primers are stable at  $-20\text{ }^{\circ}\text{C}$  for at least 1 year and can be frozen and thawed at least 10 times.

### Preparation of primers

The arbitrary primer (0.5 nmole for each primer) in each well is lyophilized. Before use, dissolve the primers in 50  $\mu\text{l}$  of TE buffer (10 mM Tris, pH 8.0, with 0.001 mM EDTA). The dissolved primer set may be divided between 2 or 3 plates. Use only one plate repeatedly until depleted, before switching to a new plate to avoid cross-contamination.

### Related Products

GeneSnare Differential Expression Kit,  
Product Code GS 0010  
GeneSnare Arbitrary Primer, Set 2,  
Product Code GS 0050  
GeneSnare Arbitrary Primer, Set 3,  
Product Code GS 0060  
GeneSnare RedAccuTaq™ RT-PCR Kit,  
Product Code GS 0030  
GeneSnare Differential Expression Primer Kit,  
Product Code GS 0020

Under License from Seegene, Inc.

Annealing Control Primer and ACP are trademarks of Seegene, Inc.  
GeneSnare is a trademark and RedAccuTaq is a registered trademark of Sigma-Aldrich Biotechnology LP

AH,PHC 09/10-1