



3050 Spruce Street
Saint Louis, Missouri 63103 USA
Telephone 800-325-5832 • (314) 771-5765
Fax (314) 286-7828
email: techserv@sial.com
sigma-aldrich.com

Product Information

ArrayHyb™ Hybridization Buffer

Product No. **A7718**
Technical Bulletin No. MB-745
May 2000
Store at room temperature

TECHNICAL BULLETIN

Product Description

ArrayHyb™ Hybridization Buffer is optimized for use with microarrays spotted on glass slides. Hybridization of fluorescently labeled nucleic acids using ArrayHyb Hybridization Buffer consistently provides superior signal with minimal background when compared to standard and other commercially available hybridization buffers. For instances in which lower hybridization temperatures (42-50°C) are desirable, Sigma recommends formamide-containing ArrayHyb™ LowTemp Hybridization Buffer (Product No. A 3095).

The ArrayHyb formulation increases the rate of hybridization and can provide results for analytical purposes in as little as 2 hours. Approximately 80% of the maximum signal is achieved within 6 hours, however, the signal will continue to increase over time with no increase in background for greater sensitivity. This flexibility allows hybridizations to be tailored to meet specific requirements.

Precautions and Disclaimer

ArrayHyb Hybridization Buffer is for laboratory use only. Not for drug, household or other uses.

Storage/Stability

Store ArrayHyb Hybridization Buffer at room temperature. ArrayHyb has a shelf life of 1 year upon receipt. If a small amount of precipitate forms, warm the solution to >30°C until dissolved, invert to mix and store at room temperature.

Reagents and Equipment Required But Not Provided
(Sigma product numbers have been given where appropriate)

<u>Product Name</u>	<u>Product No.</u>
Humidity chamber	H 6644
Wheaton 20 slide staining dish	S 6141
22 x 22 mm Coverslips	C 9802
10% Lauryl sulfate (SDS)	L 4522
95% Ethanol	E 7148
Single strand DNA for hybridization	D 7656
Poly (dA) _n	P 0887
Yeast tRNA	R 5636
20X Saline-sodium citrate (SSC)	S 6639
Orbital shaker	

Preparation Instructions

- 1% SDS for pre-treatment of slides prior to hybridization. Prepare a 1% SDS solution by diluting 100 ml of 10% SDS (Product No. L 4522) with 900 ml of distilled deionized water.
- Wash Buffer 1 (1X SSC, 0.03% SDS). Prepare Wash Buffer 1 by combining 50 ml of 20X SSC (Product No. S 6639) with 3 ml of 10% SDS. Dilute to 1 L with distilled deionized water.
- Wash Buffer 2 (0.2X SSC). Prepare Wash Buffer 2 by diluting 10 ml of 20X SSC to 1 L with distilled deionized water.
- Wash Buffer 3 (0.05X SSC). Prepare Wash Buffer 3 by diluting 2.5 ml of 20X SSC to 1 L with distilled deionized water.

Procedure

A. Pre-treatment of Slides for Hybridization

Slides must be pre-treated after spotting of targets and prior to use in hybridization. The pre-treatment procedure outlined below was developed for SigmaScreen™ Coated Slides for Microarrays, which contain aminopropyltriethoxysilane-derived amine functional groups covalently attached to the slide surface. If a different type of slide has been utilized as a printing substrate, use a pre-treatment protocol optimized for that substrate. If the printed slides have already been pre-treated by similar procedures, proceed directly to Section B, Hybridization. This protocol has been optimized for hybridizations using ArrayHyb Hybridization Buffer. This step denatures the spotted, double stranded DNA to make it available for hybridization and washes away any excess spotted nucleic acid. If slides are pre-treated using other procedures, overall signal and consistency may be compromised.

Multiple slides may be pre-treated at the same time using this procedure. Do not write on the slide at this stage, as many writing sources are soluble in ethanol and can contribute to non-specific fluorescent signal. Slides that are not used immediately for hybridizations should be stored desiccated at 2-8 °C.

1. Begin heating distilled deionized water to 95-100 °C in a clean container filled with sufficient water to cover slides in a slide staining rack.
2. Fill a humidity chamber with ~50 ml water pre-warmed to approx. 65 °C.
3. Invert arrayed slides (DNA side down) into the humidity chamber and allow spots to re-hydrate for approx. 5 seconds.
4. Snap-dry each arrayed slide (DNA side up) on a 95-100°C inverted heat block for approx. 5 seconds.
5. Place slides (DNA side up) in a second empty humidity chamber.
6. UV crosslink the DNA to the surface with 65 mJ of 254 nm UV light. Place the slides in a slide staining rack.

7. Incubate the slides for approx. 2 minutes in a slide staining dish filled with 1% SDS on an orbital shaker. This step removes the unbound nucleic acids from the arrays and helps block non-specific binding of nucleic acids.
8. Wash the slides by dipping the staining rack several times in a dish of distilled deionized water. Fresh water should be used for each set of slides. This step removes the majority of SDS from step 7.
9. Gently plunge the slide rack into the 95-100 °C distilled deionized water for 2 minutes.
10. Remove the slide rack from the water bath and rinse the slides by plunging the rack 10-20 times in 95+% ethanol.
11. Quickly transfer the slides to a centrifuge with a swinging bucket rotor for microtiter plates (place paper towels below rack to absorb liquid) and spin at 50-100 X g for 5-10 minutes.
12. Use the treated slides immediately or store in a slide box desiccated at 2-8 °C.

B. Hybridization

1. Equilibrate pre-treated slides to room temperature and label each clearly with indelible ink.
2. Create a probe mixture by combining labeled nucleic acid and blocking agent(s), if desired (see below), in a microcentrifuge tube. Ethanol precipitate or speed-vac the mixture to dryness. Resuspend the probe mixture pellet into the appropriate volume of ArrayHyb Hybridization Buffer for the size cover slips being utilized. The optimal volume has been found to be 2.5-3 µl/cm² for ArrayHyb (12.5-15 µl for 22 x 22 mm cover slips).

<u>Common Blocking Agents</u>	<u>Final Concentration</u>
Single stranded DNA	100 µg/ml
Poly (dA)	400 µg/ml
Yeast tRNA	200 µg/ml
CoT1 DNA (LTI)	400 µg/ml

3. Heat the ArrayHyb/probe mixture at 95-100 °C for 5 minutes.
4. Centrifuge the contents to the bottom of the microcentrifuge tube and store at 70 °C until needed.
5. Carefully pipette the hybridization solution onto a coverslip.
6. Slowly lower the slide (array side down) until surface tension allows the cover slip to be raised with the slide, taking care not to introduce bubbles.
7. Incubate the slides from 6 hours to overnight at 65°C in a humidity-controlled environment. This can be achieved by placing slides in a slide moat incubator; alternatively, hybridization may be carried out in one of several commercially available hybridization chambers immersed in a temperature controlled water bath.

C. Washing

Do not allow the slides to dry out at any point during this procedure. This can cause non-specifically bound probe to become permanently attached to the slide, resulting in high backgrounds.

1. Remove the slides one at a time from the water bath, immerse the slides into a clean container filled with Wash Buffer 1 (pre-warmed to 50-60 °C), and gently remove the cover slips. Place the slides into a slide rack/staining dish filled with Wash Buffer 1 and incubate for 5 minutes at room temperature with gentle mixing on an orbital shaker.
2. During this incubation, place 200 ml of Wash Buffer 2 (0.2X SSC) in a clean staining dish containing a clean slide rack.
3. Quickly transfer the slides, one at a time, to the rack in the Wash Buffer 2. Shake off excess Wash Buffer 1 from each slide as it is transferred to the rack in Wash Buffer 2.
4. Incubate the slides in Wash Buffer 2 for 5 minutes at room temperature on an orbital shaker.
5. During this incubation, place 200 ml of Wash Buffer 3 (0.05X SSC) in a clean staining dish (no slide rack required).
6. Transfer the slide rack containing the slides to the staining dish containing Wash Buffer 3 and incubate for 5 minutes at room temperature on an orbital shaker.
7. After the third washing, quickly transfer the slides to a dry slide rack and place in a centrifuge equipped with a swinging bucket rotor for microtiter plates. Centrifuge at 50-100 X g for 5-10 minutes. Immediately remove the slides from the centrifuge and store in a light-proof slide box.

Store the slides protected from light and dust. Scan as soon as possible.

Troubleshooting Guide

Problem	Cause	Solution
Low Signal	Inefficient binding of nucleic acids to the slide during printing	DNA must be cleaned properly prior to spotting. Silica matrix columns, such as Sigma's GenElute PCR DNA purification kit, are recommended for purification of the DNA.
	Hybridization temperature too high	Decrease hybridization temperature (typically by 5-10 °C) to achieve optimal hybridization signal.
	Probe was not labeled efficiently	Check the probe for labeling efficiency. If poor label incorporation is observed, remake the probe. Protect labeled probes from exposure to light.
Background Fluorescence	Drying of hybridization solution at the edges of the cover slip	Be sure to hybridize the slides in a humidified chamber. The use of humidity chambers will prevent this problem.
	Probe was allowed to dry to the slide during transfer to Wash Buffer 1	Use extra care to prevent any drying of probe solution on the slides.
	Dust has accumulated on the slide	Dust particles will show up as isolated spots of very high fluorescence. Protect slides from general lab air as much as possible. Some dust may be removed by the use of compressed air to "dust" the slide prior to scanning.
Non-specific Hybridization Signal	Hybridization temperature too low	Increase hybridization temperature (typically by 5-10 °C) to eliminate non-specific hybridization signal. Alternatively, use ArrayHyb LowTemp Hybridization Buffer.
Spots appear as streaks or comets	Recommended pre-treatment procedure was not followed	When concentrated DNA (0.5 mg/ml) is spotted on the slide, only a small percentage becomes bound to the surface. The remaining unbound DNA must be washed away. This is accomplished in the recommended pre-treatment procedure by washing the slides in 1% SDS. This step removes unbound DNA from the spots and prevents the DNA from binding to the surrounding slide surfaces.

Related Products

<u>Product Name</u>	<u>Product No.</u>
ArrayHyb™ LowTemp Hybridization Buffer	A 3095
SigmaScreen™ Coated Slides for Microarrays	S 7934
Aminoallyl cDNA Labeling Kit	AM-ALLYL
GenElute™ Mammalian Total RNA Kits	RTN-10, RTN-70 and RTN-350
GenElute™ mRNA from Total RNA Kits	MRN-10 and MRN-70
GenElute™ PCR Purification Kit	GEN-PCR
SigmaSpin™ Post-reaction Clean-up Columns	S 5059
0.1% Poly-L-Lysine	P 8920
3-Aminopropyltriethoxysilane	A 3648
SlideMoat Hybridization Incubator	Z38,067-9
BioLink BLX UV Crosslinker	Z37,537-3
Belly-Dancer Orbital Shaker	Z36,760-5

References

1. Schena, M., *et al.* Parallel human genomic analysis: microarray-based expression monitoring of 1000 genes. *Proc. Natl. Acad. Sci. USA*, **93**, 10614-10619 (1996)
2. Schena, M., *et al.* Quantitative monitoring of gene expression patterns with a complementary DNA microarray. *Science*, **270**, 467-470 (1995)

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications.

Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.

Please see reverse side of the invoice or packing slip.