

DNA BUFFERS

0.1 M MES, pH 4.5 (COUPLING BUFFER) 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
MES (2[N-Morpholino] ethanesulfonic acid)	Sigma M2933	0.1 M	4.88 g
5 N NaOH	Fisher SS256-500	-----	~ 5 drops

Filter Sterilize and store at 4°C

0.02% TWEEN-20 (WASH BUFFER I) 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
TWEEN® 20 (Polyoxyethylenesorbitan monolaurate)	Sigma P9416	0.02%	50 µL

Filter Sterilize and store at Room Temperature

0.1% SDS (WASH BUFFER II) 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
SDS (Sodium lauryl sulfate) 10% solution	Sigma L4522	0.1%	2.5 mL

Filter Sterilize and store at Room Temperature

20% Sarkosyl 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
Sarkosyl (N-Lauroylsarcosine)	Sigma L9150	20%	50g

Filter Sterilize and store at Room Temperature

TE, pH 8.0 (SAMPLE DILUENT) 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
Tris-EDTA Buffer, pH 8.0, 100X	Sigma T9285	1 X	2.5 mL

Filter Sterilize and store at Room Temperature

1.5 X TMAC Hybridization Solution (MICROSPHERE DILUENT) 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
5 M TMAC	Sigma T3411	4.5 M	225 mL
20% Sarkosyl	-----	0.15%	1.88 mL
1 M Tris-HCl, pH 8.0	Sigma T3038	75mM	18.75 mL
0.5 M EDTA, pH 8.0	Invitrogen 15575-020	6mM	3.0 mL
H ₂ O	-----	-----	1.37 mL

Store at Room Temperature

1 X TMAC Hybridization Solution (DETECTION BUFFER) 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
5 M TMAC	Sigma T3411	3 M	150 mL
20% Sarkosyl	-----	0.1%	1.25 mL
1 M Tris-HCl, pH 8.0	Sigma T3038	50mM	12.5 mL
0.5 M EDTA pH 8.0	Invitrogen 15575-020	4mM	2 mL
H ₂ O	-----	-----	84.25 mL

Store at Room Temperature

6X SSPET (STRINGENT WASH BUFFER) 250 mL

Reagent	Catalog Number	Final Concentration	Amount/ 250 mL
SSPE, 20X (Phosphate buffer, pH 7.4, sodium chloride, EDTA)	Sigma S2015	6X concentrate	75 mL
Triton [®] X-100	Sigma T9284	0.005%-0.01%	12.5-25 µL

Filter Sterilize and store at Room Temperature