

General Recommendations for DNA Electrophoresis

This protocol is for the General Recommendations for DNA Electrophoresis

- Use the same DNA loading dye (supplied with the DNA ladder/marker) for both the sample DNA and the ladder/marker DNA.
- If possible, always load equal volumes of the sample DNA and the ladder/marker DNA. The sample can be diluted with 1X DNA loading dye.
- Avoid high salt concentrations in the DNA samples as this may cause bands to shift during electrophoresis.
- Following electrophoresis, visualize DNA by staining in 0.5 µg/ml ethidium bromide solution or SYBR® Green I.
- Choose the gel percentage according to the tables below:

Table 1. Recommended Agarose Gels for Electrophoretic Separation of DNA Fragments.

Agarose gel, %	Range of effective separation, bp	Approximate positions of tracking dyes, bp*			
		Bromophenol blue		Xylene cyanol FF	
		TBE buffer	TAE buffer	TBE buffer	TAE buffer
0.5	2000-50000	750	1150	13000	16700
0.6	1000-20000	540	850	8820	11600
0.7	800-12000	410	660	6400	8500
0.8	800-10000	320	530	4830	6500
0.9	600-10000	260	440	3770	5140
1.0	400-8000	220	370	3030	4160
1.2	300-7000	160	275	2070	2890
1.5	200-3000	110	190	1300	1840
2.0	100-2000	65	120	710	1040
3.0	25-1000	30	60	300	460
4.0	10-500	18	40	170	260
5.0	10-300	12	27	105	165

Table 2. Recommended Polyacrylamide Gels for Electrophoretic Separation of DNA Fragments (1).

Polyacrylamide gel (with BIS at 1:20), % (w/v)	Range of effective separation*	Approximate positions of tracking dyes*	
		Bromophenol blue	Xylene cyanol FF
Denaturing gels			
4.0	100-500 b	50 b	230 b
5.0	70-400 b	35 b	130 b
6.0	40-300 b	26 b	105 b
8.0	30-200 b	19 b	75 b
10.0	20-100 b	12 b	55 b
15.0	10-50 b	10 b	0 b
20.0	5-30 b	8 b	28 b
30.0	1-10 b	6 b	20 b
Non-denaturing gels			
3.5	100-1000 bp	100 bp	460 bp
5.0	80-500 bp	65 bp	260 bp
8.0	60-400 bp	45 bp	160 bp
12.0	50-200 bp	20 bp	70 bp
15.0	25-150 bp	15 bp	60 bp
20.0	5-100 bp	12 bp	45 bp

Note

* Positions of the tracking dyes can only be estimated approximately because the dye front migrates as wide band. The following guidelines are recommended:

- Only high purity agarose should be used. TopVision™ Agarose was used to prepare the gels.
- Only freshly prepared electrophoresis buffers should be used. The buffers were prepared from 50X TAE Buffer and 10X TBE Buffer.
- Choose electrophoresis conditions according to the recommendations below:

Size of the DNA	Voltage	Buffer
<1 kb	5-10 V/cm	TBE
1-5 kb	4-10 V/cm	TAE or TBE
> 5 kb	1-3 V/cm	TAE
Up to 10 kb, fast electrophoresis with Express DNA ladders	up to 23 V/cm	TAE

thermoscientific.com

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

North America
Technical Services:
 techservice.genomics@thermofisher.com
Customer Services:
 customerservice.genomics@thermofisher.com
 Tel 800 235 9880
 Fax 800 292 6088

Europe and Asia
Technical Services:
 techservice.emea.genomics@thermofisher.com
Customer Services:
 customerservice.emea.genomics@thermofisher.com

Thermo
 SCIENTIFIC

Part of Thermo Fisher Scientific