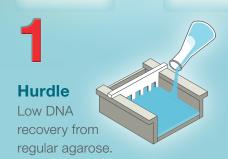
thermoscientific

Five simple wins to keep your DNA electrophoresis on track

Use these straightforward solutions to improve your results



Solution

DNA can be denatured during the melting process when using standard meltingpoint agarose since high temperatures are required to convert the agarose to a liquid. Use low-melting point (LMP) agarose to isolate DNA without the need for enzymes or electroelution. Thermo Scientific™ TopVision™ Agarose comes in standard- and low-melting temperature options.

Hurdle

Contamination or errors introduced from multiple pipetting steps.

Solution

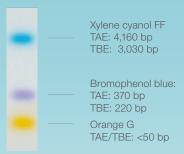
Use products for direct gel loading to avoid repetitive addition of loading buffer to samples and ladders. This can be an ideal time-saving method for high-throughput PCR.

Hurdle

Insufficient tracking of DNA migration—DNA bands run off the gel without the visual cue of lower molecular weight tracking dye.

Solution

Most loading buffer contains two tracking dyes-bromophenol blue and xylene cyanol FF. Use Thermo Scientific™ TriTrack™ DNA Loading Dye (6X) that contains a third dye, orange G, which runs around 50 bp.



Migration of dyes in TriTrack buffer.

Hurdle

Smeared DNA bands or atypical band migration.

Solution

Incorrect buffer can cause bands to smear. Use Tris-acetate-EDTA (TAE) or Tris-borate-EDTA (TBE) running buffer and follow the proper condition for voltage as shown here.

| Size of DNA | Voltage | Preferred buffer |
|-----------------|------------------|------------------|
| <1 kb | 5-10 V/cm | TBE |
| 1–5 kb | 4-10 V/cm | TAE or |
| 1–3 KD | | TBE |
| >5 kb | 1-3 V/cm | TAE |
| Up to 10 | Up to 23 V/cm | TAE |
| kb, fast | | |
| electrophoresis | | |
| with DNA | | |
| ladders | | |

Hurdle

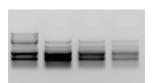
Difficulty in identifying the size of DNA fragments.

Solution

Choosing suboptimal ladders can lead to incorrect sizing of your DNA fragments. Use Thermo Scientific™ GeneRuler™ DNA ladders, which offer:

- A broad range of size options (1 kb, 100 bp, 50 bp, low range, and high range)
- Sharp and bright reference bands
- TriTrack loading buffer to easily monitor DNA migration during electrophoresis

What happened to my gel?



Diffused bands Low voltage and extended

run time



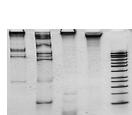
Poor separation Unsubmerged gel and/or short run time



Overloading

Unresolved bands

High salt concentration



Smeared bands

Atypical bands Incorrect loading dye, no denaturation of bound protein

