

Resolution of Linear DNA Fragments from 23 Kilobases to 6 Megabases Using Biphasic Linear Switch Time Ramping

Contributed by Donald R. VanDevanter, Tumor Institute, Swedish Hospital Medical Center, 409 Elkind Hall, 747 Summit Avenue, Seattle, WA 98104

Introduction

Analyses of uncharacterized DNA fragments using CHEF electrophoresis often require preliminary studies to determine what size ranges DNA fragments of interest fall within. It is not uncommon to run more than one gel in order to fully characterize large DNA fragments from a single cell source¹. This is because DNA fragments under 200 kilobase pairs (kbp), between 200 kbp and 2 megabase pairs (mbp), and between 2 mbp and 6 mbp are optimally separated by CHEF electrophoretic techniques using different switching frequencies, electrophoretic field strengths, angles of orientation (if using a Bio-Rad CHEF Mapper™ system), and/or agarose type and concentration². Data is presented illustrating that an initial estimate of sizes can be obtained in a single gel by using a biphasic linear switch time ramping procedure.

Results

A biphasic linear switch time ramping strategy³ with a fixed electric field strength can be used with the CHEF-DR® II system to resolve linear DNA fragments ranging from 23 kbp to 5.7 mbp in size on a single gel (Figure 1). The gel shown in Figure 1A was 0.5% Chromosomal Grade Agarose made in 0.5x TBE (45 mM Tris borate, 1.0 mM EDTA, pH 8.0), recirculated at 10 °C. The applied electric field was held constant at 1.95 V/cm (65 V), and switch times were ramped linearly from 30 seconds to 2 minutes over 33 hours, then from 2 minutes to 50 minutes over 55 hours. Under these conditions, the chromosomes of *Schizosaccharomyces pombe* (closed circles, lanes 4 and 6), *Saccharomyces cerevisiae* (open squares, lanes 3 and 5), lambda phage concatomers (open circles, lane 2), and the larger fragments of *Hind* III-digested lambda DNA (closed squares, lane 1) migrate as an approximate linear function of the log of their respective molecular weights (in mbp, Figure 1B).

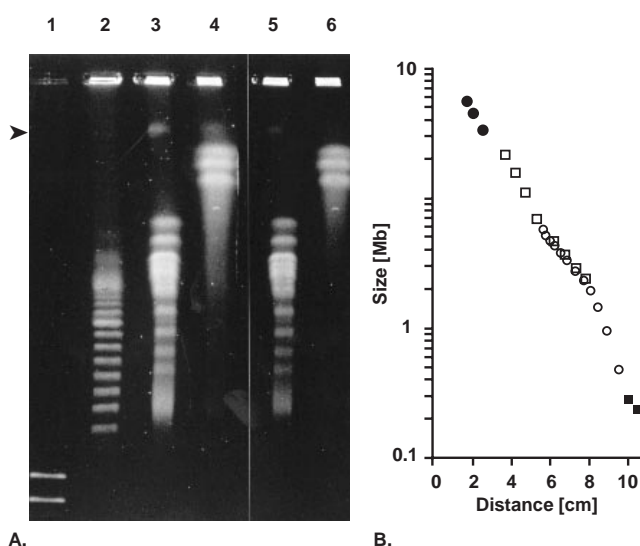


Fig. 1. Separation of DNA fragments ranging from 23 kbp to 5.7 mbp in size by biphasic linear switch time ramping. (A) Ethidium bromide stain of 27.5 kbp and 23.1 kbp lambda-*Hind* III fragments (closed squares, lane 1), lambda concatomers (open circles, lane 2), *S. cerevisiae* chromosomes (open squares, lanes 3 and 5), and *S. pombe* chromosomes (closed circles, lanes 4 and 6). The arrow indicates the electrophoretic zone of compression. Specific electrophoretic conditions are described in the text. (B) Semi-log plot of the sizes of individual fragments (in mbp) versus distance migrated from the origin (in cm). Only selected chromosomes of *S. cerevisiae* are shown.

Lowering the concentration of agarose in gels can have the effect of decreasing run times without an apparent effect on the linearity of separation (Figure 2). The migration profiles shown in Figure 2 were derived by running the *S. pombe* and *S. cerevisiae* chromosomes on 0.5%, 0.4%, 0.3% and 0.2% Chromosomal Grade Agarose gels.

Discussion

The CHEF electrophoresis strategy outlined above is not ideal for sharp resolution of bands within a particular narrow range of DNA fragment sizes, but can be extremely useful for determining what range of fragment sizes deserve closer examination. For this reason, it is a preferable approach to running multiple pulsed-field gels under different conditions when size ranges of greatest interest remain unknown.

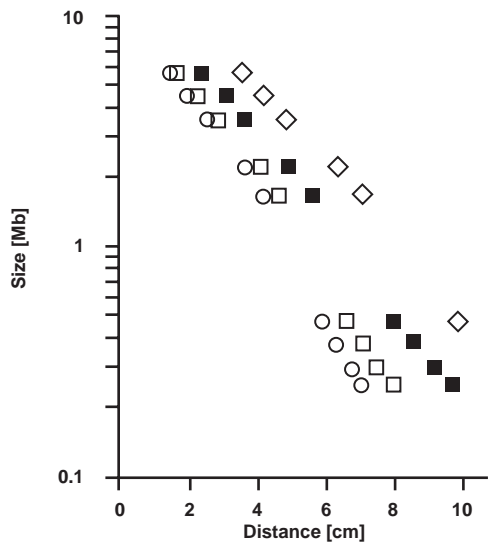


Fig. 2. Effect of agarose concentration on biphasic linear switch time ramping strategy. Semi-log plot of the migration of *S. pombe* and select *S. cerevisiae* chromosomes electrophoresed on a single 0.5% agarose gel (open circles) containing internal lanes of 0.4% (open squares), 0.3% (closed squares), and 0.2% (open diamonds) agarose. Gel parameters were identical to those in Figure 1.

References

1. VanDevanter, D.R., Piaskowski, V.D., Caspar, J.T., Douglass, E.C., and Von Hoff, D.D., *J. Natl. Cancer Inst.* **82**:1815-1821, (1990).
2. Lai, E., Birren, B.W., Clark, S.M., Simon, M.I., and Hood, L., *BioTechniques* **7**:34-42.
3. VanDevanter, D.R., *Nucleic Acids Res.* **20**:1148.

BIO-RAD

Life Science Group

2000 Alfred Nobel Drive
Hercules, California 94547
Telephone (510) 741-1000
Fax: (510) 741-1060

Eastern Regional Office, 85A Marcus Dr., Melville, New York 11747 • Phone (516) 756-2575 • Fax (516) 756-2594
European Headquarters, Bio-Rad Laboratories, Dreve du Senechal, 19, B-1180 Brussels • Phone 02 375 59 70 • Fax 02 374 61 62
Australia, Bio-Rad Laboratories Pty Limited, Unit 11, 112-118 Talavera Rd P.O. Box 371, North Ryde, N.S.W. 2113 • Phone 02-805-5000 • Fax 02-805-1920
Austria, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, A-1130 Wien • Phone 0222-877 89 01 • Fax 0222-876 56 29
Belgium, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, B-9810 Nazareth Eke • Phone 091-85 55 11 • Fax 091-85 65 54
Canada, Bio-Rad Laboratories (Canada) Ltd., 5149 Bradco Boulevard, Mississauga, Ontario L4W 2A6 • Phone (416) 624-0713 • Fax (416) 624-3019
China, Bio-Rad Laboratories, Yanshan Hotel Office Tower, #1307, A138 Haidian Road, Beijing • Phone 2563146 • Fax 2564308
France, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94203 Ivry Sur Seine Cedex • Phone 01-49 60 68 34 • Fax 01-46 71 24 67
Germany, Bio-Rad Laboratories GmbH, Heidemannstraße 164, Postfach 45 01 33, D-8000 München 45 • Phone 089-318 84-0 • Fax 089-318 84 100
Italy, Bio-Rad Laboratories S.r.l., Via Cellini, 18A, 20090 Segrate Milano • Phone 02-21609.1 • Fax 02-21609-399
Japan, Nippon Bio-Rad Laboratories, K. K., Sumitomo Seimei Kachidoki Bldg 5-3-6 Kachidoki, Chuo-Ku, Tokyo 104 • Phone 03-3534-7515 • Fax 03-3534-8027
The Netherlands, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 08385-40666 • Fax 08385-42216
New Zealand, Bio-Rad Laboratories, Pty Ltd., P. O. Box 100-051, North Shore Mail Centre, Auckland 10 • Phone 09-443 3099 • Fax 09-443 3097
Pacific, Bio-Rad Laboratories, Unit 1111, 11/F., New Kowloon Plaza, 38, Tai Kok Tsui Road, Tai Kok Tsui, Kowloon, Hong Kong • Phone 7893300 • Fax 7891257
Scandinavia, Bio-Rad Laboratories, Kanalvagen 10C, 19461 Upplands Vasby, Sweden • Phone 46 (0) 8 590-73489 • Fax 46 (0) 8 590-71781
Spain, Bio-Rad Laboratories, S. A. Avda Valdelaparra 3, Pol. Ind. Alcobendas, E-28100 Alcobendas, Madrid • Phone (91) 661 70 85 • Fax (91) 661 96 98
Switzerland, Bio-Rad Laboratories AG, Kanalstrasse, 17, 8152 Glattbrugg • Phone 01-810 16 77 • Fax 01-810 19 33
United Kingdom, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Phone 0800 181134 • Fax 0442 259118