

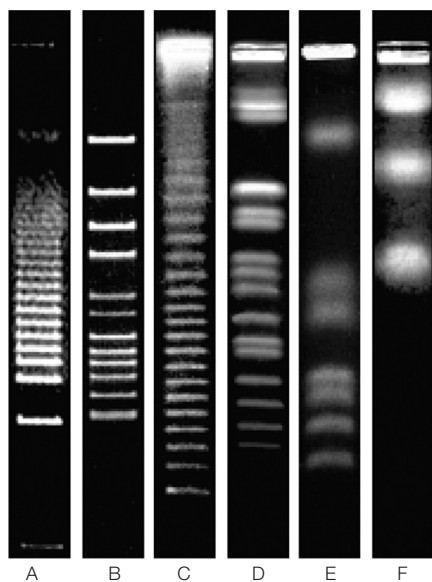
Pulsed Field Gel Electrophoresis Standards

Bio-Rad offers standards for all pulsed field gel electrophoresis (PFGE) applications, from field inversion gel electrophoresis (FIGE) separation of cosmid inserts to the largest chromosomal separations. Higher molecular weight standards are prepared in low-melt agarose blocks that can be cut to fit any well dimensions.

No more than four samples should be placed between lanes with standards. Placement of standards may be adjusted when fewer than 7 (for 10-well gels) or 11 (for 15-well gels) samples are loaded on a gel.

For 10-well gels, load standards in lanes 1, 5, and 10

For 15-well gels, load standards in lanes 1, 5, 10, and 15



Standard and markers. A, 5 kb ladder standard; B, 8.3–48.5 kb standard; C, lambda ladder standard; D, *S. cerevisiae* marker; E, *H. wingei* marker; F, *S. pombe* marker.

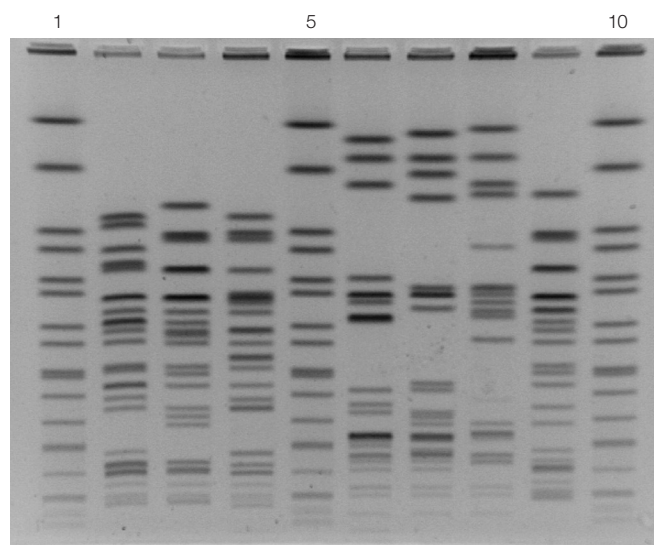
PulseNet Global Reference Standard:

Salmonella enterica, serotype Braenderup H9812

Information in this section is intended for PulseNet members.

H9812 is used as the global reference standard by PulseNet member laboratories. It was selected because when digested with XbaI, it has an optimal number of bands (18) distributed across a wide range of sizes (20.5–1,135 kb).

- H9812 is used to normalize migration variation occurring across the gel and to accurately determine sample band sizes
- H9812 is always digested with XbaI, regardless of which enzyme is used to digest the actual samples
- No more than 4 samples should be placed between H9812 lanes. Placement of standards may be adjusted when fewer than 7 (for 10-well gels) or 11 (for 15-well gels) samples are loaded on a gel:
 - For 10-well gels, load standards in lanes 1, 5, and 10
 - For 15-well gels, load standards in lanes 1, 5, 10, and 15

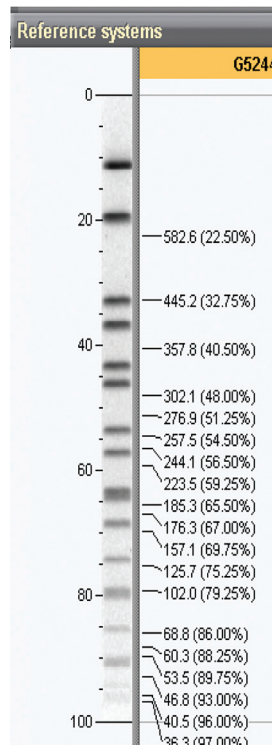


Lanes 1, 5, and 10 include the H9812 standard.

Recommendations

- Prior to using a new lot of H9812 plugs, compare it with a previously used lot to confirm that the new lot produces an acceptable PFGE pattern, that is no pattern changes, no changes in band intensity, and no DNA degradation
- Pretested H9812 standard plugs serve as a positive control for XbaI digests and can be useful in troubleshooting efforts:
 - If a problem is observed in sample lanes but not the standard lanes, it was caused by something that affected only the sample (for example, sample plug preparation, BlnI digest)
 - If a problem is observed in both sample and standard lanes, it is caused by a step involving both sample and standard plugs (for example, XbaI digest, electrophoresis)

Tip: Make an H9812 plug when preparing sample plugs to control for parameters and reagents used to make plugs on that day. If H9812 and sample plugs have deteriorated but pretested H9812 standards appear appropriately, the problem occurred during plug preparation.



S. enterica, serotype Braenderup H9812 standard.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000