

## Recipes for TAE and TBE Electrophoresis Buffers

Agarose gels are generally run two types of electrophoresis buffers. Nucleic acid agarose gel electrophoresis is usually conducted with either Tris-acetate-EDTA (TAE) buffer or Tris-borate-EDTA (TBE) buffer. While TAE buffer provides faster electrophoretic migration of linear DNA and better resolution of supercoiled DNA, TBE buffers have a stronger buffering capacity for longer or higher voltage electrophoresis runs.

**Note:** Because of higher voltages and resulting higher currents often used with Sub-Cell® Model 96 and 192 cells, we recommend that only TBE buffers be used for electrophoresis with these specific systems.

### Tris-acetate-EDTA (TAE) buffer

TAE is often prepared in concentrated stock solutions of 10× or 50× in the laboratory. A 1× working solution is prepared prior to electrophoresis.

Composition of 1× TAE buffer

- 40 mM Tris (pH 7.6)
- 20 mM acetic acid
- 1 mM EDTA

### Preparation of 50× TEA stock solution

To prepare 1 liter of 50× TAE dissolve following components in 600 ml of deionized water:

- 242 g Tris base (FW = 121)
- 57.1 ml glacial acetic acid
- 100 ml 0.5 M EDTA (pH 8.0)

Adjust the final volume to 1 liter with deionized water.

To prepare a 1× working solution from 50× stock buffer mix 50× stock buffer with DNase free deionized water at 1:4 ratio.

### Tris-borate-EDTA (TBE) buffer

TBE buffer can be made and stored in concentrated stocks of 5× or 10×.

Composition of 1× TBE buffer

- 89 mM Tris (pH 7.6),
- 89 mM boric acid,
- 2 mM EDTA

### Preparation of 10× TBE stock solution

To prepare 1 liter of 10× TBE dissolve following components in 600 ml deionized water:

- 108 g Tris base (FW = 121)
- 55 g boric acid (FW = 61.8)
- 40 ml 0.5 M EDTA (pH 8.0)

Adjust final volume to 1 liter with deionized water.

To prepare a 1× working solution from 10× stock buffer, mix 10× stock buffer with DNase free deionized water at 1:9 ratio.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

---

**Web site** [www.bio-rad.com](http://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