

## Transient Gene Expression in Adherent HEK293 Cells using PEI STAR™ Transfection Reagent

### Materials

- PEI STAR™, 1 mg/mL, pH neutralized, sterile-filtered
- HEK293 expression medium or DMEM maintained at 37°C transfection is inhibited by serum. Only use media that is reduced-serum, serum-free or chemically defined.
- HEK293 cells, ~50,000 cells/cm<sup>2</sup>, seeded the day prior to transfection
- Transfection-grade plasmid DNA (pDNA) with gene of interest, 1 µg per 10<sup>6</sup> cells to be transfected.

### General Guidance

We list recommended reagent quantities per 10<sup>6</sup> cells in the table below. If performing many very similar transfections (e.g., making many similar viruses), we would recommend co-varying on these parameters over optimization ranges first.

Parameter	Recommended	Optimization Range
<b>PEI STAR™ Concentration (per 10<sup>6</sup> cells)</b>	3.0 µg	2.0 to 4.0 µg
<b>DNA Concentration (per 10<sup>6</sup> cells)</b>	1.5 µg	1.0 to 2.0 µg
<b>PEI STAR™/DNA Complex Time</b>	10.0 min	5.0 to 15.0 min

Even if performing **co-transfection**, keep the total DNA concentration range the same (1.0 to 2.0 µg/10<sup>6</sup> cells).

If high toxicity is observed, then reduce the quantity of PEI and/or DNA first. We **do not** recommend replacing the media post-transfection to reduce toxicity except as a last resort.

### Before Transfection

Seed the cells the day before transfection to reach 50-80% confluence on the day of transfection. Seed approximately 50,000 cells per cm<sup>2</sup>.

### Transfection

1. Measure the cell density to determine the transfection parameters based on the number of viable cells.
2. Dilute 1.0 µg pDNA per 10<sup>6</sup> cells in 5% final culture volume of DMEM or media. Mix well.
3. Dilute 3.0 µg PEI STAR™ per 10<sup>6</sup> cells in 5% final culture volume of DMEM or media. Mix well.

4. Mix the solution of pDNA and PEI STAR™ together by quickly and briefly vortexing or inverting the tube several times.
5. Allow the mixture of pDNA and PEI STAR™ to rest for 10 minutes at room temperature.
6. Gradually add the pDNA and PEI STAR™ mixture to the cells dropwise while swirling the plate.
7. Incubate the cells at typical incubation conditions.

## References

**Boussif *et al*** (1995) A versatile vector for gene and oligonucleotide transfer into cells in culture and *in vivo*: polyethylenimine. *Proc.Natl.Acad.Sci.U.S.A.* **92** 7297. PMID: [7638184](#).

**Longo *et al*** (2013) Transient mammalian cell transfection with polyethylenimine (PEI). *Methods Enzymol.* **529** 227. PMID: [24011049](#).

## Support

Support is available by emailing us at [techsupport@bio-techne.com](mailto:techsupport@bio-techne.com).