

# Transfection Quality Plasmid DNA in As Little As Ten Minutes Using the PureYield™...

Transfection Quality Plasmid DNA in As Little As Ten Minutes Using the PureYield™ Plasmid Miniprep System

---

## ABSTRACT

*Here we introduce the PureYield™ Plasmid Miniprep System that allows researchers to isolate high-quality plasmid DNA suitable for a broad range of downstream applications, including transfection, in vitro expression, sequencing and cloning. The system allows direct purification of plasmid from as much as 0.6 ml of bacterial culture without cell harvesting for from up to 3 ml of culture when the cells are pelleted. The entire procedure can be completed in as little as 10 minutes.*

*Cell Notes 21, 3–5.*

**Don Smith and Eric Vincent**

*Promega Corporation*

Publication Date: 2008



## Download Article

(276 KB)

---