

# Plasmid Minipreps with a 10-year guarantee.



## Wizard® *Plus* SV Plasmid Minipreps.

**Ten years** of extracting clean plasmids.  
**Thousands** of citations.



**Promega**

## SPECIAL OFFERS NOW\*

**BUY 2 KITS GET 1 FREE**  
Wizard® *Plus* SV Minipreps,  
250 preps (A1460)

**BUY 1 KIT AT 30% OFF**  
Wizard® *Plus* SV Minipreps,  
1000 preps (A1465)

### ENQUIRIES?

Call us

6513 3450

Email us

[sg\\_custserv@promega.com](mailto:sg_custserv@promega.com)

\*Terms apply: Valid until 31 December 2012. Not valid with any other existing promotions. Terms subject to change without prior notice.

# Greater plasmid sizes at >10kb are possible.



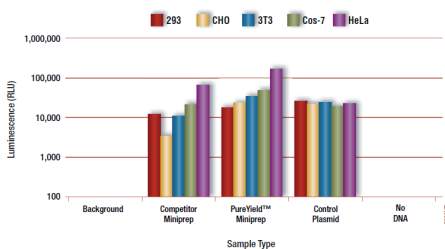
## PureYield™ Plasmid Minipreps.

for demanding applications  
such as transfection and  
cell-free expression

**Greater** plasmid purity.  
**Greater** speed - purify in 10 minutes.

### Transfection-grade Plasmid DNA

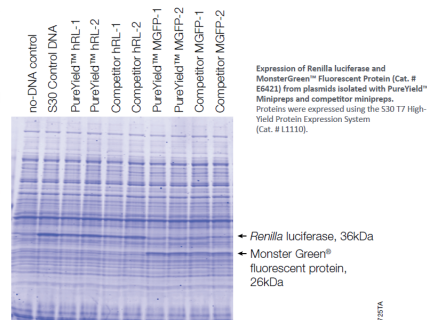
Go directly from plasmid prep to transfection



Transfection of SV40 promoter-driven firefly luciferase vector (pGL4.13 Vector) and subsequent assay in 5 different cell lines.

### Cell-Free-Expression-Grade Plasmid

Express protein in vitro directly after the prep, no further clean-up necessary



## SPECIAL OFFER\*

**BUY 2 KITS GET 1 FREE**  
PureYield™ Plasmid  
Minipreps, 250 preps  
(A1222)



ENQUIRIES? Call us 6513 3450 • Email us [sg\\_custserv@promega.com](mailto:sg_custserv@promega.com)

\*Terms apply: Valid until 31 December 2012.  
Not valid with any other existing promotions.  
Terms subject to change without prior notice.