

Fast-n-Easy Plasmid Preps



- ▶ *For PCR, sequencing, restriction digest and transformation/transfection*
- ▶ *Pure DNA within 10*/40** min*
- ▶ *Come at costs one can't ignore*

** Plasmid Mini-Prep Kit (vacuum-based protocol)*

*** Plasmid Midi-Prep Kit, Plasmid Maxi-Prep Kit (vacuum-based protocol)*

Fast-n-Easy Plasmid Preps



Plasmid Mini-Prep Kit (PP-204)



Plasmid Midi-Prep Kit (PP-211)



Plasmid Maxi-Prep Kit (PP-212)

DNA Yields	up to 20 µg	up to 200 µg	up to 1 mg
Efficiency	hands-on time ≤ 10 min	hands-on time ≤ 40 min	hands-on time ≤ 40 min
Versatility	centrifuge- and vacuum-based purification		
Quality	high-purity plasmid/cosmid DNA suitable for all downstream application		
Convenience	all steps performed at room temperature no need for phenol-chloroform-extraction or alcohol precipitation		
Economy	column refill packs available		
Customer Support	available in English, Spanish, French, Portuguese & German (info@jenabioscience.com)		

Request a sample pack

free shipping within Germany – visit www.jenabioscience.com/plasmid
or contact one of our international distributors for purchases outside of Germany.

