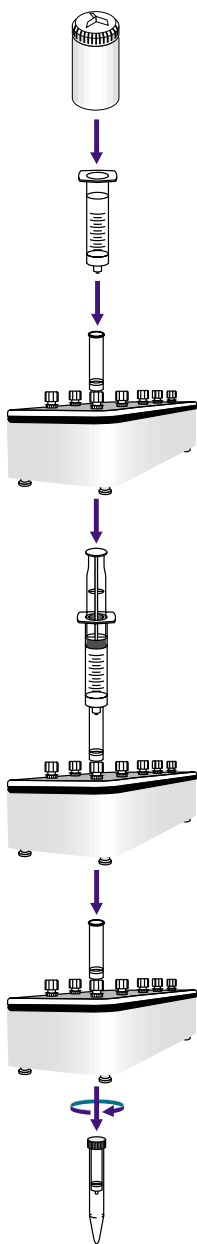


# GENELUTE HP PLASMID MIDIPREP KIT



## Vacuum Format

Preparation: See Technical Bulletin for Details

¥ Add RNase A to the Resuspension Solution

¥ Add Ethanol to the Wash Solution 2

¥ Chill the Neutralization Solution P

### 1 Harvest & Lyse Bacteria

Pellet 50 ml of an overnight culture at 5,000 x g, 10 min. Discard supernatant.

Resuspend cells in 4 ml of Resuspension Solution. Pipet up and down, or vortex.

Add 4 ml of Lysis Solution and gently invert 6-8 times to mix. Do not vortex. Allow to clear, 3-5 min.

Remove the plunger from a filter syringe and place the barrel in an upright position.

### 2 Prepare Cleared Lysate

Add 4 ml of Neutralization Solution P to the lysed cells & gently invert 6-8 times to mix.

Add 3 ml of Binding Solution G & gently invert 1-2 times to mix.

Immediately add the mix to the barrel of the filter syringe & let sit for 5 min.

### 3 Prepare Column

Place a binding column onto the vacuum manifold & apply the vacuum.

Add 4 ml of Column Preparation Solution to the column & allow it to pass through.

### 4 Bind Plasmid DNA to Column

Remove the binding column from the collection tube & place on a vacuum manifold & apply the vacuum.

Hold the filter syringe over the column & gently insert the plunger to expel the cleared lysate. Allow the lysate to pass through the column.

### 5 Wash to Remove Contaminants

Add 4 ml of Wash Solution 1 to the column & allow to pass through.

Add 4 ml of Wash Solution 2 to the column & allow to pass through.

Leave vacuum on for 10 min to dry the column.

### 6 Elute Purified Plasmid DNA

Transfer the column to a collection tube provided.

Add 1 ml of Elution Solution & spin in a swinging bucket rotor at 3,000 x g, 5 min.

See [Technical Bulletin](#) for detailed procedure