

Quantum Prep[®] Vacuum Manifold

Introduction

The Quantum Prep Vacuum Manifold is designed for use with Quantum Prep HT/96 Plasmid Miniprep Kits. This product is specifically designed to provide users with added convenience in high-throughput purification of nucleic acids.

Product Components

1 Manifold Housing	1 Hose Barb
1 Manifold Base	1 3.5 ft piece Tygon tubing
1 Bottom Gasket with pressure sensitive adhesive backing	
1 Top Gasket with pressure sensitive adhesive backing	

See Quantum Prep HT/96 instruction manuals for details on use of this manifold for vacuum filtration-mediated high-throughput plasmid preps.

Initial Assembly

1. Remove the self adhesive protection strip from the smaller (top) gasket provided in the kit, and seat this gasket on top of the Manifold Housing.
2. Remove the self adhesive protection strip from the larger (bottom) gasket provided in the kit, and seat this gasket in the groove on the Manifold Base.
3. Plug the Hose Barb into the Manifold Housing.
4. Attach vacuum tubing from the Hose Barb now plugged in to the Manifold Housing to your vacuum trap and connect the vacuum trap to either house vacuum or a vacuum pump.

Optional: A vacuum regulator may be placed between the vacuum trap and the Manifold, or between the vacuum source and the vacuum trap to monitor and regulate vacuum pressures.

Notes

Optimal performance is obtained at a vacuum pressure of 20"-25" Hg.

Product Information

732-6150	Quantum Prep HT/96 ClearVac Kit, 2x96
732-6151	Quantum Prep HT/96 ClearSpin Kit, 2x96
732-6152	Quantum Prep HT/96 ClearVac Kit, 10x96
732-6153	Quantum Prep HT/96 ClearSpin Kit, 10x96