

Product Guidelines for Site Installation

Product Name:	Thermo Scientific™ Barnstead™ MicroPure™ Water Purification Systems	Item Description:	Type 1 Lab Water System
Model #	50132373, 50132370	Manufacturer	Thermo Scientific
Order #	50132373, 50132370		

NOTE: Barnstead GenPure Pro water purification systems come complete with pressure regulator, wall mounting bracket, tubing, and all necessary consumables.

A. RECOMMENDED FEED WATER SPECIFICATIONS

Colloid Index	SDI max.	<u>< 1</u>	SDI
Resistivity	µS/cm	<u>> 0.5</u>	µS/cm
TOC Value	ppb max.	<u>< 50</u>	ppb
Free Chlorine	ppm max.	<u>< 0.05</u>	ppm
Turbidity	NTU	<u><1.0</u>	NTU
Carbon Dioxide	ppm max.	<u>< 30</u>	ppm
Silicate	ppm max.	<u>< 2</u>	ppm
Pressure	PSI	<u>1.45 - 87</u>	PSI
Feed Water Temp.	°C	<u>2 - 40 C</u>	°C
Recommended Type	ASTM 1, 2, or 3	<u>2</u>	

B. ELECTRICAL POWER REQUIREMENTS

Electrical Specifications

115/240V Volts 50/60 Hertz 2 - 1 A Amps Single # of Phases

Total number of receptacles / plugs required to operate water system. 1

Length of line cord 8 Feet 0 Inches

C. PLUMBING REQUIREMENTS

Water Inlet Line

Type of Fitting Male NPT

Size of Fitting 3/4 Inch

Manual Cutoff Valve Strongly Recommend

Waste Drain

Type of Drain Atmospheric

Max. Distance from Unit 5 Feet

D. SPATIAL REQUIREMENTS

Equipment exterior dimensions (inches): 12 Width 12 Depth 21.5 Height

Equipment Weight (lbs): 50 Lbs.

Wall mount specifications: If the system is to be mounted on the wall or on a shelf the designated area must be capable of supporting up to 200 lbs.

E. Accessories / Replacement Consumables

Accessories

09.2212 System Wall Mounting Bracket

Replacement Consumables

09.1006 Ultra pure water cartridge

09.1003 0.2µm Final Filter

09.1102 Disinfection Cartridge

CMX25 Disinfection Solution, 1 Time Use

50133981 Ultrafilter

09.1002 Replacement UV lamp