

Product Guidelines for Site Installation

Product Name:	Thermo Scientific™ Barnstead™ GenPure xCAD Plus™ Water Purification Systems		Item Description:	Type 1 Lab Water System with Bench Mounted Remote Dispenser
	50136152, 50136151, 50136153, 50136146			
Model #	50136153, 50136146		Manufacturer	Thermo Scientific
Order #	50136152, 50136151, 50136153, 50136146			

NOTE: Barnstead GenPure xCAD Plus 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	<u>115/240V</u>	Volts	<u>50/60</u>	Hertz	<u>2 - 1 A</u>	Amps	<u>Single</u>	# of Phases
Total number of receptacles / plugs required to operate water system.	<u>1</u>							
Length of line cord	<u>8</u>	Feet	<u>0</u>	Inches				

C. PLUMBING REQUIREMENTS

Water Inlet Line	
Type of Fitting	<u>Male NPT</u>
Size of Fitting	<u>3/4</u> Inch
Manual Cutoff Valve	<u>Strongly Recommend</u>
Waste Drain	
Type of Drain	<u>Atmospheric</u>
Max. Distance from Unit	<u>5</u> Feet

D. SPATIAL REQUIREMENTS

Equipment exterior dimensions (inches):	<u>14.7</u>	Width	<u>13.3</u>	Depth	<u>24.2</u>	Height
Remote Dispenser dimensions (inches):	<u>12.3</u>	Width	<u>21</u>	Depth	<u>28.5</u>	Height
Equipment Weight (lbs):	<u>55</u>	Lbs.				
Remote Dispenser Weight (lbs):	<u>26.5</u>	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.

Remote Dispenser can be mounted no more than 9 feet away from the water system. Additional remote dispensers maybe mounted up to 9 feet away from each other.

E. Accessories / Replacement Consumables

Accessories

AY1137X1 Log Printer for connection to RS232 port (110V)

Replacement Consumables

09.2005 Ultra pure water cartridge

09.1003 0.2μm Final Filter

09.2201 Disinfection Cartridge

CMX25 Disinfection Solution, 1 Time Use

50133980 Ultrafilter

09.2002 Replacement UV lamp

50136494* Additional bench mounted "client" xCAD remote dispenser

50136505* Additional wall mounted "client" xCAD remote dispenser

* Up to 2 additional "client" xCAD remote dispensers can be connected per water system in any mounting configuration (wall or bench).