

Product Guidelines for Site Installation

Product Name:	Thermo Scientific™ Barnstead™ GenPure xCAD Plus™ Water Purification Systems	Item Description:	Type 1 Lab Water System with Wall Mounted Remote Dispenser
Model #	50131317, 50131315, 50131321, 50131323	Manufacturer	Thermo Scientific
Order #	50131317, 50131315, 50131321, 50131323		

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

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