

Product Guidelines for Site Installation

Product Name:	Thermo Scientific™ Barnstead™ Smart2Pure™ 12LPH Water Purification Systems	Item Description:	Tap to Type 1 Lab Water System
Model #	50129890, 50129845	Manufacturer	Thermo Scientific
Order #	50129890, 50129845		

NOTE: Barnstead Smart2Pure water purification systems come complete with pressure regulator, tubing, and all necessary consumables. Required 1 micron pretreatment filter, 09.4003, and reservoir is not included with the system therefore must be ordered separately.

A. RECOMMENDED FEED WATER SPECIFICATIONS

Blocking Index	SDI max.	<3	SDI
Manganese	mg/l	> 0.05	mg/l
Iron	mg/l	> 0.05	mg/l
Free Chlorine	mg/l	< 0.1	mg/l
pH Range	pH	4-11	pH
Pressure	PSI	1.45 - 87	PSI
Feed Water Temp.	°C	2-35	°C
Recommended Type	ASTM 1, 2, or 3	Tap	DIN2000

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	15.7	Depth	21.5	Height
30L Reservoir Dimensions (inches):	15	Width	15	Depth	23.5	Height
60L Reservoir Dimensions (inches):	15	Width	15	Depth	35.9	Height
Pretreatment dimensions (inches):	5.1	Width	5.1	Depth	12.2	Height
Equipment Weight (lbs):	40	Lbs.				
30L Reservoir Weight (full) (lbs):	70	Lbs.				
60L Reservoir Weight (full) (lbs):	140	Lbs.				
Pretreatment Weight (lbs):	15	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.
 Wall mount specifications: If the **reservoir** 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

Required Accessories

09.4003	1 micron prefilter
06.5040	30L Storage Reservoir (must select 30L or 60L but not both)
06.5070	60L Storage Reservoir (must select 30L or 60L but not both)

Optional Accessories

09.2212	Wall mounting bracket for system
09.4001	5 micron filter and hardness stabilizer for customers with SDI >3
50138221	Hand dispenser
06.5015	30L Reservoir wall mounting bracket
06.5016	60L Reservoir wall mounting bracket

Replacement Consumables

09.1020	Ultra pure water cartridge
09.1003	0.2µm Final Filter
09.2012	RO Membrane with integrated pretreatment for 12LPH system
CMX25	Disinfection Solution, 1 Time Use
50133981	Ultrafilter
09.1002	Replacement UV lamp
06.5001	Sterile Overflow Filter
06.5002	0.2 micron filter with CO2 Adsorber