

Product Guidelines for Site Installation

Product Name:	Thermo Scientific™ Barnstead™ Pacific™ TII UV Water Purification Systems	Item Description:	Tap to Type 2 Lab Water System
Model #	50132129, 50132131, 50132132, 50131982, 50132133	Manufacturer	Thermo Scientific
Order #	50132129, 50132131, 50132132, 50131982, 50132133		

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

A. RECOMMENDED FEED WATER SPECIFICATIONS

Resistivity	µS/cm	> 0.001	µS/cm
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):	14.6	Width	13	Depth	23.7	Height
30L Reservoir Dimensions (inches):	15	Width	15	Depth	23.5	Height
60L Reservoir Dimensions (inches):	15	Width	15	Depth	35.9	Height
100L Reservoir Dimensions (inches):	15	Width	15	Depth	49.2	Height
Pretreatment dimensions (inches):	12	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.				
100L Reservoir Weight (full) (lbs):	220	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.

E. Accessories / Replacement Consumables

Required Accessories

09.4000	5 micron prefilter with activated carbon and hardness stabilizer
06.5032	30L Storage Reservoir with level display and recirculation pump (must select 30L or 60L but not both)
06.5062	60L Storage Reservoir with level display and recirculation pump (must select 30L or 60L but not both)
06.5082	100L Storage Reservoir with level display and recirculation pump (must select 30L or 60L but not both)

Optional Accessories

AY1137X1	Printer
50138221	Hand dispenser
06.5015	30L Reservoir wall mounting bracket
06.5016	60L Reservoir wall mounting bracket

Replacement Consumables

09.4011	High purity deionizer cartridge
22.0046	RO membrane with integrated pretreatment for 3, 7, 12, and 20LPH systems (order 2 for 20LPH)
22.0087	RO Membrane with integrated pretreatment for 20LPH system (order 2 for 40LPH)
CMX25	Disinfection Solution, 1 Time Use
09.4002	Replacement UV lamp
06.5001	Sterile Overflow Filter
06.5002	0.2 micron filter with CO2 Adsorber

