

Product Guidelines for Site Installation

| | | | |
|---------------|---|-------------------|---------------------------------------|
| Product Name: | Thermo Scientific™ Barnstead™ Smart2Pure™ 3LPH and 6LPH Water Purification Systems | Item Description: | Tap to Type 1 Lab Water System |
| Model # | 50129872, 50129688, 50129885, 50129887 | Manufacturer | Thermo Scientific |
| Order # | 50129872, 50129688, 50129885, 50129887 | | |

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

A. RECOMMENDED FEED WATER SPECIFICATIONS

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

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>12</u> | Width | <u>15.7</u> | Depth | <u>21.5</u> | Height |
| Pretreatment dimensions (inches): | <u>5.1</u> | Width | <u>5.1</u> | Depth | <u>12.2</u> | Height |
| Equipment Weight (lbs): | <u>40</u> | Lbs. | | | | |
| Pretreatment Weight (lbs): | <u>15</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.

E. Accessories / Replacement Consumables

Required Accessories

09.4003 1 micron prefilter

Optional Accessories

09.2212 Wall mounting bracket for system

09.4001 5 micron filter and hardness stabilizer for customers with SDI >3

Replacement Consumables

09.1020 Ultra pure water cartridge

09.1003 0.2µm Final Filter

09.2003 RO Membrane with integrated pretreatment for 3LPH system

CMX25 Disinfection Solution, 1 Time Use

09.2006 RO Membrane with integrated pretreatment for 6LPH system

50133981 Ultrafilter

09.1002 Replacement UV lamp

