



## STA 8000: Site Readiness Instructions

### a) Order Review:

Please review this order. Record any discrepancies between the PerkinElmer order and your Purchase Order, along with any agreements or commitments made by your PerkinElmer Sales Representative that are NOT listed on the order. Let your Customer Care Representative (1 800 762 4000) know about these discrepancies and/or commitments with your Site Readiness Confirmation

### b) Site Requirements:

#### Bench Space

STA 8000	450 Watts Maximum
Circulator	1440 Watts Maximum

A bench top of 100 cm (39.4in) W x 65 cm (24 in) D will accommodate an STA 8000 system with no accessories.

#### Peripherals, Accessories

Circulating chiller fan must be positioned so warm air out of the chiller is not blown onto either the instrument or autosampler (if equipped).

### c) Electrical Requirements:

#### Power Consumption

STA 8000	450 Watts Maximum
Circulator	1440 Watts Maximum

#### Power Specifications

STA 8000	100-240V/ 50/60 Hz.
Circulator	120 VAC, 12 Amps or 240 VAC, 7 Amps

### Power Outlets

Each system component (STA 8000, Computer, Printer, etc) requires a separate power outlet. All outlets should share a common earth ground.

This equipment is designed to operate within 10% of the selected line voltage. The supply must be smooth, clean and free of transient voltages over 40 volts.

Earth grounding: less than 1 ohm resistance between the grounds of any 2 components of the system

### d) Gas Requirements:

NOTE: ALL GASSES AND REGULATORS MUST BE SUPPLIED BY CUSTOMER. IF THE REQUIRED GAS QUALITY IS NOT AVAILABLE A FILTER DRIER IS REQUIRED.

IF SPLITTING A SINGLE GAS SOURCE IS NECESSARY, PLEASE CONTACT YOUR LOCAL SERVICE REPRESENTATIVE FOR INFORMATION ON THE REGULATOR INSTALL KIT PART NUMBER N519-0462 SERVICE 1 800 762 4000

A Nitrogen balance purge of ~ 40 cc/min. is required. This gas should be clean and dry having minimum purity of 99.995%. Regulator Outlet Pressure **2-3 bar (30-45 psi)**

Sample purge gases can be: dry argon, air, nitrogen or oxygen, with a minimum purity of 99.995%.

The regulator shutoff valve should have a 1/4" NPT male thread on the outlet side.

Regulator Outlet pressure **2-3 bar (30-45 psi)**

### e) Coolant requirements:

The STA 8000 **requires** a coolant. Performance will vary with the coolant temperature and stability. A controlled coolant temperature of  $\pm 0.1$  °C is required for best performance. The cooling can be one of the following types:

**Circulating water:** A source of water and a drain are required.

**Flow rate: 0.5-3.0 l/minute required (1 l/min optimum)**

**Water temperature:  $\geq 15$  °C controlled to  $\pm 0.1$  °C**

**Chiller:** A liquid circulating device such as the PolyScience Model 9102. The 9102 water circulator does not include the cooling liquid. 50/50 Ethylene Glycol (non automotive grade, Fisher P/N E1774), Distilled Water and Algaecide (N0776059 8 drops per gallon) is required as the cooling media to 15 °C. Minimum temperature control  $\pm 0.1$  °C is required.

**Coolant temperatures: 15 °C  $\geq$  35 °C (Non-condensing coolant temperature)**

**Flow rate: 0.5-3.0 l/minute required (1 l/min optimum)**

## f) Environmental Requirements:

### Laboratory Environment

Temperature range	15 to 40°C
Humidity	<80% (Non-condensing)

Clean and dust-free

Level, vibration-free work surface

### Water Requirements

The STA 8000 comes equipped with a tap water circulating system. In this configuration tap water or water circulating device (i.e. PolyScience type chiller) must be located conveniently near the STA 8000. The coolant tubes provided (2) are approximately 5 feet long (1.5 meters) having an OD of 12 mm (0.47 in) and ID of 8 mm (0.31 in). If additional tubing length is required, it is the responsibility of the customer to provide the additional tubing.

**Flow rate: 0.5-3.0 l/minute required (1 l/min optimum)**

## g) Safety Requirements:

### Gas Cylinders and Gas Delivery Lines

Lock down straps should be present on all gas cylinders.

### Ventilation

As required

## h) PC Configuration:

Due to numerous differences in PC hardware, PerkinElmer cannot guarantee that our software will run on a customer-supplied computer, even if the system meets the minimum specifications described on this page.

PerkinElmer installation of a customer-supplied computer is available for an additional fee.

PerkinElmer is not responsible for problems caused by unspecified system components, software, and/or accessories. The additional time it takes to verify this type of problem is **billable** at the current Service rate.

PerkinElmer **may not** provide maintenance service on the customer-supplied computers.

### Recommended for the PYRIS Thermal Analysis Systems

If providing a custom PC it should be equivalent to the system specified below.

- Model: Lenovo ThinkCentre M82 Tower
- Processor: Intel® Core™ i5 -3550 CPU @ 3.30 GHz, 3.30 GHz
- Factory Installed Operating System: Microsoft® Windows® 7 Ultimate 32 bit
- Memory: 4 GB
- Hard Drive: 1 TB 7200rpm SATA
- Networking: Integrated Intel Gigabit Ethernet
- External I/O Ports: 4 USB 3.0 (4 rear), 4 USB 2.0 (2 front, 2 rear), 1 Ethernet (RJ45), 2 serial (9-pin), 1 VGA out
- Expansion Slots: 2 PCI, 1 PCIe
- Removable Media: DVD Recordable

- Video: Integrated Graphics
- Mouse: Edge Mouse - USB 2-button optical mouse with scroll
- Keyboard: USB Keyboard
- Audio: Integrated Audio

### **Operating System Requirements**

Win7/32 bit

## **i) Installation Overview:**

**At receipt of the instrumentation, customer is to determine if any obvious shipping damage has occurred and is to notify the shipping company immediately.**

The customer is to unpack and verify the layout of Instruments and accessories with an eye to possible missing or damaged parts.

You should unpack the instrument and accessories. The instrument should be placed on the bench along with associated components.

### **Customer to hold Packing Material for CSE**

The STA 8000 is to be installed by a PKI Service Engineer, you should hold the instrument packing material for the Customer Service Engineer (CSE).

### **Physical Installation (Instrument Only)**

The physical installation will vary based on system configuration.

### **Physical Installation (Software)**

A PKI Service Engineer will install the instrument control and data analysis software. If the customer IT person is required to be in attendance to perform the software installation it is the responsibility of the user to make the arrangements with their IT Department prior to the installation to confirm their presence.

### **Physical Installation (Accessories)**

[As Required]

### **Installation Test Standards**

A PKI Service Engineer will test the instrument to insure that its performance meets PerkinElmer's specifications. Reference materials are supplied with the instrument to test its functionality.

### **Customer Orientation**

The customer must be available for an instrumentation orientation.

## **i) Miscellaneous:**

If IQOQ validation is required for this instrument, please contact your local Service representative at 1800 762 4000.

If a journal is required for this analyzer, please contact your local Sales /Service representative at 1 800 762 4000. (The Journal is a binder that should be located near the instrument. It provides an organized means of storing miscellaneous instrument information such as: service records, certificates tracking usage etc.)