

# Sample-Draw Transmitter

**Instruction Manual**

Part Number 71-0114

5May2010

© 2007 Thermo Fisher Scientific Inc. All rights reserved.

Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Thermo Fisher Scientific  
Air Quality Instruments  
27 Forge Parkway  
Franklin, MA 02038  
1-508-520-0430  
[www.thermo.com/aqi](http://www.thermo.com/aqi)

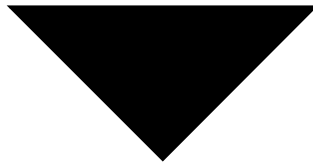
# WEEE Compliance

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the following symbol:



Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on Thermo Fisher Scientific's compliance with these Directives, the recyclers in your country, and information on Thermo Fisher Scientific products which may assist the detection of substances subject to the RoHS Directive are available at: [www.thermo.com/WEEERoHS](http://www.thermo.com/WEEERoHS).





# WARNING



THIS INSTRUMENT IS DESIGNED TO DETECT ONE OR MORE OF THE FOLLOWING:

FLAMMABLE VAPORS, OXYGEN CONTENT, AND/OR TOXIC GAS AND TO GIVE WARNING BEFORE THEY REACH HARMFUL CONDITIONS. IN ORDER TO ENSURE THAT IT WILL WARN OF DANGEROUS CONCENTRATIONS, IT IS ESSENTIAL THAT THE INSTRUCTIONS IN THIS MANUAL, PARTICULARLY THOSE CONCERNING START UP, OPERATION, CALIBRATION, AND MAINTENANCE, BE READ, UNDERSTOOD, AND FOLLOWED.

## NOTATION CONVENTIONS

Notices are used in this operator's manual to alert you to hazardous conditions to person or instrument and to notify you of additional information. This operator's manual uses the following notices.



---

### **WARNING**

Notifies you of potential danger that can result in personal injury or death.

---



---

### **CAUTION**

Notifies you of potential damage to equipment.

---



---

### **NOTE**

Notifies you of additional or critical information.

---

## **SERVICE LOCATIONS**

For additional assistance, service is available from exclusive distributors worldwide. Contact one of the phone numbers below for product support and technical information or visit us on the web at [www.thermo.com/aqi](http://www.thermo.com/aqi).

**866-282-0430 Toll Free**  
**508-520-0430 International**

# Warranty

Seller warrants that the Products will operate or perform substantially in conformance with Seller's published specifications and be free from defects in material and workmanship, when subjected to normal, proper and intended usage by properly trained personnel, for the period of time set forth in the product documentation, published specifications or package inserts. If a period of time is not specified in Seller's product documentation, published specifications or package inserts, the warranty period shall be one (1) year from the date of shipment to Buyer for equipment and ninety (90) days for all other products (the "Warranty Period"). Seller agrees during the Warranty Period, to repair or replace, at Seller's option, defective Products so as to cause the same to operate in substantial conformance with said published specifications; provided that (a) Buyer shall promptly notify Seller in writing upon the discovery of any defect, which notice shall include the product model and serial number (if applicable) and details of the warranty claim; (b) after Seller's review, Seller will provide Buyer with service data and/or a Return Material Authorization ("RMA"), which may include biohazard decontamination procedures and other product-specific handling instructions; and (c) then, if applicable, Buyer may return the defective Products to Seller with all costs prepaid by Buyer. Replacement parts may be new or refurbished, at the election of Seller. All replaced parts shall become the property of Seller. Shipment to Buyer of repaired or replacement Products shall be made in accordance with the Delivery provisions of the Seller's Terms and Conditions of Sale. Consumables, including but not limited to lamps, fuses, batteries, bulbs and other such expendable items, are expressly excluded from the warranty under this warranty.

Notwithstanding the foregoing, Products supplied by Seller that are obtained by Seller from an original manufacturer or third party supplier are not warranted by Seller, but Seller agrees to assign to Buyer any warranty rights in such Product that Seller may have from the original manufacturer or third party supplier, to the extent such assignment is allowed by such original manufacturer or third party supplier.

In no event shall Seller have any obligation to make repairs, replacements or corrections required, in whole or in part, as the result of (i) normal wear and tear, (ii) accident, disaster or event of force majeure, (iii) misuse, fault or negligence of or by Buyer, (iv) use of the Products in a manner for which they were not designed, (v) causes external to the Products such as, but not limited to, power failure or electrical power surges, (vi) improper storage and handling of the Products or (vii) use of the Products in combination with equipment or software not supplied by Seller. If Seller determines that Products for which Buyer has requested warranty services are not

covered by the warranty hereunder, Buyer shall pay or reimburse Seller for all costs of investigating and responding to such request at Seller's then prevailing time and materials rates. If Seller provides repair services or replacement parts that are not covered by the warranty provided in this warranty, Buyer shall pay Seller therefor at Seller's then prevailing time and materials rates. ANY INSTALLATION, MAINTENANCE, REPAIR, SERVICE, RELOCATION OR ALTERATION TO OR OF, OR OTHER TAMPERING WITH, THE PRODUCTS PERFORMED BY ANY PERSON OR ENTITY OTHER THAN SELLER WITHOUT SELLER'S PRIOR WRITTEN APPROVAL, OR ANY USE OF REPLACEMENT PARTS NOT SUPPLIED BY SELLER, SHALL IMMEDIATELY VOID AND CANCEL ALL WARRANTIES WITH RESPECT TO THE AFFECTED PRODUCTS.

THE OBLIGATIONS CREATED BY THIS WARRANTY STATEMENT TO REPAIR OR REPLACE A DEFECTIVE PRODUCT SHALL BE THE SOLE REMEDY OF BUYER IN THE EVENT OF A DEFECTIVE PRODUCT. EXCEPT AS EXPRESSLY PROVIDED IN THIS WARRANTY STATEMENT, SELLER DISCLAIMS ALL OTHER WARRANTIES, WHETHER EXPRESS OR IMPLIED, ORAL OR WRITTEN, WITH RESPECT TO THE PRODUCTS, INCLUDING WITHOUT LIMITATION ALL IMPLIED WARRANTIES OF MERCHANTABILITY OR FITNESS FOR ANY PARTICULAR PURPOSE. SELLER DOES NOT WARRANT THAT THE PRODUCTS ARE ERROR-FREE OR WILL ACCOMPLISH ANY PARTICULAR RESULT.





# TABLE OF CONTENTS

## Chapter 1

### Introduction

Overview .....	1
Description .....	1
Target Gases .....	2
Specifications .....	3

## Chapter 2

### Installation

Mounting the Transmitter .....	5
Connecting the Sample Lines .....	6
Wiring the Transmitter .....	7

## Chapter 3

### Start Up & Operation

Preparing for Start Up .....	9
Setting the Channel Parameters at the Controller .....	9
Setting the Fresh Air Signal .....	10

## Chapter 4

### Calibration

Preparing for calibration .....	12
Calibrating the Detector .....	12

## Chapter 5

### Maintenance

Preventive Maintenance .....	17
Troubleshooting .....	18
Replacing the Toxic or Oxygen Sensor .....	22
Replacing the Combustible Gas Sensor .....	23
Replacing the Toxic and Oxygen Amplifier .....	24
Replacing the Combustible Amplifier .....	25
Replacing the Pump .....	26

**Appendix A**


Parts List ..... 29

**Appendix B**

External Wiring Instructions ..... 33

**Appendix C**

Calibration Response Charts ..... 37



Chapter  
1

## INTRODUCTION

### Overview

The Sample-Draw Transmitter is a series of oxygen, combustible gas, and toxic gas transmitters.

### Description

The Sample-Draw Transmitter assembly is comprised of a sample-drawing apparatus, main board, gas sensor, a temperature compensation circuit, and an amplifier/transmitter electronic assembly all enclosed in a weather and dust resistant housing.

## Target Gases

Table 1-1 lists the target gas and detection ranges for the Sample-Draw Transmitter.

**Table 1-1 Target Gases and Detection Ranges**

<b>Target Gas</b>	<b>Detection Range</b>
Ammonia	0 to 100 ppm
Arsine	0 to 1 ppm
Carbon Monoxide	0 to 500 ppm
Chlorine	0 to 10 ppm
Chlorine Dioxide	0 to 2 ppm
Diborane	0 to 1 ppm
Fluorine	0 to 10 ppm
Hydrogen Chloride	0 to 30 ppm
Hydrogen Cyanide	0 to 50 ppm
Hydrogen Fluoride	0 to 10 ppm
Hydrogen Sulfide	0 to 100 ppm
Nitric Oxide	0 to 100 ppm
Nitrogen Dioxide	0 to 20 ppm
Oxygen	0-30% vol
Ozone	0 to 1 ppm
Phosphine	0 to 1 ppm
Silane	0 to 20 ppm
Sulfur Dioxide	0 to 20 ppm
Combustible gases	0 to 100% LEL <sup>1</sup>

<sup>1</sup>. Lower Explosive Limit (may also be used in ranges of 0 to 5,000 or 10,000 parts per million for some hydrocarbons).

## Specification

Table 1-2 lists the Sample-Draw Transmitter performance, electrical and environmental specifications.

**Table 1-2 Specifications**

Amplifier Outputs	4 to 20 mA analog signal (to controller) 100 to 500 mV analog test signals (at test jacks)
Fault Condition	0 mA loop current
Sampling Method	Sample-Draw
Accuracy	± 10% detection range
Repeatability	± 5% detection range
Housing	Fiberglass (with Lexan <sup>1</sup> window), NEMA 4X
Dimensions	10.0 in. H x 7.15 in. W x 4.5 in. D (25.4 cm H x 18.2 cm W x 11.4 cm D)
Weight	4 lbs (1.8 kg)
Enclosure Rating	NEMA 4X
Area Classification	Non-hazardous locations
Power Source	12 to 24V DC
Temperature Range	- 4° F to 113° F (-20° C to 45° C)
Standard Accessories	Manual In-line hydrophobic filter (some versions) <sup>2</sup>
Optional Accessories	Calibration kits, gas cylinders, gas collection bag

<sup>1</sup>. Lexan is a registered trademark of General Electric Company.

<sup>2</sup>. This filter is not furnished for use with reactive gases Cl<sub>2</sub>, ClO<sub>2</sub>, F<sub>2</sub>, HCl, HF, NH<sub>3</sub>, NO<sub>2</sub> and O<sub>3</sub>. These gases react with or are absorbed by entrapped water or other residue, and thus would be lost prior to entering the sample-drawing system.



# INSTALLATION

## Mounting the Transmitter



### WARNING

Perform all installation procedures in a fresh air environment (known to be free of combustible and toxic gases and of normal oxygen content). The transmitter is not in operation as a gas monitoring system until the start-up procedure is complete.



### CAUTION

The Sample-Draw Transmitter is not suitable for Class I hazardous areas. Mount the Sample-Draw Transmitter in a non-hazardous area

1. Select a mounting area that is indoors or sheltered from rain or snow. Make sure there is enough room to mount the housing, open the housing door, and to make wiring and sample line connections at the bottom of the housing.
2. Secure the housing to the vertical surface using bolts or screws through the mounting flanges (see Figures 2-1).

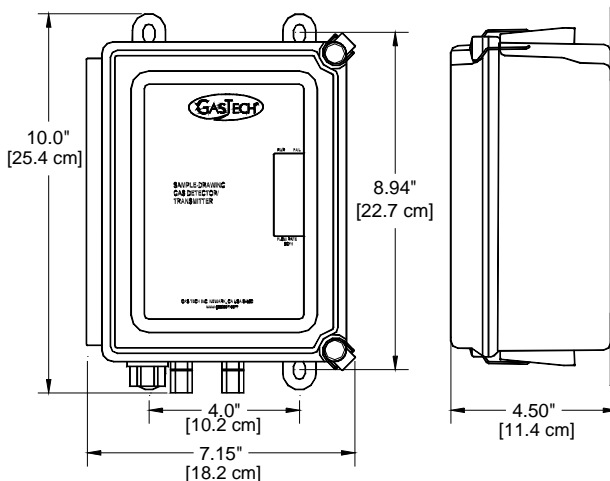
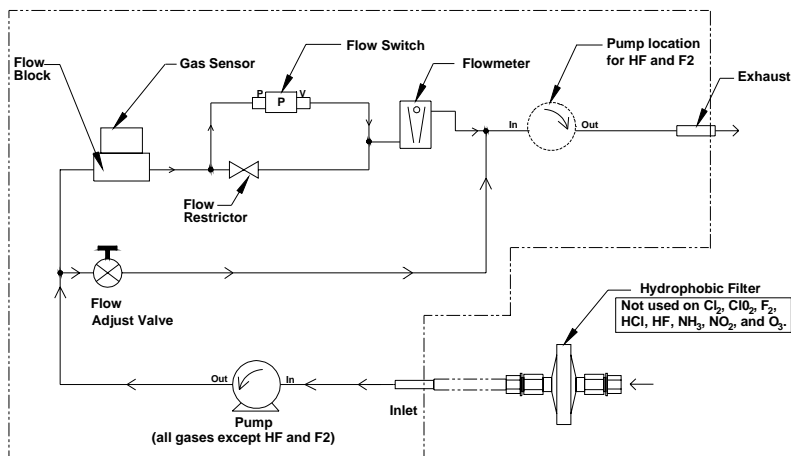


Figure 2-1 Outline and Mounting Dimensions

## Connecting the Sample Lines

Refer to Figure 2-2 Flow Diagram and Figure 2-3 for Inlet and Outlet connections.



**Figure 2-2 Flow Diagram**

1. Attach 1/4 in. OD TEFLON® tubing to the inlet fitting. The other end of the tubing should be at the monitoring location.
2. If supplied, insert the filter into the incoming sample line at a point near the inlet fitting, of the sample-draw transmitter.
3. If necessary, connect a 1/4 in. OD plastic or stainless steel tube to the EXHAUST fitting. Keep the exhaust tubing as short as possible. Route the exhaust to an open area where the sample can safely dissipate.

®DuPont registered trademark.



---

## Wiring the Transmitter

Use a three-conductor, shielded cable, or run the wiring within metal conduit to reduce Radio Frequency Interference (RFI) and Electromagnetic Interference (EMI).

Use 18 AWG wire or larger. The two-way wire and receiver resistance must not exceed 200 ohms for 12 volts systems or 600 ohms for 24 volts systems.



---

### WARNING

**Make sure all power to the controller is turned off during all wiring procedures.**

---

- The power circuit (“+12-24 volts” and “—”) supplies both the flow system (pump and signal LEDs) and the signal loop circuit.
- The signal loop circuit (amplifier/transmitter plus all external components such as controller, displays, etc.) all of which share the common 4-20 mA current loop, which begins at “+12-24 volts” and terminates back at controller common “—”.



---

### WARNING

**Do Not Overload the Transmitter Signal loop or signal limiting due to excessive loop resistance could result in understatement of hazardous conditions. From Ohm’s Law, a signal loop powered from a 12 volt source cannot deliver more than 20 mA (full scale reading) if the total loop resistance, including all wiring, this transmitter (25 ohms internal), the receiver, and all other loop device input circuits, exceed 600 ohms. Operation from 24 volts permits about twice the loop resistance. Do not operate from voltages below 12 volts.**

---



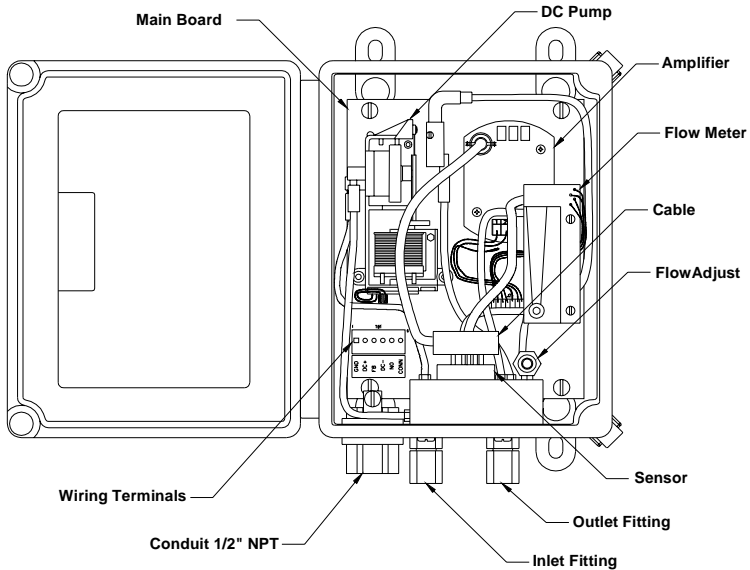
---

### CAUTION

**Do not run transmitter and AC power wiring through the same conduit.**

---

1. Connect the sample-draw transmitter to the receiver, controller or other 12-24 volt DC power source.



**Figure 2-3 Component Location**

2. Ground the shield cable or metal conduit by connecting it to the SHIELD or GND terminal at the receiver or controller. Do not ground any of the three power/signal terminals in the sample drawing transmitter to its housing.
3. Confirm that the detector is connected to the amplifier.
4. Confirm that the terminal block on the amplifier is connected to the PC board cable.
5. Turn on the power to the system at the power source.
6. Confirm that the flowmeter "PILOT" light is on and the flowmeter indicates a flow rate of approximately 0.6 SCFH. Adjust if necessary.
7. Confirm that the incoming sample line is not leaking. To test the sample line, put your thumb over the end of the incoming sample tubing. If the flowmeter ball drops to the bottom of the column and the flow fail circuit activates, the sample tubing is ready to use.



**CAUTION**

**Always ground the cable or conduit at the controller and never at the transmitter.**

---

## START UP & OPERATION

Complete the following procedure to place the controller and the transmitter into normal operation.

### Preparing for Start Up

1. Complete the mounting and wiring procedures described in the Installation chapter of this manual.
2. Connect incoming power to the controller as described in the receiver or controller manual.



---

#### NOTE

Allow the sensor to stabilize for 1 hour after power is introduced before calibrating.

---

### Setting the Channel Parameters at the Controller

The sample drawing transmitter may be used with a wide variety of receivers and controllers which provide 12 to 24 volts DC power at 400 mA and a means of accepting and interpreting a 4-20 mA signal proportional to gas concentration in the ranges shown in Table 1-1 of this manual

When you install the transmitter, set the parameters for that channel at the controller. For instructions on setting the channel parameters at the controller, see the receiver or controller manual.

## Setting the Fresh Air Signal

---



### WARNING

**If you suspect the presence of target gas during the start-up procedure, use the calibration kit described in the calibration chapter and the zero-emission air cylinder to introduce fresh air to the detector and confirm an accurate zero setting (span setting for oxygen monitoring systems).**

---

1. Confirm that the receiver, controller or DC power source are on.
2. Open the housing cover.



### NOTE

Some controllers have a feature that enables you to disable the alarm LEDs, buzzer, and relays during the calibration procedures, response tests, and so on. Refer to the controller manual for information. Make sure you re-enable full alarm functionality after the calibration procedure is complete.

---

3. Confirm that the sample line inlet is sampling from a source of clean air.
4. Select the millivolt (mV) range on the multimeter. Plug the positive multimeter lead into the white (+) test jack: plug the negative lead into the blue (-) test jack.
5. Confirm a reading of 100 mV (toxic/combustible) or 379 mV (oxygen) on the multimeter. Adjust the potentiometer on the amplifier if necessary.
  - Toxic/combustible - Adjust “ZERO” pot until the multimeter reading is 100 mV
  - Oxygen - Adjust “SPAN” pot until the multimeter reading is 379 mV
6. Remove the multimeter leads from the test jacks, and secure the cover to the housing.

The transmitter is now in operation.

## CALIBRATION

This chapter describes the optional calibration kit used for calibrating the transmitter. See Parts List, in Appendix A of this manual for ordering information.



---

### WARNING

**Accurate calibration of the transmitter is essential to ensure accurate readings of toxic gas concentrations. Incorrect calibration can impair the performance of the transmitter and place you in unnecessary danger if hazardous conditions exist.**

---

The calibration kit contains all of the equipment you need to introduce a calibration sample to the detector. It includes the following components:

- Storage Case (safely stores the components of the calibration kit).
- Cylinder (contains a known concentration of target gas).
- Regulator (controls the flow of the sample from the cylinder to the detector)
- Tubing (connects components of the calibration kit).
- Gas collecting Bag (permits matching of regulator flow rate to the sample pump flow rate). One of two types may be furnished:
  - TEDLAR™ bag may be used for all gases.
  - Vinyl bag is suitable for Nitrogen, hydrogen, propane, hexane and carbon monoxide, or Zero Air.
- Pinch Clamp and Y fitting (controls gas flow between the gas collecting bag and transmitter).

For calibration it is preferable to introduce the calibration gas directly to the inlet of the sample line. If this is not practical, temporarily remove the sample line tubing from the Inlet fitting of the sample-draw adapter, connect a short piece of tubing in its place and admit the calibration gas there. Be sure to replace the sample line when calibration is complete.

## Preparing for Calibration

This section describes how to prepare the transmitter for calibration. The procedure includes step-by-step instructions for preparing the calibration kit and the transmitter.



---

### WARNING

**Calibrate the detector in a fresh air environment (environment known to be free of toxic gases). If the in-line hydrophobic filter is used during normal operation, calibrate the detector with the filter in place.**

---

## Calibrating the Detector

This chapter describes how to prepare the gas cylinder, set the clean air base line and gas response reading for the toxic gas and oxygen transmitters.

### Calculating the Calibration Gas Response Reading

The 100 to 500 mV test signal at the test jack of the transmitter amplifier is used to calibrate the transmitter. The following formula describes how to calculate the output test signal as a function of the gas concentration:

$$\text{Test signal} = ((\text{gas concentration}/\text{full scale range}) \times 400 \text{ mV}) + (100 \text{ mV})$$

- For example, if you are using a gas cylinder of 5 ppm chlorine to calibrate a transmitter whose full-scale range is 0-10 ppm:

$$((5 \text{ {ppm conc.}})/10 \text{ {ppm full-scale conc.}}) \times 400 \text{ mV {range}} + (100 \text{ mV {offset}}).$$

$$= (5/10 \times 400 \text{ mV}) + (100 \text{ mV}) = 300 \text{ mV gas signal setting on multimeter.}$$

- For oxygen, the measurement range is 0-30% by volume and the recommended calibration gas is clean air, which contains 20.9% oxygen, so the calculation is:

$$((20.9/30) \times 400) + (100) = 378.7 \text{ (or } 379) \text{ mV.}$$

For user convenience, graphs of signal versus gas concentration are shown in Appendix C of this manual for each gas range.

---

## Preparing the Gas Cylinder

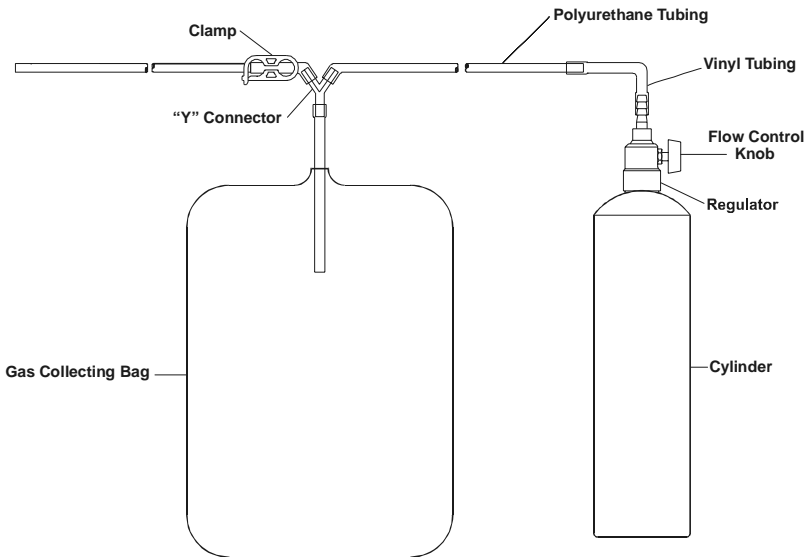
---

**NOTE**

When performing the following steps, use gas samples of known concentration. Cylinders of known concentration are available from Thermo Fisher Scientific (see Parts List in Appendix A). Specify gas and concentration when ordering gas cylinders.

---

1. Verify the regulator flow control valve is closed, then carefully screw the regulator onto the cylinder.
2. Verify that all tubing connections are tight and secure.



**Figure 4-1 Calibration Kit with Gas Cylinder and Gas Collecting Bag**

## Calibrating the Toxic Gas or Combustible Detector

1. Verify the absence of toxic/combustible gas or any atmosphere other than clean air at the calibration site, by ventilation with clean air or by testing with a portable gas analyzer.
2. Open the housing.
3. Set the multimeter to 0-500 DC mV range. Plug the positive multimeter lead into the white (+) test jack and the negative lead into the blue (-) test jack of the amplifier.
4. Confirm that the multimeter shows a reading of 100 mV. If the reading is other than 100 mV adjust the ZERO potentiometer on the amplifier so that it does so. Leave the multimeter probes connected to the test jacks for setting the toxic gas response.



---

### NOTE

For Cl<sub>2</sub>, ClO<sub>2</sub>, F<sub>2</sub>, HF, NO<sub>2</sub> and O<sub>3</sub> transmitters, turn the zero control counter-clockwise to increase reading, for all other gases turn it clockwise to increase reading.

---

## SETTING THE TOXIC OR COMBUSTIBLE GAS RESPONSE READING

1. Set the multimeter to 0-500 DC mV range. Plug the positive multimeter lead into the white (+) test jack and the negative lead into the blue (-) test jack.
2. Close the clamp, then open the flow control knob until the gas collecting bag is approximately 3/4 full.
3. Open the clamp and allow the sample-draw pump to draw gas from the gas collecting bag. Adjust the regulator so the gas collecting bag remains 3/4 full.
4. After the reading stabilizes (normally 1 to 2 minutes), confirm a reading of the SPAN pot on the amplifier until the multimeter reads the value calculated at the beginning of this section or adjust the SPAN potentiometer on the amplifier so that it does so.



---

### CAUTION

**If the reading on the multimeter has not stabilized after 2 minutes, see the Troubleshooting section of this manual for slow response.**

---

5. Remove the multimeter test probes and secure cover to housing.
6. Close the flow control knob and disconnect the calibration gas cylinder. Flatten the gas collecting bag to expel any remaining gas.

The transmitter is now in normal operation.



---

## Calibrating the Oxygen Detector

### SETTING THE OXYGEN ZERO READING

1. Set the multimeter to 0-500 DC mV range. Plug the positive (+) multimeter lead into the white (+) test jack and the negative lead into the blue (-) test jack of the amplifier.
2. Attach the regulator to the cylinder of 100% nitrogen.
3. Connect the regulator output to the sample line inlet.
4. Observe the millivolt reading on the multimeter, which should decline towards 100 mV as nitrogen enters the detector and displaces residual air.
5. When the multimeter reading has stabilized, turn the “ZERO” pot on amplifier until reading is 100 mV.
6. Leave the multimeter connected to the amplifier test jacks.
7. Close the flow control knob on the regulator and disconnect the nitrogen cylinder. Flatten the gas collecting bag to expel any remaining gas.

### SETTING THE OXYGEN RESPONSE READING

For calibration of oxygen transmitters, the recommended calibration gas is clean fresh air, if its purity can be assured by ventilation or by test with a portable oxygen indicator. If this is impractical, a cylinder of certified ZERO AIR should be used. For clean fresh air, the calibration setting should be 379 mV at the test jacks (equivalent to 15.1 mA in signal loop), to correspond with the 20.9% oxygen present in clean air. For ZERO AIR, calculate the proper span setting for the analyzed oxygen content (if other than 20.9%) from the formula or from the chart in Appendix C of this manual, then proceed as follows:

1. If using clean air for calibration, assure that the sample line inlet is relocated to a source of air known to be free of contamination, then go directly to step 5 (otherwise connect the regulator to a cylinder of certified ZERO AIR).



### NOTE

The ZERO AIR cylinder is not included in the standard calibration kit, as clean air is most often readily available to the detector by normal methods of ventilation. It may be ordered separately (81-0076) and uses the same regulator (81-1003) as the nitrogen cylinder.

---

2. Attach the flow regulator to the cylinder of zero air and the Y fitting of the gas collecting bag.
3. Close the clamp and then open the flow control valve until the gas collecting bag is approximately 3/4 full.
4. Open the clamp and allow the sample-draw pump to draw gas from the gas collecting bag. Adjust the flow regulator so the gas collecting bag remains 3/4 full.
5. After the reading stabilizes (normally 1 to 2 minutes), confirm a reading of 379 mV (or calculated value for cylinder oxygen content, if different from 20.9%) on the multimeter. If necessary, adjust the "SPAN" pot on the amplifier until the multimeter reads correctly.
6. Close the flow control knob and disconnect the ZERO AIR gas cylinder. Flatten the gas collecting bag to expel any remaining gas. If using the method of relocating sample line inlet, be sure too move the sample line inlet back to the desired monitoring location.

The transmitter is now in normal operation.

# MAINTENANCE



---

## WARNING

**Perform all maintenance activities in a non-hazardous environment.**

---

## Preventive Maintenance

This schedule describes daily, monthly, and quarterly procedures to ensure the performance and durability of the transmitter.

### Daily

1. Verify that the receiver indicates a reading near 0 ppm (toxic), 0% LEL (combustible) or near 21% volume (oxygen). Investigate significant changes in the display reading and signal output.
2. Confirm that the transmitter pilot light (above the flowmeter) is on.
3. Confirm that the flowmeter indicates a flow rate of approximately 0.6 SCFH. Adjust if necessary using the flow adjustment valve.

### Monthly



---

## NOTE

Some controllers have a feature that permits disabling of the alarm devices during maintenance and calibration procedures. Refer to the controller manual for information. Be sure to enable full alarm functionality after the calibration or response test procedure is complete.

To test any visual, audible and relay alarm indications by the controller during the response test, use a concentration of gas greater than the alarm setpoints.

If you have evacuation alarms or alarms that are forwarded to the fire department, be sure to notify the appropriate people before you test the alarms.

---

1. Confirm that the controller display reading is approximately 0 ppm (toxic) and 21% (oxygen). If not, set the fresh air signal at the amplifier to 100 mV (toxic) or 379 mV (oxygen), as described in the Start Up & Operation chapter of this manual.
2. Assemble the calibration kit and introduce the gas to the detector as described in the Calibration chapter of this manual.
3. After the reading stabilizes (normally 1 to 2 minutes), confirm that the display reading for the controller responds to the gas sample as the sample is introduced to the detector and is within  $\pm 10\%$  of the desired reading. If not, perform the calibration procedure as described in the Calibration chapter of this manual.
4. Confirm that the receiver or controller alarm functions (if any) respond appropriately.
5. Turn off the calibration gas. Disassemble and store the calibration kit as described in the Calibration chapter of this manual.

### Quarterly



---

#### **CAUTION**

**Calibrate the transmitter at least once every three months. Some applications may require a more frequent calibration schedule.**

---

Perform the calibration procedure as described in the Calibration chapter of this manual.

## Troubleshooting

The section describes symptoms, probable causes, and suggested responses for problems you may encounter with the transmitter.

### Fault Condition

#### **Symptoms:**

- FAULT or FAILURE indication.
- Negative reading at receiver or controller.
- Transmitter flow “FAIL” light is on.
- Transmitter flowmeter indicates less than 0.6 SCFH.

**Probable Causes:**

- The power supply, controller, Sample-Draw Transmitter termination PC board, amplifier or detector wiring connections are incorrect or incomplete, see Chapter 2.
- Sample-Draw Transmitter ZERO potentiometer is adjusted incorrectly, see Chapter 3.
- Low flow due to clogged filter, obstructed sample line, failed pump, etc., see Chapter 3.
- Sensor is missing, improperly installed or defective.
- Amplifier improperly set or defective.

**Suggested Response:**

- Check all components for proper connection to power and continuity of signal loop (see wiring diagram in Appendix B).
- Verify other components are properly connected and appear to be operating normally, check the setup of the transmitter. A fault problem can arise in either (or both) of two areas: the sample-draw subsystem or the electronic subsystem.

**CHECK THE SAMPLE-DRAW SUBSYSTEM AS FOLLOWS:**

1. Check that the flowmeter indicates approximately 0.6 SCFH. If this reading is low, attempt to set the correct rate using the flow adjust valve. If you cannot set the correct rate using the flow adjust valve, replace the filter in the sample line and check the sample line for obstructions or kinks.
2. If you still cannot set the correct rate using the flow adjust valve, replace the pump.
3. If the fault condition continues, contact Thermo Fisher Scientific for further instructions.

**CHECK THE ELECTRONIC SUBSYSTEM AS FOLLOWS:**

1. Confirm that the wiring to the controller terminal strip, the main board and the amplifier terminal block is complete and correct.
2. Confirm that the sensor cable is connected to the amplifier.
3. Confirm that the sensor is installed in the flow block.
4. Confirm that the sensor is plugged into the cable to the amplifier.
5. Set the fresh air signal as described in the Start Up & Operation chapter of this manual.

IF THE CONDITION CONTINUES, FURTHER ISOLATE THE PROBLEM AS FOLLOWS:

1. Disconnect the transmitter wires at the controller terminal strip and connect them to the terminals of a different channel or controller known to be operating correctly. If the fault condition clears, the controller terminal strip or main circuit board is bad. Contact Thermo Fisher Scientific for further instructions. If the fault continues, go to the next step.
2. Check the fresh air signal at the amplifier test jack as described in the Start Up & Operation chapter of this manual. If you can set the fresh air signal to 100 mV (toxic/combustible) or 379 mV (oxygen), then the detector amplifier and sample-draw subsystem are operating correctly. Contact Thermo Fisher Scientific for further instruction. If the fault condition continues, go to the next step.
3. Disconnect the detector and connect a detector known to be operating correctly. If the fault condition clears, replace the detector. You can replace the plug-in sensor as an alternate to replacing the entire detector assembly. If the fault condition continues, go to the next step.
4. Connect the detector in question to an amplifier known to be operating correctly. If the fault clears, replace the amplifier.
5. If you replaced the amplifier or detector assembly, perform the calibration procedure.
6. If the fault condition continues contact Thermo Fisher Scientific for further instruction.

---

## Difficult or Unable to Calibrate, Slow or No Response

### Symptoms:

- Unable to accurately calibrate the transmitter.
- Slow or no response to calibration gas during monthly response test.
- Transmitter requires frequent calibration.



### NOTE

Under normal conditions, the transmitter requires calibration approximately every 3 months. Some applications may require a more frequent calibration schedule.

---

### Probable Cause:

- Sample in gas cylinder is low, exhausted or out-dated.
- Sensor is outdated or is reaching the end of its useful life.
- Low flow due to clogged filter, obstructed sample line, failed pump, .

### Suggested Response:

1. Make sure the calibration cylinder has an adequate supply of fresh gas.
2. Replace the plug-in sensor if it is outdated or shows signs of leakage or liquid contamination (investigate for source of liquid, if present).
3. Check the in-line hydrophobic filter (if installed). If the filter is installed, remove it. If the gas response changes after you remove the filter, the filter is contaminated. Replace the filter.



### NOTE

Filter is not recommended for reactive gases such as Chlorine, Fluorine, etc.

---

4. Check the sample line for leaks by plugging the inlet to the sample line and assuring that the flowmeter float ball drops to the bottom of the flow column and that the flow “FAIL” light (above the flowmeter) is lit. If not, find and correct the leak. Unplug the sample line inlet and perform a calibration.
5. If calibration difficulties continue, contact Thermo Fisher Scientific for further instruction.

## Replacing the Toxic or Oxygen Sensor

1. Turn off all incoming power at the power source.
2. Open the housing.
3. Unplug the cable from the sensor, (pull straight up), and pull old sensor from the flow block with fingers (if sensor appears to be stuck, pry **gently** with a small screwdriver).
4. Remove the replacement sensor from its container and **if installed** remove the spring placed between the pins marked "R" (reference) and "S" (sensing).



### NOTE

Only certain toxic sensors are shipped with these shorting springs, others either do not benefit from shorting "Ref" to "Sens" pins, or may be degraded by it. Do not be concerned if the spring is missing on the new sensors.

---



### CAUTION

**Do not remove the spring until you are ready to perform the start-up procedure. The detector will take longer to stabilize if the spring is removed before start-up.**

---

5. Press the new sensor into the cavity until it is firmly seated.
6. Reconnect the cable to the sensor.
7. Turn the power on.
8. Perform the calibration procedure described in the Calibration chapter of this manual. You can zero the detector right away, but for best results, let the detector stabilize for 1 hour before calibrating. (Oxygen sensors may be zeroed and calibrated immediately.)



## Replacing the Combustible Sensor

1. Turn off all incoming power at the power source.
2. Open the housing, disconnect the three sensor wires (green, red, white) from the terminal block (R, A, C) on the amplifier.
3. Remove the sensor and adapter from the flow block by pulling straight up with your fingers. If the sensor appears to be stuck, pry **gently** with a small screwdriver.
4. Remove the four screws from the retaining plate, and slide the adapter and gasket from the sensor. Keep the gasket, adapter, and screws for the next step.
5. Attach the new sensor to the adapter, aligning the gasket between the adapter and retaining plate. Tighten all four screws.
6. Push the sensor adapter into the flow block until it is firmly seated.
7. Reconnect wires to the amplifier terminal strip as shown in Table 5-1.

**Table 5-1 Combustible Detector Wiring**

<b>Combustible Detector</b>	<b>Amplifier Terminal Block</b>
Green (reference)	R (terminal 4)
White (common)	C (terminal 6)
Red (active)	A (terminal 5)
No connection	O <sub>2</sub> (terminal 7), J (terminals 8 & 9)

8. Test the connection by gently pulling the leads from the terminal connections.
9. Turn the power on.
10. Allow the new detector to warm up for 15 minutes, then calibrate the detector. Perform the calibration procedure described in the Calibration chapter of this manual.

## Replacing the Toxic and Oxygen Amplifier

1. Turn off all incoming power at the power source.
2. Open the housing.
3. Unplug the cable from the amplifier.
4. Disconnect wires from the amplifier (green, violet).
5. Remove the two screws that secure the amplifier to the housing and remove amplifier.
6. Place the new amplifier into the housing so the holes in the housing line up.
7. Secure the amplifier to the housing using the same screws. Do not over-tighten.
8. Reconnect the wires from the main board to the amplifier terminal block, refer to table 5-2.

**Table 5-2 Toxic and Oxygen Amplifier Wiring Connections**

<b>Amplifier Terminal Block</b>	<b>Main Board J2 Connector</b>
+	Green
FB	Violet

9. Plug cable in to the amplifier socket.
10. Turn the power on.
11. Perform the calibration procedure described in the Calibration chapter of this manual. You can zero the detector right away, but for best results, let the detector stabilize for 1 hour before calibrating.

## Replacing the Combustible Amplifier

1. Turn off all incoming power at the power source.
2. Open the housing.
3. Disconnect sensor wires from the amplifier (green, red, white).
4. Disconnect power/signal wires from the amplifier (green, blue, violet).
5. Remove the two screws that secure the amplifier to the housing and remove amplifier.
6. Place the new amplifier into the housing so the holes in the housing line up.
7. Secure the amplifier to the housing using the same screws. Do not over-tighten.
8. Reconnect the wires from the main board to the amplifier terminal block, refer to Table 5-3.

**Table 5-3 Amplifier Wiring Connections**

3-Point Terminal Block		6-Point Terminal Block	
Amplifier Terminal	Main Board "J2" Connector	Amplifier Terminal	Sensor Wires
1 (+)	Green	4 (R)	Green
2 (-)	Blue	5 (A)	Red
3 (FB)	Violet	6 (C)	White
-----	-----	7, 8, & 9	No Connection

9. Reconnect the sensor wires to the amplifier terminal block, refer to Table 5-3.
10. Turn the power on.
11. Perform the calibration procedure described in the Calibration chapter of this manual. You can zero the detector right away, but for best results, let the detector stabilize for 1 hour before calibrating.

## Replacing the Pump

1. Turn off all incoming power at the power source.
2. Open the housing and locate the DC pump (see Figure 2-3).  
Disconnect the pump power cable by pulling the white push-on connector forward.
3. Remove the three screws that secure the pump to the short aluminum standoffs from the printed circuit board. The two lower screws are readily visible. The third screw, best removed last while holding the pump, is located inside the pump frame, just above and to the right of the rubber diaphragm. A screw-holding screwdriver is beneficial, but not necessary.
4. Examine the tubing connections from the old pump to the flow system, noting where each of the two white elbow fittings connects to other components in the flow system (see Figure 2-3), and observe that the upper fitting (outlet) is identified with a "+" mark and the lower fitting (inlet) with a "-", both molded into the pump frame. There is also a molded "/" line at an angle showing the proper alignment of each of the keyed white elbow fittings. Note that the new pump has its own new set of the white elbow connectors, but to simplify pump replacement it is recommended that these be removed from the new pump and the elbow and tubing from the old pump transferred intact to the new pump, as follows.
5. Remove each white elbow fitting from the old pump, one at a time, while holding the pump in one hand and pulling the fitting with a slight twist, and transfer it to the new pump, aligned with the "/" mark on the pump body as before. Push the elbow firmly in place (again with a slight twist) to the proper angle. Repeat this with the second fitting. Use care to keep the keyed elbow fittings properly aligned with the keyed nipples to avoid breaking the fittings.
6. Remove the three barbed rubber pump mounts from the new pump (these are furnished to accommodate mounting in an older version of sample draw unit, but not used in later versions) by pulling them out at an angle from the pump.
7. Insert one mounting screw with its lockwasher in the upper mounting hole, holding it in place with a screwdriver. Align the two lower mounting holes with their standoffs, as a guide for starting the upper screw in its standoff (do not tighten yet). Start the other two screws (with their lockwashers in place) in their respective standoffs. Tighten all three screws until snug, but do not over-torque.
8. Plug the pump cable into the connector on the new pump. The connector is polarized to ensure that it is plugged in properly.

9. Restore power to the Sample-Draw unit and allow a few minutes to warm up, then perform the system checks outlined in Chapter 5.

## **Returning for Repair**

Before you remove the transmitter from the monitoring area, first contact a Thermo Fisher Scientific representative.

The Thermo Fisher Scientific representative may guide you through certain diagnostic procedures with the transmitter in place. If you cannot correct the malfunction, the representative will assist you in returning the transmitter for repair.



Appendix  
A

## PARTS LIST

Table A-1 lists part numbers for the Sample-Draw Transmitter replacement parts and accessories.

**Table A-1 Parts List**

<b>Part No.</b>	<b>Description</b>
06-1156	Tubing, TEFLON <sup>®</sup> 1/4" OD x 1/8" ID
06-1253	Tubing, polyurethane, 1/4" OD x 3/16" ID
07-0009	Gasket, flange for Combustible Detector
07-6117	O-ring for Oxygen Sensor Adapter
10-0120	Retaining Screw for Combustible Sensor (4-required)
14-2112	Adapter for Combustible Sensor
14-2113	Adapter for Oxygen Sensor
17-0605	Connector, Y barb (gas collecting bag)
30-0011	Pump, sample-draw transmitter
33-0151	Filter, sample-draw transmitter
57-7045-01	Amplifier, Combustibles
57-7210	Amplifier, AsH <sub>3</sub> , B <sub>2</sub> H <sub>6</sub> , CO, HCN, H <sub>2</sub> S, NH <sub>3</sub> , O <sub>2</sub> , PH <sub>3</sub> , SiH <sub>4</sub> , SO <sub>2</sub>
57-7210-01	Amplifier, Cl <sub>2</sub> , ClO <sub>2</sub> , F <sub>2</sub> , HF, NO <sub>2</sub> , O <sub>3</sub>
57-7210-02	Amplifier, HCl, NO
61-0121-01	Sensor, combustible
65-1061	Sensor, oxygen (O <sub>2</sub> )
65-2417	Sensor, sulfur dioxide (SO <sub>2</sub> )
65-2425-01	Sensor, carbon monoxide (CO)
65-2425-02	Sensor, hydrogen sulfide(H <sub>2</sub> S)
65-2431-01	Sensor, chlorine (Cl <sub>2</sub> ), fluorine (F <sub>2</sub> )
65-2431-02	Sensor, hydrogen cyanide (HCN)
65-2431-03	Sensor, nitric oxide (NO)

**Table A-1 Parts List (Continued)**

<b>Part No.</b>	<b>Description</b>
65-2431-04	Sensor, hydrogen chloride (HCl)
65-2431-05	Sensor, hydrogen fluoride (HF)
65-2431-07	Sensor, ammonia (NH <sub>3</sub> )
65-2431-08	Sensor, arsine (AsH <sub>3</sub> ), diborane (B <sub>2</sub> H <sub>6</sub> ), phosphine (PH <sub>3</sub> ), silane (SiH <sub>4</sub> )
65-2431-09	Sensor, ozone (O <sub>3</sub> )
65-2431-10	Sensor, nitrogen dioxide (NO <sub>2</sub> )
65-2431-11	Sensor, chlorine dioxide (ClO <sub>2</sub> )
71-0114	Sample-Draw Transmitter Operator's Manual
81-0065	Cylinder, CO-in-air (100 ppm)
81-0076	Cylinder, Zero-emission air
81-0078	Cylinder, 100% nitrogen (N <sub>2</sub> )
81-0151	Cylinder, H <sub>2</sub> S-in-N <sub>2</sub> (25 ppm)
81-0170	Cylinder, SO <sub>2</sub> -in-N <sub>2</sub> (5 ppm)
81-0175	Cylinder, NO-in-N <sub>2</sub> (25 ppm)
81-0180	Cylinder, NO <sub>2</sub> -in-N <sub>2</sub> (5 ppm)
81-0188	Cylinder, SiH <sub>4</sub> -in-N <sub>2</sub> (5 ppm)
81-0190	Cylinder, Cl <sub>2</sub> -in-N <sub>2</sub> (5 ppm) <sup>1</sup>
81-0191	Cylinder, NH <sub>3</sub> -in-air (25 ppm)
81-0192	Cylinder, HCN-in-N <sub>2</sub> (10 ppm)
81-0193	Cylinder, PH <sub>3</sub> -in-N <sub>2</sub> (0.5 ppm) <sup>2</sup>
81-0194	Cylinder, HCl-in-N <sub>2</sub> (10 ppm)
81-1003	Regulator, CO, O <sub>2</sub> , Combustible gases
81-1051-01	Regulator, toxic gases (except Cl <sub>2</sub> & HCl)
81-1062	Regulator, Cl <sub>2</sub> , HCl
81-1131	Gas collecting bag, vinyl
81-1132	Clamp, tubing (gas collecting bag)
81-1133	Gas collecting bag, TEDLAR™, reactive gas
81-1134	TEDLAR™ gas collecting bag kit



**Table A-1 Parts List (Continued)**

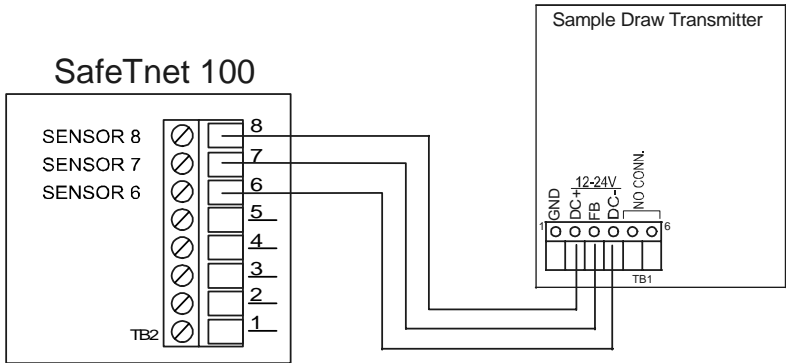
<b>Part No.</b>	<b>Description</b>
81-6414-01	Calibration kit, H <sub>2</sub> S (25 ppm)
81-6414-02	Calibration kit, Cl <sub>2</sub> (5 ppm) <sup>1</sup>
81-6414-03	Calibration kit, HCl (10 ppm)
81-6414-04	Calibration kit, HCN (10 ppm)
81-6414-05	Calibration kit, NH <sub>3</sub> (25 ppm)
81-6414-06	Calibration kit, PH <sub>3</sub> (0.5 ppm) <sup>2</sup>
81-6414-07	Calibration kit, NO (25 ppm)
81-6414-08	Calibration kit, NO <sub>2</sub> (5 ppm)
81-6414-09	Calibration kit, SO <sub>2</sub> (5 ppm)
81-6414-10	Calibration kit, SiH <sub>4</sub> (5 ppm)
81-6415-01	Calibration kit, LEL, Methane
81-6415-02	Calibration kit, LEL, Hydrogen
81-6415-03	Calibration kit, LEL, Propane
81-6415-04	Calibration kit, LEL, Hexane
81-6415-05	Calibration kit, PPM, Methane
81-6415-06	Calibration kit, PPM, Hexane
81-6415-07	Calibration kit, PPM, Hydrogen
81-6415-08	Calibration kit, O <sub>2</sub> (100% N <sub>2</sub> )
81-6415-09	Calibration kit, CO (100 ppm)
82-0005	Screwdriver, adjustment
1. Appropriate also for fluorine (F <sub>2</sub> ) sensors. 2. Appropriate also for arsine (AsH <sub>3</sub> ) and diborane (B <sub>2</sub> H <sub>6</sub> ) sensors.	



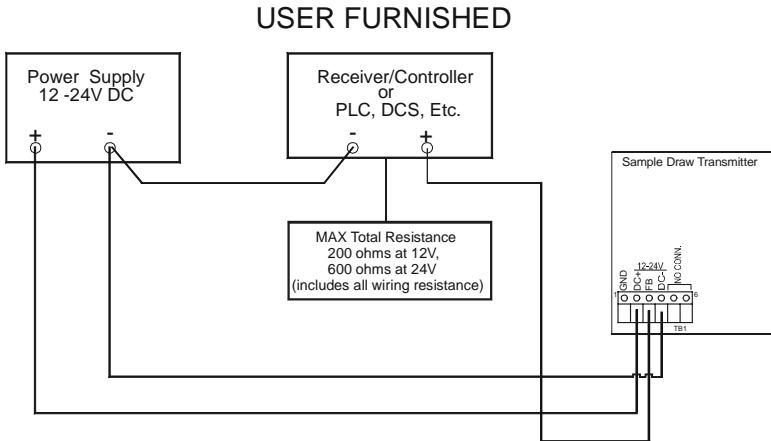
# Appendix B

## EXTERNAL WIRING INSTRUCTIONS

Appendix B shows the Sample-Draw Transmitter external wiring instructions.

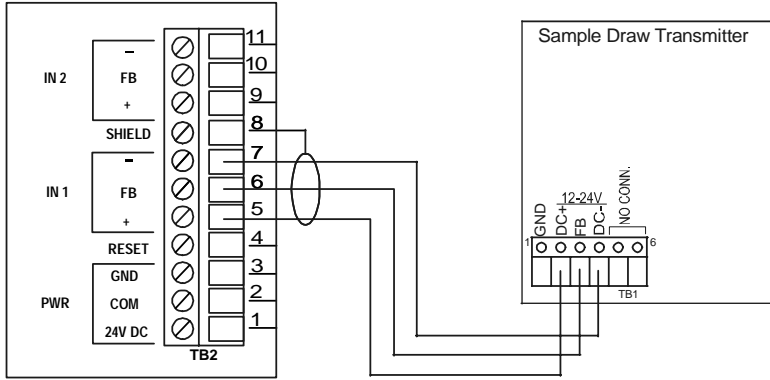


**Figure B-1 SafeTnet 100 External Wiring**

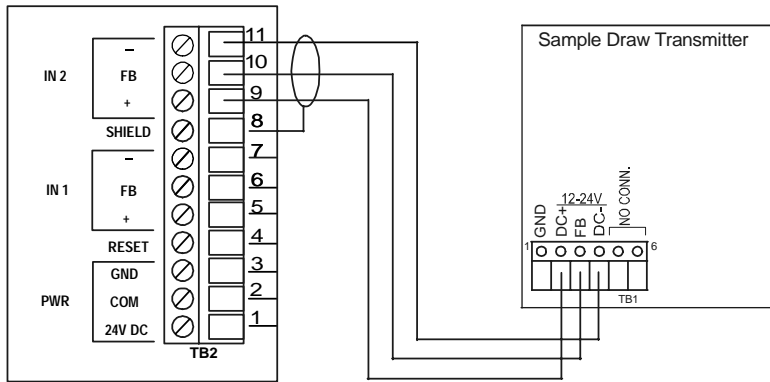


**Figure B-2 User Furnished External Wiring**

STN 2000, 210 and 410

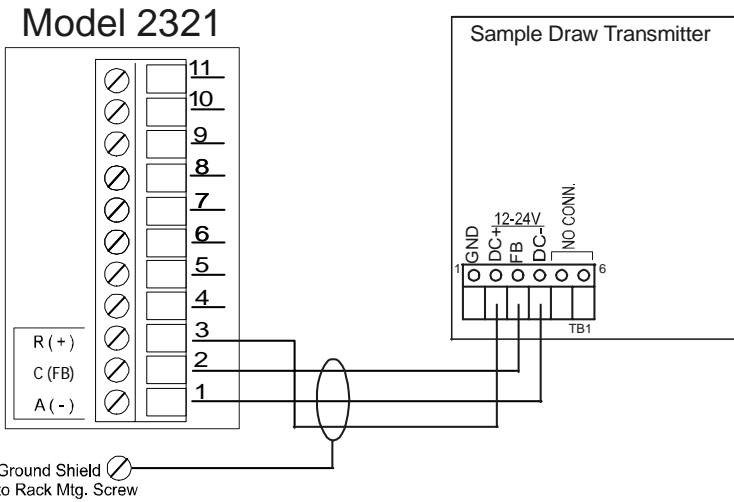


Channel 1

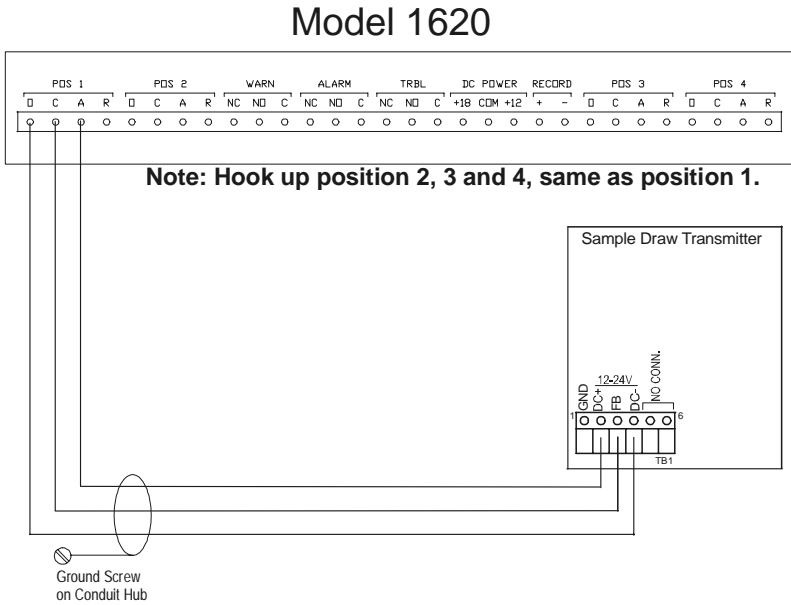


Channel 2 (optional)

Figure B-3 SafeTnet 2000, 210 and 410 External Wiring



**Figure B-4 Model 2321 External Wiring**



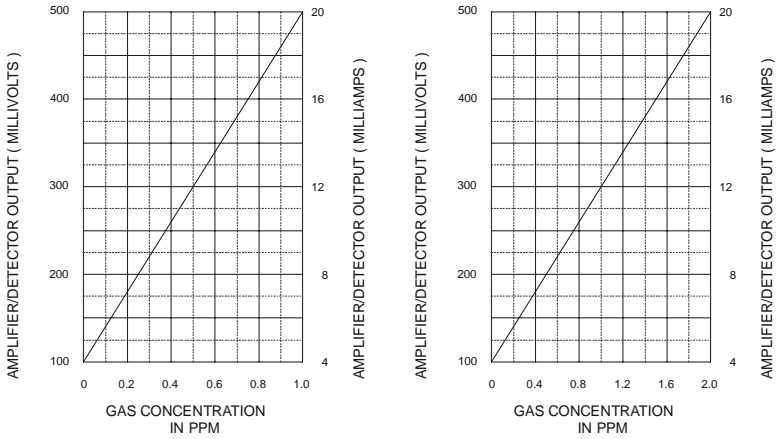
**Figure B-5 Model 1620 External Wiring**



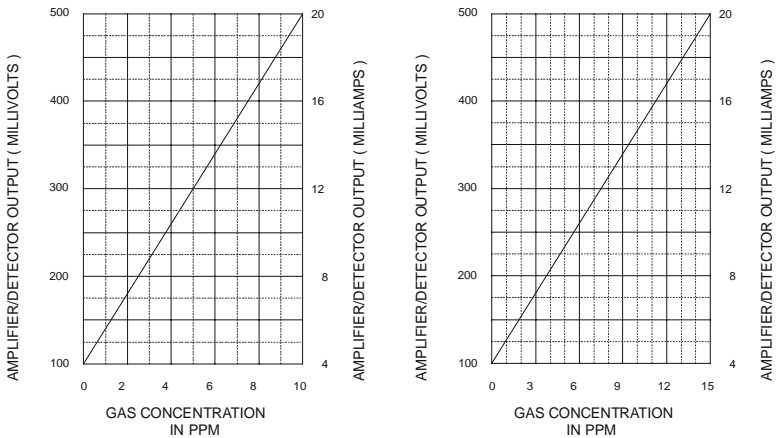
Appendix  
C

# CALIBRATION RESPONSE CHARTS

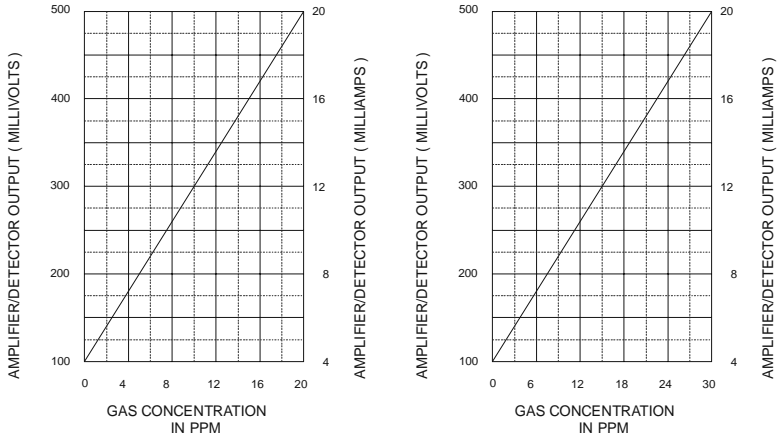
Appendix C shows the Sample-Draw Transmitter calibration response charts.



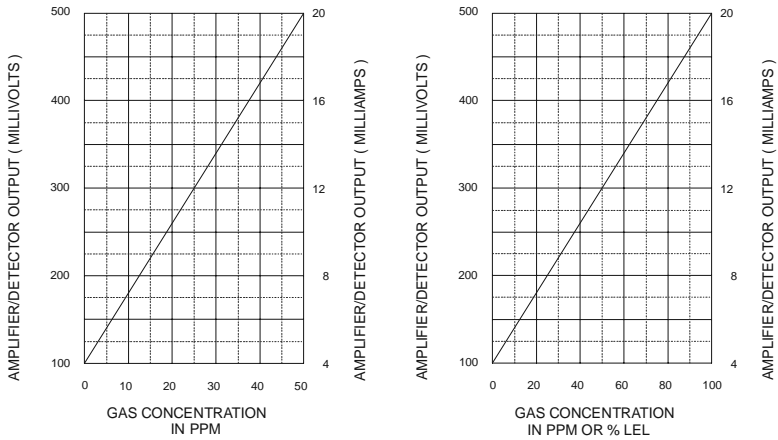
**Figure C-1 Calibration Response Charts for 0-1 and 0-2 ppm**



**Figure C-2 Calibration Response Charts for 0-10 and 0-15 ppm**

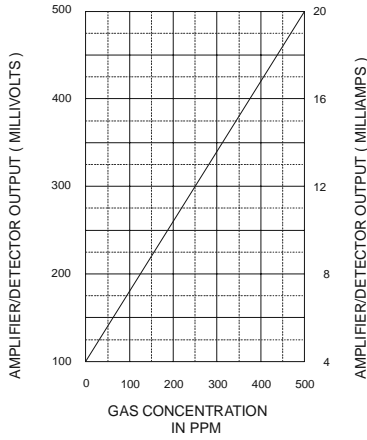


**Figure C-3 Calibration Response Charts for 0-20 and 0-30 ppm**

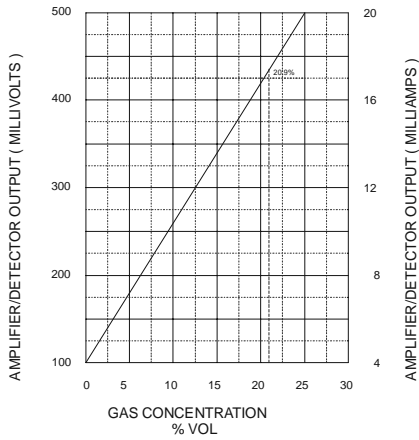


**Figure C-4 Calibration Response Charts for 0-50 ppm and 0-100 ppm or % LEL**





**Figure C-5 Calibration Response Charts for 0-500 ppm**



**Figure C-6 Calibration Response Charts for 0-30% by volume**

