

NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to [www.agilent.com/chem](http://www.agilent.com/chem).



# ***XGS-600 Unit: Tilt Stand Kit***

*FIELD INSTALLATION  
INSTRUCTIONS*

Part No. 699908401

Rev. A

August 2008

# ***XGS-600 Unit: Tilt Stand Kit***

## **Warranty**

Products manufactured by Seller are warranted against defects in materials and workmanship for twelve (12) months from date of shipment thereof to Customer, and Seller's liability under valid warranty claims is limited, at the option of Seller, to repair, replacement, or refund an equitable portion of the purchase price of the Product. Items expendable in normal use are not covered by this warranty. All warranty replacement or repair of parts shall be limited to equipment malfunctions which, in the sole opinion of Seller, are due or traceable to defects in original materials or workmanship. All obligations of Seller under this warranty shall cease in the event of abuse, accident, alteration, misuse, or neglect of the equipment. In-warranty repaired or replaced parts are warranted only for the remaining unexpired portion of the original warranty period applicable to the repaired or replaced parts. After expiration of the applicable warranty period, Customer shall be charged at the then current prices for parts, labor, and transportation.

When products are used with toxic chemicals, or in an atmosphere that is dangerous to the health of humans, or is environmentally unsafe, it will be the responsibility of the Customer to have the product cleaned by an independent agency skilled and approved in handling and cleaning contaminated materials before the product will be accepted by Varian, Inc. for repair and/or replacement.

Reasonable care must be used to avoid hazards. Seller expressly disclaims responsibility for loss or damage caused by use of its Products other than in accordance with proper operating procedures.

Except as stated herein, Seller makes no warranty, express or implied (either in fact or by operation of law), statutory or otherwise; and, except as stated herein, Seller shall have no liability under any warranty, express or implied (either in fact or by operation of law), statutory or otherwise. Statements made by any person, including representatives of Seller, which are inconsistent or in conflict with the terms of this warranty shall not be binding upon Seller unless reduced to writing and approved by an officer of Seller.

### **Warranty Replacement and Adjustment**

All claims under warranty must be made promptly after occurrence of circumstances giving rise thereto, and must be received within the applicable warranty period by Seller or its authorized representative. Such claims should include the Product serial number, the date of shipment, and a full description of the circumstances giving rise to the claim. Before any Products are returned for repair and/or adjustment, written authorization from Seller or its authorized representative for the return and instructions as to how and where these Products should be returned must be obtained. Any Product returned to Seller for examination shall be prepaid via the means of transportation indicated as acceptable by Seller. Seller reserves the right to reject any warranty claim not promptly reported and any warranty claim on any item that has been altered or has been returned by non-acceptable means of transportation. When any Product is returned for examination and inspection, or for any other reason, Customer shall be responsible for all damage resulting from improper packing or handling, and for loss in transit, notwithstanding any defect or non-conformity in the Product. In all cases, Seller has the sole responsibility for determining the cause and nature of failure, and Seller's determination with regard thereto shall be final.

If it is found that Seller's Product has been returned without cause and is still serviceable, Customer will be notified and the Product returned at its expense; in addition, a charge for testing and examination may be made on Products so returned. 3/1/00

*This page intentionally left blank.*

# Varian Field Instruction Sheet

## XGS-600S: Tilt Stand Kit

### Preface

#### **Documentation Standards**

This manual uses the following documentation standards:

#### **NOTE**

Notes contain important information.



#### **CAUTION**

Cautions appear before instructions, which if not followed, could cause damage to the equipment or data loss.



#### **WARNING**

Warnings appear for a particular procedure or practice which, if not followed correctly, could lead to serious injury or death.



#### **Kit Components**

The *Tilt Stand Kit* is comprised of:

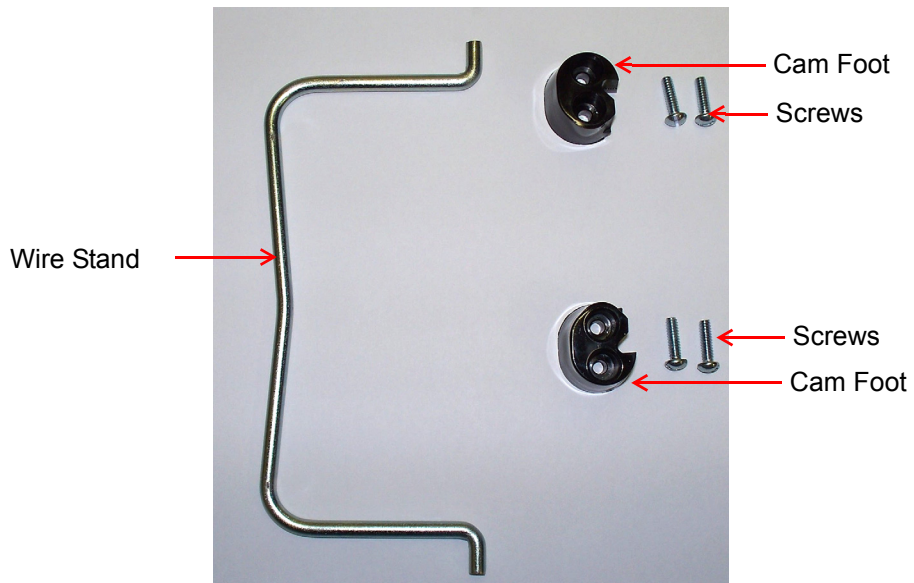
- One (1) universal 1427D5 Hammond Manufacturing Stand kit
- Two (2) 648099034 self-adhesive rubber bumper feet
- One (1) 699908401 instruction sheet

Since the Hammond Stand kit is a generic kit, some of the parts are not used during this install.

The Hammond Stand kit parts used are shown in Figure 1: Hammond Stand Kit: Parts Used:

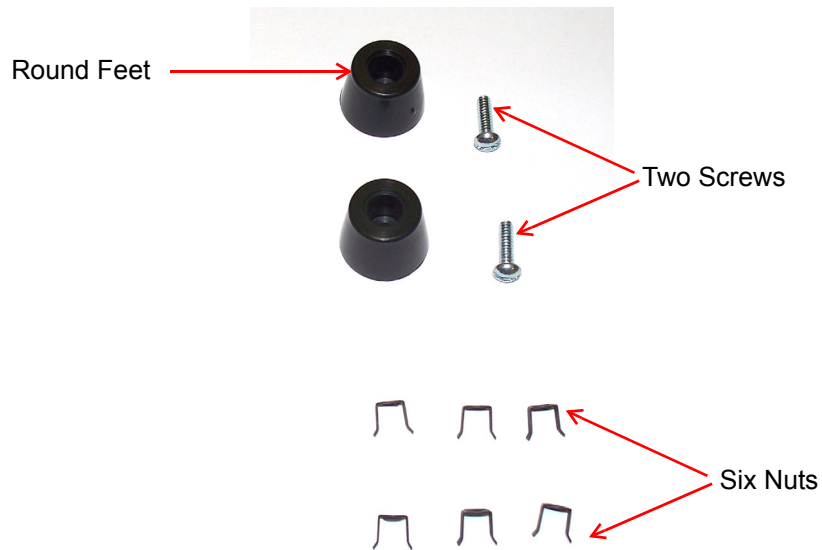
- One (1) stand wire form
- One (1) right and one (1) left cam foot
- Four (4) mounting screws





**Figure 1: Hammond Stand Kit: Parts Used**

Discarded items appear in Figure 2: Hammond Stand Kit: Discarded Parts.



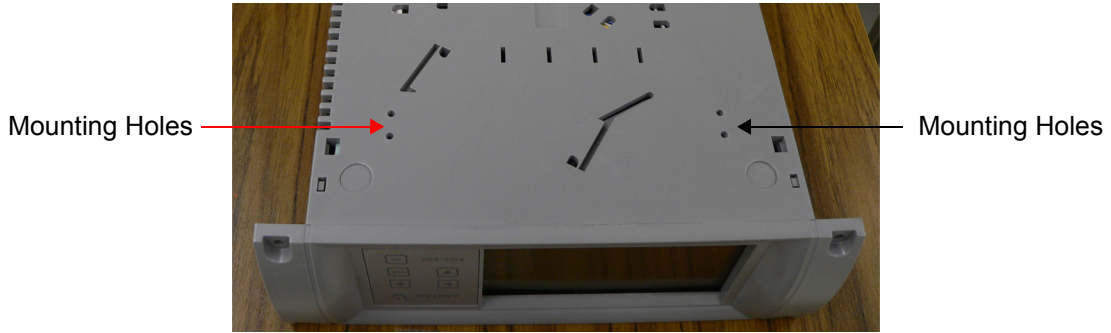
**Figure 2: Hammond Stand Kit: Discarded Parts**

**Equipment Required**

- Flat head screw driver

### Installation Procedure

1. Turn off the power to the unit, wait 30 seconds and disconnect any cables to the unit.
2. Flip the unit on its back with the Front Panel facing front and locate the four mounting holes (Figure 3: Mounting Holes).



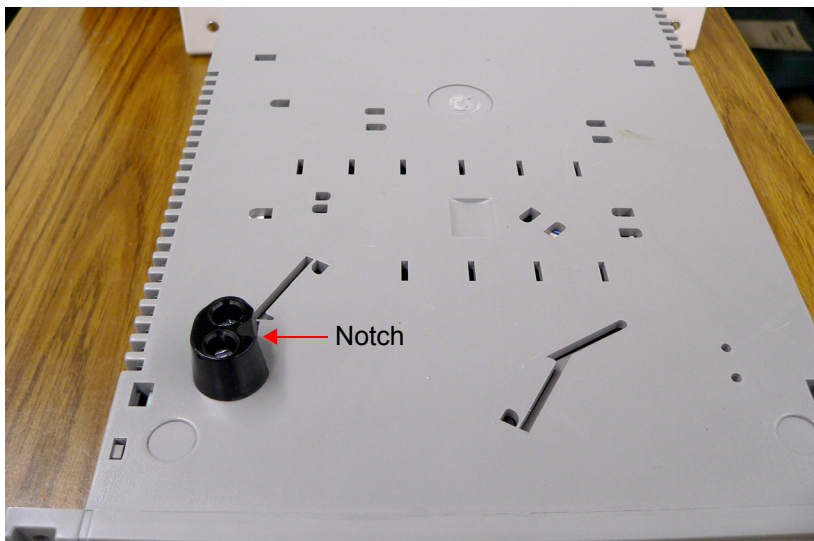
**Figure 3: Mounting Holes**

Figure 4: Cam Feet shows the cam feet. There is one left and one right foot.



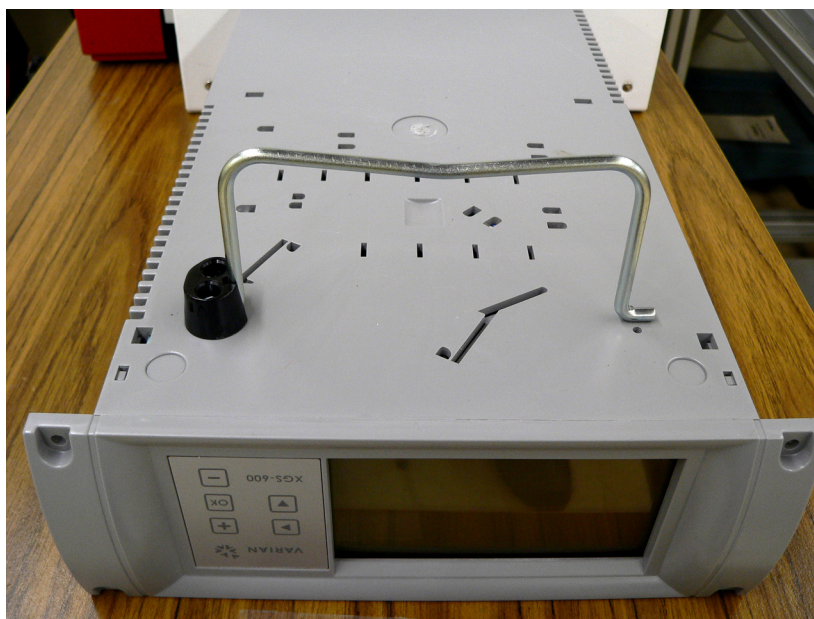
**Figure 4: Cam Feet**

3. Place one of the feet with the notch for the wire stand inward (Figure 5: One Foot Installed) and secure it with two of the screws. Tighten screws snug to the feet. **DO NOT OVER TIGHTEN.**



**Figure 5: One Foot Installed**

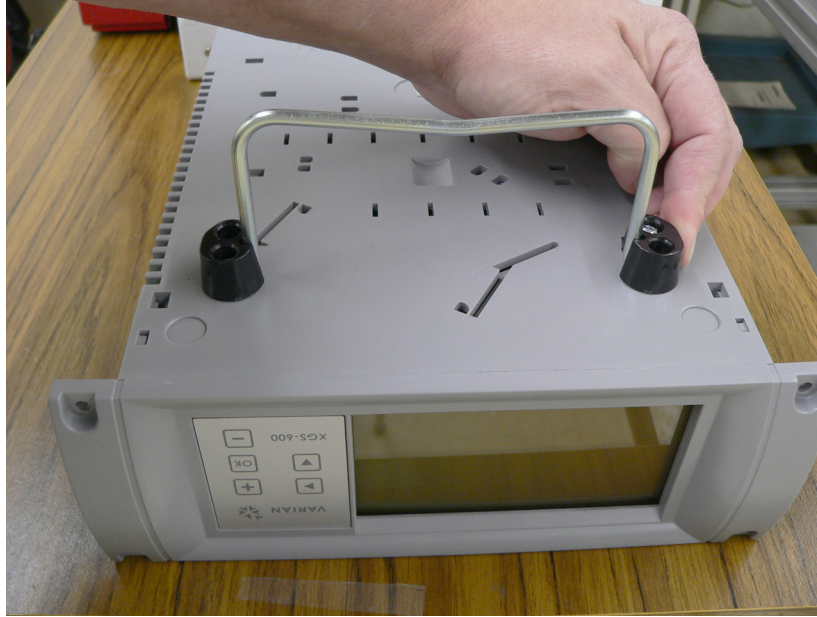
4. Slip the wire stand into the notch of the mounted foot and leave the wire stand in the upright position (Figure 6: Wire Stand with One Foot).



**Figure 6: Wire Stand with One Foot**



- Slip the other end of the wire stand into the second foot and apply pressure from the outside of the foot to align it with the second set of holes (Figure 7: Foot with Pressure Applied).



**Figure 7: Foot with Pressure Applied**

- Insert the two remaining screws and tighten down.
- Add the two self-adhesive rubber bumper feet into the round depressions on the rear of the XGS-600 enclosure, if not already installed (Figure 8: Adhesive Rear Feet Installed).



648099034 self-adhesive rubber bumper foot

**Figure 8: Adhesive Rear Feet Installed**





## Sales and Service Offices

### Canada

#### Central coordination through: Varian, Inc.

121 Hartwell Avenue  
Lexington, MA 02421  
USA  
Tel.: +1 781 861 7200  
Toll-Free: +1 800 882 7426  
Fax: +1 781 860 5437

### China

#### Varian Technologies China, Ltd.

Room 1648  
Central Tower South Wing  
Beijing Junefield Plaza  
No. 10 XuanWuMenWai Street  
Beijing 100052  
P. R. China  
Tel.: +86 (10) 6310 8550  
Toll-Free: 800 820 6556  
Fax: +86 (10) 6310 0141

### France

#### Varian s.a.

7 avenue des Tropiques  
Z.A. de Courtaboeuf - B.P. 12  
91941 Les Ulis cedex  
France  
Tel.: +33 (0) 1 69 86 38 84  
Fax: +33 (0) 1 69 86 29 88

### Benelux

#### Varian Vacuum Technologies

Herculesweg 8  
4338 PL Middelburg  
The Netherlands  
Tel.: +31 118 671570  
Fax: +31 118 671569

### Germany & Austria

#### Varian Deutschland GmbH

Alsfelder Strasse 6  
Postfach 11 14 35  
64289 Darmstadt  
Germany  
Tel.: +49 (0) 6151 703 353  
Fax: +49 (0) 6151 703 302

### India

#### Varian India Pvt. Ltd.

205-A, "A" wing of Galleria  
2nd floor, Hiranandani Gardens  
Powai, Mumbai-400 076  
India  
Tel.: +91 22 2570 8595/8597  
Fax: +91 22 2570 8599  
Mobile: +91 98 679 55969

### Italy

#### Varian, Inc.

via F.lli Varian 54  
10040 Leini, (Torino)  
Italy  
Tel.: +39 011 997 9111  
Toll-Free: 00 800 234 234 00  
Fax: +39 011 997 9350

### Japan

#### Varian Technologies Japan, Ltd.

8th Floor  
Sumitomo Shibaura Building  
4-16-36 Shibaura Minato-ku  
Tokyo 108  
Japan  
Tel.: +81 3 5232 1253  
Toll-Free: 0120 655 040  
Fax: +81 3 5232 1710

### Korea

#### Varian Technologies Korea, Ltd.

Shinsa 2nd Bldg. 2F  
966-5 Daechi-dong  
Kangnam-gu, Seoul  
Korea 135-280  
Tel.: +82 2 3452 2452  
Toll-Free: 080 222 2452  
Fax: +82 2 3452 2451

### Mexico

#### Varian, S. de R.L. de C.V.

Concepcion Beistegui No 109  
Col Del Valle  
C.P. 03100  
Mexico, D.F.  
Tel.: +52 5 523 9465  
Fax: +52 5 523 9472

### Taiwan

#### Varian Technologies Asia, Ltd.

14F-6, No. 77, Hsin Tai Wu Road, Sec. 1  
Hsi chih, Taipei Hsien  
Taiwan, R.O.C.  
Tel.: +886 2 2698 9555  
Toll Free: 0800 051 342  
Fax: +886 2 2698 96782

### UK & Ireland

#### Varian Ltd.

6 Mead Road  
Oxford Industrial Park  
Yarnton, Oxford OX5 1QU  
UK  
Tel.: +44 (0) 1865 291570  
Fax: +44 (0) 1865 291571

### United States

#### Varian, Inc

121 Hartwell Avenue  
Lexington, MA 02421  
USA  
Tel.: +1 781 861 7200  
Toll-Free: +1 800 882 7426  
Fax: +1 781 860 5437

### Other Countries

#### Varian Vacuum Technologies

via F.lli Varian 54  
10040 Leini, (Torino)  
Italy  
Tel: (39) 011 997 9 111  
Fax: (39) 011 997 9 350

### Customer Support and Service:

#### North America

Tel: 1 (800) 882-7426 (toll-free)  
vtl.technical.support@varianinc.com

#### Europe

Tel: 00 (800) 234 234 00 (toll-free)  
vtl.technical.support@varianinc.com

#### Japan

Tel: (81) 3 5232 1253 (dedicated line)  
vtj.technical.support@varianinc.com

#### Korea

Tel (82) 2 3452 2452 (dedicated line)  
vtk.technical.support@varianinc.com

#### Taiwan

Tel: 0 (800) 051 342 (toll-free)  
vtw.technical.support@varianinc.com

### Worldwide Web Site, Catalog and On-line Orders:

www.varianinc.com

Representatives in most countries



# VARIAN