

Guide to Good Laboratory Practice

Varian 400-MR

Pub. No. 01-999338-00, Rev. B1106

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**.



Guide to Good Laboratory Practice

Varian 400-MR

Pub. No. 01-999338-00, Rev. B1106

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**.



Guide to Good Laboratory Practice
Varian 400-MR
Pub. No. 01-999338-00, Rev. B1106

Revision history:
A0906 – Initial release
B1106 – Specification correction

Technical contributors: Iain Green
Technical writes: Mike Miller
Technical editor: Everett Schreiber

Copyright ©2006 by Varian, Inc.
3120 Hansen Way, Palo Alto, California 94304
1-800-356-4437
<http://www.varianinc.com>
All rights reserved. Printed in the United States.

The information in this document has been carefully checked and is believed to be entirely reliable. However, no responsibility is assumed for inaccuracies. Statements in this document are not intended to create any warranty, expressed or implied. Specifications and performance characteristics of the software described in this manual may be changed at any time without notice. Varian reserves the right to make changes in any products herein to improve reliability, function, or design. Varian does not assume any liability arising out of the application or use of any product or circuit described herein; neither does it convey any license under its patent rights nor the rights of others. Inclusion in this document does not imply that any particular feature is standard on the instrument.

Trusted VnmrJ and VnmrJ are trademarks of Varian, Inc. Other product names are trademarks or registered trademarks of their respective holders.

Table of Contents

Chapter 1. Introduction	7
1.1 21 CFR Part 58, Subpart D -- Equipment	7
1.2 Yes/No Choice	8
1.3 Glossary	8
Chapter 2. Design Qualifications	11
2.1 400-MR and Magnet Specifications	11
2.2 Standards Compliance	11
ISO Compliance Certificate	12
CE Conformity	13
2.3 Instrument Selection Criteria	14
Chapter 3. Installation Qualifications	15
3.1 Review the Installation Planning Guide	15
3.2 Site Survey	16
3.3 Review and Confirmation of Shipped Items	16
3.4 Computer Audit	16
3.5 System Installation Checklist	19
3.6 Supercon Shim Values	20
Chapter 4. Operation Qualifications	21
4.1 AutoTest	21
4.2 AutoTest Sample	21
4.3 List of Tests Performed by AutoTest	24
Chapter 5. Performance Qualifications	27
5.1 AutoTest for Maintenance Testing	27
5.2 AutoTest Sample	28
5.3 Additional Performance Qualification Tests	28
Appendix A. 400-MR System Specifications	29
Compact, high performance 2 channel NMR spectrometer	29
High band Linear Amplifier	31
Broadband Linear Amplifier	32
Additional Optional Modules	32
Varian 400 MHz Magnet	32
Appendix B. Corrective Actions	33
Site Survey Corrective Actions	33
Review and Confirmation of Shipped Items Corrective Actions	34
System Installation Corrective Actions Checklist	34
Index	35

Table of Contents

List of Figures

Figure 1. ISO Certificate of Conformance	12
Figure 2. CE Declaration of Conformity	13
Figure 3. Autotest Configuration Panel with <i>All Standard Tests</i> Selected	22
Figure 4. Autotest Test Library Showing the Full List of Available Channel 1 Tests	22
Figure 5. Autotest History Panel Showing the Graphical Output of the MOD1 Test	23
Figure 6. Autotest Test Report Panel Showing the Text of a Completed Report	23
Figure 7. AutoTest Configuration panel showing a minimized selection of tests, potentially suitable for a PQ operation	27

List of Tables

Table 1. Probe Specification List	11
Table 2. Instrumentation Selection Criteria.	14
Table 3. Additional Performance Qualification Tests	28

Chapter 1. Introduction

Sections in this chapter:

- 1.1 “21 CFR Part 58, Subpart D -- Equipment,” this page
- 1.2 “Yes/No Choice,” page 8
- 1.3 “Glossary,” page 8

Good Laboratory Practice (GLP) regulations (see FDA regulation 21 CFR part 58) dictate a system designed to assure the quality and integrity of the safety data attained from "nonclinical laboratory studies that support or are intended to support applications for research or marketing permits for products regulated by the Food and Drug Administration, including food and color additives, animal food additives, human and animal drugs, medical devices for human use, biological products, and electronic products" (Section 58.1 Scope).

Although, for example, certain preliminary pharmacological studies and pharmacokinetic studies are still exempt from FDA regulations, laboratories have extended the reach of GLP to enable greater quality management throughout their processes.

Varian, Inc. is not responsible for compliance to GLP regulations, but the information detailed under the framework of this document, may or may not, depending on any individual customer's internal procedures, be used in conjunction with various SOPs to aid in creating a compliant environment.

1.1 21 CFR Part 58, Subpart D -- Equipment

GLP has several subparts covering the total suite of requirements. Subpart D, Equipment, contains those requirements for which analytical instrument vendors can aid compliance by providing certain capabilities, information, and tools. These areas are generally contained within the 4Q model - Design Qualification, Installation Qualification, Operational Qualification and Performance Qualification.

The definitions used in this manual are as follows (excerpted from "The Development and Application of Guidance on Equipment Qualification of Analytical Instruments", Peter Bedson, Mike Sargent, *Journal of Accreditation and Quality Assurance*, 1996, 1:265-274):

Design Qualification	Defines the functional and operational specifications of the instrument and details the conscious decisions in the selection of the supplier.
Installation Qualification	Establishes that the instrument is received as designed and specified, that it is properly installed in the selected environment and that this environment is suitable for the operation of the instrument.

Operation Qualification	The process of demonstrating that an instrument will function according to a series of predefined operational specifications in the selected environment.
Performance Qualification	The process of demonstrating that an instrument performs according to a specification appropriate for its routine use.

Varian, Inc. has developed this guidance document to:

- Contain important information relating to the 4Q model in one location
- Provide an avenue for aiding the tracking of the various processes that occur during the individual qualification stages

1.2 Yes/No Choice

At several stages in this document, Yes/No choices are to be made. If the choice is **YES**, that criterion has been met. However, if the Customer or Varian Representative decide that a certain criterion has not been met and apply **NO**, then the various appendices at the end of the document allow the tracking of the corrective actions required until all the elements of that criterion have finally been concluded. This approach collects and maintains all of records pertaining to the elements in the document in one place.

1.3 Glossary

The following is a glossary of terms that are used in this manual.

21 CFR Part 11	Code of Federal Regulations, Title 21, Food and Drugs, Part 11 “Electronic Records: Electronic Signatures Final Rule”, Federal Register 62 (54), 13429-12466. www.fda.gov/ora/compliance_ref/part11/Default.htm
21 CFR Part 58	Code of Federal Regulations, Title 21, Food and Drugs, Part 58 “Good Laboratory Practices for Nonclinical Laboratory Studies”, April 1, 2003. www.fda.gov/ora/compliance_ref/glp/default.htm
administrator	The operating system login account that owns the VnmrJ system files. The name of this administrator is selected during installation of VnmrJ for Secure Environments in the User Name: entry field. By default, the name <i>vnmr1</i> is used. The password for this account should be kept secure.
Analytical system	A system that combines instrument, computer, and analytical method.
AutoTest	Automated test software used to demonstrate that a Varian NMR instrument functions according to the operation specifications. AutoTest is designed to demonstrate key console and probe performance criteria, measuring the output using NMR data derived from a single sample. In this way, a large series of tests is automatically completed; data from one system is easily compared to similar systems.
Batch	A specific quantity or lot of a test or control article that has been characterized according to Section 58.105(a). (58.3)

Control article	Any food additive, color additive, drug, biological product, electronic product, medical device for human use, or any article other than a test article, feed, or water that is administered to the test system in the course of a nonclinical laboratory study for the purpose of establishing a basis for comparison with the test article. (58.3)
FDA	United States Food and Drug Administration (www.fda.gov)
FID	An NMR term that stands for Free Induction Decay. A FID is the time-domain data acquired by the NMR spectrometer. The frequency-domain NMR spectrum is then obtained from the FID by Fourier transform.
GAP	Good Auditing Practices
GLP	Good Laboratory Practices
GMP	Good Manufacturing Practices
IQ	Installation Qualification. Applies to new and used equipment that is to be installed. See IQ, OQ, PQ .
IQ, OQ, PQ	Installation Qualification, Operational Qualification, and Performance Qualification. Verification of equipment and instruments used for FDA (and other regulatory bodies) and client audits.
NMR	Nuclear Magnetic Resonance. A scientific technique for determining molecular structure.
Nonclinical laboratory study	In vivo or in vitro experiments in which test articles are studied prospectively in test systems under laboratory conditions to determine their safety. The term does not include studies utilizing human subjects or clinical studies or field trials in animals. The term does not include basic exploratory studies carried out to determine whether a test article has any potential utility or to determine physical or chemical characteristics of a test article. (58.3)
noncontrolled	Data and records that are not controlled, monitored, or audited by a regulatory compliant system, or system administrator. See controlled .
OQ	Operational Qualification. Verification of equipment or instrument performance to prove operation before and after movement. See IQ, OQ, PQ .
Part 11	The Electronic Records; Electronic Signatures section of the Title 21-Food and Drugs regulation.
permissions	Granted in the OS by the system administrator to allow users access to files or resources.
PQ	Performance Qualification or Performance Verification. Operational verification that is performed at least annually for equipment and instruments. See IQ, OQ, PQ .
qualify	In terms of electronic records regulations, a company has the SOPs and software in place so that their records qualify as compliant.

raw data	Any laboratory worksheets, records, memoranda, notes, or exact copies thereof, that are the result of original observations and activities of a nonclinical laboratory study and are necessary for the reconstruction and evaluation of the report of that study. In the event that exact transcripts of raw data have been prepared (e.g., tapes which have been transcribed verbatim, dated, and verified accurate by signature), the exact copy or exact transcript may be substituted for the original source as raw data. Raw data may include photographs, microfilm or microfiche copies, computer printouts, magnetic media, including dictated observations, and recorded data from automated instruments. (58.3)
record	Directories where FID and processed data are saved.
root path	Directory path of a primary directory.
SOP	Standard Operating Procedures. A company or site has SOPs in place for regulatory compliance, among other things.
trusted data	Same as controlled data, which is protected, monitored, and audited under VnmrJ for Secure Environments.
VnmrJ	Data acquisition, processing, and display software for Varian, Inc. NMR spectrometer systems.
XML	The eXtensible Markup Language used for data exchange. VnmrJ uses XML to format data for the Locator database.

Chapter 2. Design Qualifications

Sections in this chapter:

- 2.1 “400-MR and Magnet Specifications,” this page
- 2.2 “Standards Compliance,” page 11
- 2.3 “Instrument Selection Criteria,” page 14

Varian, Inc. is one of the world's leading suppliers of NMR spectrometers and the major provider of research level, high magnetic field functional MRI spectrometers. Innovative design for NMR performance is the trademark of Varian's line of spectrometers. Engineering to the highest standards and using leading-edge technologies has brought to the NMR community spectrometers that support all applications without compromise.

2.1 400-MR and Magnet Specifications

A full suite of specification for the 400-MR is available in [Appendix A, “400-MR System Specifications,”](#) and additional information can be requested from Varian, Inc.

Individual probe specification sheets are available upon request and should be appended to this document and summarized in [Table 1.](#)

Table 1. Probe Specification List

<i>Probe description and P/N</i>	<i>Specification sheet attached Yes/No (see next column)</i>	<i>Reason for 'No' attachment</i>	<i>Customer Signature/Date</i>

2.2 Standards Compliance

Varian's continuing commitment to the highest standards of design, manufacture, and in-house servicing of NMR spectrometers is confirmed by compliance with:

- ISO 9001 (see [Figure 1](#))
- All applicable directives required for CE conformity (see [Figure 2](#))

ISO Compliance Certificate



Figure 1. ISO Certificate of Conformance

CE Conformity


Declaration of Conformity


VARIAN
NMR Systems

Manufacturer's Address: 3120 Hansen Way, Palo Alto, California 94304, USA
European Representative: Regina Shuck
Alsfelder Strasse 6, 64289 Darmstadt, Germany

Hereby declare that the product:

Spectrometer, Models 400-MR and 400-MRH

to which this declaration relates is in conformity with the following standards:

Low-Voltage Directive, 73/23/EEC, applicable Safety Standards:
EN 61010-1:2001

And EMC Directive, 89/336/EEC, applicable EMC Standards:
EN61326:1997/A1:1998/A2:2001 & 50011:1998, Group II, Class A

I, the undersigned, hereby declare that the equipment specified above conforms to the above Directives and Standards.

September 25, 2006
Date


Tom Ingram, Operations Director

Figure 2. CE Declaration of Conformity

2.3 Instrument Selection Criteria

In many cases, instrument selection is based upon criteria related to specific applications, data management, data processing techniques, or specific site criteria, **Table 2.** provides a place to record such criteria. This table can be copied so that additional pages can be appended to this document.

Table 2. Instrumentation Selection Criteria.

<i>Selection Criteria</i>	<i>Description of, name of document or other path that identifies additional criteria. append copies of appropriate pages as necessary.</i>	<i>Customer Name/Signature/Date</i>

Chapter 3. Installation Qualifications

Sections in this chapter:

- 3.1 “Review the Installation Planning Guide,” this page
- 3.2 “Site Survey,” page 16
- 3.3 “Review and Confirmation of Shipped Items,” page 16
- 3.4 “Computer Audit,” page 16
- 3.5 “System Installation Checklist,” page 19
- 3.6 “Supercon Shim Values,” page 20

Various factors pertain to the installation of an instrument in a selected environment.

3.1 Review the Installation Planning Guide

Reviewing the applicable *Pre-Installation Instructions* for the 400-MR will aid determination of site suitability and provide specifications for such requirements as operating temperature, humidity, provision for air and electrical supply.

<i>Pre-Installation Instructions Rev. No.</i>	<i>Reviewed Yes/No</i>	<i>Customer Name/Signature/Date</i>

Pre-installation Instructions for the 400-MR spectrometer is available from Varian Inc. Sales Representatives or the local Sales Office. A list of sales representatives and local sales offices is located at:

http://www.varianinc.com/cgi-bin/nav?products/nmr/contact_nmr/sales_offices&cid=JKQOKLIJFI.

3.2 Site Survey

A site survey can be purchased for the 400-MR if required. The site survey is performed before the installation to check the site suitability specifically for vibrations, temperature, humidity control, and external rf interference.

<i>Site Survey Required Yes/No</i>	<i>If Yes, Performed on Date</i>	<i>Corrective action? Yes/No and description. If Yes, see "Site Survey Corrective Actions," page 33</i>	<i>Varian Representative Name/Signature/Date</i>	<i>Customer Name/Signature/Date</i>

3.3 Review and Confirmation of Shipped Items

All of the sales orders should be reviewed and checked for accuracy.

<i>Sales Order No.</i>	<i>Correct? Yes/No and description. If No see "Review and Confirmation of Shipped Items Corrective Actions," page 34</i>	<i>Varian Representative Name/Signature/Date</i>	<i>Customer Name/Signature/Date</i>

3.4 Computer Audit

The computer audit documents information on each computer (please copy and append this chart for additional computers if more than three computers are required for the installation).

The Varian and customer representative's signatures in these forms certify that the information in these forms is accurate and that all computers to be used to run Varian's MR workstation software, VnmrJ, or to run other software to process data obtained on this spectrometer, have been included in the audit (including those previously registered as part of purchases of other Varian NMR spectrometers).

Information about your site (please print):

Company/University _____
 Address _____

 Principal User _____
 Phone _____ Spectrometer type _____
 Fax _____ Console S/N _____
 Sales Order No. _____ Delivery (month/day) _____

Information on each computer (additional forms are on the back of this page). As required by SOPs, include computers directly attached to the spectrometer, computers (networked or non-networked, on-site or off-site) used to process NMR data using Varian's VnmrJ software, and computers (on-site and off-site) used to process data collected on this spectrometer with software from other vendors.

Information on computer ____ of ____ (e.g., 1 of 3)	
Manufacturer _____	Model no. _____
Computer S/N _____	Purchased from _____
Memory (Gbytes) _____	Screen size (in.) _____
Peripherals: Internal hard disk (Mbytes) _____	
CD-ROM drive model _____	Serial no. _____
Printer model _____	Serial no. _____
Plotter model _____	Serial no. _____
Terminal model _____	Serial no. _____
Other peripheral _____	Serial no. _____
Computer function: NMR host _____	
Workstation running VnmrJ _____	on-site or off-site
Workstation running other NMR software _____	on-site or off-site
Workstation running VnmrJ and other NMR software _____	on-site or off-site
VnmrJ version _____	Operating system _____

The above computer audit was performed during installation of the system.

Varian Representative _____ Date _____

I certify that the information on this form is accurate and that all computers to be used to run VnmrJ software or to run other software to process data obtained on this spectrometer, have been included, as required, in the audit (including those previously registered as part of purchases of other Varian NMR spectrometers).

Customer Representative _____ Date _____

Use these forms for additional computers. If more forms are needed, copy this page. Attach all copies to the Computer Audit.

Information on computer ____ of ____ (e.g., 2 of 3)	
Manufacturer _____	Model no. _____
Computer S/N _____	Purchased from _____
Memory (Gbytes) _____	Screen size (in.) _____
Peripherals: Internal hard disk (Mbytes) _____	
CD-ROM drive model _____	Serial no. _____
Printer model _____	Serial no. _____
Plotter model _____	Serial no. _____
Terminal model _____	Serial no. _____
Other peripheral _____	Serial no. _____
Computer function: NMR host _____	
Workstation running VnmrJ _____	on-site or off-site
Workstation running other NMR software _____	on-site or off-site
Workstation running VnmrJ and other NMR software _____	on-site or off-site
VnmrJ version _____	Operating system _____

Information on computer ____ of ____ (e.g., 3 of 3)	
Manufacturer _____	Model no. _____
Computer S/N _____	Purchased from _____
Memory (Gbytes) _____	Screen size (in.) _____
Peripherals: Internal hard disk (Mbytes) _____	
CD-ROM drive model _____	Serial no. _____
Printer model _____	Serial no. _____
Plotter model _____	Serial no. _____
Terminal model _____	Serial no. _____
Other peripheral _____	Serial no. _____
Computer function: NMR host _____	
Workstation running VnmrJ _____	on-site or off-site
Workstation running other NMR software _____	on-site or off-site
Workstation running VnmrJ and other NMR software _____	on-site or off-site
VnmrJ version _____	Operating system _____

3.5 System Installation Checklist

Fill in the system installation checklist. If corrective action is required (No check box selected), use “System Installation Corrective Actions Checklist,” page 34.

Company/University _____
 Address _____

 Principal User _____
 Phone _____ Spectrometer type _____
 Fax _____ Console S/N _____
 Sales Order No. _____ Magnet S/N _____

Shipment Damage: _____

Preinstallation Preparation:

Line voltage measured (Vac): console _____ accessory _____
 Line pressure: air _____ N₂ _____
 Air conditioning: _____
 Cryogenics (liters): LHe _____ LN _____

Testing and Customer Familiarization:

1. Acceptance tests and computer audit

yes	no	Acceptance tests procedures finished
yes	no	Test results forms completed and signed
yes	no	Computer audit completed and signed

2. System documentation review

yes	no	Software Object Code License Agreement (acceptance of product constitutes acceptance of object code license regardless of whether agreement is signed or not)
yes	no	Varian and OEM manuals
yes	no	Explanation of warranty and where to telephone for information

3. Magnet demonstration

yes	no	Posting requirements for magnetic field warning signs
yes	no	Warning signs posted
yes	no	Cryogenics handling and safety
yes	no	Magnet refilling
yes	no	Flowmeters
yes	no	Homogeneity disturbances

4. Console and probe demonstration

CAUTION! Avoid possible preamp. damage; make sure probe is connected and tuned to resonance.

yes	no	Loading programs
yes	no	Experiment setup
yes	no	Basic operation to obtain typical spectra

Customer Representative: _____ **Date** _____

Varian Representative: _____ **Date** _____

3.6 Supercon Shim Values

Fill in the following information:

Magnet Frequency and Serial Number:

Magnet Frequency _____

Serial Number _____

Measurement in:

Amps _____

Measurement	1. Date:	2. Date:	3. Date:
Z0			
Z1			
Z2			
Z3			
Z4			
X			
Y			
ZX			
ZY			
XY			
X2-Y2			
Drift			
Spacers			
Main Field Current			
Customer Signature:			
Varian Representative Signature:			

Chapter 4. Operation Qualifications

Sections in this chapter:

- 4.1 “AutoTest,” this page
- 4.2 “AutoTest Sample,” page 21
- 4.3 “List of Tests Performed by AutoTest,” page 24

An NMR spectrometer is a complex instrument, available with a variety of optional items and capable of performing a variety of different experiments. AutoTest (Figure 3) is used to demonstrate that an instrument functions according to a series of operation specifications.

The customer’s representative or a Varian representative can run the Operations Qualifications test. The customer SOP dictates the operator. Operations qualification testing by a Varian Inc. representative is not included in standard installation, warranty, or service contracts. A specific quotation, purchase order, and sales order are required. The sales order number can be recorded in this chapter. Sales Order _____.

4.1 AutoTest

The AutoTest software is designed to demonstrate key console and probe performance criteria, measuring the output using NMR data derived from a single sample. In this way, a large series of tests is automatically completed; data from one system is easily compared to similar systems.

Autotest has a variety of straightforward screens that allow the basic configuration to be set up (operator name, console type etc.) (Figure 3). A library of tests is provided (Figure 4). A history log graphically represents the data over time (Figure 5) and a report page provides a textual output for each test (Figure 6). The report can be printed providing a hard copy record.

Some tests run by AutoTest have specifications attached; some use visual data inspection. For quick reference, all tests are listed in 4.3 “List of Tests Performed by AutoTest,” page 24. System specifications are listed in the *Console Acceptance Tests and Specifications* manual.

For an OQ process, it is recommended that all applicable tests for the instrument are run. This same suite can be run after service calls, probe changes, or other major operations, according to the organization’s SOPs.

4.2 AutoTest Sample

AutoTest uses 0.1% ^{13}C methanol, 0.1% acetonitrile in 1% H_2O /99.8% D_2O .

The sample is doped with gadolinium chloride at a concentration of 0.30 mg/ml, which produces a ^1H T1 relaxation time of about 50 to 75 ms. The sample supplied with every spectrometer, has an expiration date of one year and is available for order using Varian P/N 01-901855-06, 5-mm sample.

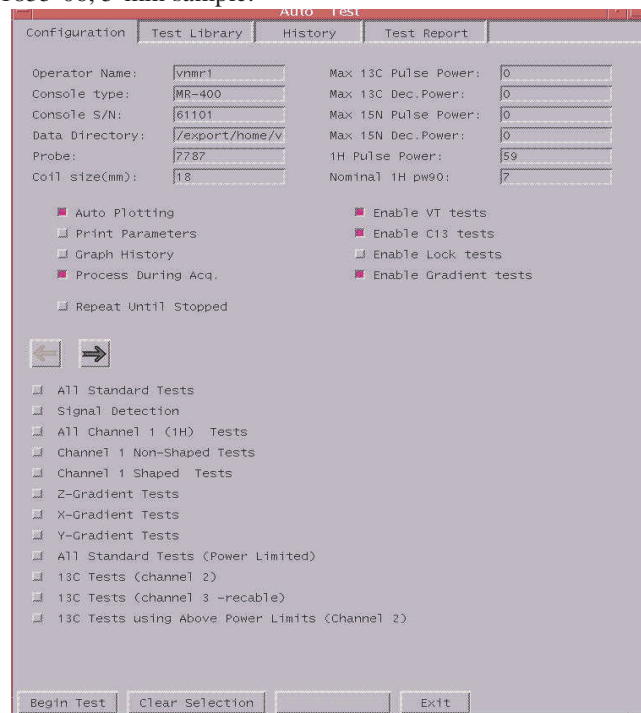


Figure 3. Autotest Configuration Panel with All Standard Tests Selected



Figure 4. Autotest Test Library Showing the Full List of Available Channel 1 Tests

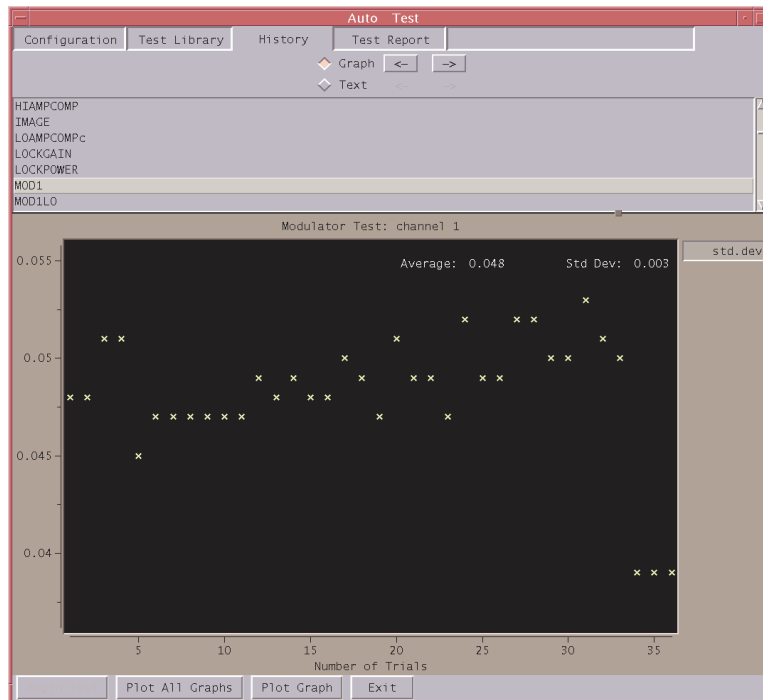


Figure 5. Autotest History Panel Showing the Graphical Output of the MOD1 Test

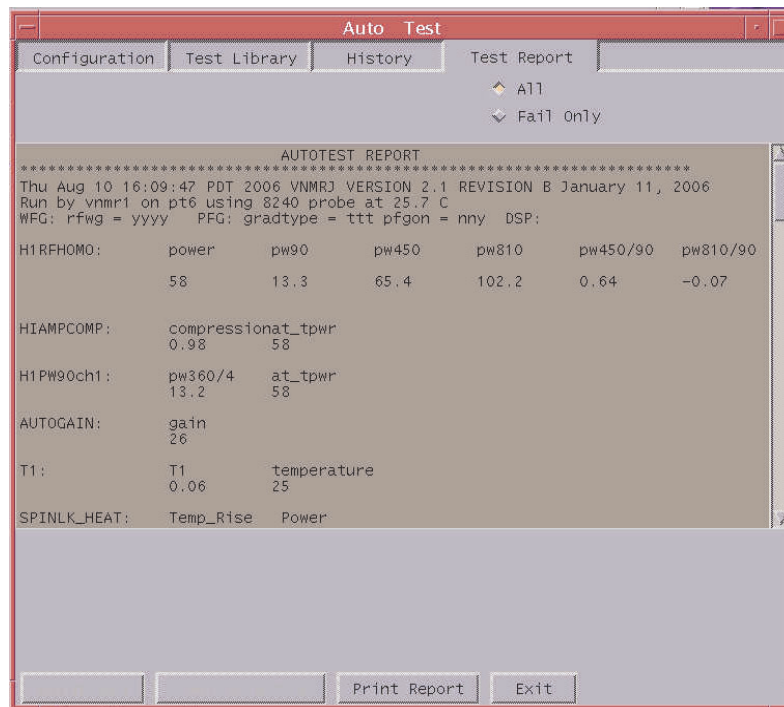


Figure 6. Autotest Test Report Panel Showing the Text of a Completed Report

4.3 List of Tests Performed by AutoTest

For more detail, AutoTest is fully documented in the *AutoTest for VnmrJ*, Pub. No. 01-999247-00.

- RF homogeneity test: ^1H
- RF homogeneity test: ^{13}C
- Water resonance frequency
- T_1
- ^1H PW90 determination
- ^{13}C PW90 determination
- High-band amplifier compression
- Low-band amplifier compression
- ^{13}C phase modulation decoupling profile
 - GARP-1 modulation
 - WALTZ-16 modulation
 - XY-32 modulation
 - MLEV-16 modulation
- ^{13}C adiabatic decoupling profiles
 - STUD modulation
 - WURST modulation
- Receiver gain tests
- Small-angle phase test
- 90° pulse stability (see Note 1)
- Phase cycle cancellation – 2 scans (see Note 2)
- Phase cycle cancellation – 4 scans (see Note 2)
- 30° amplitude stability (see Note 1)
- Pulse turnon time
- Attenuator linearity tests
 - 400 MHz systems and above
- Attenuator linearity test (reduced power)
 - 400 MHz systems and above
- Modulator linearity ($t_{\text{pwr}} = 40$)
- Modulator linearity ($t_{\text{pwr}} = -16$)
- Phase stability error
- Gradient level for 10 G/cm along Z
- X-gradient echo stability (10 G/cm)
- Z-gradient recovery stability (10 G/cm)
- Z-axis signal recovery (rect)
- Z-axis signal recovery (sine)
- Gradient phase cycle cancellation – 4 scans (see note 2)
- Gaussian 90° stability
- Gaussian 13° phase error
- GaussianSLP 13° phase error
- Temperature rise in spinlock test
- Lock power test: corr. coef.
- Lock gain test: corr. coef.

Chapter 5. Performance Qualifications

Sections in this chapter:

- 5.1 “AutoTest for Maintenance Testing,” this page
- 5.2 “AutoTest Sample,” page 28
- 5.3 “Additional Performance Qualification Tests,” page 28

Autotest can also be used for routine maintenance testing. The customer’s representative or the Varian representative can run the performance qualifications test. The customer SOP dictates the operator. Performance qualification by a Varian Inc. representative is not included in standard installation, warranty, or service contracts. A specific quotation, purchase order, and sales order are required. The sales order number can be recorded in this chapter.

5.1 AutoTest for Maintenance Testing

The AutoTest software is designed to demonstrate key console and probe performance criteria. However, for PQ, rather than run a full suite, individual tests should be chosen to best represent some of the more important criteria for the organization (see [Figure 7](#)). These tests can be run as required by the organization SOPs. The value of each selected test can be compared against earlier tests over time and against the values measured for the same test run during an OQ operation.

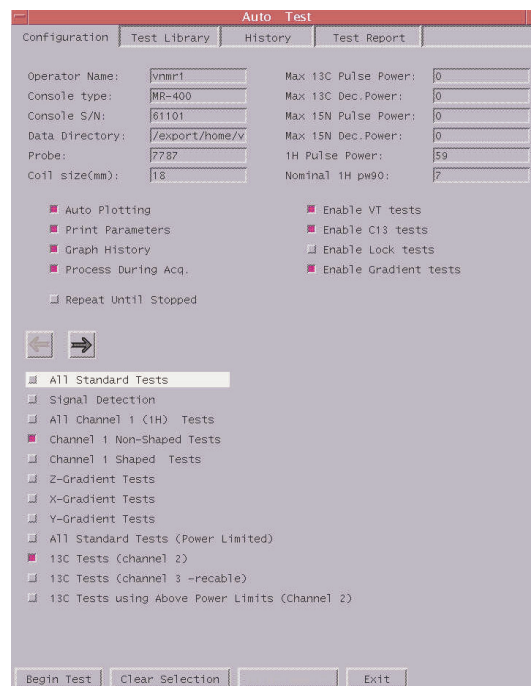


Figure 7. AutoTest Configuration panel showing a minimized selection of tests, potentially suitable for a PQ operation

5.2 AutoTest Sample

AutoTest uses 0.1% ¹³C methanol, 0.1% ¹⁵N acetonitrile in 1% H₂O/99.8% D₂O.

The sample is doped with gadolinium chloride at a concentration of 0.30 mg/ml, which produces a ¹H T1 relaxation time of about 50 to 75 ms. The sample supplied with every spectrometer, has an expiration date of one year and is available for order using Varian P/N 01-901855-06, 5-mm sample.

5.3 Additional Performance Qualification Tests

Additional tests using a particular sample with a documented result can be applied to an SOP. These tests can be documented in [Table 3](#).

Note: A Varian Inc. quotation for performance qualification services may include sample preparation and validation.:

Table 3. Additional Performance Qualification Tests

<i>Test</i>	<i>Sample</i>	<i>Result</i>	<i>Customer Representative Signature</i>	<i>Varian Sales Order (if applicable)</i>

Appendix A. **400-MR System Specifications**

Sections in this chapter:

- “Compact, high performance 2 channel NMR spectrometer,” this page
- “High band Linear Amplifier,” page 31
- “Broadband Linear Amplifier,” page 32
- “Additional Optional Modules ,” page 32
- “Varian 400 MHz Magnet,” page 32

Compact, high performance 2 channel NMR spectrometer

Frequency Sources

Frequency ranges: 18.5 - 245 MHz & 370 - 425 MHz

Frequency resolution of 0.1 Hz (not including phase continuous frequency shifting)

High Band Linear Amplifier

For ¹H & ¹⁹F, 50 watts pulsed power

Lowband, Linear Amplifier

For ¹⁰⁹Ag - ³¹P, 300 watts pulsed power, 18.5-245 MHz

RF Controller

Power PC & FPGA

64MB RAM & 32MB Flash

> 30 MB controller memory per RF channel for waveform definition:

Phase Shift Generation

50 ns minimum event time and 12.5 ns timing resolution

50 ns settling time

0.044° phase resolution

0.04° phase accuracy

Coarse amplitude control

79 dB in 1 dB steps

Fine amplitude control

60 dB in 4096 linear steps

50 ns minimum event time and 12.5ns timing resolution

50 ns settling time

Phase Coherent and Phase Continuous Frequency Shifting

50 ns minimum event

50 ns settling time

Infinitesimally fine resolution for infinitely long pulse

PFG Controller

Power PC & FPGA

64MB RAM & 32MB Flash

> 30 MB controller memory for PFG channel for waveform definition:

Gradient Shaping

12.5 ns timing resolution

Gradient amplitude control in $\pm 32,768$ linear steps

Minimum gradient pulse length 2.4 ms

Preamplifiers

High band (^{19}F and ^1H) GaAs FET preamplifier.

Active (linear) transmit/receive switch.

1 μs recovery, 1 μs switching time.

Lowband (^{109}Ag - ^{31}P) preamplifier

Active (linear) transmit/receive switch.

1 μs recovery, 1.5 μs switching time.

DDR Controller

Power PC & FPGA

64MB RAM & 32MB Flash

DirectDigital™ Receiver

Digitizer/ Max over sampling rate, 14 bit at 80 MHz, (eff. 20 bit, 10 KHz)

Maximum spectral width, 5 MHz

Maximum spectral width with no quadrature images, 5 MHz

On-the-fly Digital filtering

Lock Controller

Power PC & FPGA

64MB RAM & 32MB Flash

Adaptive 2H Lock

2H frequency +/- 1 MHz

Quadrature detection with simultaneous sampling during lock capture.

Continuous field correction for short-term field fluctuations.

Synchronization with External Events

3 TTL output line per RF channel for user programming

1 additional Ethernet connection

Broadband (^{109}Ag - ^{31}P) preamplifier

Active (linear) transmit/receive switch.

1 μs recovery, 1.5 μs switching time.

High dynamic range, low noise and fast recovery.

Note: Nuclei using frequency ranges 61 to 72 MHz and 130 to 143 MHz are not supported and cannot be viewed.

Note: Deuterium observe uses the lock channel.

High band Linear Amplifier

Nominal power	50 W
Nuclei	^1H , ^{19}F
Max pulsed power (W)	50
Max CW power (W)	15
Full power duty cycle (%)	10
Full power pulse length (ms)	250
Rise time (ns)	200
Droop (%)	5
Unblank/Blank time (μs)	1

Broadband Linear Amplifier

Nominal power	300 W
Frequency (MHz)	185 - 245
Nuclei	^{109}Ag - ^{31}P
Max pulsed power (W)	300 (full range)
Max CW power (W)	30
Full power duty cycle (%)	10
Full power pulse length (ms)	100
Rise time (ns)	200
Droop (%)	5
Unblank/Blank time (μs)	1

Additional Optional Modules

Hardware Accessories

Low Temperature Sample Management

Software Accessories

NMRAnalyst	Analyzes data from the carbon-carbon connectivity experiment, revealing carbon connectivities at previously unattainable levels of sensitivity
DOSY	Analyzes data from Diffusion Ordered Spectroscopy experiments

Varian 400 MHz Magnet

Drift	< 8 Hz/hr
Superconducting shim coils	Z1, Z2, Z3, X, Y, ZX, ZY, XY, X2-Y2
Radial 5G Stray Field from Magnet Center	55 cm/21.7 inches
Axial 5G Stray Field from Magnet Center	100 cm/39.4 inches
Axial 5 G height above floor	208 cm/81.9 inches
Axial 5 G depth below floor	N/A
System weight, operational 650 kg	650 kg/1433 lbs
Min. ceiling height	294 cm/115.7 inches
Min. ceiling height (optional transfer)	260.5 cm/102.6 inches
Liquid helium refill volume	95 L
Liquid helium hold time	< 270 days (9 months)
Liquid nitrogen refill volume	51 L
Liquid nitrogen hold time	< 14 days

Appendix B. Corrective Actions

- “Site Survey Corrective Actions,” this page
- “Review and Confirmation of Shipped Items Corrective Actions,” page 34
- “System Installation Corrective Actions Checklist,” page 34

Site Survey Corrective Actions

<i>Action Required (from description)</i>	<i>Completed Yes/No</i>	<i>Customer Name/Signature/Date</i>

Review and Confirmation of Shipped Items Corrective Actions

<i>Action Required (from description)</i>	<i>Completed Yes/No</i>	<i>Customer Name/Signature/Date</i>

System Installation Corrective Actions Checklist

<i>Action Required (from description)</i>	<i>Completed Yes/No</i>	<i>Customer Name/Signature/Date</i>

Numerics

21 CFR Part 11, 8
 21 CFR Part 58, 8
 4Q model, 7

A

administrator, 8
 AutoTest, 8, 21
 Autotest, 21
 AutoTest sample, 28

B

Batch, 8

C

CE conformity, 11
 computer audit, 16
 Control article, 9

D

Design Qualification, 7

F

FDA, 9
 FID, 9

G

GAP, 9
 glossary, 8
 GLP, 9
 GLP regulations, compliance, 7
 GMP, 9

I

Installation Qualification, 7
 IQ, 9
 ISO 9001, 11

N

NMR, 9
 Nonclinical laboratory study, 9
 noncontrolled, 9

O

Operation Qualification, 8
 OQ, 9
 OQ process, 21

P

Part 11, 9
 Performance Qualification, 8
 Performance qualification, 27
 permissions, 9
 PQ, 9
 preinstallation checklist, 19
 Pre-installation Instructions, 15

Q

qualify, 9

R

Raw data, 10
 record, 10
 root path, 10

S

shipment damage, 19
 site survey, 16
 SOP, 10
 Subpart D, Equipment, 7
 system administrator, 8
 system installation checklist, 19

T

trusted data, 10

V

VnmrJ, 10

W

web address, 15

X

XML, 10

Y

Yes/No choices, 8