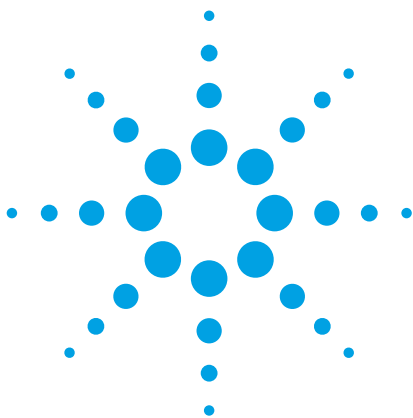




Agilent VnmrJ Software

User Rights in a multi-user environment

Technical Overview

Advantage Statement

The User Rights software tool allows an administrator to configure Agilent VnmrJ software with respect to each user, matching the varied needs within your organization. The features are extremely valuable for a multi-user system that is shared by several people or groups of people. For all users, menus and experimental protocols can be streamlined to display only the functions required or appropriate for the user's tasks, simplifying operation.

Rights Management

A pop-up window available to the VnmrJ administrator, (Figure 1) simplifies the task of creating profiles for either individuals or groups. Rights, tools, and protocols to be included in a profile are selected using checkbox and saved. One or more of up to 23 available rights and tools, and an unlimited number of protocols can be granted as appropriate to an infinite number of different profiles.



Agilent Technologies

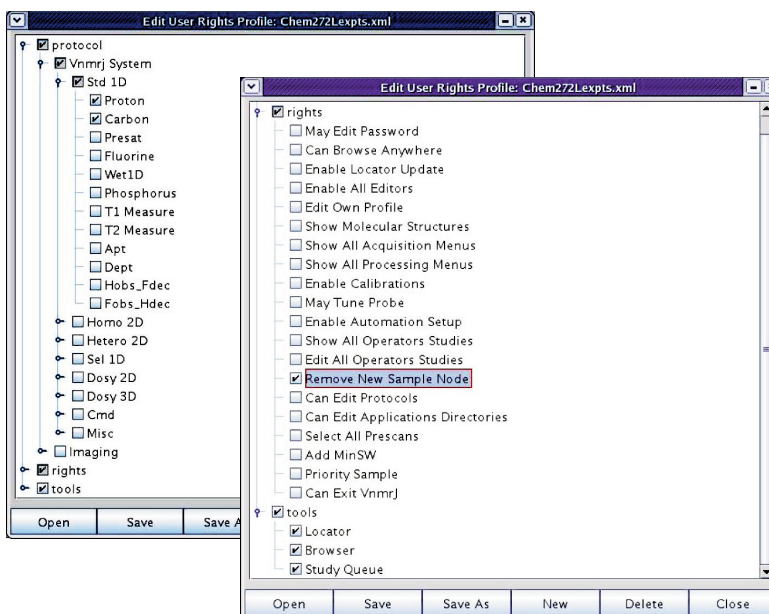


Figure 1. In this example, the profile created for Chemistry 272 laboratory (a first year organic chemistry class) allows for only a Proton and Carbon protocol to be available, as well as removal of a sample node, and access to the Locator, Browser and Study Queue tools.

Each user is assigned a single profile, as shown in Figure 2. Privileges, such as access to the command line, level of panels available, and the ability to send e-mail, are provided in the same pop-up where profiles are assigned.

Modify Operators		Delete Operator			
Operator	Email	Panel Level	Full Name	Show Command Line	Profile Name
CHEM272L		9		no	Chem272Lexpts
CHEM273L		9		no	Chem272Lexpts
CHEM333L		20		no	Chem333Lexpts
CHEM423		20		no	Chem423expts
CHEM443		20		no	Chem443expts
DavisG	davisg@chemist...	30	G. Davis	yes	AllLiquids
JohnsonM	johnson@chemi...	30	M. Johnson	yes	AllLiquids
SmithJ	smithj@chemistr...	30	J. Smith	yes	AllLiquids

Figure 2. The Chemistry 272 laboratory user, CHEM272L, has been assigned the Chem272Lexpts profile, with access to a subset of VnmrJ panels, but not to the command line or e-mail function. Research advisors Davis, Johnson, and Smith are given more access to the functions in VnmrJ.

Streamlined Menus and Experimental Protocols

Users access the VnmrJ interface through a login screen, which can be configured with unique passwords to guard the user's session and prevent its use by others. Once logged in, the VnmrJ interface is displayed with appropriate privileges. The three interfaces in Figures 3–5 reflect different profiles and privileges in a university's multi-user environment. The CHEM272L and CHEM273L users have functions specific to those course requirements, while the CHEM333L, CHEM423 and CHEM443 users have access to more advanced protocols, such as the gradient HSQC experiment. The research advisors have access to even more protocols and functions, such as the Hadamard HSQC experiment and pulse sequence panels.

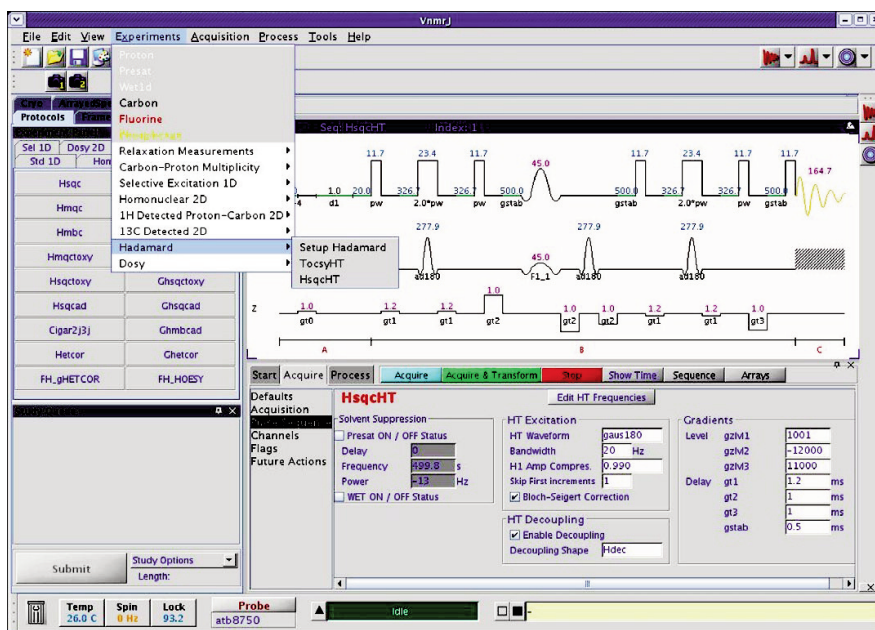


Figure 5. The research advisors have been given access to all liquid protocols, the entire set of menus (File, Edit, View, Experiments, Acquisition, Process, Tools, and Help), all panels and the command line to assist in advanced research.

www.agilent.com

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2011

Printed in the USA

March 31, 2011

SI-1661



Agilent Technologies