



# Agilent 2100 Expert Security Pack

## Backup and Migration of the Secured Area

### Introduction

With the release of 2100 Expert Software B.02.09 and higher versions, the upgrade and migration process of the *Secured Area* has slightly changed compared to previous software versions. This document provides an overview on how to perform a successful software upgrade to a higher version. In addition, detailed information are provided for backup and migration of the *Secured Area* content on the same PC or between different PCs.

This includes the migration of:

- Existing Methods and Workflows
- Existing Data
- Security Pack users and their appropriate roles

Please note that there are differences for the backup process comparing 2100 Expert Version B.02.08 and B.02.09 (or higher).

#### NOTE

After the migration is completed Agilent recommends performing an IQ-service. Performing the *Installation Verification* from the “Validation” context can be an alternative approach, if this fulfills your internal software update policy. Please be aware and consider before performing the *Secured Area* migration.

### About the tool

The “ExtractUsersAndRoles.exe” is an external tool designed to assist the migration of the *Secured Area* for 2100 Expert version B.02.08 as shown below.

	B.02.08 → B.02.08	B.02.08 → B.02.09	B.02.09 → B.02.09	B.02.09 → B.02.10
Migration on same PC	-	X	-	-
Migration between PCs	X	X	-	-



## 2100 Expert B.02.08 - Successful backup of the Secured Area and uninstallation

### Initial preparation

It is highly recommended to create a general backup of your system and data before proceeding.

Before the update or migration can start, you need to perform the following tasks while logged on to Windows as an Administrator:

- Archive your existing method and data files, using the feature within the 2100 Expert Software (restricted to users with either the Administrator or Backup operator roles).
- Clarify how the current generic “2100System” account is setup (local or network/domain). It is important to note the user name if you are working within a company network/domain.
- Uninstall only the 2100 Expert Software from the current PC using the standard operating system process depending on your Microsoft Windows version.
- Using an Administrator user account, copy the “ExtractUsersAndRoles.exe” to the <\System> directory of the *Secured Area*.

### Creating the *Secured Area* backup directory

- 1 Execute the “ExtractUsersAndRoles.exe” tool while logged on to Windows as an Administrator.
- 2 The “UsersAndRoles.xml” file will be created in the <\System> directory of the *Secured Area*. (This will fail if you did not uninstall the 2100 Expert Software).
- 3 Uninstall the 2100 Expert Secured Area from the current PC using the standard Windows process.
- 4 The remaining *Secured Area* directory can now be used as your backup directory. Copy the complete <\Secured Area> directory to the desktop, or an external backup drive.

For Windows 7 the *Secured Area* can typically be found at:

[C:\Program files \(x86\)\Agilent\2100 bioanalyzer\2100 expert\Secured Area](C:\Program files (x86)\Agilent\2100 bioanalyzer\2100 expert\Secured Area).



## 2100 Expert B.02.09 or higher - Backup of the Secured Area during uninstallation

It is highly recommended to create a general backup of your system and data before proceeding.

Before the migration can start, you need to perform the following tasks while logged on to Windows as an Administrator:

- Clarify how the current generic “2100System” account is set-up (local or network/domain). It is important to note the user name if you are working within a company network/domain.
- Uninstall only the 2100 Expert Software from the current PC using the standard operating system process depending on your Microsoft Windows version.
- A backup will be created automatically in the root directory of your previous installation. The default path would be “C:\SecuredArea.bak”.

The backup folder contains all information required for the migration to a different PC. You can copy it to an external hard drive to use it during upcoming installations. The external tool is not required for version B.02.09 or higher.

### NOTE

If you want to install 2100 Expert with Security Pack on a new PC from an existing backup without uninstallation on your current PC, make sure that your backup contains the “Users & Roles” information that was generated with the external tool for the previous software revision B.02.08. In case you have any doubts, contact your Agilent Support Team for clarification upfront.

---



## Migrating on the same PC

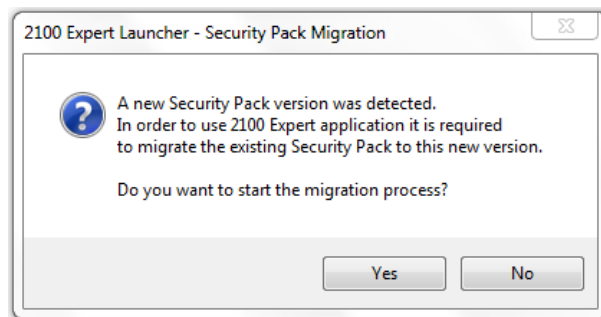
The following section describes the standard *Secured Area* migration on the same PC (in case of a software upgrade).

### **Important:**

If you want to **upgrade from 2100 Expert Software B.02.08 to B.02.09 or higher**, you need to use the external tool and completely uninstall the software as described in the section “*2100 Expert B.02.08 - Successful backup of the Secured Area and uninstallation*”. Then follow the instructions for migration between PCs in the next chapter.

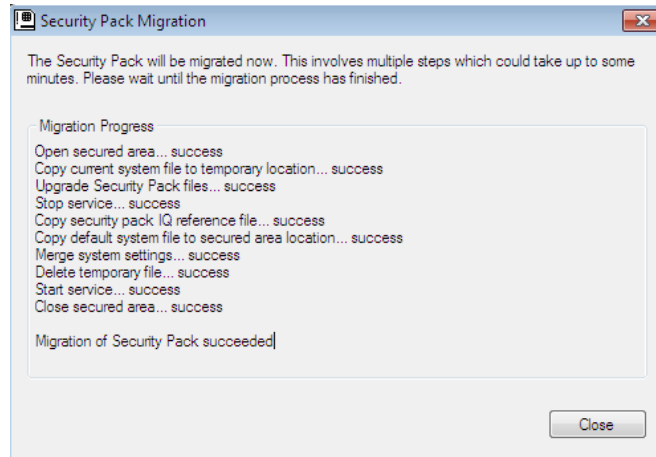
When installing the same 2100 Expert Software version on the same PC, the external tool is ***not*** required. This also applies to an upgrade from B.02.08 to B.02.08 (SR1, SR2 or SR3).

- 1 Install the 2100 Expert Software and follow the on-screen instructions.
- 2 If upgrading to a higher version of the 2100 Expert Software, the launcher recognizes the upgrade. Specific system settings as well as files need to be migrated from the old version. Click “Yes” to start the migration process (“No” would cancel the software launch).





- 3 A window will automatically open and show the progress of the migration.

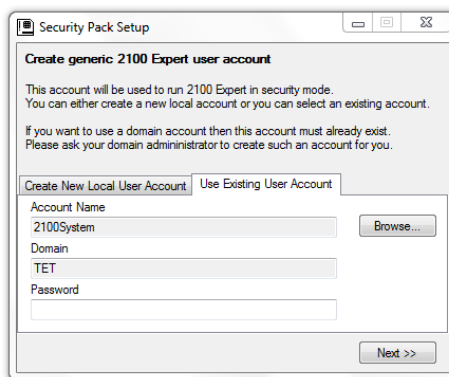


- 4 Upon successful migration of the *Secured Area*, click “Close” to launch the 2100 Expert Software.

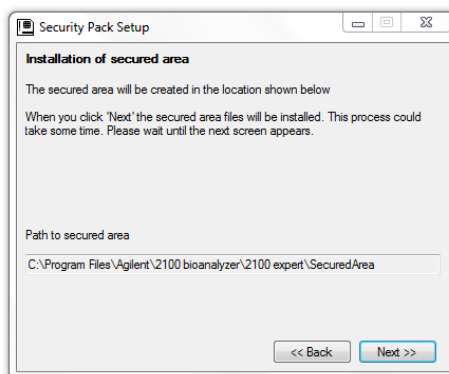
## Migration between different PCs

For the migration process between PCs, you first need to create a backup of your *Secured Area* as described in the previous chapters for 2100 Expert B.02.08 and B.02.09 or higher.

- 1 Install the 2100 Expert Software (version B.02.09 or higher) on the new PC.
- 2 Enter all the licenses including the Security Pack license and start the Security Pack setup.
- 3 Please use the same account names for the generic “2100System” account and the initial administrator account as used on the previous PC or for your previous 2100 Expert installation.

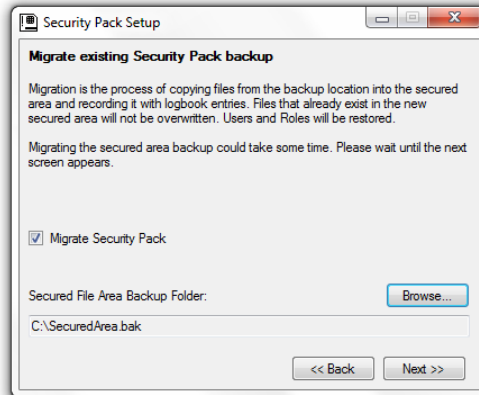


- 4 A new *Secured Area* is prepared on the PC in the default location (this can take a couple of minutes).

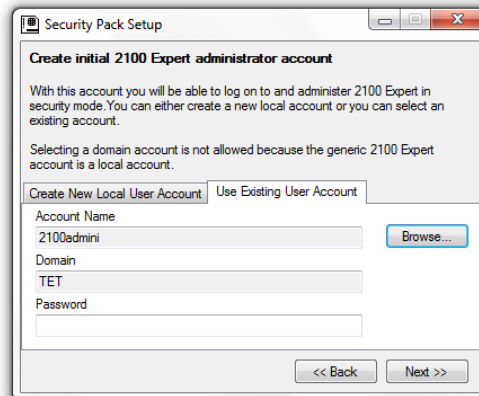


## Migration from the backup of the previous *Secured Area*

- 5 When prompted for the migration of the *Secured Area*, check the “Migrate Security Pack” box and select the directory of your *Secured Area* backup.



- 6 Specify the initial 2100 Expert administrator user account. You can create a new local Windows administrator account or choose an existing account from a domain.



- 7 Finish the Security Pack setup following the on-screen instructions. (When upgrading from B.02.08 to B.02.09, a previous Security Pack installation might be detected and you must confirm migration of system files).
- 8 Eventually, restore your files by de-archiving via the 2100 Expert Software.

## General Notes on Migration of the Secured Area between PCs

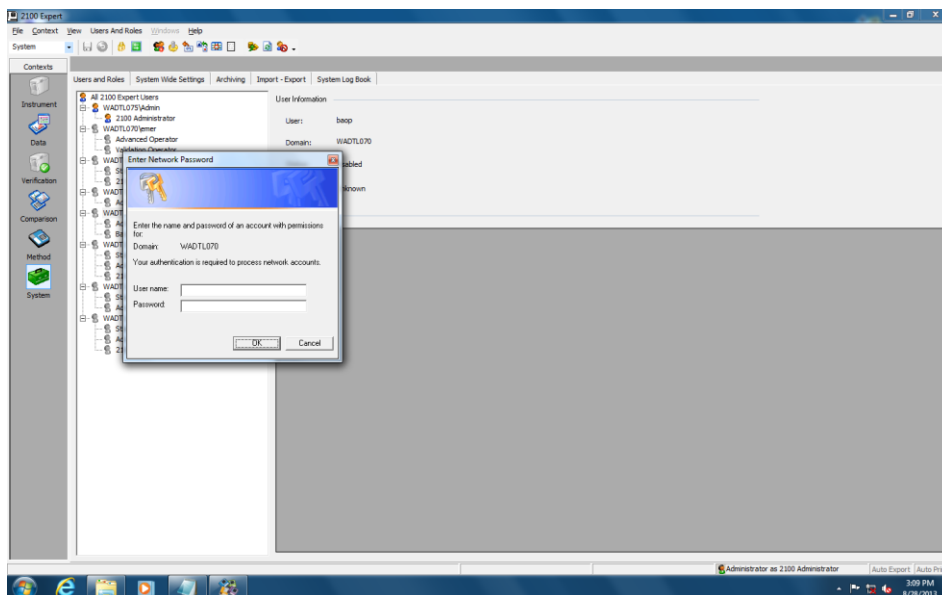
### NOTE

For domain accounts, the users are automatically migrated and do not need to be set up in the Security Pack software again. For local accounts the migrated accounts can only be used as a reference for previously acquired data. The users need to be set up on the PC and in the Security Pack software again.

## NOTE

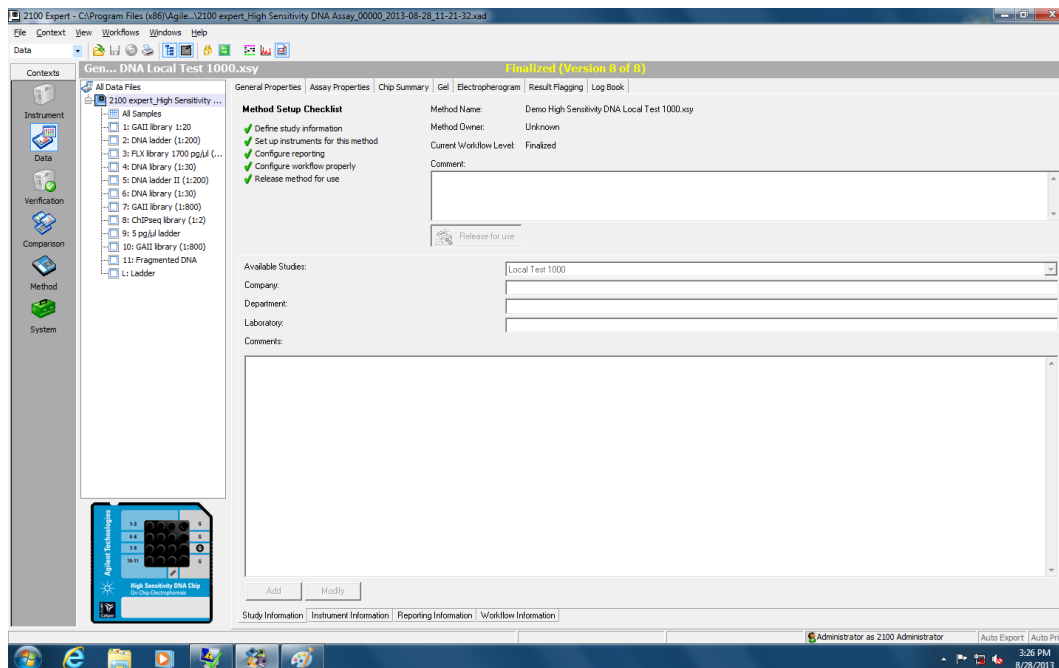
The data and methods from a previous PC can be opened for review on the new PC. However, for local user accounts, they can no longer be edited and old methods cannot be used unless imported by a new local user, who will become the new owner of the method.

In the “System” context tab named “Users and Roles”, the old users will be greyed out. This means they were migrated to the new PC without functionality. When trying to enable the user, there will be a prompt for the user’s logon credentials (i.e. user ID that was used on the old PC).





If a migrated method which was created by a local account is opened, the method owner will be changed to “unknown” as the local users from the old PC are unknown to the 2100 Expert Software installation on the new PC.



## NOTE

Important information: If the “ExtractUsersAndRoles.exe” is executed again on the new PC where the data was migrated to, it will overwrite any user information from the old PC and create a new UsersAndRoles.xml with the current user information.

**For Research Use Only**  
Not for use in diagnostic procedures

**Page 9 of 9**



G2938-90161 Rev.B

Part Number: G2938-90161 © Agilent Technologies, Inc. 2013, 2017, 2018

Edition 05/18  
Printed in Germany

Agilent Technologies  
Hewlett-Packard-Strasse 8  
76337 Waldbronn, Germany