

AriaMx & AriaDx – Steps to Activate HRM and ET Software Licenses

Introduction

This document provides a step-by-step guide for customers to activate their HRM (High Resolution Melt) Analysis and ET (Electronic Tracking) SW licenses. Customers who purchase an AriaMx Real-Time PCR instrument have the option of the HRM Analysis and ET software packages both at the time of instrument purchase and after initial instrument purchase. The HRM license is not offered for the AriaDx.

Aria SW add-on options	Part Number
High Resolution Melt (HRM) Analysis for AriaMx	G5381AA
High Resolution Melt (HRM) Analysis for AriaDx	N/A
Electronic Tracking (ET) for AriaMx	G5380AA
Electronic Tracking (ET) for AriaDx	K5822AA

Steps

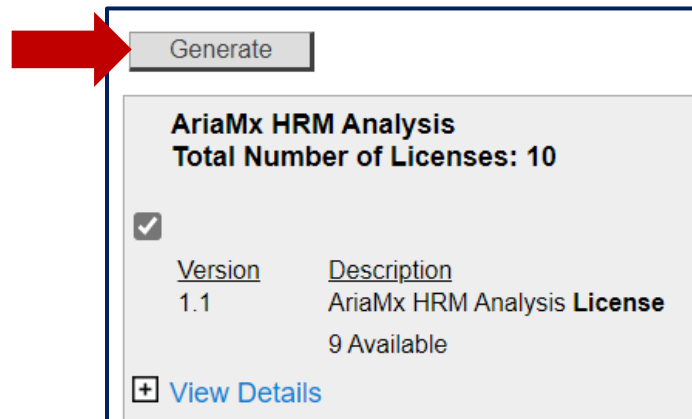
1. With the purchase of the HRM or ET software, a physical certificate with an Authorization Code is shipped. See [Appendix 1 – Example of SW Certificate with Authorization Code](#) for an example. This paper includes instructions for creating a profile in SubscribeNet and for configuring the license code. Authorization Codes are specific to the account for the same Agilent Genomics product.
2. Navigate to <https://agilent.subscribenet.com/>, follow the instructions to create a profile, and then login to the account.
3. Once logged in, use the Authorization Code on the certificate to register for a license. Click “Register Authorization Code” on the left tab.



- Once the Authorization Code has been accepted, navigate to “Generate or View Licenses” on the left tab.

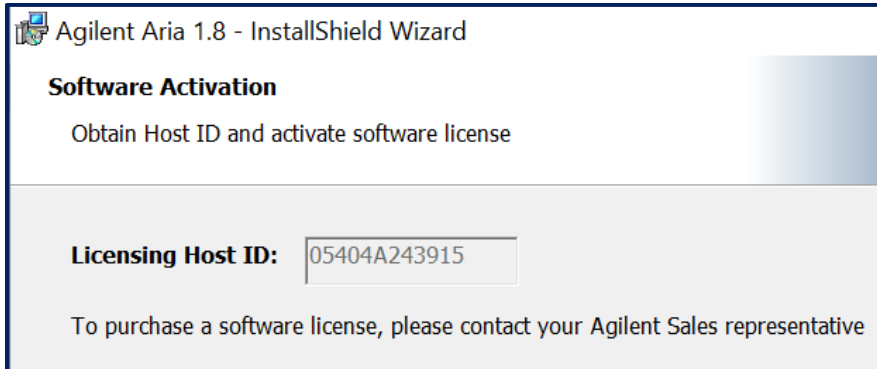


- Select the license type and click “Generate.”



6. Enter the Host ID number, which is specific to the customer’s PC. Enter “VHD=” followed by the Host ID number, which is the twelve-digit hexadecimal number displayed in the 'Software Activation' dialog box of your product software. This dialog box is found during the installation of the AriaMx software (with ET or HRM features enabled) or AriaDx software (with ET features enabled) on the customer’s PC. *

*For customers who already have downloaded the Aria software on their PC and would like to activate add-on licenses, you must first uninstall and reinstall the Aria software to identify the host ID and activate the license.



Click “Generate”.

Electronic Software and License Delivery

Generate Licenses

For each available license, enter 'VDH=' followed by the Host ID and click the 'Generate' button.

For AriaMX ET licenses, the Host ID is the twelve-digit hexadecimal number displayed in the 'Software Activation' dialog box of your product software.

After generation, the number of available licenses will be decremented by one.

		Number of Licenses Available
<u>Version</u>	<u>Description</u>	
1.1	AriaMx HRM Analysis License	9

Host ID

Generate

- To generate more licenses, click “Generate or View Licenses” in the left bar and repeat Steps #5 and #6. For HRM, the system will allow the user to generate a total of 10 licenses per authorization code.
For ET, each authorization code is good for 1 license.
- If you would like to switch or upgrade a PC, there is an option to retire an old HRM license file without losing the number of licenses in your pool. To return a license back to your License Pool, navigate to “Generate or View Licenses” and then click on the “View Details” link. Click on the “Return” button. The return increments the available quantity in the License Pool for the related license product.

AriaMx HRM Analysis
Total Number of Licenses: 10

<u>Version</u>	<u>Description</u>
1.1	AriaMx HRM Analysis License
	9 Available

[View Details](#)

License Quantity: 1 Expiration Date: permanent

Server Host ID: VDH=000000000000 **Return**

Generated By: on Jun 14, 2022

[About Return](#)
[About Upgrade](#)

NOTE: You only have two returns for each license product quantity in your License Pool. For example, for a license product with a quantity of two, you would have four returns. When you have used all your returns for a given license product, the 'Return' button will no longer appear. If you need additional returns, please contact [Agilent SubscribeNet Support](#).

Appendix 1 – Example of SW Certificate with Authorization Code



Agilent Technologies

Software Entitlement Certificate

This certificate evidences Agilent Technologies' grant to customer of licenses for the Agilent Genomics software product purchase as provided per the Agilent Genomics Authorization Code. The Agilent Genomics Authorization Code is specific to the software product purchased.

Please Note: Your Authorization Code is provided in the lavender license envelope in form of a separate label. Please keep this filed for future reference. Please place your label on the space below. The following represents an **EXAMPLE** only:

Feature Extraction 1-year Upgrades CML

Product: G4460AA Qty: 1

Authorization Code: D1A 021 68C3



Place your label here

To activate your license, please visit the following URL:

<https://agilent.subscribenet.com>

Each Agilent Genomics Product requires an individual account in SubscribeNet. Authorization Codes must never be registered to an account with any other Agilent product.

If you do not have a login yet, navigate to the registration page and provide your profile information including the authorization code from above. In turn you will receive an e-mail with your password.

If you have your login for the same Agilent Genomics product, go into SubscribeNet and register the additional authorization code for this product using "Register Authorization Codes". After that you can configure your license using "Generate and View Licenses" on the left navigation bar, and download the license file to your computer.

To obtain support, please contact your local Agilent Account Manager or Agilent Technical Support.

Rights and restrictions on the use, transfer, and the copying of software are set forth in Agilent Technologies' Software License agreement.



G2566-90080

Part Number: **G2566-90080**

Edition 06/14
Printed in USA

© Agilent Technologies, Inc. 2014

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051
USA