

Agilent Genomic Workbench ライセンスの取得方法

必要なもの

1) Order Number

Genomic Workbenchに添付のライセンス権利証明書 の
右上に記載されています

2) Certificate Number

Genomic Workbenchに添付のライセンス権利証明書 の
右上に記載されています

3) Host Name

下ページをご参考ください

4) メールアドレス

Agilent Technologies

Order Number: [Redacted]
Certificate Number: [Redacted]

Software Entitlement Certificate

This certificate evidences Agilent Technologies' grant to customer of licenses for the following software product purchase:

Product	Description	Quantity
G4147AA-001_V81	Feature Extraction Upgrade for Scanner Version 8.1	1

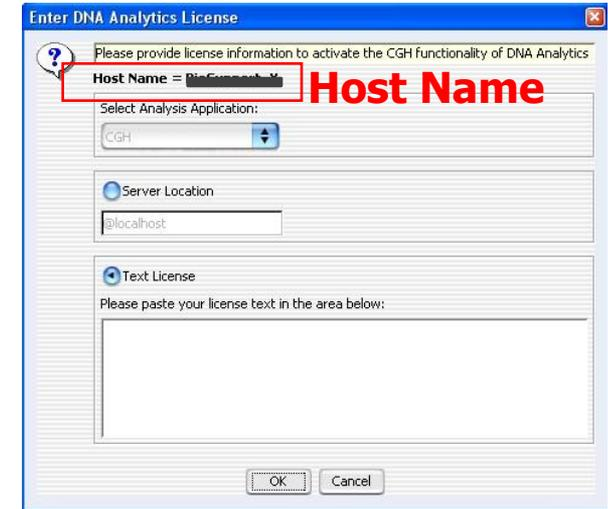
NOTICE:
Agilent Feature Extraction software allows you to visualize IFF images without a license. To enable data processing you must obtain a license by visiting our website at:
<http://www.agilent.com/ind/softwarelicense>
You will need to provide both the Order Number and the Certificate Number printed above, as well as the Host ID of the computer that will have the software installed.
To find the Host ID for the computer:
1. Install and Run the Feature Extraction software.
2. At the top of the screen, click on: Help
3. Select: About
The Host ID will be listed in the resulting screen.
The license file will be sent to you via email when the redemption process has been completed.

Rights and restrictions on the use, transfer, and copying of this software are set forth in Agilent Technologies' Software License agreement.
CE 106 Page 1 of 1

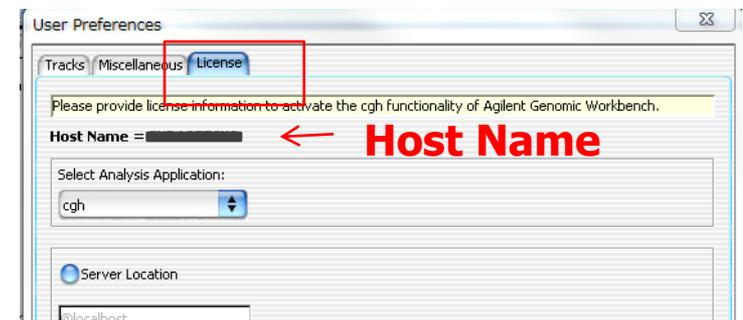
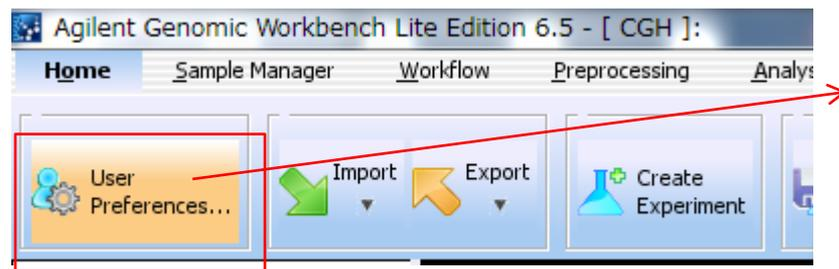
Host Name 確認方法

Host Name は各PC固有の文字列です。下記のいずれかの方法でご確認ください。

1) ライセンスが切れている・もしくはライセンスが入っていないAgilent Genomic Workbenchを立ち上げた際に
“Invalid License”のメッセージの後に表示されるライセンスボックス内



2) Agilent Genomic WorkbenchのHomeメニュー>User Preferences で表示されるボックスの中のLicense タブ内



1. 下記のAgilentウェブを開いてください。

<https://software.business.agilent.com>

2. 下記のページの“Click here to start software license redemption”をクリックしてください

The image shows a screenshot of the Agilent Software License Redemption web application. At the top left is the Agilent Technologies logo. To the right are navigation links: Search, Site Map, Login, Contact Us, and Home. Below these is a search bar with the text "Quick Search:" and a "GO" button. A secondary navigation bar contains links for About Agilent, Products & Services, Industries, International, and Online Stores. The main content area has a heading "Welcome to Agilent Software License Redemption." followed by two paragraphs of text explaining the application's purpose. A blue arrow points to the underlined link "Click here to start software license redemption." at the bottom of the main content area. The footer contains copyright information "© 2000 - 2002 Agilent Technologies", a link to "Terms of Use", and a link to "Privacy Statement".

3. [Login]のページにて Order NumberとCertificate Numberを入力してください

 **Agilent Technologies** **Close**

Help

**Agilent Software Licensing
Step 1 -- Login**

1 Login **2** Redemption Options **3** Enter Request **4** Review Request **5** Complete Request

Enter your Agilent Order number

Enter your Agilent Certificate number

I do not have a certificate number

Clear **Next**

4. [Redemption Options]ページのProduct 名を確認後、クリックしてください

Agilent Software Licensing
Step 2 -- Redemption Options

1 Login2 Redemption Options3 Enter Request4 Review Request5 Complete Request

Our Records Indicate You Have the Following for Order **Internal Use Only** CGH1005

Product	Description	Quantity Ordered	Quantity Available
G4175AA	CGH Analytics Node-Locked Annual License	120	28

|< << >> >| Page : 1 of 1

Request License(s) for:

- One or more products on a single instrument or host computer ?
- A single product on one or more instruments or host computers ?

[View Previously Issued Licenses For This Order](#)
[To Upload a License Input File For This Order](#)

BackNext

5. [Enter Request]

- ① ご購入いただいたライセンスをプルダウンからご覧いただけます。ライセンス発行するものを選択してください。(FeatureExtractionやChIPモジュールなど、一度に数種類のソフトウェアをご注文いただいた場合、ご注意ください。)
- ② “Add”ボタンをクリックしてください。
- ③ Host IDを入力してください
- ④ Nextボタンを押してください

Agilent Software Licensing
Step 3 - Request License(s)

1 Login 2 Redemption Options 3 Enter Request 4 Review Request 5 Complete Request

Please Select Products for Order Internal Use Only CGH1005

Please select one ①

Quantity Available Add

Enter Desired Quantity

License Requests (Product, Qty):

Back Next

Agilent Software Licensing
Step 3 - Request License(s)

1 Login 2 Redemption Options 3 Enter Request 4 Review Request 5 Complete Request

Please Select Products for Order Internal Use Only CGH1005

G4175AA - CGH Analytics Node-Locked Annual License (28 available)

Quantity Available 28 ② Add

Enter Desired Quantity 1

License Requests (Product, Qty):

Back Next

Agilent Software Licensing
Step 3 - Request License(s)

1 Login 2 Redemption Options 3 Enter Request 4 Review Request 5 Complete Request

Please Select Products for Order Internal Use Only CGH1005

G4175AA - CGH Analytics Node-Locked Annual License (28 available)

Quantity Available 28 Add

Enter Desired Quantity 1 Remove

License Requests (Product, Qty):
G4175AA, 1

Enter Node ID ③

Back Next ④

8. [Complete Request] 画面にて情報を確認後、画面を閉じてください

Agilent Software Licensing (ASL)
Step 5 -- Confirmation

1 Login2 Redemption Options3 Enter Request4 Review Request5 Complete Request

Nov 10, 2006 2:7:2

Dear Agilent Technologies customer:

Your request for license was submitted **successfully**.

The following is a summary of your request:

- Submitted request for host=BioSupport_Y product=G4175AA quantity= 1

In a few moments you should receive your requested licenses via email. Please remember to keep your **entitlement certificate** in a safe place. You can obtain order license status, view previously redeemed licenses, and request additional licenses from Agilent Software Licensing by using your entitlement certificate.

We thank you for your business.

Sincerely,

Agilent Software Licensing Team

P.S. To print this page for your records, click on the **Print** button below.

Please click [here](#) to request more licenses.

Print

9. ライセンスファイルが添付されたメールが指定アドレスに届きます。
*メールアドレスに届くファイルは".lic"という拡張子をもっています。
このファイルをソフトウェアがインストールされたPCに保管してください

10.

- ① Agilent Genomic Workbench を開いてください
- ② Switch Moduleメニューから、ご購入いただいたモジュールを選択してください
- ③ "Invalid License"のメッセージが出たらOKボタンを押してください。



- ④ ライセンスファイル”をノートパッドやテキストエディターなどのソフトウェアで開き、全文をコピーして
Text License ボックスにキーボードのコントロールキーとVを同時に
押してペーストしてください。

