

## 下記の項目をチェックください。

- インストールするバージョンの System Requirement および Software Compatibility をご確認ください。
   参照; <u>http://www.genomics.agilent.com/article.jsp?pageId=2579</u>
   なお、English OS でのみ検証されており、日本語 OS の場合予期せぬエラーや不具合が起こり得ます。
   何卒ご了承ください。
- 2. インストールする PC/Mac のログインアカウントに Administrator 権限があることを確認ください。
- 3. Widnows7の場合、アカウント権限の設定(p.2)を行ってください。
- 4. Common Storage フォルダの設定(p.3)の設定を行ってください。
- 5. 上記の設定を行った後、インストーラーを右クリックし'Run As Administrator'からインストールを実行してください

ご不明な点は、カストマコンタクトセンタまで(電話受付 9:00~12:00、13:00~18:00 土、日、祝日
 電話 0120-477-111
 FAX 0120-565-154
 E-mail: email\_japan@agilent.com/che



## ■アカウント権限の設定: Windows7 にインストールしている場合下記の2つの設定を行ってください。

1. Windows ログインアカウントのセッティングについて

1)WindowsのStartメニューより

> Control Panel > User Accounts > User Accounts を選択ください。

2)現在ログインで使っているアカウント名を選択後、Change User Account

Control Settings をクリックください

3)画面にスライダーが表示されますので、一番下の"Never notify me"まで下げ

てください。

2.プログラムの設定について

1) desktop の Cytogenomics のショートカット、

もしくは Start>すべてのプログラム> Agilent Cytogenomics

ver···> Agilent Cytogenomcis を選択し、

右クリック> Properties を選択ください

2)表示されたボックスの互換性 (Compatibility) タブを選択くださ

い。

3) 一番下にある項目 「特権レベリレ(Privilege Level)」にて、

管理者としてこのプログラムを実行する(Run this program as an

0120-477-111

0120-565-154

administrator)、にチェックを入れてください。

電話で

FAXOO





ご不明な点は、カストマコンタクトセンタまで(電話受付 9:00~12:00、13:00~18:00 土、日、祝日

E-mail: email\_japan@agilent.com http://www.agilent.com/che



## ■Common Storage フォルダの設定

- 1) C:¥Agilent¥AgilentCytoGenomicsEdition<version名>¥CytoGenomics\_Server"フォルダを開いてください。
- 2) CommonStorage\_XXXXX"という名称のフォルダがいくつかあります。

全てのフォルダについて 2)-3)のステップを繰り返してください。

GytoGenomics_Server		
マイル(日) 編集(日) 表示(以) お死に入り(日)	ツールロ ヘルプ国	
) #5 · 🕤 🤺 🖉 ## 🌔	7#119	
FER 🖗 😋 C#Agilen/#Agilen/CytoGenomics	Edition2.0.6.0#CytoGenomics_Server	
771#278#50920	7 6H	CommonStorage_201.0_13600
● 新しいフォルズを打ちまる ● このフォルズを Web に公開する ■ このフォルズを共有する	CommonStorage,2010,14206	GommonStorage 2010_14673
7018 ×	data (	Docs
AcientOvtoOxnomicsEdition		

3) フォルダを右クリック>共有 (Share with) > 特定のユーザー(Specific people…)を選択してください。表

 第 ファイルの共有

 ポットワーク上の共有相手となるユーザーを選択してください

 名柄を入力して[追加]をクリックするか、または、矢印をクリックして相手を除来してください。

 ・
 道加(A)

 名柄
 アクセス計可のレベル

 夏とveryone
 夏み取り/書き込み ・

 土丸の問題の詳細
 共有(H)

示されたボックスの中で、"Everyone"を追加ください。

4) 「アクセス許可のレベル(Permission Level)」にて、読み書き/書き込み(Read/Write) を選択してくださ

い。

ご不明な点は、カストマコンタクトセンタまで(電話受付 9:00~12:00、13:00~18:00 土、日、祝日

電話 20120-477-111 FAX 20120-565-154 E-mail: email\_japan@agilent.com http://www.agilent.com/che