## グリッドモードによる Agilentマイクロアレイの数値化 FeatureExtraction 11.x以降

#### Spot Finding of the Four Corners of the Array



#### 数値化後にグリッドがず れていた場合、マニュア ル操作で数値化を実行で きます。



## 1. Project 完了後にグリッドに関するWarningが出るの でYesを押してください>> Gridが保存されます



もしグリッドがずれてしまったスライドが1枚の場合、ステップ3に進んでください。

グリッドずれが起こったスライドが複数枚の場合、

ステップ2に進んでいただき、1スライドごとに数値化してください。



## 1. (別法)Project Explorerをセットします

#### File> New> Standard Projectを選択してください。 Project ExplorerにTifイメージ、Grid Name,Protocol Nameを1組だけ(1スライ ド分だけ) セットします。

FE_Project1 [New] [Standard - Config]							
Project Explorer P ×	😭 Project Properties 📲 Extraction Set Configuration						
- 🔂 FE_Project1	Extraction Set Name $\nabla$	Grid Name	Protocol Name	Output Name			
US82800151_251486829638_	US82800151_25148682	_F_20100118	GE1_107_Sep -	US82800151_25			
014868_D_F_20100118							
<b>€</b> GE1_107_Sep09							







#### FE Projectのツリー上で右クリックし、 Open Current Extraction in Grid Modeを選 択します



## 3. Mainグリッドの調整

#### 3-1)Adjust Main Grid ボタンを押した後全 体のMain Gridをあわせます





4. Subグリッドの調整

**4-1) Adjust SubGrid** ボタン を選択します。

4-2) SubGridCornerViewにてグリッドの位置 調整を行います



Select Corners and automatically fit Gridボタンが

・ONになっていた場合、コーナースポットをク リックすると自動的にサブグリッドがその位置に きます。

・OFFになっていた場合マウスを当てると Grid ↓ カーソルが表示されます。右クリックを押しなが らグリッドを移動し位置を調整します









## 5. グリッドファイルの保存 すべてのサブグリッドをあわせた後

#### 画面右上のGrid画面のClosexボタンを押します(Windows7の場合、 File>Save Gridを押すとエラーを起こすので必ずこの方法で閉じてください)

Agilent Feature Extracti	ion - [[Grid Mode] C:\Documents and Settings\chiho yoshida\Deskt	ppDemoDataVGE_TR_FE1010VJS11063876_25 🗐 🗖 🔀
Ele yew Edit Color Io	ools Feature Egtraction Window Help	- <del>0</del> ×
🔝 🗈 Green Channel	• 🗊 🔍 🔍 🔩 Auto Color Scaling 🛛 📲 1 🛛 🥊	. 🕫 🗶 🖂 🥻
15 15 10 10 .		
2000000	版 ● ■ Q .	
Grid Adjustment 4	9 ×	
Hain grid adjustment	0	
Array label 0 ⊐ Too left corner		
Left 961		
Top 469	×	
Main oxid adjustment		
riani gria adjusciniciik		CONCERNMENT OF CONTRACTORS

# Gridファイルを上書きするかどうか尋ねるボックスが表示されますので、Yesボタンを選択します





## 6. Feature Extractionの実行 すでに保存したグリッドがGrid欄に入っています。

### 6-1) Project Propertiesタブの Overwrite Previous ResultsをTrueに変更してください。

A discar Francisco Francisco FFF Designational February Academy Academi	🔮 Project Properties 📲 Extraction Set Configuration   🕴 4 🕨		
Ne Aglient Feature Extraction - [FE_Frojecti [New] [Standard - Config]]	Output and Data Transfer		A
The Mount File Distance Tests (Distance Dist	Outputs		
<u>File view Ealt Project Tools Window H</u> elp	MAGE	None	
	JPEG	None	
	TEXT	Local file only	
	Output Package	Compact	
	Zip Text File	False	
	Visual Results	Local file only	
	Grid	None	
Project Explorer 9 × 1 🚓 Project Property 📫 Extraction Set Configuration	QC Report	Local PDF file only	
	FTP Send Tiff File	False	
Erre Projecti Eutropian Set Name T Grid Name Protocol Name Output Name Service	Crop multipack image to individual arrays	False	
Extraction Set Name V Child Name Protocol Name Output Name Scan	<ul> <li>Local File Folder</li> </ul>		
■ ■ Get 2000101_201400020000 ULC020001E1 25140002 C01 gid opp CE1 107 Cop ▼ ULC020001E1 25 ULC020	Same As Image	True	=
JS82800151_2514868296: US62600151_25146662_301_gild.csv de1_107_3ep US62600131_23 US626	Results Folder		
	FTP Setting		
0002000101_2014000290	Post Processing Plug Ins		
G GE1 107 Sep09	Automatic Protocol Assignment		
	Highest Priority Default Protocol	Grid Template Default	
	Project Default Protocol		
	Automatic Grid Template Assignment		
	Use Grid file if available	False	
	External DyeNorm List File		
	Overwrite Previous Results	Irue	<b>_</b>
			•

#### 6-2) Projectのスタートボタンを押してください

