スキャナーマニュアル (v.7.0対応, 簡易版)



- Step 1—マイクロアレイスキャナの起動
- Step 2—スライドをスライドフォルダに取付ける
- Step 3—スライドフォルダをカローセルに入れる
- Step 4—スキャナコントロールソフトウェアを起動する
- Step 5—ユーザーの名前の入力とスロット範囲を設定する
- Step 6—カローセルの空のスロットをチェックする
- Step 7—スキャンの設定を変更する
- Step 8—カローセルランを開始する
- Step 9—アウトプットファイルを選択する
- Step 10—システムを終了する

Step 1—マイクロアレイスキャナの起動

- 1. マイクロアレイスキャナの電源スイッチを入れます。スイッチはスキャナ前面の左下にあります。
- 2. PCの電源スイッチを押してください。Windowsのユーザー名とパスワードを入力し、ログオンしま す。ログオンの初期設定は以下の通りです。
 - ユーザーアカウント: Administrator
 - パスワード: 3000hanover

あるいは、

- ユーザーアカウント: User
- パスワード: なし
- 3. 以上の順番でスイッチをONにすると以下の動作が行われます。
 - a. 換気がスタートします(背面の冷却ファンが動き出します)。
 - b. スキャナのセルフチェック(Power on Self Test)が始まります。通常約2分程度かかります。
 - c. セルフチェックが終了しPCとの通信接続が確立すると、黄色と緑色のインディケーターランプ が点灯した状態になります。この後、レーザーのウォームアップに数分~20分程かかります。
 - d. スキャナの準備が完了すると、黄色のランプが消え、緑色のランプのみが点灯した状態になり ます。



Figure 1 スキャナ電源スイッチとインディケータランプの位置

- ※ ご使用 20 分前にはスキャナとPC の電源を入れてください。スキャナの電源だけを入れても、レー ザーのウォームアップが終了しません(黄色のランプが消えない)のでご注意ください。
- ※ 必ずスキャナの電源を入れてから、PC の電源を入れてください。順序が逆になるとPC がスキャ ナを認識しないことがあります。

Step 2—スライドをスライドフォルダに取付ける

スライドはスキャン前にスライドフォルダに差し込んでおく必要があります。これにより、スキャン中ス ライドを保護することになります。

指紋等の汚れは、蛍光となってイメージに現れます。スライドを扱うときは手袋をし、スライドの端以 外は触らないようにします。また、バーコード以外のシール類はスライドに貼らないようにして下さい。

1. スライドフォルダをアジレントのロゴがある方を上にして持ちます。親指をカバーの中央に置き、押 しながら上部にスライドしてカバーを外します。(Figure 2参照)



Figure 2 スライドフォルダのカバーを開ける

- 2. スライドを取付ける前に、以下のことに注意してください。
 - バーコードがない端を奥に入れます。
 - アレイ面がカバーの方になるようにスライドを置きます。アジレントのマイクロアレイスキャナは
 アレイの裏側からイメージを読み取ります。

アレイによっては、バーコードが片面あるいは両面に貼り付けられている場合があります。どちらの場合も、常にアレイ面をカバー側にして、スライドを置くようにして下さい。

詳細は英文マニュアル123ページの "Barcode specifications"をご参照ください。

3. Figure 3のように、スライドをファルダに置き、カバーを倒して押しながら手前にスライドさせて、カ バーを固定します。



Figure 3 スライドをフォルダに取付ける



Figure 4 スライドオリエンテーション

Step 3—スライドフォルダをカローセルに入れる

スライドをスライドフォルダに差し込んだ後は、スライドフォルダをカローセルにセットします。カローセ ルはスライドフォルダが常に正しい位置に入るようにデザインされています。

この項ではカローセルがスキャナに設置されている場合のステップを説明します。カローセルをスキャナから取り出し、別の位置でスライドフォルダを入れる場合は、英文マニュアル86ページの"Mount or remove the carousel" をご参照ください。

- スライドフォルダをカローセルのスロット1にいれます。スロットの順番はホームスロット(ホームポ ジション)から右回りとなります。
 スライドフォルダをホームスロットに入れないように注意してください。(ホームスロットは他のスロットに比べて幅がやや広く、またHomeという表示がついています。)
- バーコードラベル側がカローセルの中心部に向くように、スライドフォルダを入れてください。スライ ドガラスは手前にくるようにします。
- 3. スライドフォルダがきちんとカローセルの底まで位置したことを確認します。



Figure 5 カローセル

(オプション) カローセルの蓋を取付けます。タブを持ち上げ、きちんと蓋がカローセルにはまるようにします。



Figure 6 カローセルカバー

5. スキャナの蓋を閉めます。

以下の動作はスキャナの蓋を閉めないと行えません。

- スキャナのイニシャライゼーション(スキャンコントロールソフトウェアの起動)
- カローセル内の空のスロットの確認(Check Carousel)
- バーコードの確認(Read Barcode from Carousel)
- カローセルランの開始(Scan Slot)

Step 4—スキャナコントロールのソフトウェアを起動する

PCワークステーションにあるスキャナコントロールソフトウェアを使って、スキャンを行います。スキャンの設定、スタート、停止(アボート)はこのソフトウェア画面から操作します。

1. デスクトップ上のアジレントスキャナのアイコンをダブルクリックして、ソフトウェアをスタートします。



Figure 7 スキャンコントロール アイコン

Agilent Scan Controlのウィンドウが表示され、スキャナのイニシャライズが開始されます。

Agilent Scan Control									
jie Settings Help									
Current c	Current carousel settings								
	e <u>×tended Dynamic Range Scan Mode</u>								
<u>O</u> perato	Operator: XDR Red PMT (%)								
			- 1						
St <u>a</u> rt slo	t: 1 <u>▼</u> E <u>n</u> d slo	t: 48 - Check Larouse							
Slot #	Slide ID/Barcode	Scan Region(mm)	Red PMT(%)	Green PMT(%)	Scan Resolution(µm)	Dye Channel	Output Path		
1		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
2		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:A		
3		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:V		
4		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
5		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
6		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
7		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
8		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
9		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
10		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
11		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:A		
12		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
13		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\		
Scan re	gion option: Use selected :	scan region for the slides in the carouse	l	5µm scanning m	ode: Single Pass	Hes	Edit Slot Values >>		
_									
icanner sta	atus: Lasers are warming				<u>^</u>	<u>S</u> car	n Slot 1-48 Abort		
					7				

Figure 8 スキャンコントロールメイン画面

イニシャライズ後、レーザーのウォームアップがまだ終了していない場合は、Scanner statusメッセ ージ欄に"Lasers are warming…"と表示されます。

レーザーがウォームアップ中も、次のステップ5の作業は可能です。レーザーのウォームアップには数分~20分程度かかります。

レーザーのウォームアップが完了しないうちは、カローセルランを開始することができません。ウォ ームアップが完了すると、メイン画面のScanner statusメッセージに"Scanner ready"と表示されま す。(併せて、スキャナ本体前面のLEDを確認して下さい。緑のLEDだけが点灯している状態が、 準備完了です。緑と黄色、両方のLEDが点灯している場合は、レーザーのウォームアップが完了 していません)

×	Agilent !	Scan Control							
Ei	le S <u>e</u> tting	gs <u>H</u> elp							
Γ	Current c	arousel settings				tended Dunamic Bange	Scan Mode		
	Onerator Onerator								
	Hit 100 Y Lo: 10 Y Lo: 10 Y								
	Start slot: 1 💌 End slot: 48 💌 Check Carousel								
	Slot #	Slide ID/Barcode	Scan Region(mm)	Red PMT(%)	Green PMT(%)	Scan Resolution(µm)	Dye Channel	Output Path	-
	1		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:V	
	2		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:V	
	3		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	4		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	5		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	6		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	_
	7		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:V	
	8		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	9		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:N	_
	10		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	11		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:A	_
	12		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	13		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
		1		1400	1400	140			
	Scan rej	gion option: Use selecte	ed scan region for the slides in the carous	el	5µm scanning m	node: Single Pass	Res	e <u>t</u> Selection [Edit Slot Va	ues <u>>></u>
							C	0	
	canner st	atus: Scanner ready.)			~ 	<u> </u>	1 510t 1-48 Ag	ion
_			t						
		,	\mathbf{N}						
			\mathbf{N}						
		r	1						
			Scanner Ready	/					

Figure 9 スキャンコントロールメイン画面(レーザーのウォームアップが終了している場合)

Step 5—ユーザーの名前の入力とスロット範囲の設定

初めに:「スキャン」とはスライド1枚あたりのイメージの読み取りを指します。「カローセルラン」はカローセルに ある、全てのスロットのスキャンを意味します。

カローセルの全48スロットがスライド(フォルダ)で埋まっていない場合は、スライドを入れたスロット の場所(番号)を特定することができます。これにより、スキャン作動時間の節約ができます。

カローセルランを開始すると、ユーザーが入力したスタートスロットからスキャンを開始し、エンドスロットでスキャンを終了します。

- 1. Operator のところにユーザー名を入力します。
- カローセルの48スロットを全て使用する場合(48枚スライドがある場合)は、以下の項目は無視してステップ7にお進みください。
- 3. Start slot: ボックスのドロップダウン↓をクリックします。
- 4. 最初にスキャンするスロットの番号を選択します。



Figure 10 スタートスロット番号の選択

- 5. End Slot: ボックスのドロップダウン↓をクリックします。
- 6. 最後にスキャンするスロットの番号を選択します。
 ここで入力したスロットの番号に基づき、以下のスキャンコントロールのメイン画面でテーブル表
 - 示が変更されます。
 - ・ 指定したスロット範囲のみ、Check Carousel機能が使えます。
 - 指定したスロット範囲のみ、スキャンの設定変更ができます。



Figure 11 スロット範囲の指定

Step 6---カローセルの空のスロットをチェックする

指定スロット範囲の中に、空のスロットがないかどうか調べます。これによりカローセルラン(スキャン)の時間が節約できます。

このステップを行わないと、ステップ5で指定した範囲内の全てのスロットに対して、スキャンをしようとします(スロットにスライドフォルダがあるかないかを調べます)。

1. スキャンしたいスロットを入力します(前項ステップ5をご参照ください)。

2. Check Carouselをクリックします。

×	Agilent S	ican Control							×
Eil	e S <u>e</u> tting	s <u>H</u> elp							
Г	Current carousel settings								
	Derator: XDR Red PMT (%) XDR Green PMT (%)								
	St <u>a</u> rt slot	::1 ▼ E <u>n</u> d	slot: 🚺 🧲 Check Carouse	el)					
								T	
	Slot #	Slide ID/Barcode	Scan Region(mm)	Red PMT(%)	Green PMT(%)	Scan Resolution(µm)	Dye Channel	Output Path	
	1		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:V	
	2		Scan Area (61 x /1.6mm)	100	100	10	Red&Green	D:V	
	3		Scan Area (61 21.6mm)	100	100	10	Red&Green	D:\	
	4		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	5		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:A	
	6		Scan Area 61 x 21.6mm)	100	100	10	Red&Green	D:A	
		_	/						
			Check Carousel						
			Oneek Garouser						
								Þ	
	Scan region option: Use selected scan region for the slides in the carousel 5µm scanning mode: Single Pass Reset Selection Edit Slot Values 2>								
s	canner sta	itus: Scanner ready.				A V	<u>S</u> car	n Slot 1-6 Abort	

Figure 12 Check Carouselの位置

カローセルチェック後のメイン画面は以下のようになります。

Current carousel settings							
Slot #	Slide ID/Barcode	Scan Region(mm)	Red PMT(%)	Green PMT(%)	Scan Resolution(µm)	Dye Channel	Output Path
1	No chin						
2	No chip						
3	•	Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\
4	No chip						
5	No chip						
6	No chip						
Scan region option: Use selected scan region for the slides in the carousel 5µm scanning mode: Single Pass Reset Selection Edit Slot Values >>							

Figure 13 カローセルチェック後のスキャン画面の表示例

カローセルチェック後に、スライドフォルダを入れ替えた場合(例:空のスロットにスライドフォルダを 入れたり、スライドフォルダを取り出してスロットを空にした場合)は、全てのスロット行を選択して、 Reset Selectionをクリックします。この操作により、初期設定の表示に戻り、48スロットが表示されま す。もう一度設定を変更したい場合は、Check Carouselを再度クリックします。

Step 7—スキャンの設定を変更する

アジレントマイクロアレイをお使いの場合:

スキャンコントロールの初期設定はアジレントのマイクロアレイ(高密度アレイを除く)に適した値となっています。

- 設定を変更する必要のない場合は、ステータス表示が"Scanner Ready" となっていることを確認し、ステップ8にお進みください。
- 設定を変更する場合、バーコード情報を入力する場合は以下のステップを行います。

設定を変更する場合は、Edit Slot Values>>をクリックして、メイン画面を拡張します。

Agilent Scan Control								
le S <u>e</u> tting	gs <u>H</u> elp							
- Current ca	Current carousel settings							
Operator	Potential of narrige scan mode VDD Course DMT (%) VDD Course DMT (%)							
<u>o</u> perator								
	Hit 100 💌 Lo: 10 💌 Hit 100 💌 Lo: 10 💌							
St <u>a</u> rt slot	Start slot: 1 End slot: 6 Check Carousel							
Slot #	Slide ID/Barcode	Scan Region(mm)	Red PMT(%)	Green PMT(%)	Scan Resolution(µm)	Dye Channel	Output Path	
1	No chip							
2	No chip	Corrections (C1 in C1 Correct)	100	100	10	De de Casara	D.I.	
3	No chin	Scan Area (61 x 21.6mm)	100	100	10	ned&ureen		
5	No chip							
6	No chip							
-	Edit Slot Values>>							
Scan reg	gion option: Use selected s	can region for the slides in the carousel		5μm scanning m	node: Single Pass	Res	et Selection Edit Slot Values 2>	
uarmer sta	atus. Jocariner ready.				- -	<u>5</u> car	Agon Agon	

Figure 14 Edit Slot Values>>の位置

plent Scan Control gettings Settings tells genator Page alors Start side: I I Scan Region(rm) Red PMT(2) Scan Resolution(im) Duput Path Duput Path I No chip Scan Area (SI x 21 6mm) 100 I No chip Scan region option: Use selected scan region for the slides in the carouse! Single Pasc Reset/Selection Hide Editing C Scan Area (SI x 21 6mm) Single Pasc Reset/Selection Hide Editing C Scan Area (SI x 21 6mm) Single Pasc Reset/Selection Hide Editing C Scan Area (SI x 21 6mm) Scan Region (gmm) PMT(12) Reset/Selection Hide Editing C Scan Area (SI x 21 6mm) ToO mode Scan Area (SI x 21 6mm) ToO mode Description: Read Barcode from Carouse! No chip						eXtend	led Dyna	mic Sca	an Mode
sptms Bels gendor									
		c				/			
	S <u>e</u> ttin	Ings Help		_					-
	Current c	carousel settings							
Hit 100 K Lot 10 K Hit 100 K Lot 10 K Start stor: 1 K Epd stor. 5 K Check Garousel Start stor: 1 K Epd stor. 5 K Check Garousel Start stor: 1 K Epd stor. 5 K Check Garousel Start stor: 1 K Epd stor. 5 K Check Garousel Start stor: 1 K Epd stor. 5 K Check Garousel 1 No chip 1 K Chip 1 K Chep 2 No chip 5 Can Area (51 x 21.6mm) 100 10 4 No chip 5 Can Area (51 x 21.6mm) 100 10 Rest&Green 5 No chip 5 Can Area (51 x 21.6mm) 100 10 Rest&Green D/A 6 No chip 5 Can region option Use selected scan region for the slides in the carousel Sum scanning mode: Single Pass Rest Selection Hide Editing S dt slot values Bed Green Scan Area (51 x 21.6mm) 100 K No K Single Pass Rest Selection Hide Editing S Read Bacode from Carousel Brogree Set Yalues Set Yalues Set Yalues Nutput path: D:A Brogree Set Yalues Set Yalues Set Yalues <td>Operato</td> <td>or:</td> <td></td> <td></td> <td></td> <td>ended Dynamic Range Bed Phil (%)</td> <td>Scan Mode</td> <td>'NB Green PMT</td> <td>(%)</td>	Operato	or:				ended Dynamic Range Bed Phil (%)	Scan Mode	'NB Green PMT	(%)
Start slot: 1 Egd slot: 6 Check Carouzel Start slot: 1 Egd slot: 6 Check Carouzel Start Skide ID/Barcode Scan Region(mm) Red PMT(2) Green PMT(2) Green PMT(2) Green PMT(2) Red&Green DX Red&Green DX Red&Green DX Red&Green DX Red&Green Reset Selection Hide Editing Green Scan PMT(2) Red&Green Reset Selection Hide Editing Green Scan PMT(2) Red&Green Scan PMT(2) PMT		1			Hi:	100 V Lo: 1	10 🖵	Hi: 100 🔽	Lo: 10
Start stol: 1 Egd stol: 6 Check Carousel Start stol: 1 No chip Free PMT(%) Scan Resolution(um) Dye Channel Duput Path 1 No chip 5 Scan Area (61 x 21.6mm) 100 10 Red PMT(%) Scan Resolution(um) Dye Channel Duput Path 1 No chip 5 Scan Area (61 x 21.6mm) 100 10 Reset Selection Hide Editing 4 1 Scan Region option: Use selected scan region for the sides in the carousel Sym scanning mode: Single Pass Reset Selection Hide Editing 4 Start start Scan Area (61 x 21.6mm) PMT(%) PMT(%) resolution(um) Dye channel Set Yalues Read Barcode from Carousel Scan Area (61 x 21.6mm) 100 10 Red Barcode from Carousel Description: Scan Stol 1-6 Abort Nuput path D:1 Brogree Scan Stol 1-6 Abort Scan Stol 1-6 Abort								, _	, _
Side II //Barcode Scan Region(mm) Red PMT(2) Green PMT(2) Scan Resolution(µm) Dype Channel Dutput Path 1 No chip Scan Area (61 x 21 6mm) 100 10 Red&Green D/ 3 Scan Area (61 x 21 6mm) 100 10 Red&Green D/ 4 No chip Scan Area (61 x 21 6mm) 100 10 Red&Green D/ 5 No chip Scan Area (61 x 21 6mm) 100 10 Red&Green D/ 5 No chip Scan Area (61 x 21 6mm) 100 10 Red&Green D/ 5 No chip Scan region option: Use selected scan region for the sides in the carousel Sym scanning mode: Single Pass Reset Selection Hide Editing set of the sides in the carousel 5 No chip Scan region (gm) PMT(2) resolution(µm) Que channel Set Yalues 6 Scan Area (61 x 21.6mm) 100 100 10 Red&Breen Set Yalues 7 Scan Slot 1.6 Scan Slot 1.6 Abort Set Yalues Set Yalues	St <u>a</u> rt slo	ot:: 1 💌 E <u>n</u> d slo	ot: 6 💌 Cł	neck <u>C</u> arousel					
1 No chip Image: Scan Area (61 x 21.6nm) 100 100 10 Red&Green D/A 4 No chip Image: Scan Area (61 x 21.6nm) 100 100 10 Red&Green D/A 5 No chip Image: Scan Area (61 x 21.6nm) 100 100 10 Rest Selection Hide Editing S ati slot values Image: Scan region (gm) Eled Green Scan Scan region (gm) Dye channel 3 Scan Area (61 x 21.6nm) 100 100 I00 I00 Rest Selection Hide Editing S ati slot values Brogyse Description: Set Values Set Values Tread Barcode from Carouse] Immer status: Scanner ready. Scan Stot 1-6 Abort August path. D/A Brogyse Scan Stot 1-6 Abort	Slot #	Slide ID/Barcode	Scan Region(mm)	Red PMT(%) Green PMT(%)	Scan Resolution(µm)	Dye Channel	Output Path	
2 No chip Scan Area (61 x 21 6mm) 100 10 Red&Green 01 4 No chip Image: Scan Area (61 x 21 6mm) 100 10 Red&Green 01 5 No chip Image: Scan Area (61 x 21 6mm) 100 10 Red&Green 01 4 No chip Image: Scan Area (61 x 21 6mm) 100 10 Rest Selection Hide Editing Scan region option: 5 No chip Image: Scan region (for the slides in the carouse! Sum scanning mode: Single Pass Rest Selection Hide Editing Scan Area (61 x 21.6mm) 5 Scan Area (61 x 21.6mm) Y 100 Y 100 Rest/Side (01 P Area Scan region (for the slides in the carouse! Set Yalues 7 Read Bacode from Carouse! PMT(%) PMT(%) Y 100 Y 100 Set Yalues 7 Read Bacode from Carouse! Perception: Set Yalues Set Yalues 9 Scan Slot 1-6 Abort Set Yalues Set Yalues	1	No chip							
4 No chip No chip 5 No chip Image: Second Sec	3	No chip	Scan Area (61 x 21 6mr	n) 100	100	10	Bed&Green	D:\	
5 No chip 6 No chip 6 No chip 4 Scan region option: Use selected scan region for the slides in the carousel 5µm scanning mode: Single Pass Scan region option: Use selected scan region for the slides in the carousel 5µm scanning mode: Single Pass Reset Selection dit slot values: Bed Green Scan slot values: Bed Green Scan ilot Slide [D/Barcode Scan region (mm) PMT(%) PMT(%) PMT(%) 3 Scan Area (61 x 21.6mm) 100 100 Read Barcode from Carouse] Description: Pread Barcode from Carouse] Browse Description: Scan Slot 1-6 Abort mmer status: Scanner ready. Scan Slot 1-6 Abort	4	No chip		ny 100	100		1100acaroon		
6 No chip ▲ Scan region option: Use selected scan region for the slides in the carousel Sµm scanning mode: Single Pass Reset Selection Hide Editing ≤ dit slot values Bide [D/Barcode Scan region (mm) PMT(X) Perception Set Values 3 Scan Area (61 x 21.6mm) 100 100 PMT(2) Description: Read Barcode from Carouse! Browse Description: Scan Slot 1-6 Abort Inner status: Scanner ready. Scan Slot 1-6 Abort	5	No chip							
Scan region option: Use selected scan region for the slides in the carousel Sum scanning mode: Single Pass Reset Selection Hide Editing d the slot values Bed Scan Region (mm) PMT(%) PMT(%) PMT(%) Resolution(µm) Dye channel Set Values Read Barcode from Carousel Dutput path D:\ Read Barcode from Carousel Scanner ready. Scan Stot 1-6 Abor: Set Values Set Values	6	No chip							
dit slot values Stot Slide [D/Barcode Scan region (mm) PMT(2) PMT(2) resolution(µm) Dye channel Scan Area (61 x 21.6mm) 100 100 10 Red&Green Description: Read Barcode from Carouse! Dutput path: D:\ Inner status: Scanner ready. Scan Slot 1-6 Abort. Set Values									
Side [D/Barcode Scan region (mm) PMT(% PMT(%) resolution(µm) Dye channel Scan Area (61 x 21.6mm) 100 100 10 Pectigrion: Read Barcode from Carouse Dutput path: D:\ Inner status: Scanner ready. Scan Slot 1-6 Abort Set Values Set Values	▲ Scan re	egion option: Use selected	scan region for the slides i	n the carousel	5µm scanning n	node: Single Pass	Res	et Selection	Hide Editing <
s Scan Area (61 x 21.6mm) 100 100 I00 IRed&Green Set Values Read Barcode from Carouse] Description: Description: Description: Set Values Inner status: Scanner ready. Scan Slot 1-6 Abort.	Scan re Edit slot v	egion option: Use selected values	scan region for the slides i	n the carousel	5µm scanning n	node: Single Pass	Res	set Selection	Hide Editing ≤
Read Barcode from Carouse! Description: Dutput path: D:\ inner status: Scanner ready. Scan Slot 1-6 Set Values	Scan re Edit slot v Slot	egion option: Use selected values Slide JD/Barcode	scan region for the slides i Scan region (<u>m</u> m)	n the carousel Red PMT(%) P	5μm scanning n reen Scan MT(%) resoluti	node: Single Pass cn(μm) Dye channel	Res	set Selection	Hide Editing <
Dutput path: D:\ Immer status: Scanner ready. Scan Slot 1-6 Abort Set Values	Scan re Edit slot v Slot	agion option: Use selected values Slide [D/Barcode	scan region for the slides i Scan region (<u>m</u> m) [Scan Area (61 × 21.6mm	n the carousel Red PMT(%) P	5µm scanning n reen Scan MT(%) resoluti 00 ▼ 10	node: Single Pass on(μπ) Dye channel	Res	tet Selection	Hide Editing <u><</u> Set <u>V</u> alues
Inner status: Scanner ready.	Scan re Edit slot • Slot 3 Read	egion option: Use selected values Slide [D/Barcode d Barcode from Carouse]	scan region for the slides i Scan region (<u>m</u> m) Scan Area (61 x 21.6mm	n the carousel Red PMT(왕) P a)	5µm scanning n reen Scan MT(%) resolyti 00 ▼ 10	ode: Single Pass an(µm) Dye channel V Red&Green Description:	Res		Hide Editing ⊴ Set ⊻alues
Inner status: Scanner ready.	Scan re Edit slot • Slot Reac Output p	agion option: Use selected values Side [D/Barcode d Barcode from Carouse] path: [D:\	scan region for the slides i Scan region (<u>m</u> m) [Scan Area (61 × 21.6mm	n the carousel PMT(%) P a)	5µm scanning n reen Scan MT(%) resoluti 00 ▼ 10 Browse	node: Single Pass on(µm) Dye channel ▼ Red&Green Description:	Res		Hide Editing ≤ Set ⊻alues
Set Values	Scan re Edit slot v Slot 3 Read	agion option: Use selected values Slide [D/Barcode d Barcode from Carouse] path: D:\	scan region for the slides i Scan region (<u>m</u> m) [Scan Area (61 x 21.6mm	n the carousel Bed PMT(%) P 100 ¥ 1	5µm scanning n reen Scan MT(%) resoluti 00 ▼ 10 Browse	ode: Single Pass on(μm) Dye channel ▼ Redt&Green Description:	Res		Hide Editing ⊴ Set ⊻alues
Set Values	Scan re Edit slot v Slot Read Output p	egion option: Use selected values Slide ID/Barcode d Barcode from CarouseI path: D:\ tatus: Scanner ready.	scan region for the slides i Scan region (<u>m</u> m) Scan Area (61 × 21.6mm	n the carousel <u>B</u> ed PMT(%) P) ▼ 100 ▼ 1	5µm scanning n reen Scan MT(≋) resolyti 00 ▼ 10 Browse	node: Single Pass on(µm) Dye channel ▼ Red&Green Description:	Res	n Slot 1-6	Hide Editing ≤ Set ⊻alues Abort
	Con re Scan re Slot Slot Reac Output p anner st	egion option: Use selected values Slide JD/Barcode d Barcode from Carouse] path: D:\ tatus: Scanner ready.	scan region for the slides i Scan region (<u>m</u> m) Scan Area (61 x 21.6mm	n the carousel Bed PMT(%) P 100 V [5µm scanning n reen Scan MT(%) resoluti 00 ▼ 10 Browse	ode: Single Pass	▼ <u>S</u> ca	n Slot 1-6	Hide Editing ≤ Set ⊻alues

任意のスキャン設定項目を変更後、右端のSet Valuesをクリックします。設定変更後、Set Values を押さないとCurrent slot settings内のテーブルに反映されません。

<注意>

アジレントの高密度アレイ 4 x 44K フォーマットをスキャンする場合は、必ず eXtended Dynamic Scan Mode にチェックを入れ、Scan resolution を 5 μ m に設定します。

244K, 2x105K フォーマットのアレイをスキャンする場合は、チェックを入れず、Scan resolution のみ 5 μ mに設定します。

XDR(e<u>X</u>tended <u>Dynamic Range</u>)機能の詳細については、英文マニュアルの 20, 62, 65 ページを 参照してください。

初期設定を変更する場合は、	Settings>Modify Default Settingsをクリックして、Defaul	t
Settings画面を表示します。		

Default Settings	
Scan configuration	
Scan region	Dye <u>c</u> hannel: Red&Green 💌
Regio <u>n</u> (mm): Scan Area (61 x 21.6mm)	Scan resolution (μm): 10 💌
Options	PMT sensitivity level
Attempt to retrieve from XML (GEML) files	🔲 eXtended Dynamic Range scan mode
Set XML (GEML) File Path	Red (%) Green (%)
	Standa <u>r</u> d: 100 💌 100 💌
	XDR Hi 100 🔽 100 💌
5μm scanning mod <u>e</u> : Single Pass 💌	XDR Lo 10 - 10 -
Description:	<u>^</u>
	v
Scan image file handling	
Output path: JD:N	<u>B</u> rowse
Automatic file naming (the format is: Prefix1_Prefix2_scanne	umber.TIF)
Prefix2	
Slot Number Customized:	
Split and rotate TIFF image Compress TI	FF image
Help	OK Cancel

Figure 16 Default Settings画面

スキャン設定についての詳細は英文マニュアルのChapter 3 How Do I...?をご参照ください。

スキャン前の設定チェックリスト(アジレント、アジレント以外のスライドをお使いの場合)

カローセルランを開始する前に設定を変更する場合は、以下のリストをご参考ください。それぞれの 設定項目については、英文マニュアルの3 How Do I...?.に記載してあります。

設定チェックリスト	英文マニュアル
	参照ページ
スキャナで読み取れるバーコードについて	56ページ
バーコードのマニュアル入力	58ページ
スキャン設定の初期設定の変更	60ページ
表設定の変更	65ページ
アジレントスライドのスキャン範囲の設定	67ページ
non-アジレントスライドのスキャン範囲の設定	69ページ
non-アジレントスライド用のスキャン範囲の新規設定	71ページ
感度、分解能、レーザーの設定	75ページ
ファイル保存先の指定、ドライブのスペース確認	76ページ
自動ファイル名記入の設定変更	78ページ
スキャンイメージの回転と1色イメージへの変換	79ページ
スキャンイメージの圧縮	80ページ

Table 1 スキャン前の設定チェックリスト

Step 8—カローセルランを開始する

開始する前に:

マイクロアレイスキャナのウォーミングアップが終了していることを確認します。(Statusタブが Scanner Ready と表示されているかご確認ください。)

カローセルランの開始

1. メイン画面のScan Slot m-nをクリックします。

ж	🔀 Agilent Scan Control								
Eile	S <u>e</u> tting	ıs <u>H</u> elp							
Г	Current ca	arousel settings							
					□ e⊻	ended Dynamic Range	Scan Mode		
	Operator.	r.				Red PMT (%)	×	DR Green PMT (%)	
					Hi:	100 V Lo:	10 🖵 🛛 H	Hi: 100 🔻 🛛 Lo: 10	
	Charles and Friday and Charles Caround								
	Start slot								
	Slot #	Slide ID/Barcode	Scan Region(mm)	Red PMT(%)	Green PMT(%)	Scan Resolution(µm)	Dye Channel	Output Path	
	1	No chip							
	2	No chip							
	3		Scan Area (61 x 21.6mm)	100	100	10	Red&Green	D:\	
	4	No chip No chip							
	6	No chip							
		No chip							
	Soonroo	tion option: Use colocted a	approximation for the elides in the entry we		Euro coopping p	odo: Single Page	Bese	et Selection Edit Slot Valu	10533
	scarrieg	gion option. To se selected s	can region for the sides in the carouser		opun scanning n	ioue. Joingle Fass			<u>403 Z</u> /
S	canner sta	atus: Scanner ready.				÷	Scar	n Slot 1-6	ort
							200		
							1		
							/		
	Fi	iqure 17 Scar	n Slot の位置				2		
		.g				Scan Slot	n-m		

m には最初のスライドが位置するスロット、n には最後のスライドが位置するスロットのそれぞれの番号が表記されます。

スキャンを開始すると、以下の操作が自動的に行われます。

- スキャナのドアにロックがかかります。
- Scan Progressスクリーンが現れ、スキャン動作のステータスを表示します。

Scan Progress	
Current scan info	mation
File name:	D:\\US45102908_SLOT03_S01.tif
Barcode:	16011535030005
Slot number:	3
- Status	
Scanning	×
Progress	
Scan progress	20 %
Carousel progre:	ss 0 % (0 of 1 slides completed)
🔲 Ium laser p	ower off after the current carousel run completed
View <u>B</u> ep	ort Close

Figure 18 Scan Progress スクリーン

Scan Progressスクリーンを見ることで、スライド1枚あたりどれだけスキャンが終了しているか (Slide Progress)、カローセルのなかで何枚のスライドがスキャン済みか(Carousel Progress) を確認できます。

61 x 21.6 mm スキャン範囲のスキャン時間は10ミクロンの分解能で約8分、5ミクロンの分解能で Double Passの場合は、約15分かかります。各スキャンが終了すると、データは自動的に保存先へ 送信されます。(詳細は 英文マニュアルの76ページ"Change the storage directory and check drive capacity" をご参照ください。

スキャン時間はスキャナがスライドフォルダを取り込み、スライドをスキャンし、フォルダをカローセルに返すま での一連の作業を意味します。

2. (オプション) データが指定のディレクトリに保存されると、Agilent Feature Extraction Software あるいはその他のプログラムでスキャンイメージを開くことができます。

3. スキャンが終わると自動的にCarousel Report Log が表示されます。

Carousel Report Log にはカローセルラン中のスキャン情報とエラーメッセージが表記されます。



Figure 19 Carousel Report Log

<注意>

スキャナが作動中は、スキャンデータファイルをコピーしたり、移動したりすることは**絶対に**しないで 下さい。この操作により、スキャンデータが消失してしまいます。

カローセルランが終了すると、Statusの項目に "Carousel run completed." と表示されます。この 表示が出た後は、カローセルやスライドフォルダを移動できます。また、

C:¥Program Files¥Agilent¥Microarray¥CarouselRunLog.htm

にCarousel Report Logが自動的に保存されます。このLogファイルはカローセルランごとに上書きされていきます(ファイルには常に1つのLogファイルしか保存されていない状態です)。

スキャンを中止する

カローセルラン中に、スキャンを中止することができます。

• Scan Progressスクリーンの Abort ボタンをクリックします。以下のメッセージが表示されます。

Carousel	Run Abort			×					
?	Do you want to Click Yes to ab Click No to abo Click Cancel to	o abort the caro ort the run now, rt the run after cancel your abo	usel run now? ; you may lose s the current sca ort command.	scan data. n is done.					
<u>Yes</u> <u>N</u> o Cancel									

Figure 20 Carousel Run Abort メッセージ

スライドフォルダをカローセルに戻すため、スキャン終了には数分かかります。

スキャン中止前のデー	-タは指定ディレクトリ	に保存されています。
------------	-------------	------------

Scan Progress	
Current scan info	rmation
File name:	D:\\US45102908_SLOT03_S02.tif
Barcode:	16011535030005
Slot number:	3
Status	
Carousel run at	ported. Operation aborted by user.
Progress	
Scan progress	17%
Carousel progre	ss 100 % (1 of 1 slides completed)
—]
L <u>T</u> urn laser p	ower off after the current carousel run completed
View <u>R</u> ep	ortClose

Figure 21 スキャン中止後のScan Progress スクリーンの例

Step 9—ファイルの移動とデータの保管

スキャンが終了すると、ファイルを他のディレクトリに移動することができます。 ファイルは以下の4箇所に保存できます。

- アウトプットで指定したローカルドライブの指定ディレクトリ
- ローカルドライブのほかのディレクトリ
- ネットワーク上のディレクトリ
- CDまたは、DVD

アジレントではまずローカルドライブにファイルを保存することをお勧めしています。これはネットワークに問題があった時のファイルデータ消失を防ぐためです。

データを定期的に移し変えないと、ハードディスクの残り容量が少なくなります。ディスクがいっぱいになるとスキャンは中止され、エラーメッセージが表示されます。

Step 10—システムを終了する

レーザーの耐用期間を有効に使うため、4時間以上スキャナを使用しない場合は、レーザーをオフに するか、スキャナの電源を切ってください。

- 1. スキャナコントロールのソフトウェアを終了します。
- 2. マイクロアレイスキャナの電源を切ります。
- 3. 最後にPCの電源を切ります。

付録— TiffSplitterの使い方

アジレントのスキャナを用いてアレイのイメージを読み取ると、基本的に16bitのtifファイルが作成され ます。このtifファイルは2色合成イメージであるために、アジレント社以外のソフトウェアで読み取ること が出来ません。アジレント社以外のDNAマイクロアレイを読み取ったイメージを、他社のソフトウェアで 読み込み、数値化を行うためには、2色のイメージを1色のイメージに分割する必要があります。 アジレントのスキャナには、イメージを分割するためのソフトウェア(TiffSplitter)が標準装備されていま す。これを用いると、2色のイメージを1色のイメージに分割するだけではなく、イメージの反転、回転も 同時に行なうことが出来ます。

1. TiffSplitterの起動

Windows画面左下のStartボタンをクリックし、Programs→Agilent-Life Sciences→TiffSplitterを 選択します。画面上にTiffSplitterのアイコンがある場合は、アイコンをクリックして起動することもで きます。



2. Browseボタンを押して、分割したいtifイメージファイルを選択します。

fSplitter: Agilent's Tiff Image	Channel Splitter Utility (A.6.1.)	1)
- Agilent Two Channel Tiff / Tiffs		
		Browse
Impachianin ulation		
magemanipulation	C Eliz aluna main diagonal	
🔲 Flip/Rotate Image	 Filp along main diagonal 	₩¥
	C Flip along opposite diagonal	
Output Directory		
🔲 Write Split tiffs back to input	directory	
		Browse
Split Tiff	Cancel	

3. 分割したイメージを裏返してかつ反転させたいときには、ImageManipulationの Flip/Rotate Imageの項目をチェックし、かつ反転の向きを選択します。

アジレントのスキャナは、アレイを裏側から、バーコードを左においた状態でスキャンをするので、 読み取ったイメージも、この向きになっています。例えば、このイメージを、アレイを表側にしてバー コードが下側になる向きに反転/回転させたい場合は、Flip along opposite diagonal の方をチェ ックします。

- 4. 分割したファイルをOutputするディレクトリを選択します。もとのイメージファイルと同じ場所に保存 したい場合は、Write Split tiffs back to input directoryの項目をチェックします。
- 5. Split Tiffボタンをクリックします。1色づつに分割されたイメージファイルが、指定されたディレクトリ に保存されます。ファイル名は自動的に付けられます。分割する前のファイル名に、それぞれ、緑 色のチャンネルのデータはGreen、赤色のチャンネルのデータにはRedが接頭語として付加されま す。