



本ユーザーズガイドは2100 エキスパート ソフトウェアの簡易版取り扱い説明書です。 詳細機能については **"Agilent 2100 Bioanalyzer** 2100 expert Guide(PN ; G2946-90004)"を ご覧ください。2100 expert ソフトウェアのHelpメニューの Contexts and Index からご覧いただけます。

ラボチップの調製については、各キットに添付のReagent kit をご覧ください。

## Ver.08.16.13







はじめに~2100 エキスパート ソフトウェアの構成 3
1. バイオアナライザの準備
2.分析を実行する
3. データを表示する 8
・ゲルイメージ
・エレクトロフェログラム10
4. データを解析する 12
・サブタブの紹介12
・Peak Table サブタブ13
・Region サブタブ16
・Results サブタブ18
・Fragment サブタブ20
5. 分析設定を変更する
・セットポイントエクスプローラー機能21
・マニュアルインテグレーション機能23
・任意のレーンだけ抽出する(SaveSelectedSample)
6. データのアウトプット
・印刷機能
・画面のコピー27
・転送(Export)
7. チップ間比較(Comparison コンテキスト)
8. ハードウェア診断
付録1.フラグ機能
付録 2 . Support Package





## はじめに~ **2100エキスパートソフトウェアの画面表示**

タイトルバー												
メニューバー	2 <sup>9</sup> 2100 expert	Country Party (NDW) (MR 1580	Gene DNV	7500 Curl	ani,2100 asp	nt_DNR.7500_	0000_2005-0	6-19_14-0	1-Staad			-181
	Into Central	- 🙆 Li O Se 🔚 🖬 🤞			0000	0 20	<b>1</b>					
ツールハー	Carbodz	GelCR. Mix 3a (Not R	eviewed	)		And	ynen		888888888		88888	
情報バー	(Contractor)	M si véc. S C 2100 expert_044.750_00 - R Samples	General Peo	La. P.	P. So. So.	Chir Sermary 6	Electropheni Sol - Sol - Sel	ignen Res. Sali Sali	R Magging Log Br	tecel Gtb	a	
コンテキストバ		KCR Mis Sk     KCR Mis Sk     Sangle 3     Sangle 4     Sangle 5     Sangle 6								Parenal Entergration da Entergration of Stope Thready Constitution for	at the (c)	Collepte 24 54 1,5
ツリー表示	refician Sector	- Sangle 7 - Sangle 7 - Sangle 8 - Sangle 9 - Sangle 9 - Sangle 13	10390 5000 2000 2000	=			=		- 10390 - 7000 - 2000 - 1500	Head (filter sad	Ad Pul In [c]	L.
タブ		Sample 12	2000			=	=	=-	- 1000 - 700			
サブ-タブ	Sinten		300					==	- 300			
			50	=					- 50			
Lower Panel			**	Size (ba) 90 74 11.2 195	Canc. [19]44 8,30 8,40 1,92 0,95	Polarty [coul/] 201.0 129,1 25,1 0,5	Obervations Long Policy	Circle Coport Config Coport	an Cokens	0440		
ステイタスバー			6 7 5 Tesubs	215 281 310 412 724 book Toble	2,56 1,75 1,98 2,94 7.44 Erres	30/0 9/0 9/1 20,1 20,1	9 7 7	Marya Marya Panas Product	ly Set Lover Marke ly set Upper Marke r Pods	r •	<u>.</u>	a Help
	-		-					<b>餐</b> 种	edvanced as Advan	and Openator	[1 001	DOOT BAD NO
					t	ヱットポ	イント	・エク	スプロ・	ーラー		

## 重要 2100 エキスパートソフトウェア の構成

2100 エキスパートソフトウェアは次の6つのコンテキストから構成されます。



- ・*Instrument* コンテキスト -- 分析スタート、ストップなどの本体制御画面
- ・Data コンテキスト -- データを見る、ピーク設定を変更するなどの分析画面
- ・Verification コンテキスト -- 稼動性能適格性確認試験(OQ)のための
- ドキュメント作成画面 ・ Comparison コンテキスト -- 複数にわたるデータを比較分析する画面
- ・Assay コンテキスト 解析設定をカスタマイズしたassayファイルの作成画面
- ・Systemコンテキスト -- ファイル名やデータ保存先のなどの設定変更画面

コンテキストの切り替えは、コンテキスト バー(下図参照)のいずれかの絵をクリックすることで行えます。



2100Expertガイド

# 2100 exp

# 1. バイオアナライザの準備

 (1) デスクトップのアイコンをダブルクリックし、2100 エキス パートソフトウェアを立ち上げてください。下記の画面 (Instrument コンテキスト) が表示されます。



(2) コンテキストバーから"Instrument" コンテキストを選択してください。

(3) 複数台バイオアナライザを接続し ている場合、ツリー表示から装置を選 択 してください。



(4) ソフトウェアがバイオアナライザ本体を正常に認識しているかどうか、画面 中央のアイコンで確認してください。



本体を正常に認識していま す。蓋が開いている状態で す。



本体を正常に認識して います。



本体を認識していません。下図をご参考ください。

トラブルシュート1 バイオアナライザ本体が接続されていない場合

- ・バイオアナライザ本体の電源が入っているかどうか確認してください。
- ・PCと本体のケーブルが正しく接続されているかどうか確認してください。
- ・下図に従い、COMポートの設定を行ってください。



COM PORT欄のプルダウンか ら、数字(1,2,もしくはその他の 数字)を選択してください。 (Demoを選択するとオフライン 用となり、測定できません。)





## 2. 分析を実行する

(1) InstrumentタブのAssay ボタンから 実行したいAssayを選択します。



ご使用のキットやサンプルタイプに応じて、適切なアッセイを選択してから次のステップに進んでください。

#### DNAアッセイ 5種類

フォルダ	アッセイ名	使用キット	対象サンプル	備考
dsDNA	DNA 1000 Series II	DNA 1000 キット	dsDNA 25-1000 bp	
	DNA 7500 Series II	DNA 7500 キット	dsDNA 100-7500 bp	
	DNA 12000 Series II	DNA 12000 キット	dsDNA 100-12000bp	
	DNA 12000 Laddering Series II	DNA 12000 キット	dsDNA 100-12000bp アポトーシス細胞からの DNAサンプル	ノイズピーク除去アル ゴリズムを含みます
	High Sensitivity DNA	High Sensitivity DNA キット	dsDNA 50-7500 bp	

#### RNAアッセイ 9種類

フォルダ	アッセイ名	使用キット	サンプルタイプ	備考
RNA	Eukaryote Total RNA Nano Series II	RNA 6000 Nano キット	真核生物由来 totalRNA	RIN計算
	Prokaryote Total RNA Nano Series II	RNA 6000 Nano キット	原核生物由来 totalRNA	RIN計算
	Plant RNA Nano	RNA 6000 Nano キット	植物由来 totalRNA	RIN計算
	mRNA Nano Series II	RNA 6000 Nano キット	mRNA, cRNA	(RIN計算できません)
	Eukaryote Total RNA Pico Series II	RNA 6000 Pico キット	真核生物由来 totalRNA	RIN計算
	Prokaryote Total RNA Pico Series II	RNA 6000 Pico キット	原核生物由来 totalRNA	RIN計算
	Plant RNA Pico	RNA 6000 Pico キット	植物由来 totalRNA	RIN計算
	mRNA Pico Series II	RNA 6000 Pico キット	mRNA, cRNA	(RIN計算できません)
	Small RNA Series II	Small RNA キット	totalRNA, smallRNA抽出物	( <sub>RIN</sub> 計算できません)

#### Proteinアッセイ 3種類

フォルダ	アッセイ名	使用キット	サンプルタイプ	備考
Protein	Protein 80 Series II Protein 230 Series II	Protein 80 キット Protein 230 キット	5 - 80 KDa 14 - 230 KDa	
	High Sensitivity Protein 250	High Sensitivity Protein 250 キット	10 - 250 kDa	





(2) バイオアナライザの蓋を開け、正 しく調製されたラボチップを台座に乗 せます。





(3) 蓋を閉めた後、ソフトウェア上のアイコンがラボチップの 絵に変わることを確認して下さい。選択したアッセイによってラボチップの色は変わり ます





トラブル ; 調製済みのラボチップを入れたにも 関わらず、アイコンがラボチップの 絵に変わらない場合

ラボチップの調製に問題があります。 (液量が少ない、気泡があるなど) 再度調製を確認してください。

#### (4)オプション ; データ保存場所、ファイル 名、サンプル数、サンプル名を確認・記入し ます。

Γ	instrument Diagnostics Instru	ment Logbook			
		Name:		COM Port	Demo 💌
	13 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0	Serial#:		Assay Sel	ection: À Assavs
		Vendor:		Start/Stop	p Run: C Start
	DNA Chip On-Chip Dechaptoresis	Product ID:			Juit
		Firmware:	Simulation Mode	Assay File	:: <u>C:¥nt¥2100 bioa</u>
				Data File:	
10	Destination and Data Acquisitio	n Parametere			
K	Default     C:\Pr	ogram Files (x86)*	Agilent\2100 bioanalyzer\2100 exp	ert\Data\2013-07	<b>a</b>
	C Custom C:\Pr	ogram Files (x86) <sup>s</sup>	Agilent\2100 bioanalyzer\2100 exp	ert\Data\2013-07	-26
	File Prefix 2100	expert	(max 16 characters)		
	Run sample	1 to   1	2 렆		
	Chip Summary				
	Sample	e Name	Sample Comment	Rest. Digest	Observation
	Ladder 1			V	
u	Ladder 2			V	
	Ladder 3			V	
	DNA 1000 ladder			V	
u	Ladder 5			V	
	Ladder 6			V	
	Ladder 1			V	
Ш	Ladder 2			V	
	Ladder 3			V	
K	DNA 1000 ladder				
10	🔷 10 H			- ex 🏊	

-----

(5) Startボタンを押してください。





## **Agilent Technologies**

6



(6) 分析が始まります。スタート開始後、数分は電圧チェック、温調、 フォーカシングなどの初期化を行います。

#### 注意;

分析中は装置に振動を与えないでください。 装置のランプが点滅している間は蓋を開けないでください。



(7) 分析が終わると、自動的に初期画面にもどります。ラボチップをバイオア ナライザから取り出し、クリーニングチップで電極洗浄を行ってください。 (電極洗浄については各Reagent kit guideを 参照ください。キットごとに洗浄方法が異なります。)



Agilent Technologies

7





(1) コンテキストバーから"Data"コンテキストをクリック してください。



(2) Fileメニュー> Open(あるいはツー ルバーのOpen;右図)を選択し、フォル ダの中からファイルを選びます。(測定 中および測定終了後は、この作業を行わ なくても測定データが開かれていま す。)



#### **ノート** 2100 エキスパートソフトウェアによる分析データの拡張子は"xad" です。

(3) ファイルが開きます。



データ表示は選択タブで様々に 切り替えることができます。







## 3-1 ゲルイメージ表示機能

"Data" コンテキストにてGelタブを押すと、ゲルイメージが表示されます。



Global Scale; チップのすべてのウェルのなかで最も濃いバンドを基準にして、全ウェル を同じスケールにします。

Individual Scale;個々のウェルの最適スケールにします。従って、同じチップ内でもそれぞれのウェルのスケールは異なります。

Selected Scale; 選択したウェルが最適表示となるように、全てのウェルのスケールを合せます。

The Measure of Confidence





## 3-2 エレクトロフェログラム表示機能

"Data" コンテキストにてElectropherogramタブを押すと、エレクトロフェログラムが表示されます。



## 3-2-1 データ重ね描き表示

任意のエレクトロフェログラムを重ねて表示することができます。





# 2100 expert

## 3-2-2 エレクトロフェログラムのスケール機能

エレクトロフェログラム (シングルウェル表示の時) には2つのスケール 調整モードがあります。

(1) ツールバーの中にあるScaling Mode(下図)の下向き矢印をクリックします。

망 -100 M |鎌| 😰 🎜 Overlaid Samples

(2) 2種類のうち、いずれかをクリックしてください。



Individual Scale; 個々のウェルの夫々を最適スケールにします。従って、各ウェルのスケール は異なります。

Selected Scale; 選択したウェルが最適表示となるスケールに、全てのウェルのスケールを合せます。

## 3-2-3 その他エレクトロフェログラム画面の機能





**Agilent Technologies** 

11

## 4. データを解析する



解析用途に応じて、様々なサブタブをご用意しております。





ロブロードな形状のピーク濃度・サイズを解析する場合 ・次世代シークエンスサンプル ・smallRNAサンプル ・たんぱく租精製サンプル

⇒4-2 Region Table サブタブへ



□RNAサンプルのクオリティーチェック ・totalRNA サンプル ・mRNAサンプル ・smallRNAサンプル ⇒4-3 Results サブタブへ 200 500 1000 2000 400

□RNAサンプルのrRNA解析

- ・totalRNA サンプル
- ・mRNAサンプル

⇒4-4 Fragment Tableサブタブへ





# 2100 expert

## 4-1 Peak Tableサブタブ

PCR産物などのシャープな形状のピーク濃度・サイズを解析する場合に適しています。



Lower Marker; ピークの上に緑の数字 Upper Marker; ピークの上に紫の数字

ピークが割れている例

各ピークの境目は赤い点線で示されていま す。

正しくシングルピークとして認識している ことを確認してください。特に、Upper Markerのエリアは濃度補正に使うため、正 しく認識されている必要があります。



・各ピークの積分線を変更したい場合

・目的のピークが表示されていない場合

⇒ p14 ピークがスプリットしている場合

⇒ p15 Upper Markerピークを正しく捉えていない場合





## ◆ピークがスプリットしている場合

#### a)グラフ上で右クリック> Manual Integration選択



b)各ピークの下にドット付の青いベースライン が表示されます。 削除したいベースラインのドットをクリックし ます。(選択されたドットがグリーンに変わり ます)



c)右クリックから Remove Peakにてピーク削除します。

c) ドットをドラッグで動かします。

The Measure of Confidence







#### ◆Upper Markerピークを正しく捉えていない場合



異なるピークをLower/Upper Markerと認識している場合

正しいピークを選択し、右クリック > Manually Set Upper Markerを選 択して下さい。





## 4-2 Region Tableサブタブ

(1) Electropherogramタブを開き、 Region Tableサブタブを選択します。

◆Regionタブが表示されていない場合、

b) [Global]タブを選択します。

Analysis]にチェックを入れます。

c)[Advanced] モードを選択します。

a) ViewメニューからSetpointsを選択し、セッ

トポイントエキスプローラーを表示します。

d) [Smear Analysis]の [Perform Smear





150 200 400 500 600

(3) Region位置をドラッグで調整

(4)Tableにサイズ、濃度、Molarityな どの情報が表示されます。





◆全レーン同一のRegionを設定する方法



(1) Assay Properties タブを選択

(2) Globalタブの中のSmear Analysisの下のTableをダブルクリック



(3) 表示されたボックスにRegionをAdd ボタンで追加、From,Toに任意の 数字を入力してOKボタンを押しま す。



ポイント

Region Tableで計算している濃度 は、緑の点線で示される始点・終 点を結んだ実線から上のエリアを 使用します。 右図のようにずれている場合、始 点・終点をずらして下さい。







[FU]

#### 2100Expertガイド



## 4-3 Results サブタブ

Electropherogramタブ
 を開き、Resultsサブタ
 ブを選択します。



#### □totalRNA用のアッセイでは

- ・RNA Area;面積値
- ・RNA Concentration;濃度
- rRNA Ratio;

Eukaryote totalRNAアッセイでは28S/18S Prokaryote totalRNAアッセイでは23S/16S Plant totalRNAアッセイでは25S/18S

・RNA Integrity Number (RIN); 分解の度合いを1-10の数字で表示

トラブルシュート RNA Integrity Number(RIN)がNAと表示される場合、次ページをご覧ください。

#### □mRNA用のアッセイでは

- ・RNA Area;面積値
- ・RNA Concentration;濃度
- ・rRNA Contamination; rRNAピークが占める割合(%)

#### □smallRNA用のアッセイでは

- ・Small RNA Concentration; Integration StartからEnd(緑点縦破線)までの濃度
- ・miRNA Concentration; 10-40 nt の濃度
- ・miRNA/Small RNA Ratio(%); 上記の比率

The Measure of Confidence



#### ◆RIN =NAと表示される場合



RINはHuman, Mouse, Ratのtotal RNAの泳動パターンを基に設計されており、 典型的なパターンと異なる場合警告の意味でフラグを立て、RINがN/Aとなり非表示となります。 フラグを解除していただくとRINが再表示されます。

1)フラグのたっているレーンを開き、グ ラフの下にある"Error"タブ①を開きま す。 ここにフラグ情報が記載されています。

2)セットポイントエクスプローラーを開 くために、②のボタンを左側にドラッグ します(左クリックを押したまま左側に 移動させる)

3) "Normal"モードから"Advance" モードに切り替えます

③ (タブから"Advance"を選択する)

4) スクロールを一番下まで下げます④ 5)RNA Integrity Numberの項目の中 で、フラグのたっている項目の数値を下 図のように変更します。





Error(Errorタブで表示されるDescriptionカラム)	Setpoint Explorerの項目	数値
Unexpected signal in pre-region	Pre Region Anomaly ···	1以上に変更
Unexpected signal in 5S-region	5S Region Anomaly…	1以上に変更
Unexpected signal in fast-region	Fast Region Anomaly…	1以上に変更
Unexpected signal in inter-region	Inter Region Anomaly…	1以上に変更
Unexpected signal in precursor-region	Precursor Region Anomaly…	1以上に変更
Unexpected signal in post-region	Post Region Anomaly…	1以上に変更
Unexpected baseline signal	Baseline Anomaly····	1以上に変更
Unexpected ribosomal ratio	Ribosomal Ratio Anomaly…	1以上に変更
Unexpected sample type	Unknown Sample Type…	1以上に変更
	Threshold Prerequisite Concentration	
Total RNA concentration too low		0に変更





## 4-4 Fragment Tableサブタブ



## Electropherogramタブを開き、 Fragmentサブタブを選択しま す。

各rRNAピークについて下記情報がテーブ ルで表示されます。

- Start Size
- End Size
- Area
- ・% of total RNA
- ドット付のベースラインは各rRNAの区 切り線を示しており、ドラッグで 動かすことができます。







画面上で右クリックからAdd Fragmentを選択することで、rRNA ピークを増やすことができます。





## 5.分析設定を変更する



## 5-1 セットポイントエクスプローラー機能

ピーク認識の設定値(ピーク幅、ピーク高)は、セットポイントエクス プローラーで変更することができます。(測定後のデータに関しても、設 定値を変更することができます。)

(1) コンテキストsから"Data"コンテキストをクリックしてください。

(2) 下記のうち、いずれかのタブを選択してください。

- Assay Properties Tab
- Electropherogram Tab (Single/Grid 表示)

• Gel Tab

(3) Gel タブ、Electropherogram タブを選択している場合は、 画面右端のバー(下図) をクリックしてください。 (Assay Properties タブを 選択している場合は画面右端にすでにセットポイントエクスプローラーが表示されています。)



(4) 下記のセットポイントエクスプローラーで、様々な設定値が変更できます。







#### (5) 変更したい項目に数値を入力した後、キーボードのenterキーをクリックして ください。





## 5-2 マニュアルインテグレーション機能



2100エキスパートソフトウェアの電気泳動分析では、マニュアルインテグレーションを行うことができますマニュアルインテグレーションにより、 ピークのベースラインを移動させたり、加えたり、削除したりすることができます。

## (1) ベースラインを移動する

**1** [Data and Assay]コンテキスト中の[Electropherogram]タブを選択し、エレクトロフェログラムを拡大し、目的のピークを大きく表示します。

- 2 ツールバー中の[Manual Integration]ボタン をクリックします。
- 3 ベースラインポイントを適切な場所に設定します。



ヒント

垂線に沿ってピークベースラインポイントを移動させるためには、CTRLキーお よび左のマウスボタンを押してドラッグしてください。エレクトロフェログラ ムに沿ってピークベースラインポイントを移動させるためには、左のマウスボ タンのみを押してドラッグしてください。





ß



#### (2)ピークの追加

エレクトロフェログラム上で右クリックし、表示メニューより[Add Peak]を選択 します。



## (2) ピークの削除

ベースラインポイントを右クリックし、表示メニューから[Remove Peak]を選択します。2つのベースラインポイントとそれを結ぶ線が消えます。





2100Expertガイド

ШШ

2100 Expert - C:¥Program Files (x86)¥Agilent¥2100 b

Windows

Help

Ц

II.

٦

Ctrl+O

Ctrl+S

All Data Files Demo DNA 1000 Series II.. All Samples 1: Ladder 1 2: Ladder 2

3: Ladder 3
4: DNA 1000 ladder

- 5: Ladder 5 - 6: Ladder 6 - 7: Ladder 1 - 8: Ladder 2 - 9: Ladder 3 - 10: DNA 1000 ladder - 11: Ladder 5 - 12: Ladder 6 - 12: Ladder

File Context View Electropherogram

Save Selected Sample...

## 5-3 任意のレーンだけ抽出する(SaveSelectedSample)

Instrumer Data Comparison Assay

À Open...

Close Close All

E Save

Revert
Save As...

1880 E

Import...



分析した全レーンの中から、あるレーンだけを抽出して保存したい、 余分なレーンを外したい、という場合下記の方法でデータを抽出し て別名保存することができます。

- (1) Dataコンテキストで該当のデータ を開きます。
- (2)File メニューから Save Selected Sample を選択します。

- (3) データ内のサンプルが表示されま す。抽出したいサンプルにチェックを 入れ、Applyボタンを押します。
- Save Selected Samples ... Data File: Demo DNA 1000 Series II2.xad Location: C:¥Program Files (x86)¥Agilent¥2100 bioanalyzer¥2100 expert¥data Catego ID Sample Name Com Ladder 1 V 1 Sample 2 Ladder 2 Sample 5 3 4 5 Ladder 3 Sample 5 DNA 1000 ladder Sample Sample Ladder 5 Ladder 6 Sample 6 7 Ladder 1 Sample 8 Ladder 2 Sample Г 9 Ladder 3 Sample 10 DNA 1000 ladder Sample 11 Ladder 5 Sample 12 Ladder 6 Sample V 13 Ladder Ladder Cancel Help Apply

Save selected samples

(4) 名前を付けて任意の場所に保存します。





## **Agilent Technologies**

25

2100Expertガイド



## 6. データをアウトプットする

#### 6-1 印刷機能

(1) コンテキストバーから"Data " コンテキス トをクリックしてください。

(2) 複数のデータが開かれている場合、 印刷したいデータを ツリー 表示 パネルで選択してくださ い。



- (3) FileメニューからPrintを選択してください。
- (4) 下記のダイアログが現れます。印刷したいアイテムにチェックを入れ、 Printボタンをクリックしてください。

Print Item	🗖 Electropherogram	🗖 Calibration Curve		印刷内容を選択
🗖 Assay Details	🗖 Gel Like	🗖 Standard Curve		
	🗖 Result Tables			
-Wells		Options		
🖸 Ali Wells		1 per page		
O Wells		Exclude Marker		
Enter sample separated by	number and/or sample ranges, commas. Example: 1,2,3-6	🗖 Include Ladder		
- Save To File				
🗖 PDF 🛛 File Path:	C:¥izing_DNA-7500_00000	0_2003-01-07_17-54-12.r		出力形式を選択 (PDF/HTML)
🗖 HTTML 🛛 File Path:	C:¥zing_DNA-7500_00000	_2003-01-07_17-54-12.F	1	
Page Setup Printer.	. Preview Cancel	Help Print	1	
eviewをクリック イアウトを予め	クし 確認			







測定データ画面を簡単にコピーすることができます。

6-2画面のコピー

Electropherogram > Copy Electropherogram を選択し、図を貼り付ける先のアプリ ケーション上にペーストします(Controlを押しながらVを押す)



Gel > Copy Gel を選択し、図を貼り付ける先のアプリケーション上にペーストします(Controlを押しながらVを押す)









## 6-3 データを転送(Export)する

- (1) コンテキストバーから"Data and Assay" をクリックしてください。
- (2) 複数のデータが開かれている場合、転送したいデータを ツリー 表示 パネルで選択してください。



- (3) FileメニューからExportを選択してください。
- (4)下記のダイアログが現れます。転送したいアイテムと形式に チェックを入れ、保存先を選択した後、Exportボタンをクリックしてください。







# 7. チップ間比較(Comparisonコンテキスト)

Comparisonコンテキスト画面では、複数のデータから任意のウェルを選択し、比較表示したり、ファイルに保存することができます。

(1) コンテキストバーから" Comparison"コンテキストをク リックしてください。



(2) Fileメニュー> Open(あるいはツールバーのOpen; 下図)を選択してください。 表れたダイアログボックスにて、フォルダの中から比較したいデータファイル(.xad) をいくつか選び、Openボタンをクリックしてください。Dataコンテキストで開いてい るデータは再度開く必要はありません。



(3) 開いたデータファイルは、画面中央 下の"Select Data Files"にリストされ ます。





## **Agilent Technologies**

29





、、、 を選び、**右クリック**を押してください。

(6) "Add Sample to New Comparison File" をクリックします。

🏹 🖴 🛛 Select Data Files		а 
Demo DNA 1000 Serie	es II 👻	
Ladder 1		
: Ladder 2		
3: Ladder 3		
4: DNA 1000 ladder		
i: Ladder 5	Ade	d Sample to New Comparison File
i: Ladder 6		
7: Ladder 1		
3: Ladder 2		
9: Ladder 3		



新しくComparison ファイルが作 成され、選択されたウェルがツ リー表示に表示されます。



30



(5) Select Data Filesリストから、比較したいウェルを次々に足していきます。

追加する場合は" <b>Add</b>	Select Dat	aFiles	
Sample to	Demo DNA 100	) Series II 🔻	
Comparison File"をク	1: Ladder 1		
リックします。(新しく Comparisonファイルを	2: Ladder 2 3: Ladder 3		
作成したい場合のみ"Add	4: DNA 1000 ladder	r 🛛	
Sample to <b>New</b>	5: Ladder 5		
Comparison File"を選択	6: Ladder 6		Comparison Comments :
します)	8: Ladder 2		
	9: Ladder 3		
	10: DNA 1000 ladd	Add Caraala	ta Campanisan Fila
	11: Ladder 5	Add Sample	to comparison File
	12: Ladder 6	Add Sample	to New Comparison File

L: Ladder

#### ノート

違うアッセイで得られたデータを比較することはできません。(例えば DNA7500アッセイで得られたデータと、DNA1000アッセイで得られたものを 比較することはできません。)違うアッセイで得られたデータを選択した場合、 下記のメッセージが表示されます。

-	2100 Expe	rt	
-	▲	Assay class does not match. Please choose a sample class DNA 1000 Series II	with assay
l			ОК

(6) Comparison ファイルからウェルを削除したい場合、ウェルを選択して右クリック後、Delete Sample From Comparison Fileを選択してください。





(7) Comparison ファイル内のエレクトロフェログラムを重ね描きする場合は、下記 を実行してください。

①Electropherogramタブを選択します。

②ツールバーのOverlaid Samplesから、重ね描きしたいウェルを選択します。



(8) 作成したComparisonファイルを保存する場合、FileメニューからSaveを選択してください。

ノート

Comparisonファイルの拡張子は"xac" です。 デフォルト名は分析の種類に由来します。(「ComparisonFileX[分析の種 類].xac」、Xは自動的に与えられた番号です。) 例: 「ComparisonFile 0 Protein 200.xac」







## 8. ハードウェア診断

Agilent 2100 バイオアナライザシステムには、ソフトウェア中にハードウェアの診断ツールが用意されています。 この診断ツールによりユーザーご自身でバイオアナライザ本体装置の状態についてのチェックを 行うことが可能です。

診断ツールのテスト結果は、" passed "もしくは" failed "で表示されます。" failed "は、不完全なハード ウエアコンポーネンツの存在を示しております。この結果が出た場合は、弊社サポートまでお問い合わせく ださい

#### 【ご用意いただくもの】

#### ① 未使用のラボチップ 1個

RNA, DNA, Protein ラボチップのいずれか。埃の入らない環境で保管いただければ、次回の測定に使用できます。

#### ② テストチップセット

型番;G2938-68100もしくはG2938-68300 1セットは装置に付属しております Expire Date(使用期限)が銀色の袋のシールとテストチップに記 載されています。期限内であることをご確認ください。



## ハードウェア診断の操作手順

(1) コンテキストバーから"Instrument"コンテキストを選択してください。

(2) 複数台バイオアナライザを接続している場合、ツリー表示から診断したい装置を選択してください。

(3) Diagnosticsタブを選択してください。



ノート

Diagnosticsタブは装置とソフトウェアが正常に通信されていない場合、 選択できません。事前に装置電源が入っているか、接続ケーブルが適切 につながっているかどうかを確認してください。

ノート

2100エキスパートソフトウェアが測定を行っている間は、ハードウェア 診断を行うことはできません。





#### (4) Diagnosticsタブにて、診断したい項目のApplyボックスにチェックを 入れて下さい。

Instru	ment	Diagnostics				
lame :	DEL17	00058		Firmware : C.01.03F		
ierial 4	¢: DEL1	700058		Product ID : G2938B		
Aval	lable T	ests:				
$\square$	Apply	Name	Description	Stabus	-	Start
	×	Communication Test	Tests if communication with instrument is funct	🐺 Selected		JACAN
	×	Electronics Test	Tests instrument electronics.	Selected		Select All
	×	Fan Test	Tests if instrument fan is working.	Selected		
	×	Ud Senfor Test	Tests if the lid sensors are working.	Selected		Unselect Al
	×	Temperature Test	Checks if the temperature sensors and heater	😴 Selected		E-law.
	×	Stepper Motor Test	Tests if horizontal and vertical motors are wor	Selected		14899
	X	Electrope / Diode Test	Tests conductivity of channels (pin to pin).	Selected		
		Utab Valence Shability Test	Tests bisk rollings, service rand stability	Salastad	-	

	Name	主なチェック項目
1	Electronics test	電源ボードのチェック
2	Fan test	ファン機能チェック
3	Lid sensor test	Lid(蓋) センサーチェック
4	Stepper motor test	Stepper Motor稼動チェック
5	<i>Temperature test</i>	チップ台座の温調機能チェック
6	HV Stability and Accuracy Test	16個の高圧電源の精度と安定性をチェック
7	HV accuracy test (on-load)	リファレンスチャネルを使用した高圧電源コ ントローラーチェック
8	Short circuit test	漏れ電流チェック(電極に水分がついている 場合や、温度25℃相対湿度60%以上の部屋の 場合、このテストがfailします)*
9	Electrode diode test	電極ピン間の伝導度チェック
10	Optics test	LEDとレーザーのDark Current値をチェック
11	Electrophoresis autofocus test	レーザーのフォーカスと強度チェック
12	Laser stability test	レーザーの安定性チェック

注)湿度が高い部屋では、漏れ電流値は高くなります。適切な結果を得るために、室温25℃ 相対湿度60%以下で実施ください

(5) Startボタンを押して下さい。

(6) 診断が始まります。表示されるダイアログボックスの指示に従い、各ハードウェア診断項目を 進めてください。

2122100 exper					
File Context	<u>View Assers Tools Windows</u>	Help			
Instrument	• 👌 🛐				
Contexts	DE11700058 - mRN/	\ Pico			
	🕼 Al Instruments	Instrument	Diagnostics		
Instrument	DELL/20058	Name : DEL170	0158		
14	Bioanalyzer Diagnose	Sena 4: DCI1)	00056	×	1
Data and Assay	Please inset	pl	Des		
<b>1</b>	Close the lid		ts if communication		
	<b>V</b>				ts instrument electr
					ts it instrument tan
100					cks E the terrorat
				Lancel	ts if horizontal and
Comparison		×	Electrode / Diode	Test Ti	ests conductivity of d
		Test Proper	loak och car erd ties	/ Ann 7	ente konk ondrene ente
		1D:	4		
		Barner	Terroration Te	**	







## 各項目における手順

1.Electoronic Test 2 Fan Test	Notes that the second s	hardware test Please insert cartridge! Close the lid and press OK!	×.
上記の画面が現れます。 (1) 装置本体の蓋を閉めてく (2) 画面の"OK"ボタンを押 3 Lid Sensor Test	ください。 してくださ		OK Cancel
	👈 Bioanalyzer	hardware test	x
	2	Please open the Lid!	

上記の画面が現れます。

- (1) 装置本体の蓋を開けてください。
- (2) 画面の"OK"ボタンを押してください

#### **4 Stepper Motor Test**

この項目の操作は不要です。 自動的ソフトウェアが診断を進行します。

#### **5** Temperature Test



ΟK

Cancel

#### 上記の画面が現れます。

- (1) 空のラボチップをバイオアナライザにセットしてください
- (2) 装置本体の蓋を閉めてください。
- (3) 画面の"OK"ボタンを押してください

The Measure of Confidence





## 6 HV Stability and Accuracy Test

この項目の操作は不要です。 自動的ソフトウェアが診断を進行します。

8 Short Circuit Test



lectrode diode test	🔧 Bioanalyzer	hardware test		×I
Optics Test	2	Please insert Electrode Cartridge and the Electrode/Diode Test Chip! Close the lid!		
		1	OK Cancel	
<ul> <li>(1)Electrode/Diode testチップをバイ</li> <li>(2)装置本体の蓋を閉めてください。</li> <li>(3) 画面の"OK"ボタンを押してください</li> </ul>	オアナライ ハ	イザにセット	トしてください	$\left \right $
名前を確認ください ―――				
期限内であることを確認くだ			TestChip Electrode/Diode	
Electrophoresis autofocus te	est			
Laser stability test		Bioanalyzer har	rdware test sase insert Electrode Cartridge and Autofocus Test C d enter chip values! sse the lid!	<b>≥</b> hip
(1)ト記のダイマログボックフロー Auto	focus			

11 Electrophoresis autofo

- 12 Laser stability test
  - (1)上記のダイアログボックスに、Autofocus testチップ情報を入力してください。各 Autofocus testチップにはそれぞれ Offset値とIntensity値が書いてありま す。その値を画面に入力してください。 (2)Autofocus testチップをバイオアナライザに セットしてください
  - (3) 装置本体の蓋を閉めてください。
  - (4) 画面の"OK"ボタンを押してください

名前を確認ください 期限内であることを確認ください

000 OOC**Test Chip** Autofocus Exp.xxxxxx Lot xxxxxx Offset\*\*\*\* Intensity\*\*

Offset: 13

Intensity:

620

Cance

Offset値とIntensity値 は各チップ固有の値です



#### (7) 各診断項目のStatus欄には、テスト結果が表示されます。



/		
Description 🚺 Status 🍾	<ul> <li>Stop</li> </ul>	- Executing (谁行中)
mnunication with instrument is fun <mark>it 🥑</mark> Executed, passed		
ument electronics. 🚽 🚽 Executed, passed	Select	- Execution pending(ペンディング状態)
trument fan is working. 🔰 🚽 Executed, passed 🚽		
elid sensors are working. 🚽 Executed, passed 💡	Unselec	- Executed, passed (診断に八人した状況)
he temperature sensors and heater 🔨 Executing	Dilour	
rizontal and vertical motors are wo <mark>r</mark> 🚱Execution pending	19397	- EXecuted, Talled (診断に異常値か見つかつに状況)
uctivity of channels (pin to pin). 📩 🍋 Execution pending 🛛 🧎	_	
sultan anna an an an taitiste 🥄 🖓 🖓 🖓 🖓 🖓 🖓 🖉	-	

- (8) "Failed"と表示された項目に関しては、再度診断を行ってください。
- (9) 再度 "Failed"と表示される項目が残っている場合、下記のファイルをメールに添付の上、

#### email\_japan@agilent.com

にお送りください。

ハードウェア診断ファイル;拡張子.xdy files 場所; Local Drive内の ¥Program Files¥Agilent¥2100 bioanalyzer ¥2100 expert¥diagnosis

t¥2100 bioanalyzer¥2100 expert¥diagnosis					
	名前 △	サイズ	種類	更新日時	
	🖻 Diagnosis_28-07-2004_14-00-06.xdy	71 KB	XDY ファイル	2004/07/29 17:40	
	Diagnosis_28-07-2004_13-00-31.xdy	86 KB	XDY ファイル	2004/07/29 17:40	
diagnosis	Diagnosis_28-07-2004_13-02-37.xdy	57 KB	XDY ファイル	2004/07/29 17:40	
	🖻 Diagnosis_28-07-2004_13-08-09.xdy 👘	98 KB	XDY ファイル	2004/07/29 17:40	
オブジェクトを選択すると、その説明が表 示されます。	Diagnosis_28-07-2004_12-56-09.xdy 🖻	62 KB	XDY ファイル	2004/07/29 17:40	
関連項目: マイドキュメント マイネットワーク					
<u> マイコンピュータ</u>					







## 付録1.フラグ機能

say Properties	Chip Sun	mary Gel	Electrop	herogram	🖉 Re	sult Flaggi	ng Log	Book		_	_		
[bp] La	ider Ladde	r 1 Ladder 2	Ladder 3	DNA 1	Ladder S	Ladder 6 I	adder 1	Ladder 21	adder 3	DNA 1	Ladder 5	Ladder 6	[bp]
				-	-		-	-	-	-	-	-	
1500 -				_		_							- 1500
1000 - 700 -	==		—	_	_	_	_		_	=	_		- 1000
500 -						_						_	- 500
400 -	_	_											- 400
300	_	_								_			- 300
200 -		_											- 200
150 -	-												- 150
100 - 50 - 15 -													- 100 - 50 - 15
	L 1	2	3	4	5	6	7	8	9	10	11	12	

#### フラグ機能では

試料に対してユーザー定義の色分け設定を行うことがで きます。この機能を使えば、特定の特徴をもつ試料を簡 単に同定することができます。

フラグ則は、「Result Flagging ]タブ上で定義でき ます。 このタブは[Data and Assay]コンテキスト上で電 気泳動測定データファイル(.xad.)またはアッセイ ファイル(.xsy)を選択しているときに使用できま す。



下記の2つのモードがあります。 このモードでは、あらかじめいくつかのフラ グ条件がある程度設定されております。 あとは基準値を入力するだけです







Editor Mode このモードでは、よりフレキシブルにフラグの定義を行うことができます。複 雑な演算をおこなう場合には このモードを使用します。



一度設定したフラグ条件は、

- ・Save Resultsボタンから名前を付けて保存できます。保存形式は.xmlです。
- ・Import Rules from Fileから 取り込むことができます。



# 2100 expert

## 例1)totalRNAアッセイでRINの値によって色分けする

totalRNAアッセイでは、Form Mode の'Search RIN'が適応されています。	Image: Select Form       Image: Se
デフォルトでは、 ・RIN1以上のレーン ; 薄いブルー という条件のみが適応されています。	1. Specify minimal       Search Pattern         Set primary threshol       Search Fragments in Range         Search Fragment with Concentration       ✓         Search RIN       ✓
Assay Properties       Chip Summary       Gel	3. Specify results Samples with RIN equal to or h Shall be colored Change Shall be labelled RIN:
例えば ・RINが9以上は青 ・RIN 8-9 は緑 ・上記以外はオレンジ と表示するには、下記のように設定します。	L 1 2 3 4 5 6 7 8 9 10 11 12
<ol> <li>1) 1.Specify minimal RIN value Set primary threshold の設定値を9</li> <li>2) 2.Specify options Set secondary threshold をONにし、設定値を8</li> </ol>	Set primary threshold       9.0       (1.0 <= Value <= 10.0)
3) 3. Specify resultsのShall be coloredの項目からそれぞれ上記の色 を決定	Shall be tabeled All Other Samples
<ul> <li>4)実行ボタンを押します。</li> <li>結果がゲルイメージ上に表示されます</li> </ul>	
	L 1 2 3 4 5 6 7 8 9 10 11 12







## 例2)DNA・Proteinアッセイで目的のサイズの濃度によって色分けする

例えば、DNAアッセイにて目的の400-500bpのピークが

- ・検出されていれば緑
- ・10ng/ul以上あればオレンジ
- ・それ以外であれば赤
- と表示するには下記のように設定します。
  - 1) Form Modeを選択します。
  - 2) Select Formから'Search Fragment in Range'を選 択します。

Select Form Search Fragment Search Fragment with Purity Search Fragment with Purity Search Fragments in Range

- 1) 1.Specify search rangeにて、
  Set lower limit; 400 bp(任意の値)
  Set upper limit; 500 bp (任意の値)
  2) 2.Specify options
  Set secondary threshold をONにし、
  設定値を10 ng/ulに設定(任意の
  値)
  3) 3.Specify resultsのShall be coloredの項目からそれぞれ上記の
- 色を決定
- 4)実行ボタンを押します。





結果がゲルイメージ上に表示されます





# 付録2. Support Package

ソフトウェア機能に異常がある、動作が不安定、という症状の場合、テクニカルサ ポートスタッフが状況を把握するため、Support Packageファイルを要求すること があります。

## Support Package 作成手順

- 1) 2100 expertソフトウェアのHelpメ ニューから Create Support Packageを 選択します。
- もしくは、WindowsのStartメニューから All programs>Agilent 2100 Bioanalyzer>Utilities>Create Support Packageを選択してください。
- 2) 以下のダイアログが表示されますので、全 ての項目を選択し、 "Collect"をクリックして ください。



💱 File Collect		×
- Task		
<ul> <li>Diagnose Hang Mode</li> </ul>		
- Task		
<ul> <li>Trace Files</li> </ul>		
- Task		
<ul> <li>Application Minidumps</li> </ul>		
- Task		
<ul> <li>Instrument Diagnostics</li> </ul>		<b>I</b>
– Task		
Instrument Packets		S .
- Task		
<ul> <li>System Information</li> </ul>		S .
– Task		
<ul> <li>License Information</li> </ul>		Sector 1
– Task		
System File		S .
- Task		
<ul> <li>List Installation Files</li> </ul>		
- Task		
<ul> <li>Registry Information</li> </ul>		<b>I</b>
- Task		
System Event Logs		<b>I</b>
- Task		
<ul> <li>Logged on User Information</li> </ul>		
– Task		
<ul> <li>Security Service Configuration</li> </ul>		<b>I</b>
	Collect Open	Exit

3) デスクトップ上に""Expert"から 始まる名称のzipファイルが作成され ます。

4)ファイルをメールに添付の上、

#### email\_japan@agilent.com

にお送りください。



