# Expertソフトウェア ver01.02版 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

【ご用意していだたくもの】 ① 未使用のラボチップ 1個 (DNA用 RNA用 Protein用のいずれでも良い)

 ② テストチップセット(G2938-68100もしくは G2938-68300;1セットは装置に付属しておりま す)



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。

nstru	ment	Diagnostics				
ame	: DEL17	00058		Firmware : C.D1.03F		
erial	¢: DEL1	700058		Product ID : G2938B		
Ava	i <del>lable</del> T	ests:				
i	Apply	Name	Description	Status	-	Start
11	×	communication Test	Tests if communication with instrument is funct	🗟 Selected		
1	×	Electronics Test	Tests instrument electronics.	Selected		Select All
1	×	Fan Test	Tests if instrument fan is working.	Selected		
1	×	Ud Sensor Test	Tests if the lid sensors are working.	Selected		Unselect All
1	×	Temperature Test	Checks if the temperature sensors and heater	😇 Selected		Dila.uu
1	×	Stepper Motor Test	Tests if horizontal and vertical motors are wor	🐷 Selected		TABIN
1	×			ected		
1		₩→ 診断県	日はハイオアナライサに、	∧ ortent	-	
les	Proper		ライヤンスの種類に依左			
i.			ノービンハッ川主族に以行			

選択いただく項目;下記の二重線以外の項目



二重線で引いた項目は行わないでください。

(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。



### 各項目における手順=

## **1.Electoronic Test**

#### 2 Fan Test

🔧 Bioanalyze	r hardware test		×
2	Please insert cartridge! Close the lid and press (	)K!	
		ОК	Cancel

上記の画面が現れます。 (1)装置本体の蓋を閉めてください。 (2)画面の"OK"ボタンを押してください



## **3 Lid Sensor Test**

🔧 Bioanalyze	r hardware test		×
2	Please open the Lid!		
		ОК	Cancel

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 Temperature Test**



#### 注意!

デフォルトではCancelボタンが選択されていま す。(3)のステップでは必ず"OK"ボタンを選択 して下さい。キャンセルボタンを選択すると、こ のテスト項目は"Failed"となります.)

### **5 Stepper Motor Test**

この項目の操作は不要です。自動的にソフトウェアが診断を進行します。

## **6 HV Stability Test**

🔧 Bioanalyze	r hardware test	x
2	Please insert Electrode Cartridge and Empty Chip! Close the lid!	_
	OK Cancel	

#### 前項で(1)と(2)を行っている場合、そのまま(3)を行ってください。



(3) 画面の"OK"ボタンを押してください

### 9 Short Circuit Test



(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。前項で用いている ラボチップが古い場合や汚れている場合は、新しいチップに変えてくだ さい。



(2) 装置本体の蓋を閉めてください。

# **11 Optics Test**



(1) Electrode/Diode test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

#### **12 Electrophoresis autofocus test**

🍾 Bioanalyz	er hardware te	st		×
2	Please insert E and enter chip Close the lid!	lectrode Cartridge and values!	Autofocus Test Chip	
		Offset:	18	
		Intensity:	620	
Edit		ОК	Cancel	

- 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



## 13 Laser stability test

🔧 Bioanalyzer	hardware test	×	
2	Please insert Electrode Cartridge and Autofocus Test Chip and enter chip values! Close the lid!		
	Offset: 13		
	Intensity: 620		
Edit	OK Cancel		

000

 $\bigcirc \bigcirc \bigcirc$ 

OOC

OO

**Test Chip** 

Autofocus Offset \*\*

Intensity \*\*\*\*\*

- 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

### 14 Electrode diode test

🔧 Bioanalyze	r hardware test	×
2	Please insert Electrode Cartridge and the Electrode/Diode Test Chip! Close the lid!	_
	OK Cancel	

(1) Electrode/Diode test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

(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 I	-	Stop
mnunication with instrument is funct	🚽 Executed, passed		
ument electronics.	🚽 Executed, passed		Select
trument fah is working.	🚽 Executed, passed		
ild sensors are working.	🚽 Executed, passed		Lineler
he temperature sensors and heater	CExecuting		E-lau
izontal and vertical motors are wor	Execution pending		1939
uctivity of channels (pin to pin).	Execution pending		
coltrast solutions and stability	The Execution panding	-	

- Executing
- Execution pending
- Executed, passed
- Executed, failed

(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、弊社に 下記のファイルのご提供をお送りください。

> 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

送付先; email\_japan@agilent.com

	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
iagnosis	🖻 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
ブジェクトを選択すると、その説明が表 だれます。	Diagnosis_28-07-2004_12-56-09.xdy 🖻	62 KB	XDY ファイル	2004/07/29 17:40
連項目:				
<u>1    + + + + + + + + + + + + + + + + + +</u>				
<u>イネットワーク</u>				
<u>1 コンピュータ</u>				