Agilent 2100 バイオアナライザ







Contents

* Agine facturation	1 City	バイオアナライザ	豆知
102			

- ・サイズの算出方法
- ・濃度の算出方法 DNA
 - RNA Nano/Pico small RNA Protein Protein2
- ・RINの計算フロー

識

З

4 5

6 7 9

10

▶ よく報告されるトラブル・ご質問

- **Expert**編 ・ピークが認識されない 13
- ・間違ったピークがマーカーと認識される 14 15
 - ・RINが計算されない
 - 16 ・rRNAのピークを調整したい
 - ・秒表示からサイズ表示に切り替わらない 18
- 実験操作編

20



便利な機能のご紹介

・スメアassay	
DNA	10
DNAUS	11
· Comparison機能	12

Comparison饿肥





^{豆知識} サイズの算出方法

例) DNA7500 assay

ヾイオアナライザ



- ①ラダは既知のサイズを用いているため
 ソフトウェアにはあらかじめ
 ピークサイズがインプットされています
- ②ラダレーンのピークのMigration timeを もとに、Standard Curveを描きます
- ③それぞれのピークのMigration timeから Standard Curveをもとにサイズを算出 します

Ladder size はAssay Propertyから Standard Curve:はChip Summaryから確認可能





Image: Image: Image: State S





Overlaid Samples

 Markerに入っているUpper Markerは 既知濃度です

②各レーンのUpper Markerのエリアを基準に インジェクション量の補正を行います

③各ピークがもつサイズをもとに、correction factorをかけます correction factor: Agilentがもつ経験的な値。 インジェクション時のサイズ によるバイアスを補正する ためのもの

④サイズに応じてモル濃度計算をおこないます

Peak Conc.= (The Area)* (UpperMarker known conc./UpperMarker Area) * (injectoin bias factor)



Image: Image: Image: Selection of the selection of the



①Ladderに入っているRNAは既知濃度です

Nano: 150ng/ul Pico : 1000pg/ul

②Ladderのエリアを基準に濃度を算出します

Ladder Area : Ladder Conc. = Sample Area : Sample Conc.

Nano RNA Conc. (ng/ul)= (The Area) *(Ladder Area) / (150 * Sample Area)

Pico RNA Conc. (pg/ul)= (The Area) *(Ladder Area)/(1000 * Sample Area)

Page 5



Image: Instance of the second secon

Ladder



①Ladderに入っている各ピークは既知濃度です

②Ladderの各ピークのエリアを基準に サイズごとの補正値を計算します

③各レーンのLower Markerは既知濃度 (500pg/ul)です



④ ②と③で求めたfactorでmiRNA、small RNAと
 想定される領域の濃度を計算します









■ バイオアナライザ 豆知識 濃度の算出方法 - High Sensitivity Protein 250 assay-



Ladder のtime corrected areaと濃度(4,167 pg/ µL*)を基に Sampleのtime corrected areaから算出

*通常プロトコルで調製時のSample Bufferと混合するときのladder濃度 (ladder のもとの濃度= 1 mg/mL)



Image: Instance of the second secon

染色効率はタンパク質の種類により異なります Upper Markerのみでの補正(前頁)の他、標品による濃度補正も可能です

Assay Pro	operties	Chip Summary	Gel	Electroph	erogram	Result Flagging	Log Book
	_		Dal	ta File :	Demo Pr	rotein 80 Series II	.xad
logie	1-3		Loc	tation :	C:¥…t¥	2100 bioanalyzer ^y	∉2100 expert¥dat
schno	4-6	000	Cre	eated :	July 11,	2008 21:45:00	
ent Te	7-9		Мо	dified :	June 03	, 2009 1:46:21	
Agil		# G	Sol	ftware :	Created	by version B.02.0	07.SI437
*	Pro 0n-02	ptein Chip ip-Electrophoresis	Ass	say :	Protein	Analysis 5 - 80 kD	a Diagnostics, v4.
2							

	Sample Name	Sample	Com	Use For Calibration	Conc.[ng/ <i>f</i> Êl]	Statu	
	IgG non-red.				0	~	
2	IgG 1:5 (red.)			V	200	~	
3	IgG 1:20 (red.)				50	~	
4	IgG 1:10 (red.)				100	~	
5	IgG 1:2 (red.)				500	~	
6	IgG (pur, red.)				1000	~	
7	IgG non-red.	_			0	~	
8	Low Range ladder			Calk Y = 1.01	ration Curve *X R*2 = 0.986		
9	Low Range ladder		1000				
10	PBS blank		800-E				
		(E) (E)	800				
	Chip Lot # Reagen	t Kit Lot # 🖁	500-E				
		Relativo	400				
			300-				
			100	< ·			
			100	200 300 400 50 Stary	0 600 700 800 Sard Conc. (ng)(伊)	900	1000

- ①10ウェルのうち、何ウェルかに
 濃度をふった標品、残りにサンプルを
 入れて測定します
- ②データファイルを開き、 Chip Summaryに標品の情報を入力 します
- ③ソフトウェアが標品のレーンの メインピークをもとにStandard Curveをかきます

```
④③をもとに濃度を補正し算出します
```







① エレクトロフェログラムの region分けをおこないます



- Anomaly thresholdによる 異常データの検出をおこないます
- ●フラグ 赤字→Critical
 ⇒RIN値が計算されず
 ●フラグ 黒字→Criticalではない
 ⇒RIN値は計算される

Unexpected baseline signal Unexpected signal in pre-region Unexpected signal in 5S-region Unexpected signal in fast-region Unexpected signal in inter-region Unexpected signal in precursor-region Unexpected signal in post-region Unexpected ribosomal ratio Unexpected sample type Unexpected lower marker











Page 11

-🔅 💮 Agilent Technologies

便利な機能のご紹介 スメアアッセイ – DNA assay以外-

DNA assay以外の場合、スメアアッセイはデフォルト設定ではありません。

(Region Tableタブがありません)







異なるチップからのデータを選択抽出して、ひとつのファイルに統合できます

最大48データまで可能です

便利な機能のご紹介





よく報告されるトラブル・ご質問 (Expert編) ピークが認識されない

設定されている閾値以下のピークは認識されません

(例:シグナルが弱い、ピークがブロード)

認識させるためにSetpointの値を変更する必要があります





Lage Content of the second content of

マーカーと近い泳動度を持ったピークがある場合、

正しくマーカーが認識されない場合があります







Logic Laboration Laboration

プログラム上の想定範囲外のデータの場合、エラーとなりRINが表示されません ⇒p.9「RIN計算のフロー参照」

RINを表示させるため、フラグのたった項目の設定を変更する必要があります







Page 17

- Agilent Technologies

「 よく報告されるトラブル・ご質問 (Expert編) rRNAのピークを調整したい

設定したピークに名前(真核生物の場合、18S/28S)を設定します







Let a c a b b b c matrix of the constraint of the constraint

以下の原因が考えられます

- ・マーカーが認識されていない場合
 →マーカーのピークが低い場合、ピークとして認識されません
 ピークを認識させ、マーカーを設定してください
 (p.13 の「ピークが認識されない」参照)
- ・分析途中でストップボタンを押した場合
 →分析途中のデータがあるとサイズ表示されません (次ページ参照)
- Comparison機能の場合

 →Comparisonモードではサイズ表示ができません
 Dataモードでは同ーチップ上のラダを基にサイズを計算しています
 このため、複数のチップ上のデータを比較するComparisonモードでは
 計算上、サイズ表示が難しく秒表示になります
- ・上記以外 → 弊社サポート窓口へお問い合わせください





しょく報告されるトラブル・ご質問 (Expert編) 砂表示からサイズ表示に切り替わらない

・分析途中でストップボタンを押した場合 →分析途中のデータがあるとサイズ表示されません 最後のデータを除く必要があります





実験操作編 -チェックポイント-

データに問題がある場合は以下の点についてチェックしてください

【試薬】・Expire Dateを過ぎた試薬・チップを使用していませんか?

*チップにも使用期限があります。古いチップを使用すると

泳動パターンが乱れる可能性がございますのでご注意ください

・Gel-dye Mixの調製時に

試薬は常温に戻し、よくMixしてから使用していますか?

試薬の量、遠心速度、時間は正確ですか?

・チップにアプライする時に





- よく報告されるトラブル・ご質問

実験操作編 -チェックポイント-

データに問題がある場合は以下の点についてチェックしてください

【サンプル】

・スペックよりも高濃度ではありませんか?

Kitにより適正濃度が異なります

- ・高分子Genomic DNAが高濃度に入っていませんか? DNase処理、希釈等を行ってください
- ・サンプルの塩濃度は適正ですか?

DNA HS, RNA Pico, Small RNAの場合、ご注意ください





よく報告されるトラブル・ご質問

実験操作編 -チェックポイント-

データに問題がある場合は以下の点についてチェックしてください

【プライミングステーション、ゲル充填時】

・プライミングステーションをメンテナンスしていますか?

ガスケットに汚れ、亀裂はありませんか?

シリンジは定期的に交換していますか?(シリンジはKitについています)

シリンジはスタンドにきちんとセットされていますか?

・充填時のストッパーの位置、充填時間は正しいですか?

・充填後、流路に気泡が入っていませんか?(気泡が入った場合は調製し直してください)





よく報告されるトラブル・ご質問

実験操作編 -チェックポイント-

データに問題がある場合は以下の点についてチェックしてください

【電極】

- ・分析後の電極の洗浄を適切に行っていますか?
 泳動終了後、直ちに洗浄を行ってください
 クリーニングチップでの洗浄は液量 350 ul を守っていますか?
 洗浄液がクリーニングチップ内で偏っていませんか?
- ・数か月に一度、超音波による洗浄を行ってください









Agilent 2100 バイオアナライザ

日本アジレントゲノミクスウェブサイト http://Agilentgenomics.jp

製品に関するお問い合わせ先

Phone: 0120-477-111

- Fax: 0120-565-154
- Mail: email_japan@agilent.com

		75004	こついて「お聞きを生」ご	28 70-AL+AL				
● 長をすービス	+£→ <u>\$</u>		K	ME				
		~	Hong	5 84784 ED X 5 878 8				
ゲノミウス								
			~					
		7		A State of the				
		~						
14:0-IA-2	JAME			A STATE OF A STATE OF A STATE				
e-crice-meter	ヒルデノムの構成によって、科学 実験内に存在する全ての研想の	には新しい時代が別本しました の最佳平均環時間まする技術は	、被補加生体を表現的に明らかに 今日の研究療法においては、会	することが可能になったのです とり飲のものとなっています。				
NUESEE(月)勝定で用 8日初にた肥料後たSナー								
lat 10 にて [アジレント企業論 創がございます] テーマ『マーカ	デジレントは、温祉干売用の分布 解決する和手他いをいたします		法戦争として、世界の主要な群落	活躍の視覚する思想に同想を				
- としての MRILA の可能性」								
CESCTL-C ERR	ナビゲーション かめ	の内容ジョンプするナドドニシュ	177.					
マ: マ:	◆ 然長・レリューション							
	 ● 施文·技術資料・アプリテー5 ● デデート(通知二)(1)/2 							
7レイ スキャナ キャンペーン (2010.10.29主で)		A CONTRACTOR OF THE OWNER						
-Engli	豊島 : アジレント ぎノムク	スートーオルンラューション						
- スモバシーテンシング : Aglast 2100 パイナアナライデカ合発	/(c+7++c+	01147400764	武泉(発現)	試験(デンム)				
30425-90500'940'90 5 485	-	ink-jet technology	A CONTRACTOR	Carlotter State				
•56299-2				and a second				
■ アジビジャーテクノロジー、ヒト デノムの金エクソン構築を与って、		I STATION						
PやするSureSelectターデット SILK 産業与いたを発表(10/15)		Construction of the	the pass	Sec. 2				
750.546 JD-JSR84	14++++++	豊臣平便県アレイ	遺伝干発現用:ラベル化	OCH/CN/R:54LE				
を定義のた業業是代マイクロアレ の数単分のの1	CRASH CCD Not South to Clin	NERILATE-C	Low light Galok Ang /	(発生生亡 /				
TRANK STATES	4-05- 014-016	CHY7L-F	2-54242(18/28)	ハイプロダイゼーション				
きる部分のみを経済可能にする	タンパク戦分野	CMP-m-a-shp7t-4 Op074/95/67t-4	いんでもあったーション (10回応 / 大力量)	(25日応 / 大安量) 法承し(10万元)				
どの産効素化を完成するシステ	神秘分析	2231.76-((AMB))	信奉(いっつす)	ChiP-on-a-chip/ST/L-E				
1011日11日11日	CALCULAR PROVINCE AND		(法半注量体干発病と同に)	きべれた(発生の意)				
 CIN間道:11の疾病の遺活子 約第四の結婚に向けたいTOCOT 	143/1-1/21/017			(活動)HOOH/ON//と間に)				
osiedweitijelle	A479-34-85	于-26时	1994-595	#72.4×79×				
04-09-0		And Address of the Party of the						
- スルモクジーン構築		19- × 10						
アジレントアレイ発展ページ	*			SumSoloot				
・ AAME/ログイン ・ アジレントデノSクスなる構象				Surescielu				
+#-+A-00540	24++	Feature Extraction v10	リアルタイムPCR装置	Stare Select 25				
0.309/12936	オーラン ティンパイティンパ道院長	CED Aglent Generates	DECOUP / NECCOSP) OPCR & CRT-PCRJ2R	Hansan All Econ 4-ct-운 문				
1440 <i>7</i>	1	workbanch		X装色体やいくれたナペオ エンペSCLDフラグメール				
01014740076-45476				¢.				
				へそれ民計会 47-1連続手号 10ストロを使利金付け				
	第二: 臣司司書: 7:157-052章							
	1 : Jacobie (Open Gen	ordex)						
2: 武府県計・7277-1922								
Part of state and an over	3: 237 - 2 2 3 3 5 5 5	and the second sec						



