

<u>Agilent 2200 TapeStation Analysis Table 機能についての紹介</u>

Agilent2200 TapeStation Analysisソフトウェアでは、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。



1 Peak Table機能

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。

<u>開き方</u>

1) Peak Tableを開くために、該当データ を開き、Viewメニューで Electropherogramを選択してください。





<u>設定方法</u>

✓ 各ピークの境目は青い実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、 積分線をドラッグして左右に動かすこ とで変更できます。



✓目的のピークが表示されていない場合、右クリック>Addで積分線を追加できます。



2 Region Table機能

Home

Electropherogram

View

Sample Table

N

To [bp]

🗐 🗸 File

222

===

Gel

Region Table

From [bp]

ブロード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。 任意の領域 (Region)を指定して解析します。

<u>開き方</u>

1) 該当データを開き、Viewメニューで Electropherogramを選択してください。

2) 画面下のタブでRegion Tableを選択し てください。

<u>見方</u>

設定した領域の

- ・From(任意のスタート位置)
- ・To(任意のエンド位置)
- Average Size (平均サイズ)
- ・Conc (濃度)
- ・Region Molarity (モル濃度)
- ・% of Total (領域の全体に対する割合)
- Region Comment

が表示されます。

- つ割合) Region Table Sample Table Tom [tabl Sample Table] Tom [tabl Sample Table]

任意の領域の設定方法

(1) Region Settingsを選択します。

	🔤) 👝 🚍 🖘)				0 TapeStation Analysis S	oftware A.01.04 - 2014-	08-04-02.D1000	
	II - File Home							
	Gel Electropherogram Region	Aligned Scale to Co Sample Sna	Gel Contrast Show All Lanes	EPG Reset Snapshot Zoom	Assign Insert Remove Settings	Restore Default Settings Region	Region ns Settings	
	View	Display	Gel	EPG		Analysis		
							/	
(2) 表示されたボ FromとToの欄にご カします。	ックスで任意 スタートとス	の領域を設 トップのサ	設定します。 ⁺イズを入へ	Assay Options - D1000 Regions	This page allows you to set applied to newly analyzed fi If you want to apply these r Va From [bp]	up default regions for files of tr les, or files which have been 'R regions to the current file, use t To [bp]	his protocol. These regions evert[ed] to Originai'. he 'Apply to File' button. Name	x will only be
(3)全レーンに同 to Fileボタンを押	同じ領域を設定 します。	ミする場合	、Apply		 ✓ 200 ✓ 100 	500 1000		
(4) Applyボタンな す。	をクリックして	てボックス	を閉じま				A	pply to File
			-				OK Cancel	Apply
								~

3 Sample Table 機能

totalRNA サンプルのRINeの値やrRNA比を見る場合に使います。

<u>開き方</u>

1) 該当データを開いた後、Viewメニューの設定はGel,Electropherogram、Region いずれでもSample Tableは表示されます。

2) 画面下のタブでSample Tableを選択 してください。



A0

<u>見方</u>

□ RNA用のアッセイでは下記が表示されます。

- ・ウェル情報
- ・RINeの値;分解の度合いを1-10の数字で表示
- 28S/18S (Height)
- 28S/18S (Area)
- ・濃度
- Sample Description
- Alert
- Observations
- Total RNA Area
- rRNA Area

Well	RINe	28S/18S (Height)	28S/18S (Area)	Conc. [ng/µl]	Sample Description	Alert	Observations	Total RNA Area	rRNA Area
C2	10.0	1.9	3.0	139				5.93	2.30
D2	9.9	2.3	3.3	400				17.07	6.31

- □ DNA, Proteinのアッセイでは下記が表示されます。
 - ・ウェル情報
 - ・濃度 (Peak Tableの総濃度)
 - Sample Description
 - Alert
 - Observations

電話 🔼 0120-477-111

Well	Conc. [ng/µl]	Sample Description	Alert	Observations
A0	18.0	D1000 Ladder		Ladder
A1	7.11	1		
B1	6.90	2		

