





SNaPshot[®] Kit SNP 解析 Getting Started Guide

ご使用になる前に

1

2

3

リファレンス データを使用した Panel および Bin Set の作成

Primer Focus[®] Kit データを 使用した Panel および Bin Set の作成

SNaPshot[®] Kit 解析の設定

5

4

結果の解析と検証

6

結果の印刷とエクスポート

© Copyright 2005, Applied Biosystems.All rights reserved.

研究用にのみ使用できます。診断目的およびその手続き上での使用はできません。

本マニュアルに記載されている情報は、予告なく変更されることがあります。Applied Biosystems は、本マニュアルのあらゆる誤謬に対して、一切の責任を負わ ないものとします。本マニュアルの情報は、発行の時点においては、完全かつ正確であるものとみなします。Applied Biosystems は、本マニュアルと関連、もし くは本マニュアルから生じる、偶発的、特別、複合的、派生的損害に対して、いかなる場合も責任を負わないものとします。

ご購入者への告知:ライセンス拒否

本ソフトウェアの製品のみの購入では、明示されているか否かに関わらず、または禁反言によるか否かに関わらず、Applera Corporation が所有または管理する 特許権利に基づくあらゆるプロセス、機器またはその他の装置、システム、混合物、試薬、キットの権利のいかなる権利(ライセンス)の所有も意味しません。 GeneMapper Software has not undergone specific developmental validation for human identification applications. Human identification laboratories analyzing single-source or parentage samples which choose to use GeneMapper Software for data analysis should perform their own developmental validation studies.

商標:

ABI PRISM, Applied Biosystems, GeneMapper, LIZ, Primer Focus, and SNaPshot are registered trademarks, and the AB Design, Applera, GeneMapper and SNPlex are trademarks of Applera Corporation or its subsidiaries in the U.S. and/or certain other countries.

This product includes software developed by the Apache Software Foundation

(http://www.apache.org/). Copyright © 1999-2000 The Apache Software Foundation. All rights reserved.

This product includes software developed by the ExoLab Project (http://www.exolab.org/). Copyright 2000 © Intalio Inc. Allrights reserved.

JNIRegistry is Copyright © 1997 Timothy Gerard Endres, ICE Engineering, Inc., http://www.trustice.com.

AFLP is a registered trademark of Keygene N.V.

Microsoft is registered trademark of Microsoft Corporation.

Oracle is a registered trademark of Oracle Corporation.

All other trademarks are the sole property of their respective owners.

Applera Corporation is committed to providing the world's leading technology and information for life scientists. Applera Corporation consists of the Applied Biosystems and Celera Genomics businesses.

Part Number 4363078 Rev. B 06/2005

目 次

	まえがき v
	このガイドの使い方
	詳細情報の入手方法
	サポートの入手方法 vii
第1章	ご使用になる前に 1
	SNaPshot [®] Kit 解析について
	データ例について
	SNaPshot [®] Kit 解析のワークフロー6
	GeneMapper [®] ソフトウェア の用語
	ソフトウェアの起動とログイン
	ユーザ独自のサンプル ファイルを使う場合のこのガイドの使用
	GeneMapper ソフトウェア v3.7 を使う場合のこのガイドの使用
	このガイドにおける手順のその他の方法8
第2章	リファレンス データを使用した Panel および Bin Set の作成 9
	新規プロジェクトの作成とリファレンス サンプル ファイルの追加
	アナリシス パラメータと Table Settings の設定12
	プロジェクトでのサイジングのみの実行14
	Kit、Panel、Marker、Bin Set、Bin の作成19
第3章	Primer Focus [®] Kit データを使用した Panel および Bin Set の作成 31
	新規プロジェクトの作成と Primer Focus [®] Kit サンプル ファイルの追加
	アナリシス パラメータと Table Settings の設定34
	プロジェクトでのサイジングのみの実行
	Kit、Panel、Marker、Bin Set、Bin の作成41
第4章	SNaPshot [®] Kit 解析の設定 49
	新規プロジェクトの作成とサンプル ファイルの追加
	プロジェクト用のアナリシス パラメータと Table Settings の設定
第5章	結果の解析と検証 57
	プロジェクトの解析
	結果の検証

第6章	結果の印刷とエクスポート	63
	結果の印刷	64
	結果のエクスポート	65

索引

まえがき

このガイドの使い方

このマニュアルの GeneMapper[®] Software Version 4.0 SNaPshot Kit SNP 解析 Getting Started Guide では、 互換性のある任意の Applied Biosystems 社製シーケンサ、および Data Collection ソフト ウェアを使用して SNaPshot Kit データをサイジングおよびジェノタイピングする手順を、段 階を追って簡潔に説明しています。また、トラブルシューティング、データの印刷とエクス ポート、レポート作成の方法についても記載しています。このマニュアルは、GeneMapper ソフトウェアの基本機能が素早く理解できるような構成になっています。

- 対象読者 このマニュアルは、GeneMapper ソフトウェアの初級ユーザを対象としています。
- 前提条件 このマニュアルは、対象読者が次の条件を満たしていることを前提としています。
 - 『GeneMapper[®] Software v4.0 Quick Installation and Administration Guide』 (PN 4363080)の説明に従って、GeneMapper ソフトウェア v4.0 がインストールされて いること。
 - Microsoft[®] Winows[®] XP オペレーティング システムに関する実践的な知識を備えている こと。
 - 表記法 このマニュアルでは、次の表記法を使用しています。
 - 太字は、ユーザの行う操作を示します。たとえば、次のとおりです。
 残りの各フィールドについて、0を入力し、[Enter] キーを押します。
 - ゴシック体で表記されている言葉は、初出または重要な語を示し、強調しています。たとえば、次のとおりです。
 解析の前に、必ず新しいマトリックスを調製してください。
 - ▶は、ドロップダウンメニューまたはショートカットメニューで連続して選択するコマンドを区切って示しています。たとえば、次のとおりです。
 - 「File」 ▶ 「Open」 ▶ 「Spot Set」を選択します。 サンプル列を右クリックし、「View Filter」 ▶ 「View All Runs」を選択します。
- **ユーザへの注意事項** Applied Biosystems のユーザマニュアルには、2 種類の注意事項が記載されています。各注 意事項は、次のような特定のレベルの注意と対応が必要なことを示しています。

注:製品を使用する上で役に立ちますが、必須ではない情報を記述しています。

重要! 装置の適切な操作、ケミストリキットの正しい使用法、化学物質の安全な使用法に関する必要な情報を記述しています。

ユーザへの注意事項の例を次に示します。

注: カラムのサイズは、ランタイムに影響します。

注:キャリブレーション機能は、Control Console でも使用できます。

重要! クライアントがデータベースに接続していることを確認するには、有効な Oracle ユー ザ ID とパスワードが必要です。

重要! 1 枚の **96** ウェル プレートに対して、それぞれ **1** つの **Sample Entry** スプレッドシートを作成する必要があります。

安全に関する
 このユーザマニュアルには、安全に関する注意事項も記載されています。詳細については、
 注意事項
 『GeneMapper[®] Software v4.0 Installation and Administration Guide』(PN 4363080)を参照してください。

詳細情報の入手方法

- 安全に関する情報 安全に関する情報の詳細については、『GeneMapper[®] Software v4.0 Installation and Administration Guide』(PN 4363080)を参照してください。
- ソフトウェアの保証
 すべての保証とライセンスに関する情報の詳細については、『GeneMapper[®] Software v 4.0
 とライセンス
 Installation and Administration Guide』(PN 4363080) を参照してください。
 - 関連資料 このソフトウェアには、次の関連資料が同梱されています。
 - GeneMapper[®] Software v4.0 Installation and Administration Guide GeneMapper ソフトウェア v4.0 のインストール、セキュリティ、メンテナンスの手順について記載しています。
 - GeneMapper[®] Software v4.0 Getting Started Guide GeneMapper ソフトウェアで 提供されるアプリケーションに特化したデータ例を解析する方法を記載したガイド 5 冊。このガイドでは、互換性のある Applied Biosystems 社製のシーケンサおよび Data Collection ソフトウェアで作成されたマイクロサテライト、LOH、AFLP[®] システム、 SNaPshot[®] キット、および SNPlex システム データを解析する手順について、段階を 追って簡潔に説明しています。このガイドは、GeneMapper ソフトウェアの基本機能が 素早く理解できるような構成になっています。
 - GeneMapper[®] Software v4.0 Online Help GeneMapper ソフトウェアに関する説明 と、一般的なタスクの手順を提供します。オンライン ヘルプにアクセスするには、[F1] キーを押す、「Help」 → 「Contents and Index」の順に選択する、「GeneMapper」ウィ ンドウのツールバーで ? たクリックする、という 3 つの方法があります。
 - GeneMapper[®] Software v4.0 Quick Reference Guide 解析のワークフローをタイプ別に記載すると共に、GeneMapper ソフトウェアと互換性のある装置、ソフトウェア、解析アプリケーションのリストを提供します。
 - GeneMapper[®] Software v4.0 Reference and Troubleshooting Guide 動作理論な どの参照情報とトラブルシューティング情報を提供します。

このマニュアルおよび前述の資料の PDF (ポータブル ドキュメント フォーマット) 版は、 GeneMapper ソフトウェア v4.0 Documentation CD に収録されています。

注:詳細については、vii ページの「サポートの入手方法」を参照してください。

オンライン ヘルプ GeneMapper ソフトウェアは、ユーザ インタフェースの各機能を使用する方法について説明 したオンライン ヘルプ システムを備えています。オンライン ヘルプにアクセスするには、 [F1] キーを押す、「Help」→「Contents and Index」の順に選択する、「GeneMapper」ウィ ンドウのツールバーで 2 をクリックする、という 3 つの方法があります。

連絡先 Applied Biosystems では、弊社のユーザマニュアルをより使いやすいものにするため、お客様からのご意見、ご要望をお待ちしております。次の電子メール アドレス宛にお送りください。

jptechsupport@appliedbiosystems.com

サポートの入手方法

すべての地域における最新サービスおよびサポート情報を入手するには、 http://www.appliedbiosystems.co.jp にアクセスし、「Customer Support」のリンクをク リックしてください。

「カスタマーサポート」ページでは、次のことが可能です。

- よくある質問 (FAQ) を検索する
- テクニカル サポートに質問を直接送信する
- Applied Biosystems ユーザマニュアル、MSDS、分析証明書(Certification of Analysis)、およびその他の関連資料を注文する
- PDF 文書をダウンロードする
- カスタマートレーニングに関する情報を入手する
- ソフトウェア アップデートおよびパッチをダウンロードする

さらに、「テクニカルサポート」ページには、世界各地の Applied Biosystems テクニカル サービスおよびサポート機関の連絡先の電話番号やファックス番号も掲載されています。

まえがき サポートの入手方法

ご使用になる前に





SNaPshot[®] Kit 解析について

SNP マーカ

SNP (Single Nucleotide Polymorphism) マーカは、既知の DNA シーケンス内で相違する一 塩基から成り、最大4種類のマーカアレルすなわち変異を生み出します。



G-C および A-T 塩基対が、この SNP サイトで 存在する可能性のある 2 つのアレル

SNaPshot[®] Kit 解析

SNaPshot[®] Muliplex Kit

SNaPshot Multiplex Kit では、SNP を含む領域を PCR 増幅後、蛍光標識されていないプラ イマーを一塩基伸長することにより、最大 10 ローカスまでの SNP マーカを同時に調査する ことができます。このプライマーは、SNP サイトの直前にアニーリングするよう設計されて います。いったんプライマーがアニーリングすると、蛍光標識された相補 ddNTP (ダイ ター ミネータ)をアニーリングしたプライマーに追加することで、一塩基伸長反応が起こります。 4 つの ddNTP は、それぞれ異なる色の蛍光色素によって蛍光標識されています。この結果、 異なる SNP アレルに対するマーカフラグメントは、塩基数は同じで、色が異なります。電気 泳動と蛍光検出の後では、ひとつのマーカのアレルは、ほぼ同じサイズの異なる色のピークと してエレクトロフェログラム プロットに表示されます。異なるアレル ピークのサイズは、蛍 光色素の分子量によってわずかに異なります。これで GeneMapper ソフトウェアを使用して、 データのサイジングおよびジェノタイピングを行うことができます。(図 1-1 を参照)

SNaPshot Multiplex Kit では、異なる長さのプライマーを使用することにより、最大 10 ローカスまでの SNP マーカを同時に調査することができます。場合によっては、アニーリングしないテール配列をプライマーに追加して、プライマー長を他のプライマーと十分区別できるようにする必要があります。この手順により、SNP マーカのオーバーラップを回避できます(図 1-2 を参照)。



図 1-1 SNaPshot[®] Muliplex Kit ワークフローの概要



図 1-2 テンプレートへアニーリングし伸長されたテール付きプライマー

Primer Focus® Kit

オプションで Primer Focus Kit を使用すると、目的の SNP マーカで可能な 4 つのアレルを すべて作成できます。このキットに含まれている試薬により、4 つすべての潜在的 ddNTP を、 蛍光標識されていない SNaPshot プライマーの 3' 末端に、テンプレートを使用することなく 追加できます。 これで GeneMapper ソフトウェアは Primer Focus Kit サンプルのサンプル ファイルを使用 して、次を実行できます。

- 任意のテール付き伸長産物の移動度を評価し、SNP マーカがオーバーラップしていない ことを検証
- Auto Panel 機能を使用した、各 SNP マーカアレルに対する Bin の自動作成



SNaPshot[®] Kit Applied Biosystems では、次の Kit を提供しています。

- SNaPshot Multiplex Kit (PN: 4323159, 4323161, 4323163)
- Primer Focus Kit (PN: 4329538)

互換性のある装置 SNaPshot 解析と互換性のある Applied Biosystems 社製シーケンサの詳細については、 『GeneMapper[®] Software v4.0 Quick Reference Guide』(PN 4362816) を参照してください。

データ例について

サンプル ファイル このガイドの「ご使用になる前に」に記載されている演習を実行するには、コンピュータ ハー の場所 ドドライブの次の場所に配置されている 3 つのサンプル ファイル (.fsa)、およびオプション で 6 つの Primer Focus Kit サンプル ファイルを使用します。

<drive>:\AppliedBiosystems\GeneMapper\Example Data\SNaPshot

注:上記の場所は、GeneMapper[®] ソフトウェアがインストールされているドライブによっ て異なります。デフォルトでは、D ドライブにインストールされます。

装置とサイズ サンプル ファイルは、SNaPshot Multiplex Kit のサンプルを GeneScan[™] 120 LIZ[®] サイズ スタンダード スタンダードとともに ABI PRISM[®] 3100 ジェネティック アナライザで泳動したものです。また、各マーカのすべての潜在的アレルを含むサンプル ファイルは、Primer Focus Kit のサン プルを GeneScan[™] 120 LIZ[®] サイズ スタンダードとともに ABI PRISM[®] 3100 ジェネティック アナライザ上で泳動したものです。

マーカ情報 SNaPshot Kit のデータ例には、6つのマーカが含まれています。

第2章では、リファレンス データを使用して、独自のプロジェクト用の Marker と Bin (アレル定義)を手動で作成する方法について説明します。また、この章では、サンプル データのマーカ情報およびアレル情報についても説明します。

第3章では、GeneMapper ソフトウェアの Primer Focus Kit データと Auto Panel 機能を使 用して、独自のプロジェクト用の Marker と Bin (アレル定義)を自動的に作成する方法に ついて説明します。



SNaPshot[®] Kit 解析のワークフロー

次のフローチャートは、GeneMapper[®] ソフトウェアを使用して SNaPshot 解析を実行する 手順を要約したものです。

リファレンス データを使用した Panel および Bin Set の作成 (第2章)

- 1. 新規プロジェクトを作成し、リファレンス サンプル ファイルを追加します。
- 2. プロジェクト用のアナリシス パラメータと Table Settings を設定します。
- 3. プロジェクトのサイジングのみを実行します。
- 4. プロジェクト用の Kit、Panel、Marker、Bin Set、 Bin を作成します。

Primer Focus[®] Kit データを使用した Panel および Bin Set の作成 (第3章)

- 1. 新規プロジェクトを作成し、Primer Focus Kit のサ ンプル ファイルを追加します。
- プロジェクト用のアナリシス パラメータと Table Settings を設定します。
- 3. プロジェクトのサイジングのみを実行します。
- 4. プロジェクト用の Kit、Panel、Marker、Bin Set、 Bin を作成します。



GeneMapper[®] ソフトウェア の用語

用語	定義
アナリシス パラメータ	ユーザが定義する設定の集合(Analysis Method、Size Standard、Panel を含む)。プロジェクトに使用するすべてのサンプルの解析用に GeneMapper ソフトウェアで使用されるサイジング アルゴリズムとジェノタイピング アルゴリズムを決定します。
Bin/ビン	Marker 内のアレルを定義するフラグメント サイズ(bp)と蛍光色素。Bin は、 Marker と関連付けられる可能性のある各アレルに対して作成します。
Bin Set/ ビン セット	Bin の集合(アレルの定義)。通常は、一連の実験条件に対して固有の Bin Set 作成します。
Marker/ マーカ	SNP Marker は、名前とフラグメント サイズ レンジ(bp)によって定義されます。
Panel/パネル	Marker のグループ。GeneMapper ソフトウェアでは、Panel を Bin Set と関 連付けることにより、Marker の Bin 定義を設定します。
Kit/ キット	パネルのグループ。

ソフトウェアの起動とログイン

GeneMapper ソフトウェアを起動してログインするには、次の手順を実行します。

- 1. 「Start」→「All Program」→「Applied Biosystems」→「GeneMapper」→「GeneMapper 4.0」の順に選択します。
- 2. 「Login to GeneMapper」ダイアログボックスで、次の手順を実行します。
 - a. システム管理者によって割り当てられたユーザ名とパスワードを入力します。
 - **b**. 「**OK**」をクリックします。

ユーザ独自のサンプル ファイルを使う場合のこのガイドの使用

このガイドを使用して、ソフトウェアで供給されるデータ例を解析するだけでなく、ユーザ独 自のサンプルファイルを解析する際の一般的な SNaPshot Kit 解析のワークフローを段階的 に実行できます。ソフトウェアの高度な機能に関する詳細については、GeneMapper[®] Software Online Help を参照してください。

GeneMapper ソフトウェア v3.7 を使う場合のこのガイドの使用

このガイドに記載されているワークフローおよび手順は、GeneMapper ソフトウェア v3.7 で も有効です。

重要! GeneMapper ソフトウェア v3.7 には、このガイドの演習で使用する正確なデータ例 が含まれていません。このため、GeneMapper ソフトウェア v3.7 を使用する場合、参照され るデータ例の解析に、このガイドは使用しないでください。ただし、ユーザ独自のサンプル ファイルを解析する際の SNaPshot Kit 解析ワークフローにはこのガイドを使用できます。



このガイドにおける手順のその他の方法

概要

このガイドでは、GeneMapper[®] ソフトウェアを使用して SNaPshot Kit データを解析できる いくつかの解決策を記載しています。このガイドの演習を完了後、ご自身の研究室の要件に特 化してプロセスをカスタマイズする場合に備え、この節では、いくつかの代替方法の要約と詳 細情報の参照先を提供します。

 自動解析を使用した プロジェクトの設定
 GeneMapper ソフトウェアには自動解析機能が備わっています。これを利用すれば SNaPshot Kit 解析プロジェクトに関連するタスクの大部分を排除できます。第2章、第3章、 第4章のほとんどが、SNaPshot Kit プロジェクトで使用するプロジェクトを手動で作成し、 サンプルを追加し、解析する方法の説明です。自動解析を設定すると、GeneMapper ソフト ウェアは、Data Collection ソフトウェアと連携して、自動的に該当のタスクを実行します。 自動解析機能を使用して SNaPshot Kit プロジェクトを設定する方法の詳細については、 『GeneMapper[®] Software v4.0 Quick Installation and Administration Guide』(PN 4363080) を参照してください。

コマンド ライン イGeneMapper ソフトウェアでは、コマンド ライン インタフェースを介して、ソフトウェアの
主要機能を大部分実行することができます。コマンド ライン インタフェースは SNaPshot
したプロジェクトの
設定設定Kit プロジェクトを解析する際の便利なツールとして使用でき、第2章、第3章、第4章で説
明されているタスクの多くを自動化します。コマンド ライン インタフェース、およびこれを
使用して GeneMapper ツフトウェアの機能を自動化する方法の詳細については、
『GeneMapper® Software v4.0 Quick Installation and Administration Guide』(PN 4363080)
を参照してください。





新規プロジェクトの作成とリファレンス サンプル ファイルの追加

概要 「GeneMapper」ウィンドウ内でプロジェクトを作成し、プロジェクトにサンプルを追加します。

プロジェクトの 新規作成とサンプル ファイルの追加 プロジェクトを新規作成してサンプル ファイルを追加するには、次の手順を実行します。 1. ご をクリックします (「File」 ▶ 「New Project」)。

Project Type	
🔘 Generic	
🔘 Microsatellite	
💿 SNaPshot®	
🔘 OLA Analysis	
O SNPlex™	
O AFLP	

- 2. 「New Project」ダイアログ ボックスで、SNaPshot を選択し、「OK」をクリックします。 GeneMapper v3.7 ではこのダイアログボックスは表示されません。そのままステップ 3 へ進みます。
- 3. 「GeneMapper」ウィンドウで、 🔂 をクリックします(「File」 ▶ 「Add Samples to Project」)。
- 4. 「Add Samples to Project」ダイアログボックスの「Files」タブで、次のパスを確認します。 <drive>:\AppliedBiosystems\GeneMapper\Example Data\SNaPshot

注:上記のパスは、GeneMapper[®] ソフトウェアがインストールされているドライブに よって異なります。デフォルトでは、D ドライブにインストールされます。



5. SNaPshot フォルダを選択し、「Add to List」、「Add」の順にクリックします。

注: このガイドでは、SNaPshot フォルダ内の3つのサンプルファイルをすべて追加しましたが、フォルダ内のファイルの一部(1つまたは2つ)を追加することも可能です。これを行うには、左側のペインでフォルダを展開し、[Ctrl] キーを押したまま、ファイルを個々に選択して、「Add To List」をクリックします。

📾 Add Samples to Project	X
Edit View	
Files GM Database	Samples To Add:
Wr Computer Wr Computer	 ■ SNaPshot ◆ 2_B03_03.fsa ◆ 5_E03_09.fsa ◆ 6_F03_11.fsa
Add To List >>	Clear Add Add & Analyze Eance

SNaPshot フォルダ内の 3 つのサンプル ファイルが、互換性のある Applied Biosystems 社製シーケンサの Data Collection ソフトウェアに入力した情報と共に、「Samples」タ ブ上のテーブルに表示されます。

🗇 GeneMapper - *	Untitle	d - gm l	ls Logged In Da	tabase FRMINF	ODEVD01					
File Edit Analysis View Tools Help										
😂 😂 🔒 🛛 😼	2	III 🖸	I 💷 📖 🕴 🖬		🕨 💣 🕴 Tal	ole Setting: S	NaPshot Defa	ault 🔽 🛙	🗏 🔎 🦫	, 🛛 🕐
😑 🚠 Project	Sample	es Gen	otypes							
🗄 🧰 SNaPshot		Status	Sample File	Sample Name	Comments	Sample Type	Analysis Me	ethod Panel	Size S	Standard
	1	\$	2_803_03.fsa		None	Sample	None	None	None	
	2	N	5_E03_09.fsa	5	None	Sample	None	None	None	
	3	N	6_F03_11.fsa	6	None	Sample	None	None	None	
		<)				>
Progress Status								0	%	Stop



12ページの説明に従って、プロジェクト用のアナリシスパラメータと表示設定を作成します。



概要

アナリシス パラメータと Table Settings の設定

「GeneMapper」ウィンドウ内で、プロジェクト用のアナリシス パラメータと表示設定を確認 します。

アナリシス パラメータには、次の項目が含まれます。

- Analysis Method (Bin Set を含む)
- Size Standard
- Panel (Marker の集合)

プロジェクトにおけるすべてのサンプル ファイルを解析するために GeneMapper[®] ソフト ウェアが使用する、ピーク検出アルゴリズム、サイジング アルゴリズム、ジェノタイピング アルゴリズムを決定するアナリシス パラメータを設定します。

表示設定には、Table Settings と Plot Settings が含まれます。

アナリシス パラ メータの設定 プロジェクト用のアナリシス パラメータを設定するには、次の手順を実行します。

- 1. 「GeneMapper」ウィンドウで「Samples」タブを選択します。
 - 2. 「Analysis Method」カラムで最初の列をクリックし、ドロップダウン リストから SNaPshot Default を選択します。



3. 「Panel」カラムの設定を None のままにします。

注: パネルの選択は、オプションです。Panel を選択しない場合、GeneMapper ソフト ウェアでは、データをサイジングしますが、ジェノタイピングは行いません。このプロ ジェクトを作成している目的は、リファレンス データのサイジング専用解析を実行して、 Panel と Bin Set の作成に使用するためです。従って、アナリシス パラメータの一部と して Panel を選択することはありません。

- 4. 「Size Standard」カラムで最初の列を選択し、ドロップダウン リストから GS120LIZ を選択します(これは、サンプルでランされたサイズ スタンダードです)。
- 5. 「Samples」タブのすべてのサンプル列に、次のとおり選択をフィル ダウンします。
 - a. 「Analysis Method」、「Panel」、「Size Standard」の各カラム ヘッダにまたがって クリッしてクドラッグし、3 つのカラムのすべての列をハイライト表示します。

Analysis Method	Panel	Size Standard
SNaPshot Default	None	GS120LIZ
None	None	None
None	None	None

b. 「Edit」 → 「Fill Down」の順に選択します(または、[Ctrl] キーと [D] キーを同時 に押します)。

Table Settings の
選択「GeneMapper」ウィンドウの最上部で、「Table Settings」ドロップダウン リストから
SNaPshot Default を選択します。

Microsatellite Default	~
New	
AFLP Default	
Microsatellite Default	
SNPlex_v3	
SNaPshot Default	

Table Settings は、解析後に「**Samples**」タブおよび「**Genotypes**」タブで表示される情報を 制御します。**SNaPshot Default** は、**GeneMapper** ソフトウェアで提供されるデフォルトの **Table Settings** の **1** つです。

GeneMapper Manager でカスタマイズした Table Settings を編集および作成することもで きます。詳細については、GeneMapper[®] Software Online Help を参照してください。

プロジェクトの保存 プロジェクトを保存するには、次の手順を実行します。

- 1. $\boxed{1}$ $\boxed{1}$
- **2.** 「Save Project」ダイアログ ボックスで、「SNaPshot Ref Data Project」と入力し、「OK」 をクリックします。

Save Pr		
٩	Project name: SNaPshot Ref Data Project OK Cancel	

「GeneMapper」ウィンドウのタイトル バーに、SNaPshot Ref Data Project が表示され ます。

次の手順

14 ページ の説明に従って、プロジェクトのサイジングのみを実行します。



プロジェクトでのサイジングのみの実行

概要

プロジェクト用のサンプル ファイルの追加と、アナリシス パラメータの設定が完了したら、 次に、サイジングのみを実行します。これにより、サンプル ファイルをリファレンス データ として利用し、**Bin**(アレル定義)を作成できるようになります。

サイジングのみを実行するには、次の手順を行います。

- プロジェクトの解析
- SQ および関連する PQV の閲覧
- サイズスタンダードの検証
- サンプル情報の確認(Raw Data を含む)
- サンプル プロットの確認

プロジェクトの解析 > をクリックします (「Analysis」 **>** 「Analyze」)。

GeneMapper[®] ソフトウェアは、プロジェクト内の各サンプルを解析し、「GeneMapper」ウィンドウの左下にあるステータス バーに進捗状況を表示します。

SQ と PQV の閲覧 Size Quality (SQ) および関連する Process Quality Value (PQV) を閲覧するには、次の手順を実行します。

- 1. 「GeneMapper」ウィンドウの左下にあるステータス バーに、「Analysis Completed」と 表示されていることを確認します。
- 2. 「Samples」タブの右側をスクロールして、SQを閲覧します。

😁 GeneMapper - St	NaPsho	t Ref Data Proj	ject - gm Is I	Logged In Datab	ase FRMINFO	DEVD01								
File Edit Analysis View Tools Help														
😂 🖨 📓 🍒	2	M 📰 🖻 🏨	🖾 🏛	🛅 🕨 💣	Table Setting	g: SNa	Pshot (Default		~ [P	b	🔤 🕐
😑 🚠 Project	Sample	s Genotypes												
🕀 🧰 SNaPshot		Matrix	Run Name	Instrument Type	Run Date & Tim	ne	REF	SQI	SFNF	MNF	SNF	OS	SQ	UD1
	1		SNaPshot	ABI3100	2001-07-31 16	:50:38.0								
	2		SNaPshot	ABI3100	2001-07-31 16	:50:38.0								
	3		SNaPshot	ABI3100	2001-07-31 16	:50:38.0								i
		<							11					>
									_					
Analysis Completed.														Stop

スクロール バーを右方向へドラッグして「SQ」カラムを表示

手順に従ってこのガイドに示されるデータ例を使用した場合、各サンプルに対する **SQ** は **■** (Pass) になります。一方、**SQ** (SFNF、MNF、SNF、OS) の生成に関連する **PQV** も **■** になります。



黄色の 🝐 と赤い 🔵 SQ の原因調査 **重要!** ユーザ独自のデータを解析する際、SQ が ▲ (Check) または ● (Low Quality) とな り、関連付けられた PQV (SFNF、MNF、SNF、OS) が ▲ になることがあります。これ は、サイズ スタンダード、データ、またはアナリシス パラメータに問題があることを示して います。これらの問題の原因を調査して修正するには、15 ページの「サイズ スタンダードの 検証」を参照してください。

注: 💣 をクリックすると、SQ スコアに基づいて、サンプルをソートすることができます。 「Samples」タブの一番上に、🔵 SQ の付いたサンプルがリスト表示されます。

サイズ スタンダード の検証

サイズ スタンダードを検証するには、次の手順を実行します。

- 1. 「Edit」 ▶ 「Select All」の順に選択して、「Samples」タブ内のすべてのサンプルを選択 します。
- 2. <u>Ш</u>をクリックして、Size Match Editor を開きます(「Analysis」 → 「Size Match Editor」)。



Size Quality (SQ) スコア

- 図 2-1 Size Match Editor 「Size Matches」タブ
 - 3. 「Size Matches」タブをクリックして、選択されたサンプルに対する次の項目を表示し ます。
 - Size Quality (SQ) スコア
 - サイズ スタンダード ピーク
 - サイズ スタンダード ピーク ラベル
 - 4. サンプルの Sizing Quality スコア(図 2-1)を確認します。このスコアは、サイズ スタ ンダードからのデータが、ソフトウェアで選択されたサイズ スタンダードにどの程度適 合しているかを示します。このスコアによって、SQ が ■ (Pass)、 (Check)、 (Low Quality) のいずれを表示するかが決まります。

このガイドの指示に従った場合、Sizing Quality は > 0.75 となり、SQ は 📄 (Pass) を 表示します。

一方、独自のデータを解析すると、Sizing Quality は低下し、SQ は ▲ (Check)、また
 は ● (Low Quality) を表示します。トラブルシューティングについては、16 ページの
 表 2-1 を参照してください。

5. サイズスタンダード内のすべてのピークが存在し、正確に標識されているかを判定します。 このガイドの指示に従った場合、図 2-1 に示すとおり、すべてのピークが存在し、正確 に標識されます。

一方、独自のデータを解析すると、一部のサイズスタンダード ピークは正確に標識されないか、失われてしまいます。トラブルシューティングについては、16ページの表 2-1 を参照してください。



6. 「Size Calling Curve」タブをクリックして、選択したサンプルのサイズ スタンダード曲 線を表示します。サイズ スタンダードのフラグメントを表す赤いデータ ポイントと、黒 いベストフィット曲線が表示されます。



- 7. Size Match Editor の左側ペインで別のサンプルを選択し、手順 3~6を繰り返します。
- 8.「OK」をクリックして、Size Match Editor を閉じます。
- 表 2-1 サイズ スタンダードに関するトラブルシューティング

問題	操作
Sizing Quality スコアが低下し、 SQ は 📥 (Check) または 🌍 (Low Quality)を表示するが、すべ てのサイズスタンダード ピークが 存在し、正確に標識されている。	「Size Matches」タブの最上部で「Override SQ」をクリック して、Sizing Quality を上書きします (図 2-1)。 Sizing Quality スコアの変更を 1.0 に上書きします。これは、ユーザがサイズ スタンダードを確認したことを示します。
いくつかのサイズ スタンダード ピークが正確に標識されない。	「Size Matches」タブで、サイズ ラベルを編集、削除、追加します。次に「Apply」をクリックし、更新されたサイジング情報でデータを再解析します。詳細については、GeneMapper [®] Software Online Help を参照してください。
いくつかのサイズ スタンダード ピークが存在しない。	ソフトウェアで、カスタム サイズ スタンダードを作成します。 詳細については、GeneMapper [®] Software Online Help を参 照してください。

サイジングに関するトラブルシューティングやその他の情報は、『GeneMapper[®] Software Reference and Troubleshooting Guide』を参照してください。

サンプル情報の表示 個々のサンプル ファイルに関連付けられた情報や Raw Data を表示するには、左側のナビ ゲーション ペインでサンプル ファイルを選択してから、「Info」タブ、または「Raw Data」 タブを選択します。

	\varTheta GeneMapper - SNaPsho	t Ref Data Project - gm is Logged in Database FRMINFODEVD01							
	File Edit Analysis View Tools Help								
	😂 😂 📗 🛛 🍒 🚰	🔟 📰 🔛 📗 🛅 🕨 💕 Table Setting: SNaPshot Default 🛛 🖬 🔎	🎍 🛛 🙆						
サンプル ファイル ― を選択し、 サンプル情報と Raw Data を表示	Approject SNAPshot 2,803,03 faa 5,503,09 faa 5,503,09 faa 5,503,09 faa 5,503,11 faa	Mto Rawy Data EPT Data Sample Information Sample File : 6_F03_11.fsa Sample Origin Path : C:\AppliedBiosystems\GeneMapper\Example Data SNaPhot\SNaPshot\G_F03_11.fsa Status Message : Good Firer Message : Disk media Error Message : None Current Settings							
	Analysis Completed.		Stop						

00177

サンプル プロット の表示

- サンプルのプロットを表示するには、次の手順を実行します。
 - 1. 「View」 ▶ 「Samples」の順に選択し、「Samples」タブを表示します。
 - 2. 「Samples」タブでサンプル (列) を選択します。複数のサンプルを選択するには、[Shift] または [Ctrl] キーを押したままにします。すべてのサンプルを選択するには、「Edit」 → 「Select All」の順に選択します。
- 3. **IIII** をクリックします (「Analysis」 ▶ 「Display Plots」)。

「Samples Plot」ウィンドウに、選択された各サンプルに対するエレクトロフェログラム が表示されます。



4. Plot Setting には SNaPshot Default を選択します。

注: Plot Settings は、解析後、「Samples Plot」ウィンドウに表示される情報を制御し ます。SNaPshot Default は、GeneMapper ソフトウェアで提供されるデフォルトの Plot Settings の 1 つです。GeneMapper Manager で Plot Settings をカスタマイズし、編集 および作成も可能です。詳細については、GeneMapper[®] Software Online Help を参照 してください。

5. Samples Plot 内の X 軸および Y 軸上で拡大表示する方法は次のとおりです。

目的	操作
X 軸の特定の領域を 拡大表示	すべてのプロットを拡大するには、上部 X 軸上にカーソルを置き、 を 左右ヘクリックしてドラッグします。選択したプロットのみを拡大する には、[Shift] キーを押したまま、クリックしてドラッグします。
	または
	上部 X 軸で右クリックし、「Zoom To」を選択して、範囲を入力し、「OK」 をクリックします。
Y 軸の特定の領域 を拡大表示	左側の Y 軸上にカーソルを置き、🔍 を上下にクリックしてドラッグします。
	または
	左側の Y 軸を右クリックし、「Zoom To」を選択します。最大値を入力 し、オプションで「Apply to all electropherograms」を選択して、「OK」 をクリックします。
縮小表示	X 軸または Y 軸をダブルクリックします。
	または
	X 軸または Y 軸を右クリックし、「Full View」を選択します。



「Samples Plot」 ウィンドウでの データの検証 「Samples Plot」ウィンドウでは、その他にも次のようなタスクを実行できます。

- X 軸(塩基対またはデータポイント)のスケール調整
- Y 軸 (個々のサンプルの最大値、すべてのサンプルの最大値、特定の値)のスケール調整
- 特定の蛍光色素色のピークの表示と非表示
- 個々のピークに対するステータス ラインの表示
- 検出された各ピークに対するサイジング情報の列を表示する Sizing テーブルの表示
- 検出された各ピークに対するジェノタイピング情報の列を表示する Genotypes テーブ ルの表示
- ピークの選択と、Sizing テーブル内での対応するデータ列のハイライト

18ページの図 2-2 では、上記の機能の一部を図解して示しています。

上記の機能の使用に関する詳細については、[F1] キーを押し、GeneMapper[®] Software Online Help から該当の項目を選択してください。

注:「Samples Plot」ウィンドウを使用して、異なるサンプル ファイルのアレル ピークを容易に比較し、リファレンス データとして使用するサンプル ファイル、および Bin に対する最小フラグメント長と最大フラグメント長を決定することができます。この情報は、リファレンス データを Kit に追加する際 (21 ページ)、および Marker や Bin を作成する際 (24 ページ) に有用です。

Samples Plot の表示を確認したら、区をクリックして、ウィンドウを閉じます。

次の手順

19ページの説明に従って、Kit、Panel、Marker、Bin Set、Bin を作成します。







Panel Manager 内に次の階層的オブジェクトを作成します。

- Kit Panel のグループ
- Panel Marker のグループ
- Marker フラグメント サイズ レンジ (bp)

注: このガイドでは、Panel と Marker の作成方法について説明します。ただし、Marker 情報を含む Panel (テキスト ファイル) をインポートすることもできます。Panel のインポート に関する詳細については、GeneMapper[®] Software Online Help を参照してください。

また、Panel Manager を使用して、次を作成します。

- Bin Set Bin の集合
- Bin アレル定義。すなわちフラグメント サイズ(bp) と蛍光色素色

Bin Set を作成する前に、Kit を選択する必要があります。SNP Kit には、Bin Set を 1 つだ け作成できます。作成された Bin Set は、その SNP Kit 内の任意の Panel と関連付けること ができます。

Bin を作成する前に、Panel および Bin Set を選択する必要があります。Bin は、選択された Panel 内の Marker と関連付けられ、選択された Bin Set 内に保存されます。

注: この章では、リファレンス データを使用して Bin を手動で作成する方法について説明し ます。ただし、Primer Focus Kit データと Auto Panel 機能を使用して、Bin を自動的に作成 するか、または Bin 情報を含む Bin Set (テキスト ファイル) をインポートすることもできま す。Bin の自動作成に関する詳細については、第3章「Primer Focus® Kit データを使用し た Panel および Bin Set の作成」を参照してください。Bin Set のインポートに関する詳細に ついては、GeneMapper[®] Software Online Help を参照してください。

Kit と Panel の 作成

概要

- Kit と Panel を作成するには、次の手順を実行します。
 - 1. III をクリックして Panel Manager を開きます(「Tools」) 「Panel Manager」)。
 - 左側のナビゲーション ペイン最上部で Panel Manager を選択し、 で をクリックしま す(「File」) 「New Kit」)。





3. 「New Kit」ダイアログ ボックスで、「Kit Name」に「SNaPshot Kit」と入力し、「Kit Type」では「SNP」を選択して、「OK」をクリックします。

New Kit	
Kit name	SNaPshot Kit
Kit type:	SNP
	OK Cancel

左側のナビゲーション ペインに SNaPshot Kit が表示されます。

🗢 Panel Manager
File Edit Bins View
👑 🗙 📑 🖬 🖬 🗰 🛤 Bin Set:
 ■ 品 Panel Manager ● ☐ Microsatellite Kit ● ☐ LOH Kit ● SNaPshot Kit

- 4. ナビゲーション ペインで SNaPshot Kit を選択し、 「New Panel」)。
- 5. Panel Manager の右側のペインで「New Panel」を選択し、「Panel Name」に「SNaPshot 6plex Panel」と入力してから、[Enter] キーを押します。

	Panel Name	Comment
1	SNaPshot 6plex Panel	none

左側のナビゲーション ペインで、SNaPshot Kit の下に SNaPshot 6plex Panel が表示さ れます。





Bin Set の作成 Bin Set を作成するには、次の手順を実行します。

- 1. 左側のナビゲーション ペインで、19 ページ で作成した SNaPshot Kit を選択します。
- 2. III をクリックします (「Bins」 ▶ 「New Bin Set」)。
- 3. 「New Bin Set」ダイアログ ボックスで、「Bin Set Name」に「SNaPshot Bin Set」と 入力し、「OK」をクリックします。

New Bin Set	X
Please enter name of new Bin Set. SNaPshot Bin Set	

Panel Manager 最上部の「Bin Set」ドロップダウン リストに、SNaPshot Bin Set が追加されます。これで、SNaPshot Bin Set を SNaPshot Panel(または SNaPshot Kit に追加された他の任意の Panel)と関連付けることができます。

リファレンス データ
 注: プロジェクト内のすべての、または一部のサンプル ファイルをリファレンス データとし
 て追加できます。サンプル ファイルの数が少ないため、すべてのサンプル ファイルをリファレンス データとして使用します。サンプル ファイルの数が多い場合は、サンプル ファイル内
 に存在するすべてのアレルが含まれるサンプル ファイルに絞って選択します。

SNaPshot Kit にリファレンス データを追加するには、次の手順を実行します。

- 日本
 日
- 1. 左側のナビゲーション ペインで、19ページ で作成した SNaPshot Kit を選択します。

2. $\boxed{\mathbf{k}}$ \mathcal{E} \mathcal{E} 2. \mathbf{k} \mathcal{E} 2. \mathbf{k} \mathcal{E} 2. \mathcal{E} 2. \mathbf{k} \mathcal{E} 2.

「Add SNP Reference Data」ダイアログ ボックスが開きます。



3. 「Search」をクリックします。

左下のペインに、GeneMapper データベース内に保存されている、サイジング済みのす べてのプロジェクトが表示されます。

4. SNaPshot Ref Data Project を展開し、SNaPshot フォルダを選択して、「Add to List」、 「Add」の順にクリックします。

SNaPshot Data フォルダ内の 3 つのサンプル ファイルがすべて、リファレンス サンプ ルとして SNaPshot Kit に追加され、Panel Manager のナビゲーション ペインの下半分 に表示されます。

😔 Panel Manager 🛛 🔀					
File Edit Bins View					
	Bin Set: SNaPshot Bin Set	✓ IIÎî	🈼 🔳 🖹 🕌		
😑 🚠 Panel Manager	Panel Name	Comment.			
🕀 🚞 Microsatellite Kit	1 SNaPshot 6plex Panel	none			
Reference Samples					
2 B03 03.fsa					
5_E03_09.fsa					
6_F03_11.fsa					
OK Cancel Apply					



リファレンス データ の Panel への追加 リファレンス データを Kit に追加した後は、Kit 内の個々の Panel にサンプル ファイルを追 加できます。これは、リファレンス データのパネル化とも呼ばれます。

SNaPshot 6plex Panel にリファレンス データを追加するには、次の手順を実行します。

- **1.** 上部のナビゲーション ペインで、**19** ページ で作成した **SNaPshot 6plex Panel** を選択 します。
- 2. 「Bins」 → 「Panel Reference Data」の順に選択します。

SNP Reference Data		
SNP Kit reference data:		SNP Panel reference data:
SNaPshot Kit ● 2_B03_03.tsa ● 5_E03_09.tsa ● 6_F03_11.tsa	>	 SNaPshot 6plex Panel
Remove		Remove
0	КЦС	Cancel

- 3. 「SNP Reference Data」ダイアログ ボックスで、SNaPshot Kit フォルダを選択し、 [Shift] キーを押したまま、3 つのサンプル ファイルをすべて選択します。
- 4. をクリックすると、SNaPshot Kit フォルダの 3 つのサンプル ファイルが SNaPshot 6plex Panel に追加されます。
- 5. 「OK」をクリックします。



リファレンス データ 3 つのリファレンス サンプル ファイルから Marker と Bin を手動で作成するには、次の手順 を実行します。
 Bin の手動による 作成 1.「Panel Manager」ウィンドウで、「Plot」タブを選択します。
 2. ナビゲーションペインで、次を行います。

a. 上部のナビゲーションペインで、SNaPshot 6plex Panel を選択します。

b. 下部のナビゲーションペインで、5_E03_09.fsa サンプルファイルを選択します。

5_E03_09.fsa サンプル ファイルのエレクトロフェログラムが、「Plot」タブに表示されます。



- 3. 次のさまざまな方法で、実際にデータを表示します。
 - •「View」メニューからコマンドを選択するか、X 軸および Y 軸のラベルを右クリックして、X 軸および Y 軸のスケールを変更します。
 - ツールバーのカラーボタンをクリックして、特定の蛍光色素ピークを表示または非 表示にします。

	1000		
		_	_

 エレクトロフェログラム内で、カーソルを X 軸上に置き、Q を左右にクリックしてド ラッグすることで、最初の緑のピーク(ほぼ 25 塩基対の位置に存在するホモ接合体アレ ル)を拡大表示します。



5. 「File」 → 「New SNP Marker」を選択します。

🛛 Add Marker 🛛 🗙						
Marker Name: 25mer						
Allele	Name	Min	Max			
Blue	G					
Green	A	24.5	26.0			
Yellow	c					
Red	т					
<u></u>	OK Cancel					

- 6. 「In the Add Marker」ダイアログボックスで、次の手順を実行します。
 - a. 「Marker Name」に「25mer」と入力します。
 - b. Green Allele の「Min」に 24.5、「Max」に 26.0 とそれぞれ入力します。
 - c.「Name」については、デフォルト設定である ddNTP 塩基の名前(G、A、C、T) をそのまま使用します。これらのアレル名は、Primer Focus Kit データから作成し た Bin に名前をつける際にも使用されています。
 - d.「OK」をクリックして、Marker 情報と Bin 情報を SNaPshot 6plex Panel に保存 し、エレクトロフェログラムに Marker と Bin を表示します。



7. 縮小して他のアレル ピークを表示するには、 U をクリックします(「View」 ト「Full View」)。



8. 次の SNP マーカである青および緑のピーク(ほぼ 28 塩基対の位置に存在するヘテロ接合体アレル)を拡大表示します。



9. 手順 5 ~ 6 を繰り返し、次に示すように別の Marker を作成します。

🖷 Add Marker 🛛 🔀						
Marker Name: 28mer						
Allele	Name	Min	Max			
Blue	G	27.3	28.9			
Green	А	28.3	29.8			
Yellow	c					
Red	т					
OK Cancel						

作成された Marker と Bin は、SNaPshot 6plex Panel に保存され、下の図のように表示 されます。



注: GeneMapper ソフトウェアでは、ユーザが Bin に対して定義するサイズ レンジに 基づいて、Marker のサイズ レンジを自動的に算出します。


10. 下の表に示す Bin の色、名前、レンジを使用して上記の手順を繰り返し、残り 4 つの SNP マーカ (39mer、48mer、54mer、62mer) について Marker と Bin を作成します。

Marker Name	Allele Color	Allele Name	Min	Мах
39mer	Blue	G	38.4	39.9
48mer	Red	Т	47.2	48.6
54mer	Yellow	С	53.3	54.5
	Red	Т	53.8	55.2
62mer	Yellow	С	61.5	62.8

注: 原則として、ビン レンジは 1.5 bp 前後の設定が適しています。



ここまでの手順を完了すると、プロットは下の図のようになります。

 下部のナビゲーション ペインで、SNaPshot 6plex Panel に追加した他の 2 つのリファ レンス サンプル ファイル (2_B03_03.fsa と 6_F03_11.fsa) を選択します。「Plot」タブ で、各サンプル ファイル内のピークが、5_E03_09.fsa リファレンス サンプル ファイル から作成した Bin にどの程度フィットしているかを確認します。



重要! 通常、各リファレンス サンプル ファイルには、先述のサンプル ファイルから作 成された Bin に該当しないアレルピークが含まれます。手順 4 ~ 7 を繰り返して、こう したピークに Bin を作成する必要があります。



Marker と Bin の 閲覧

Marker 情報および Bin 情報を閲覧するには、次の手順を実行します。

 Panel Manager で、上部ナビゲーションペインの SNaPshot 6plex Panel を選択し、下 部ナビゲーションペインでは Reference Samples を選択してから、「Plot」タブを選択 します。

Panel には、下部 X 軸に表される Marker と、上部 X 軸に表される Bin が表示されます。



Panel Manager」ウィンドウで、「Table」タブを選択します。
 Marker 情報および Bin 情報が、テーブル形式で表示されます。

Panel Manager									×
File Edit Bins View									
	B	n Set: SNaPsho	t Bin Set	•	· III 5.				
😑 🚠 Panel Manager	Tal	ole Plot							
GMicrosatelite Kit		Marker Name	Marker Min	Marker Max	Bin 1 Name	Bin 1 Min	Bin 1 Max	Bin 1 Dye	
E CONADebat Kit	1	25mer	24.50	26.00	A	24.50	26.00	Green	^
SNaPshot 6plex Panel	2	28mer	27.30	29.80	G	27.30	28.90	Blue	
	3	39mer	38.40	39.90	G	38.40	39.90	Blue	
	4	48mer	47.20	48.60	T	47.20	48.60	Red	
	5	54mer	53.30	55.20	C	53.30	54.50	Yellow	
	6	62mer	61.50	62.80	с	61.50	62.80	Yellow	
And Andrews Samples									×
	<							,	
			OK Cano	el Apply					

3. 「Apply」ボタンをクリックして、作成した Marker、Bin 情報を保存します。

Bin Set の採用 新規 Bin Set を採用して、Panel Manager を閉じるには、「OK」をクリックします。

Bin と Marker の
編集(オプション)このガイドの実験を完了するのに、Marker や Bin を追加、編集、削除する必要はありませ
ん。ただし、Panel Manager を開き、SNaPshot Kit や SNaPshot 6plex Panel を選択して、
これらの機能をテストすることはできます。

重要! 29 ページからの説明に従い、SNaPshot 6plex Panel の Marker や Bin を編集または 削除した場合は必ず、Panel Manager の下部にある「Cancel」をクリックしてください。 「OK」または「Apply」をクリックすると、解析の結果に悪影響が出る場合があります。

Marker に Bin を追加するには

- 1. 「Plot」タブ(下部 X 軸)または「Table」タブ(列)で、Marker を選択します。
- 2. **i** をクリックします (「Bins」 ▶ 「Add Bin」)。
- **3**. 「Edit SNP Marker」ダイアログ ボックスで、Bin の「Name」、「Min」、「Max」を入力 し、「OK」をクリックします。

Bin を編集するには

- **1**. 「**Plot**」タブで、**Bin**(上部 **X** 軸)を選択します。
- 2. をクリックするか(「Bins」 → 「Edit Bin」)、または Bin を右クリックして、「Edit SNP Marker」を選択します。
- **3**. 「Edit SNP Marker」ダイアログボックスで、Bin の「Name」、「Min」、「Max」を編集 し、「OK」をクリックします。

または

- 1. 「Table」タブを選択します。
- 2. 次の情報を編集します。
 - Bin Name
 - Bin Min
 - Bin Max

Bin を視覚的に編集するには

- **1**. 「**Plot**」タブで、**Bin**(上部 **X** 軸)を選択します。
- 2. 青いセンター ラインをクリックしてドラッグし、Bin の位置を決定します。
- 3. 右ハンドルの左側をクリックしてドラッグし、Bin のオフセット(レンジ)を決定します。

```
注:変更した内容を直前の状態に戻すには、「Edit」→「Undo」を選択します。
```





Bin を削除するには

Bin を Marker から削除するには、次の手順を実行します。

- 「Plot」タブまたは「Table」タブで Bin を選択し、
 「Delete Bin」)。
 または
- 「Plot」タブで、Bin(上部 X 軸)を選択し、Bin を右クリックして、「Delete Bin」を 選択します。

Marker を編集するには

- 1. 「Plot」タブで、Marker (下部 X 軸)を選択します。
- 2. ▲ をクリックするか(「Bins」 → 「Edit SNP Marker」)、または Marker を右クリックして、「Edit SNP Marker」を選択します。
- **3.**「Edit SNP Marker」ダイアログ ボックスで、「Marker Name」を編集し、「OK」をクリックします。

または

「Table」タブで、「Marker Name」を編集します。

Panel から Marker を削除するには

Panel から **Marker** を削除するには、「**Plot**」タブで **Marker**(下部 **X** 軸)を選択し、**Marker** を右クリックして、「**Delete Marker**」を選択します。

次の手順 第4章の説明に従って、SNaPshot 解析を設定します。





新規プロジェクトの作成と Primer Focus[®] Kit サンプル ファイル の追加

概要

「GeneMapper」ウィンドウ内でプロジェクトを作成し、プロジェクトにサンプルを追加します。

プロジェクトの 新規作成とサンプル ファイルの追加 プロジェクトを新規作成してサンプル ファイルを追加するには、次の手順を実行します。

1. 👸 をクリックします (「File」 ▶ 「New Project」)。

Project Type	
🔘 Generic	
O Microsatellite	
SNaPshot®	
🔘 OLA Analysis	
O SNPlex™	
○ AFLP	

- 2. 「New Project」ダイアログ ボックスで、SNaPshot を選択し、「OK」をクリックします。 GeneMapper v3.7 ではこのダイアログボックスは表示されません。そのままステップ 3 へ進みます。
- 3. 「GeneMapper」ウィンドウで、 [] をクリックします (「File」 ▶ 「Add Samples to Project」)。
- 4. 「Add Samples to Project」ダイアログボックスの「Files」タブで、次のパスを確認します。 *<drive>*:\AppliedBiosystems\GeneMapper\Example Data\SNaPshot

注:上記の場所は、GeneMapper[®] ソフトウェアがインストールされているドライブに よって異なります。デフォルトでは、Dドライブにインストールされます。



5. Primer Focus フォルダを選択し、「Add to List」、「Add」の順にクリックします。

注: このガイドでは、**Primer Focus** フォルダ内の 6 つのサンプル ファイルをすべて追加しましたが、フォルダ内のファイルの一部(1 つまたは 2 つ)を追加することも可能です。これを行うには、左側のペインでフォルダを展開し、[**Ctrl**] キーを押したまま、ファイルを個々に選択して、「**Add To List**」をクリックします。

🗢 Add Samples to Project	
Edit View	
Files GM Database My Computer My Computer Mig 35 Roppy (A:) CD-RW Drive (E:) Contempleter Contempleter Contempleter Contempleter	Samples To Add: Primer Focus 2 20Sample1 fsa 3 28Sample2 fsa 4 4Sample4 fsa 5 22Sample5 fsa 6 60Sample6 fsa
Add To List >> Options	ear Add Add & Analyze Cancel

Primer Focus フォルダ内の 6 つのサンプル ファイルが、互換性のある **Applied Biosystems** 社製シーケンサの **Data Collection** ソフトウェアに入力した情報と 共に、「**Samples**」タブ上のテーブルに表示されます。

🗢 GeneMapper - "Untitled - gm Is Logged In Database FRMINFODEVD01									
File Edit Analysis Viev	File Edit Analysis View Tools Help								
😂 😂 🛃 😼 🖻	😂 🔁 📊 🍹 🚰 📗 🔛 🛄 🔛 🖽 🔚 🔚 🕨 🕨 Table Setting: Microsatelite Default 🔽 🖬 👂 🌭 🖉 🥝								
😑 🚠 Project	B 器Project Samples Genotypes								
🗄 🗀 Primer Focus	[[Status	Sample File	Sample Name	Comments	Sample Type	Analysis Method	Panel	Size Standard
	1	N	20Sample1.fsa	20	None	Sample	None	None	None
	2	<u> </u>	28Sample2.fsa	28	None	Sample	None	None	None
	3	1	36Sample3.fsa 36		None	lone Sample None		None	None
	4	1	44Sample4.fsa	44	None	Sample	None	None	None
	5	1	52Sample5.fsa	52	None	Sample	None	None	None
	6	8	60Sample6.fsa	60	None	Sample	None	None	None
	<	<	1				1	1	>
< >									
Progress Status								0%	Stop



34ページの説明に従って、プロジェクト用のアナリシスパラメータと表示設定を作成します。



概要

アナリシス パラメータと Table Settings の設定

「GeneMapper」ウィンドウ内で、プロジェクト用のアナリシス パラメータと表示設定を確認 します。

アナリシスパラメータには、次の項目が含まれます。

- Sample Type
- Analysis Method (Bin Set を含む)
- Size Standard
- Panel (Marker の集合)

プロジェクトにおけるすべてのサンプル ファイルを解析するために GeneMapper[®] ソフト ウェアが使用する、ピーク検出アルゴリズム、サイジング アルゴリズム、ジェノタイピング アルゴリズムを決定するアナリシス パラメータを設定します。

表示設定には、Table Settings と Plot Settings が含まれます。

アナリシス パラ プロジェクト用のアナリシス パラメータを設定するには、次の手順を実行します。

メータの設定

- 1. 「GeneMapper」ウィンドウで「Samples」タブを選択します。
- 2. 「Sample Type」カラムで最初の列をクリックし、ドロップダウン リストから Primer Focus を選択します。



- 3. 「Analysis Method」カラムで最初の列をクリックし、ドロップダウン リストから SNaPshot Default を選択します。
- 4. 「Panel」カラムの設定を None のままにします。

注: Panel の選択は、オプションです。Panel を選択しない場合、GeneMapper ソフト ウェアでは、データをサイジングしますが、ジェノタイピングは行いません。このプロ ジェクト作成の目的は、リファレンス データのサイジング専用解析を実行して、Panel と Bin Set の作成に使用することです。従って、アナリシス パラメータの一部として Panel を選択することはありません。

5.「Size Standard」カラムで最初の列を選択し、ドロップダウン リストから GS120LIZ を選択します(これは、サンプルでランされたサイズ スタンダードです)。



- 6. 「Samples」タブのすべてのサンプル列に、次のとおり選択をフィルダウンします。
 - a. 「Analysis Method」、「Panel」、「Size Standard」の各カラム ヘッダにまたがって クリックしてドラッグし、3 つのカラムのすべての列をハイライト表示します。

Sample Type	Analysis Method	Panel	Size Standard —
Primer Focus	SNaPshot Default	None	GS120LIZ
Sample	None	None	None
Sample	None	None	None
Sample	None	None	None

b.「Edit」→「Fill Down」の順に選択します(または、[Ctrl] キーと [D] キーを同時 に押します)。

Table Settingsの「GeneMapper」ウィンドウの最上部で、「Table Settings」ドロップダウン リストから選択SNaPshot Default を選択します。

Microsatellite Default	*
New	
AFLP Default	
Microsatellite Default	
SNPlex_v3	
SNaPshot Default	

Table Settings は、解析後に「**Samples**」タブおよび「**Genotypes**」タブで表示される情報を 制御します。**SNaPshot Default** は、**GeneMapper** ソフトウェアで提供されるデフォルトの **Table Settings** の1つです。

GeneMapper Manager でカスタマイズした Table Settings を編集および作成することもで きます。詳細については、GeneMapper[®] Software Online Help を参照してください。

プロジェクトの保存 プロジェクトを保存するには、次の手順を実行します。

- 1. $\boxed{1}$ $\boxed{5}$ $\boxed{5}$
- 2. 「Save Project」ダイアログ ボックスで、「SNaPshot Primer Focus Data Project」と 入力し、「OK」をクリックします。

Save Pro	oject	\mathbf{X}
٩	Project name: SNaPshot Primer Focus Data Project	t

「GeneMapper」ウィンドウのタイトル バーに、SNaPshot Primer Focus Data Project が表示されます。

次の手順

36ページの説明に従って、プロジェクトのサイジングのみを実行します。



プロジェクトでのサイジングのみの実行

概要

プロジェクト用のサンプル ファイルの追加と、アナリシス パラメータの設定が完了したら、 次に、サイジングのみを実行します。これにより、サンプル ファイルをリファレンス データ として利用し、**Bin**(アレル定義)を作成できるようになります。

サイジングのみを実行するには、次の手順を行います。

- プロジェクトの解析
- SQ および関連する PQV の閲覧
- サイズスタンダードの検証
- サンプル情報の確認(Raw Data を含む)
- サンプル プロットの確認

プロジェクトの解析 > をクリックします (「Analysis」 **>** 「Analyze」)。

GeneMapper[®] ソフトウェアは、プロジェクト内の各サンプルを解析し、「GeneMapper」ウィンドウの左下にあるステータスバーに進捗状況を表示します。

SQ と PQV の閲覧 Size Quality (SQ) および関連する Process Quality Value (PQV) を閲覧するには、次の手順を実行します。

- 1. 「GeneMapper」ウィンドウの左下にあるステータス バーに、「Analysis Completed」と 表示されていることを確認します。
- 2. 「Samples」タブの右側をスクロールして、SQを閲覧します。

🕏 GeneMapper - SNaPshot Primer Focus Data Project - gm Is Logged In Database FRMINFODEVD01														
File Edit Analysis View Tools Help														
😂 🗁 📓 🍢 🖻	۳ ۳	Ш 🗉 Ш	🕅 🌐 🗎	🕨 🍯	Table Setting:	SN	aPshot	Default		~ [P		a 🕐
😑 🚠 Project	Sample	es Genotypes												
🕀 🧀 Primer Focus		Run Name	Instrument Type	Run Date	& Time	REF	SQI	SFNF	MNF	SNF	OS	SQ	UD1	UD2
	1	Primer Focus	ABI310	2001-03-	08 17:09:11.0						<u> </u>			
	2	Primer Focus	ABI310	2001-03-	08 17:09:11.0									
	3	Primer Focus	ABI310	2001-03-0	08 17:09:11.0									
	4	Primer Focus	ABI310	2001-03-	08 17:09:11.0									
	5	Primer Focus	ABI310	2001-03-0	08 17:09:11.0									
	6	Primer Focus	ABI310	2001-03-0	08 17:09:11.0	İ	İ –						1	
		<								ų.			<u> </u>	>
< <u> </u>														
Analysis Completed.														Stop

| スクロール バーを右方向へドラッグして「SQ」カラムを表示

手順に従ってこのガイドに示されるデータ例を使用した場合、各サンプルに対する SQ は (Pass) になります。一方、SQ (SFNF、MNF、SNF、OS) の生成に関連する PQV もほとんど (になります。



黄色の 🝐 と赤い 🔵 SQ の原因調査 **重要!** ユーザ独自のデータを解析する際、SQ が ▲ (Check) または ● (Low Quality) とな り、関連付けられた PQV (SFNF、MNF、SNF、OS) が ▲ になることがあります。これ は、サイズ スタンダード、データ、またはアナリシス パラメータに問題があることを示して います。これらの問題の原因を調査して修正するには、37 ページの「サイズ スタンダードの 検証」を参照してください。

注: 💕 をクリックすると、SQ スコアに基づいて、サンプルをソートすることができます。 「Samples」タブの一番上に、🔵 SQ の付いたサンプルがリスト表示されます。

サイズ スタンダード の検証

サイズ スタンダードを検証するには、次の手順を実行します。

- 1. 「Edit」 ▶ 「Select All」の順に選択して、「Samples」タブ内のすべてのサンプルを選択 します。
- 2. <u>Ⅲ</u> をクリックして、Size Match Editor を開きます(「Analysis」 **▶** 「Size Match Editor」)。



- 図 3-1 Size Match Editor 「Size Matches」タブ
 - 3. 「Size Matches」タブをクリックして、選択されたサンプルに対する次の項目を表示します。
 - Size Quality (SQ) スコア
 - サイズ スタンダード ピーク
 - サイズ スタンダード ピーク ラベル
 - 4. サンプルの Sizing Quality スコア (図 3-1) を確認します。このスコアは、サイズ スタ ンダードからのデータが、ソフトウェアで選択されたサイズ スタンダードにどの程度適 合しているかを示します。このスコアによって、SQ が ■ (Pass)、 (Check)、 (Low Quality) のいずれを表示するかが決まります。

このガイドの指示に従った場合、Sizing Quality は > 0.75 となり、SQ は 📄 (Pass) を 表示します。

一方、独自のデータを解析すると、Sizing Quality は低下し、SQ は ▲ (Check)、また
 は ● (Low Quality) を表示します。トラブルシューティングについては、38 ページの
 表 3-1 を参照してください。

5. サイズ スタンダード内のすべてのピークが存在し、正確に標識されているかを判定します。

このガイドの指示に従った場合、図 3-1 に示すとおり、すべてのピークが存在し、正確に標識されます。

一方、独自のデータを解析すると、一部のサイズスタンダード ピークは正確に標識されないか、失われていることがあります。トラブルシューティングについては、38ページの表 3-1 を参照してください。



6. 「Size Calling Curve」タブをクリックして、選択したサンプルのサイズ スタンダード曲 線を表示します。サイズ スタンダードのフラグメントを表す赤いデータ ポイントと、黒 いベストフィット曲線が表示されます。



- 7. Size Match Editor の左側ペインで別のサンプルを選択し、手順 3~6を繰り返します。
- 8. 「OK」をクリックして、Size Match Editor を閉じます。

問題	操作
Sizing Quality スコアが低下し、SQ は 🍐 (Check) または 🍎 (Low Quality) を表示するが、すべてのサ イズ スタンダード ピークが存在し、 正確に標識されている。	「Size Matches」タブの最上部で「Override SQ」をク リックして、Sizing Quality を上書きします(図 3-1)。 Sizing Quality スコアの変更を 1.0 に上書きします。こ れは、ユーザがサイズ スタンダードを確認したことを示 します。
いくつかのサイズ スタンダード ピークが正確に標識されない。	「Size Matches」タブで、サイズ ラベルを編集、削除、 追加します。次に「Apply」をクリックし、更新された サイジング情報でデータを再解析します。詳細について は、GeneMapper [®] Software Online Help を参照して ください。
いくつかのサイズ スタンダード ピークが存在しない。	ソフトウェアで、カスタム サイズ スタンダードを作成 します。詳細については、GeneMapper [®] Software Online Help を参照してください。

サイジングに関するトラブルシューティングやその他の情報は、『GeneMapper[®] Software Reference and Troubleshooting Guide』を参照してください。

サンプル情報の表示 個々のサンプル ファイルに関連付けられた情報や Raw Data を表示するには、左側のナビ ゲーション ペインでサンプル ファイルを選択してから、「Info」タブ、または「Raw Data」 タブを選択します。





サンプル プロット の表示

- サンプルのプロットを表示するには、次の手順を実行します。
 - 1. 「View」 ▶ 「Samples」の順に選択し、「Samples」タブを表示します。
- 2. 「Samples」タブでサンプル (列) を選択します。複数のサンプルを選択するには、[Shift] または [Ctrl] キーを押したままにします。すべてのサンプルを選択するには、「Edit」 → 「Select All」の順に選択します。
- 3. **I** をクリックします (「Analysis」 ▶ 「Display Plots」)。

「Samples Plot」ウィンドウに、選択された各サンプルに対するエレクトロフェログラム が表示されます。



4. Plot Setting には SNaPshot Default を選択します。

注: Plot Settings は、解析後、「Samples Plot」ウィンドウに表示される情報を制御しま す。SNaPshot Default は、GeneMapper ソフトウェアで提供されるデフォルトの Plot Settings の 1 つです。GeneMapper Manager で Plot Settings をカスタマイズし、編集 および作成も可能です。詳細については、GeneMapper[®] Software Online Help を参照 してください。

5. Samples Plot 内の X 軸および Y 軸上で拡大表示する方法は次のとおりです。

目的	操作
X 軸の特定の領域を 拡大表示	すべてのプロットを拡大するには、上部 X 軸上にカーソルを置き、 を左右ヘクリックしてドラッグします。選択したプロットのみを拡大 するには、[Shift] キーを押したまま、クリックしてドラッグします。
	または
	上部 X 軸で右クリックし、「Zoom To」を選択して、範囲を入力し、 「OK」をクリックします。
Y 軸の特定の領域を 拡大表示	左側の Y 軸上にカーソルを置き、Q を上下にクリックしてドラッグ します。
	または
	左側の Y 軸を右クリックし、「Zoom To」を選択します。最大値を入 カし、オプションで「Apply to all electropherograms」を選択して、 「OK」をクリックします。
縮小表示	X 軸または Y 軸をダブルクリックします。
	または
	X 軸または Y 軸を右クリックし、「Full View」を選択します。



「Samples Plot」 ウィンドウでの データの検証 「Samples Plot」ウィンドウでは、その他にも次のようなタスクを実行できます。

- X 軸(塩基対またはデータポイント)のスケール調整
- Y 軸 (個々のサンプルの最大値、すべてのサンプルの最大値、特定の値)のスケール調整
- 特定の蛍光色素色のピークの表示と非表示
- 個々のピークに対するステータス ラインの表示
- 検出された各ピークに対するサイジング情報の列を表示する Sizing テーブルの表示
- 検出された各ピークに対するジェノタイピング情報の列を表示する Genotypes テーブ ルの表示
- ピークの選択と、Sizing テーブル内での対応するデータ列のハイライト

40ページの図 3-2 では、上記の機能の一部を図解して示しています。

上記の機能の使用に関する詳細については、[F1] キーを押し、GeneMapper[®] Software Online Help から該当の項目を選択してください。

Samples Plot の表示を確認したら、区をクリックして、ウィンドウを閉じます。

次の手順

41ページの説明に従って、Kit、Panel、Marker、Bin Set、Bin を作成します。



図 3-2 Samples Plot 内の異なるサンプル ファイルからのデータの検証と比較



Kit、Panel、Marker、Bin Set、Bin の作成

概要

Panel Manager 内に次の階層的なオブジェクトを作成します。

- Kit Panel のグループ
- Panel Marker のグループ
- Marker フラグメント サイズ レンジ (bp)

注: このガイドでは、Panel と Marker の作成方法について説明します。ただし、Marker 情報を含む Panel (テキスト ファイル) をインポートすることもできます。Panel のインポート に関する詳細については、GeneMapper[®] Software Online Help を参照してください。

また、Panel Manager を使用して、次を作成します。

- Bin Set Bin の集合
- Bin アレル定義。すなわちフラグメント サイズ(bp)と蛍光色素色

Bin Set を作成する前に、Kit を選択する必要があります。SNP Kit には、Bin Set を 1 つだ け作成できます。作成された Bin Set は、その SNP Kit 内の任意の Panel と関連付けること ができます。

Bin を作成する前に、Panel および Bin Set を選択する必要があります。Bin は、選択された Panel 内の Marker と関連付けられ、選択された Bin Set 内に保存されます。

注: この章では、Primer Focus Kit データ、および Auto Panel 機能を使用して Bin を自動 的に作成する方法について説明します。ただし、Bin を手動で作成するか、または Bin 情報を 含む Bin Set (テキスト ファイル) をインポートすることもできます。Bin の手動作成に関す る詳細については、第2章「リファレンス データを使用した Panel および Bin Set の作成」 を参照してください。Bin Set のインポートまたは Bin の手動作成に関する詳細については、 GeneMapper[®] Software Online Help を参照してください。

Kit の作成

Kit を作成するには、次の手順を実行します。

- 1. III をクリックして Panel Manager を開きます(「Tools」) 「Panel Manager」)。
- 2. 左側のナビゲーション ペイン最上部で Panel Manager を選択し、 す (「File」 → 「New Kit」)。

Panel Manager								
File Ed	it Bins	View						
📑 ×				. E	Bin Set:			
	'anel Mana	iger Q						



3. 「New Kit」ダイアログ ボックスで、「Kit Name」に「Primer Focus Kit」と入力し、 「Kit Type」では「SNP」を選択して、「OK」をクリックします。

New Kit	
kît neme	Primer Focus Kit
rut nome	
Kit type:	SNP 💌
	OK Cancel

左側のナビゲーション ペインに Primer Focus Kit が表示されます。

Panel Manager	
File Edit Bins View	
📽 🗙 📲 🖩 🗎 🗎	📕 🛛 Bin Set:
 Anal Manager Microsatellite Kit CH Kit SNaPshot Kit Primer Focus Kit 	

Bin Set の作成

Bin Set を作成するには、次の手順を実行します。

- 1. 左側のナビゲーションペインで、41ページで作成した Primer Focus Kit を選択します。
- 2. $\parallel \hat{l} l$ prime pr
- 3. 「New Bin Set」ダイアログ ボックスで、「Bin Set Name」に「Primer Focus Bin Set」 と入力し、「OK」をクリックします。

New Bin	Set	×				
•	Please enter name of new Bin Set. Primer Focus Bin Set					
	OK Cancel					

Panel Manager 最上部の「**Bin Set**」ドロップダウン リストに、**Primer Focus Bin Set** が追加されます。これで、**Primer Focus Bin Set** を、**Primer Focus Kit** に追加された任意のパネルと関連付けることができます。

3

Primer Focus Kit データの Kit への 追加

- Primer Focus Kit リファレンス データを Primer Focus Kit に追加するには、次の手順を実 行します。
 - 1. 左側のナビゲーションペインで、41ページで作成した Primer Focus Kit を選択します。



2. $\boxed{1}$ $\boxed{5}$ 「Add SNP Reference」ダイアログ ボックスが開き、左下のナビゲーション ペインに、 Primer Focus Kit サンプルを含むプロジェクトがすべて表示されます。



3. SNaPshot Primer Focus Data Project を展開し、Primer Focus フォルダを選択して、 「Add to List」、「Add」の順にクリックします。



SNaPshot Primer Focus Data Project 内の 6 つの Primer Focus Kit サンプル ファイル がすべて、リファレンス データとして Primer Focus Kit に追加され、Panel Manager のナビゲーション ペインの下半分に表示されます。



3

Auto Panel 機能を 使用した Panel、 Marker、Bin の作成

Auto Panel 機能を使用して Panel、Marker、Bin を作成するには、次の手順を実行します。

1. 左側のナビゲーションペインで、41ページで作成した Primer Focus Kit を選択します。

2. 📓 をクリックします (「Bins」 ▶ 「Add Panel」)。

Save Existing	Panels	
Allele Names		٦.
Elue G	- Yellow C	
Green A	Red T	
Marker multiplex		5
Minimum multiplex	5	
Maximum multiplex	20	
Panel Names		-
Base Panel name	Primer Focus Panel	
Marker Overlap		_
Maximum o	verlan	
O No overlap	To the	
Number of b	basepairs between markers 5	

- 3. 「Auto Panel」ダイアログボックスで、次のセクションを編集します。
 - Allele Names デフォルト設定である ddNTP 塩基の名前(G、A、C、T)をそのまま使用します。これらのアレル名は、Primer Focus データから作成した Bin に名前をつける際にも使用します。

注: Marker の名前は、Primer Focus Kit サンプル ファイル名に基づいています。

- Marker Multiplex デフォルト設定である 5 および 20 をそのまま使用します。これらは、Auto Panel 機能によって Panel ごとに作成される Marker の最小数および最大数です。
- Panel Names Primer Focus Panel と入力します。これは、Panel の名前を自動で割り振る場合の基準になる名前です。Marker Multiplex 設定に基づいて、ソフトウェアで複数の Panel を作成する必要がある場合、Panel 名は、Primer Focus Panel_1、Primer Focus Panel_2 などとなります。
- Marker Overlap Maximum Overlap を選択します。Maximum Overlap を選択 すると、同じ Panel 内の Marker をオーバーラップさせることが可能になります。 ただし、同じ色の Bin 同士がオーバーラップしないことが前提になります。
- Save Existing Panels このチェックボックスは選択しないでください。このオ プションを選択すると、あらかじめ存在する Panel に Marker を追加できるように なります。
- 4. Primer Focus Kit データに Auto Panel 機能を適用するには、「OK」をクリックします。

Marker と Bin の Auto Panel 機能で作成された Panel を閲覧するには、次の手順を実行します。

- 閲覧
- 1. ナビゲーション ペインで、Primer Focus Kit を展開します。
- 2. Primer Focus Panel-1 を選択します。



3. 「Plot」タブを選択します。

Panel には、下部 X 軸に表される Marker と、上部 X 軸に表される Bin が表示されま す。下部のナビゲーション ペインでは、Panel の作成に使用された Primer Focus Kit リ ファレンス サンプルの隣に緑のチェックが表示されます。



4. 「Table」タブを選択します。

Marker 情報および Bin 情報が、テーブル形式で表示されます。

🗭 Panel Manager											×
File Edit Bins View											
🚅 🗙 📑 🖀 🖬 🔳 🗄		Bin Set: Primer	Focus Bin Set			- III 4					
😑 🚠 Panel Manager	Ta	ble Plot									
🗄 🚞 Microsatelite Kit		Marker Name	Marker Min	Marker Max	Bin 1 Name	Bin 1 Min	Bin 1 Max	Bin 1 Dye	Bin 2 Name	Bin 2 Min	_
	1	20	22.99	27.50	G	22.99	24.59	Blue	A	24.80	^
SNaPshu ru	2	28	27.74	32.00	G	27.74	29.34	Blue	A	28.92	
Primer Focus Panel-1	3	36	38.77	42.45	G	38.77	40.37	Blue	A	40.42	
🗉 🚞 Microsatelite2 Kit	4	44	46.20	49.47	G	46.20	47.80	Blue	A	47.21	
🗷 🗀 Microsatellite3 Kit	5	52	52.87	55.83	G	52.87	54.47	Blue	A	53.83	
	6	60	60.79	64.05	G	60.79	62.39	Blue	A	62.09	
Reference Samples											
											~
	<										>
				OK Can	cel Appl	y J					

Panel と Bin Set の 新規 Panel と Bin Set を採用して、Panel Manager を閉じるには、「OK」をクリックします。 採用

Bin と Marker の 編集(オプション) このガイドの実験を完了するのに、**Marker** や **Bin** を追加、編集、削除する**必要はありません**。ただし、**Panel Manager** を開き、**Primer Focus Kit** や **Primer Focus Panel-1** を選択して、これらの機能をテストすることはできます。

重要! Marker や Bin を編集または削除した場合は必ず、Panel Manager の下部にある 「Cancel」をクリックしてください。「OK」または「Apply」をクリックすると、解析の結果 に悪影響が出る場合があります。

Marker に Bin を追加するには

- 1. 「Plot」タブ(下部 X 軸)または「Table」タブ(列)で、Marker を選択します。
- 2. \mathbf{i} \mathbf{k} \mathbf{k}
- **3**. 「Edit SNP Marker」ダイアログ ボックスで、Bin の「Name」、「Min」、「Max」を入力 し、「OK」をクリックします。



Bin を編集するには

- **1**. 「**Plot**」タブで、**Bin**(上部 **X** 軸)を選択します。
- 2. をクリックするか(「Bins」 ▶ 「Edit Bin」)、またはビンを右クリックして、「Edit SNP Marker」を選択します。
- **3.**「Edit SNP Marker」ダイアログ ボックスで、Bin の「Name」、「Min」、「Max」を編集 し、「OK」をクリックします。

または

- 1. 「Table」タブを選択します。
- 2. 次の情報を編集します。
 - Bin Name
 - Bin Min
 - Bin Max

Bin を視覚的に編集するには

- **1**. 「**Plot**」タブで、**Bin**(上部 **X** 軸)を選択します。
- 2. 青いセンター ラインをクリックしてドラッグし、Bin の位置を決定します。
- 3. 右ハンドルの左側をクリックしてドラッグし、Bin のオフセット(レンジ)を決定します。





Bin を削除するには

Bin を Marker から削除するには、次の手順を実行します。

• 「Plot」タブまたは「Table」タブで Bin を選択し、 ▲ をクリックします(「Bins」 → 「Delete Bin」)。

または

• 「Plot」タブで、Bin(上部 X 軸)を選択し、Bin を右クリックして、「Delete Bin」を 選択します。



Marker を編集するには

- 1. 「Plot」タブで、Marker (下部 X 軸)を選択します。
- 2. *▲* をクリックするか(「Bins」 → 「Edit SNP Marker」)、または Marker を右クリックして、「Edit SNP Marker」を選択します。
- **3.** 「Edit SNP Marker」ダイアログ ボックスで、Marker Name を編集し、「OK」をクリックします。

または

「Table」タブで、「Marker Name」を編集します。

Panel から Marker を削除するには

Panel から **Marker** を削除するには、「**Plot**」タブで **Marker**(下部 **X** 軸)を選択し、**Marker** を右クリックして、「**Delete Marker**」を選択します。

次の手順 第4章の説明に従って、SNaPshot Kit 解析を設定します。





新規プロジェクトの作成とサンプル ファイルの追加

新規プロジェクトの プロジェクトを新規作成してサンプル ファイルを追加するには、次の手順を実行します。 作成とサンプル 1. 20 をクリックします(「File」▶「New Project」)。 ファイルの追加 1. 20 をクリックします(「File」▶「New Project」)。

🖷 New Project 🛛 🔀
Project Type Generic Microsatellite SNaPshote OLA Analysis SNPIex ⁷⁴ AFLP
OK Cancel

- New Project」ダイアログボックスで、「SNaPshot」を選択し、「OK」をクリックします。
 GeneMapper v3.7 ではこのダイアログボックスは表示されません。そのままステップ3 へ進みます。
- 3. 「GeneMapper」ウィンドウで、 🕵 をクリックします(「File」 ▶ 「Add Samples to Project」)。
- 4. 「Add Samples to Project」ダイアログボックスの「Files」タブで、次のパスを確認します。 *<drive>*:\AppliedBiosystems\GeneMapper\Example Data\SNaPshot

注:上記の場所は、GeneMapper[®] ソフトウェアがインストールされているドライブに よって異なります。デフォルトでは、Dドライブにインストールされます。



5. SNaPshot フォルダを選択し、「Add to List」、「Add」の順にクリックします。

注: このガイドでは、SNaPshot フォルダ内の3つのサンプルファイルをすべて追加しましたが、フォルダ内のファイルの一部(1つまたは2つ)を追加することも可能です。 これを行うには、左側のペインでフォルダを展開し、[Ctrl] キーを押したまま、ファイルを個々に選択して、「Add To List」をクリックします。

Files GM Database	Samples To Add:
 Wy Computer Wy Computer W (A) (A) CD-RW Drive (E) Local Disk (D) Database Defaults Database Defaults Docs Database Defaults Docs Example Data Example Data Example Data M (Croastellite NAPShot NAPShot Stree Standards Stree Standards Config Service 	SNaPshot 2_603_03.fsa 6_F03_01.fsa 6_F03_11.fsa

SNaPshot Data フォルダ内の 3 つのサンプル ファイルが、互換性のある Applied Biosystems 社製シーケンサの Data Collection ソフトウェアに入力した情報と

其に、「Samples」タブ上のテーブルに表示されます。

🖶 GeneMapper - *L	Untitled	- gm l	s Logged In Da	tabase FRMINF	ODEVD01				
File Edit Analysis V	file Edit Analysis View Tools Help								
😂 😂 🛃 🛛 🍒	2		🖪 Ш 🛛 🛛	1 🖩 🗐 🗍	🕨 💣 🕴 Tak	ole Setting: Si	NaPshot Defa	utt 🔽 🕅 📗	۶ 🗟 🔤 🕐
😑 🚠 Project	□ ♣Project Samples Genotypes								
🗄 🦲 SNaPshot		Status	Sample File	Sample Name	Comments	Sample Type	Analysis Me	thod Panel	Size Standard
	1	S	2_803_03.fsa	2	None	Sample	None	None	None
	2	1	5_E03_09.fsa	5	None	Sample	None	None	None
	3	<u> (</u>	6_F03_11.fsa	6	None	Sample	None	None	None
		<)			>
Progress Status								0%	Stop



52ページの説明に従って、プロジェクト用のアナリシスパラメータと表示設定を作成します。



概要

プロジェクト用のアナリシス パラメータと Table Settings の設定

「GeneMapper」ウィンドウ内で、プロジェクト用のアナリシスパラメータと表示設定を確認 します。

アナリシスパラメータには、次の項目が含まれます。

- Analysis Method (Bin Set を含む)
- Panel (Marker の集合)
- Size Standard

プロジェクトにおけるすべてのサンプル ファイルを解析するために GeneMapper[®] ソフト ウェアが使用する、ピーク検出アルゴリズム、サイジング アルゴリズム、ジェノタイピング アルゴリズムを決定するアナリシス パラメータを設定します。

表示設定には、Table Settings と Plot Settings が含まれます。

アナリシス パラ メータの設定

プロジェクト用のアナリシス パラメータを設定するには、次の手順を実行します。

- 1. 「GeneMapper」ウィンドウで「Samples」タブを選択します。
- 2. 「Analysis Method」カラムで最初の列をクリックし、ドロップダウン リストから「New Analysis Method」を選択します。



注:GeneMapper Manager の「**Analysis Method**」タブから **Analysis Method** を新規 作成することもできます。

3. 「New Analysis Method」ダイアログ ボックスで、Analysis Type に「SNaPshot」を 選択し、「OK」をクリックします。

New Analysis Method			
Select analysis type:			
🔘 Microsatellite	🔘 OLA Analysis	🚫 AFLP	
💿 SNaPshot®	O SNPlex™		
	OK Cancel	l	



- 4. 「Analysis Method Editor」ダイアログボックスで、5 つのタブを選択して編集します。
 - General このタブには、Analysis method に関する参照情報が表示されます。
 「Name」に「SNaPshot Analysis Method」と入力します。オプションで、説明
 や、データを作成した装置名を入力できます。

Analysis Method	Editor - SNaPShot®
General Allele Pe	ak Detector Peak Quality Quality Flags
Analysis Method De	escription
Name:	SNaPshot Analysis Method
Description:	
Instrument:	
Analysis Type:	SNaPShot®
	OK Cancel

- Allele このタブには、アレル コーリングを決定する設定が表示されます。Bin Set には、次のうち1つを選択します。
 - SNaPshot Bin Set 第2章の説明に従い、SNaPshot リファレンス データを 使用してこの Bin Set を作成した場合に選択します。
 - Primer Focus Bin Set 第3章の説明に従い、Primer Focus Kit リファレンス データを使用してこの Bin Set を作成した場合に選択します。

その他の設定は、すべてデフォルト値のままにします。

Analysis Method Editor - SNaPShot®	×
General Allele Peak Detector Peak Quality Quality Flags	
Bin Set: SNaPshot Bin Set None None Cut-off va LOH Bin Set SNaPshot Bin Set Primer Focus Bin Set	
Range Filter Factory Defaults	
ОК Саг	ncel



 Peak Detector - このタブには、ピーク検出およびピーク サイジングを決定する 設定が表示されます。「Peak Detection Algorithm」に Basic を選択します。その 他の設定は、すべてデフォルト値のままにします。

Analys	sis Method	Editor - SNaP	Shot®			X
Genera	al Allele Pe	ak Detector Pe	ak Quality 🛛 Qu	ality Flags		
Peak D	Detection Algo	orithm: Basic		~		
Minimu	um Peak Heigh	nt				
	Automs	atic				
	O User s	becified (rfu):				
	Plue	Green	Vellow	Dod	Orange	
	50	50	50	50	50	
					Contant Data Ma	
					Pactory Defaults	
					OK Cancel]

 Peak Quality – このタブには、どんな場合に特定の Process Quality Values (PQV) が緑 ■ (Pass) になるか、もしくは黄色のフラグ ▲ (Check) が示されるかを決定 する設定が表示されます。

「Min peak height ratio」に 0.3 と入力します。その他の設定は、すべてデフォルト値のままにします。

Analysis Method Editor - SNaPShot®	X
General Allele Peak Detector Peak Quality	Quality Flags
Signal level	
Homozygous min peak height	200.0
Heterozygous min peak height	100.0
-Haterozvante halance	
Min peak height ratio	0.3
Peak morphology	
Max peak width (basepairs)	1.5
-Dullum neak	
Pull-up ratio	0.1
Pull-up scan	1
Allele number	
Max expected alleles	2
SNP	
Double peak	0.5
	Factory Defaults
	OK Cancel



- Quality Flags このタブには、次の設定が表示されます。
 - 個別にフラグが示される Process Quality Values (PQV) の Genotype Quality (GQ) 全体に対する重要度を決定する設定。各 PQV は、0~1まで重み付けで きます。0 は最も重要度が低く、1 は最も重要度が高くなります。
 - Sizing Quality (SQ) および GQ に Pass
 Check
 Low Quality
 の 各フラグが示されるしきい値を決定する設定。SQ および GQ の初期スコアは 1 です。何らかのフラグが示された PQV の値を 1 から差し引き、最終的な SQ お よび GQ のスコアが生成されます。

Analysis Method Editor - SNaPSho	nt⊗		
General Allele Peak Detector	Peak Quality Qua	ality Flags	
Quality weights are between 0 ┌Quality Flag Settings) and 1.		
Spectral Pull-up (SPU) Broad Peak (BD) Narrow Bin (NB) Double Peak (DP)	0.5 0.5 0.5 0.5	Control Concordance ((.ow Peak Height (LPH) Off-scale (OS) Peak Height Ratio (PHR)	0.5 0.5 0.5 0.5
PQV Thresholds Pa Sizing Quality: From Genotype Quality: From	ss Range: 0.75 to 1.0 0.75 to 1.0	Low Quality From 0.0 to From 0.0 to <u>Facto</u>	Range: 0.25 0.25
		Ōĸ	<u>Cancel</u>

Analysis Method パラメータの詳細については、GeneMapper[®] Software Online Help を参照してください。

- **5**. 「**OK**」をクリックして **Analysis Method** を保存し、「**Analysis Method Editor**」ダイア ログ ボックスを閉じます。
- 6. 「Panel」カラムの最初の列を選択します。「Select a Panel」ダイアログボックスで、次のうちいずれか1つを実行します。
 - SNaPshot Kit を展開し、SNaPshot 6plex Panel をダブルクリックします(第2章の説明に従い、SNaPshot リファレンス データを使用してこの Panel を作成した場合)。
 - Primer Focus Kit を展開し、Primer Focus Panel-1 をダブルクリックします(第 3章の説明に従い、Primer Focus Kit リファレンス データを使用してこの Panel を 作成した場合)。





- 7. 「Size Standard」カラムで最初の列を選択し、ドロップダウン リストから GS120LIZ を選択します(これは、サンプルでランされたサイズ スタンダードです)。
- 8. 「Samples」タブのすべてのサンプル列に、次のとおり選択をフィル ダウンします。
 - a. 「Analysis Method」、「Panel」、「Size Standard」の各カラム ヘッダにまたがって クリックしてドラッグし、3 つのカラムのすべての列をハイライト表示します。

Analysis Method	Panel	Size Standard
SNaPshot Analysis Method	SNaPshot 6plex Panel	GS120LIZ
None	None	None
None	None	None

b.「Edit」 → 「Fill Down」の順に選択します(または、[Ctrl] キーと [D] キーを同時 に押します)。

Table Settingsの
選択「GeneMapper」ウィンドウの最上部で、「Table Settings」ドロップダウン リストから
SNaPshot Default を選択します。

Microsatellite Default	¥
New	
AFLP Default	
Microsatellite Default	
SNPlex_v3	
SNaPshot Default	

Table Settings は、解析後に「**Samples**」タブおよび「**Genotypes**」タブで表示される情報を 制御します。**SNaPshot Default** は、**GeneMapper** ソフトウェアで提供されるデフォルトの **Table Settings** の **1** つです。

GeneMapper Manager でカスタマイズした Table Settings を編集および作成することもで きます。詳細については、GeneMapper[®] Software Online Help を参照してください。

プロジェクトの保存 プロジェクトを保存するには、次の手順を実行します。

- 1. $\boxed{1}$ $\boxed{5}$ $\boxed{5}$
- **2.**「Save Project」ダイアログ ボックスで、SNaPshot Project と入力し、「OK」をクリックします。

Save Pr	oject	
٩	Project name: SNaPshot Project OK Cancel	

「GeneMapper」ウィンドウのタイトル バーに、SNaPshot Project が表示されます。

次の手順

第5章の説明に従って、SNaPshot プロジェクトのデータを解析および検証します。





プロジェクトの解析

概要

Bin Set を指定する Panel と Analysis Method の両方を含むアナリシス パラメータを備えた プロジェクトを作成後、サンプル ファイルを解析すると、GeneMapper[®] ソフトウェアは、 データのサイジングとジェノタイピングを行います。

「GeneMapper」ウィンドウの「Samples」タブで、次を確認します。

「Status」カラムに 🦉 アイコンが表示されていること。これは、サンプルの解析準備は できているが、「Samples」タブで選択された現在の解析パラメータでまだ解析されてい ないことを示します。

解析 をクリックします(「Analysis」)、「Analyze」)。

GeneMapper ソフトウェアは、プロジェクト内の各サンプルを解析し、「GeneMapper」ウィンドウの左下にあるステータスバーに進捗状況を表示します。

次の手順 下記の説明に従って、結果を検証します。

結果の検証

概要

サイジングおよびジェノタイピング結果の検証は、次の順序で行います。

- SQ、関連する PQV、Size Standard、サンプル情報、サンプル プロットの閲覧(14~ 18ページ、または 36~40ページ)
- GQ および関連する PQV の閲覧(59 ページ)
- 各サンプルに対するアレル コールの閲覧(60ページ)
- Genotype Plot の表示 (60 ページ)
- •「Genotypes Plot」ウィンドウでのデータの検証(61ページ)

GQとPQVの閲覧 データの Genotype Quality (GQ) を閲覧するには、「Genotypes」タブを選択し、テーブル を右へロールします。

GeneMapper - SNaPshot	Project	🔵 - gm ls Lo	gged In Da	tabase FRM	INFO	DEVDO	1								(5
le Edit Analysis View Tools	; Help																
9 🖎 🖩 🍢 🗗 🚺		🖪 Ш 🛛	🛯 🎞 🔳	🔰 🕨 💣	Tab	le Settir	ng:	SNaPs	hot De	fault		*		P	2	AB	1
Panels	Sample	Genotypes															Ī
🗄 📋 SNaPshot 6plex Panel		Peak Area 2	Data Point 1	Data Point 2	ADO	AE	OS	BIN	PHR	LPH	SPU	AN	BD	DP	NB	GQ	
	1	18240.0	2944	2944													
	2	11469.0	2985	3000				NA									
	3	23080.0	3138	3138				NA	NA								
	4	15073.0	3261	3261				NA	NA								
	5	10127.0	3347	3358				NA									
	6	18166.0	3468	3468				NA	NA								
	7	15421.0	2901	2901				NA	NA								
	8	9204.0	2942	2957		<u> </u>		NA									
	9	19070.0	3095	3095		<u> </u>		NA	NA								
	10	12303.0	3214	3214		<u> </u>		NA	NA								
							-								-	_	Î

スクロール バーを右方向へクリックしてドラッグし、「GQ」カラムを表示

手順に従ってこのガイドに示されるデータ例を使用した場合、各サンプルに対する GQ は (Pass) になります。GQ (AN、BD、BIN、DP、LPH、NB、OS、PHR、SPU)の生成に 関連する PQV も になります。

黄色の<u></u>と赤い 🔴 GQ の原因調査

重要! ユーザ独自のデータを解析する場合、GQ が ▲ (Check) または ● (Low Quality) に なり、GQ の算出に関連する PQV (AN、BD、BIN、DP、LPH、NB、OS、PHR、SPU) が ▲ になることがあります。これは、データ、Marker、または Bin の定義、アナリシス パ ラメータに問題があることを示しています。これらの問題の原因を調査して修正するには、 『GeneMapper[®] Software Reference and Troubleshooting Guide』を参照してください。

注: 💣 をクリックすると、GQ スコアに基づいて、サンプルをソートすることができます。 「Genotypes」タブの一番上に、赤い 😑 GQ の付いたサンプルがリスト表示されます。



アレル コールの 閲覧

各サンプル内の各 Marker に対するアレル コールを閲覧するには、「Genotypes」タブを選択して、「Allele 1」カラムと「Allele 2」カラムを表示します。

各	Marker	に対す	るア	レル	$\square - $	ルを	表示
			1				

Sample	es Genotypes						Т		
	Sample File	Sample Name	Panel	Marker	Allele 1	Allele 2	Size 1	Size 2	Height 1
5	2_B03_03.fsa	2	SNaPshot 6plex Panel	54mer	С	Т	53.89	54.63	1343
6	2_B03_03.fsa	2	SNaPshot 6plex Panel	62mer	С	С	62.07	62.07	2202
7	5_E03_09.fsa	5	SNaPshot 6plex Panel	25mer	A	A	25.2	25.2	1554
8	5_E03_09.fsa	5	SNaPshot 6plex Panel	28mer	G	A	28.05	29.11	3220
9	5_E03_09.fsa	5	SNaPshot 6plex Panel	39mer	G	G	39.26	39.26	2259
10	5_E03_09.fsa	5	SNaPshot 6plex Panel	48mer	T	Т	47.91	47.91	1545
11	5_E03_09.fsa	5	SNaPshot 6plex Panel	54mer	С	Т	53.82	54.58	1096
12	5_E03_09.fsa	5	SNaPshot 6plex Panel	62mer	С	С	62.07	62.07	1854
13	6_F03_11.fsa	6	SNaPshot 6plex Panel	25mer	A	A	25.21	25.21	1705
14	6_F03_11.fsa	6	SNaPshot 6plex Panel	28mer	A	A	29.23	29.23	2548
9 10 11 12 13 14	5_E03_09.fsa 5_E03_09.fsa 5_E03_09.fsa 5_E03_09.fsa 6_F03_11.fsa 6_F03_11.fsa	5 5 5 5 6 6	SNaPshot 6plex Panel SNaPshot 6plex Panel SNaPshot 6plex Panel SNaPshot 6plex Panel SNaPshot 6plex Panel SNaPshot 6plex Panel	39mer 48mer 54mer 62mer 25mer 28mer	G T C C A A	G T C A A	39.26 47.91 53.82 62.07 25.21 29.23	39.26 47.91 54.58 62.07 25.21 29.23	2: 1: 1: 1: 1: 2

Genotype Plot の 表示

サンプルの Genotypes Plot を表示するには、次の手順を実行します。

- 1. 「Genotypes」タブでサンプルと Marker (列) を選択します。複数の Marker を選択す るには、[Shift] キーまたは [Ctrl] キーを押したままにします。すべての Marker を選択 するには、「Edit」 → 「Select All」の順に選択します。

「Genotypes Plot」ウィンドウに、選択された各 Marker に対するエレクトロフェログラ ムが表示されます。



3. Plot Setting には SNaPshot Default を選択します。

注: Plot Settings は解析後、「Genotypes Plot」ウィンドウに表示される情報を制御し ます。SNaPshot Default は、GeneMapper ソフトウェアで提供されるデフォルトの Plot Settings の1つです。GeneMapper Manager で Plot Settings をカスタマイズし、編集 および作成も可能です。詳細については、GeneMapper[®] Software Online Help を参照 してください。 4. Genotypes Plot 内の X 軸および Y 軸上で拡大表示する方法は次のとおりです。

目的	操作
X 軸の特定の領域を 拡大表示	目的のプロットを拡大表示するには、上部 X 軸上にカーソルを置き、 🔍 を左右へクリックしてドラッグします。
	または
	上部 X 軸で右クリックし、「Zoom To」を選択して、範囲を入力し、 「OK」をクリックします。
Y 軸の特定の領域を 拡大表示	左側の Y 軸にカーソルを置き、🔍 を上下にクリックしてドラッグします。
	または
	左側の Y 軸を右クリックし、「Zoom To」を選択します。最大値を入力 し、オプションで「Apply to all electropherograms」を選択して、「OK」 をクリックします。
縮小表示	X 軸または Y 軸をダブルクリックします。
	または
	X 軸または Y 軸を右クリックし、「Full View」を選択します。

「Genotypes Plot」 ウィンドウでの データの検証

- 「Genotypes Plot」ウィンドウでは、その他にも次のようなタスクを実行できます。
 - X 軸(塩基対またはデータ ポイント)のスケール調整
 - Y 軸 (個々のサンプルの最大値、すべてのサンプルの最大値、特定の値)のスケール調整
 - 特定の蛍光色素色のピークの表示と非表示
 - 個々のピークに対するステータス ラインの表示
 - アレルコールの追加、名前変更、削除
 - Marker および Bin の編集と削除

上記の機能の使用に関する詳細については、[F1] キーを押し、GeneMapper[®] Software Online Help から該当の項目を選択してください。

Genotypes Plot の表示を確認したら、区をクリックして、ウィンドウを閉じます。



第5章 結果の解析と検証 結果の検証




結果の印刷

「File」→「Print」の順に選択すると、次のウィンドウおよびタブでの結果を印刷することが できます。

ウィンドウ / タブ	「GeneMapper Project」ウィンドウからのアクセス
「GeneMapper」ウィンドウ – 「Samples」タブ	「View」 ▶ 「Samples」
「GeneMapper」ウィンドウ – 「Genotypes」タブ	「View」 ▶ 「Genotypes」
「GeneMapper」ウィンドウ – 「Info」タブ	「View」 ▶ 「Sample Info」
「GeneMapper」ウィンドウ – 「Raw Data」タブ	「View」 ▶ 「Raw Data」
「GeneMapper」ウィンドウ – 「EPT Data」タブ	「View」 ▶ 「EPT Data」
「Samples Plot」ウィンドウ	「Samples」タブで、「Analysis」)「Display Plots」
「Genotypes Plot」ウィンドウ	「Genotypes」タブで、「Analysis」 → 「Display Plots」
「Report Manager」ウィンドウ	「Analysis」 ▶ 「Report Manager」

注: レポートを印刷することもできます。**Report Settings** の設定とレポート作成の詳細については、**GeneMapper[®] Software Online Help** を参照してください。

結果のエクスポート

「Samples」タブお 「GeneMapper」ウィンドウの「Samples」タブおよび「Genotypes」タブに表示された結 果をエクスポートするには、次の手順を実行します。 よび「Genotypes」 タブのエクスポート 1. エクスポートするデータの内容と形式を次のとおり準備します。 a. 「GeneMapper」ウィンドウ最上部のドロップダウン リストから、目的の Table Setting を選択します。Table Setting では、表示するカラムと、サンプルのソート 順序を制御します。 b. 必要に応じてデータをソートし、サンプルを表示する順序を決定します。「Edit」▶ 「Sort」の順に選択するか、「Samples」タブまたは「Genotypes」タブで、[Shift] キーを押しながらカラム ヘッダをクリックします。また、 💕 をクリック (「Analysis」 → 「Low Quality on Top」) すると、GQ スコアに基づいてサンプル をソートすることができます。 注: Table Settings の編集と作成、およびデータのソートに関する詳細については、 GeneMapper[®] Software Online Help を参照してください。 2. 次のコマンドの1つを選択します。 •「File」→「Export Table」-選択したタブに表示された情報をエクスポートします。 • 「File」 ▶ 「Export Combined Table」 – 両方のタブに表示される情報をエクスポート します(このコマンドは、「Samples」タブが選択されている場合のみ使用できます)。 Kit のエクスポート Panel Manager のすべての Kit をエクスポートするには、次の手順を実行します。 1. 🔠 をクリックして Panel Manager を開きます(「Tools」)、「Panel Manager」)。 2. 「File」 → 「Export All Kits」の順に選択します。 Panel の Kit 内のすべての Panel をエクスポートするには、次の手順を実行します。 エクスポート 1. 🔢 をクリックして Panel Manager を開きます(「Tools」) 「Panel Manager」)。 2. 左側のナビゲーションペインで Kit を選択します。 3. 「File」 ▶ 「Export Panels」を選択します。 Bin Set の Bin Set をエクスポートするには、次の手順を実行します。 エクスポート 1. 🔠 をクリックして Panel Manager を開きます(「Tools」)、「Panel Manager」)。 2. ナビゲーションペインで、Bin Set が関連付けられた Kit を選択します。 3. 「Bin Set」ドロップダウン リストで、Bin Set を選択します。 4. 「File」 → 「Export Bin Set」を選択します。



プロジェクト、メ ソッド、Settings、 Size Standard の エクスポート プロジェクト、Analysis Method、Table Settings、Plot Settings、Reports Settings、Size Standard をエクスポートするには、次の手順を実行します。

- 1. **[** をクリックして GeneMapper Manager を開きます(「Tools」) 「GeneMapper Manager」)。
- 2. 次のタブの1つを選択します。
 - Projects
 - Analysis Methods
 - Table Settings
 - Plot Settings
 - Report Settings
 - Size Standards
- 3. エクスポートするオブジェクトを選択します(複数可)。複数のオブジェクトを選択する には、[Shift] キーまたは [Ctrl] キーを押したままにします。
- 4. 「Export」をクリックします。

レポートの レポートをエクスポートすることもできます。Report Settings の設定とレポート作成の詳細 については、GeneMapper[®] Software Online Help を参照してください。

索引

Α

Advanced ピーク検出アルゴリズム 54 Analysis Method Bin Set の選択 53 エクスポート 66 作成 12, 34, 52 保存 55 Analysis Method Editor 「Allele」タブ 53 「General」タブ 53 「Peak Detector」タブ 54 「Peak Quality」タブ 54 「Quality Flags」タブ 55 Analysis Method Editor \mathcal{O} [Allele] \mathcal{P} 53 Analysis Method Editor \mathcal{O} [General] \mathcal{P} 53 Analysis Method Editor の「Peak Detector」タブ 54 Analysis Method Editor の「Peak Quality」タブ 54 Analysis Method Editor \mathcal{O} [Quality Flags] \mathcal{P} 55 **Applied Biosystems** テクニカル サポート vii ユーザマニュアルに関する顧客フィードバック vii 連絡先 vii Auto Panel 45

В

```
Bin

Marker に追加 29, 46

閲覧 28, 45

削除 30, 47

自動作成 45

手動による作成 24

定義 7, 19, 41

編集 29, 47

Bin Set

Analysis Method での選択 53

インポート 19, 41

エクスポート 65

採用 28, 46

作成 21, 42

定義 7, 19, 41
```

F

Fill Down 12, 35, 56

G

Genotype Quality GQ を参照 「Genotypes Plot」ウィンドウ 印刷 64 エクスポート 65 拡大表示 61 データの検証 61 表示 60 「Genotypes」タブ 59 Genotypes テーブル 18,40 GQ 閲覧 59 原因調査 59

K

Kit Primer Focus 3 Primer Focus Kit データへ追加 43 SNaPshot Multiplex 2 エクスポート 65 作成 19,41 種類 20,42 定義 7,19,41 リファレンス データの追加 21

Μ

```
Marker
Bin に追加 29, 46
Panel からの削除 30, 48
閲覧 28, 45
作成 24, 45
定義 2, 7, 19, 41
データ例の 5, 24, 25, 26, 27, 45, 46
プロット、表示 60
プロットの表示 60
編集 30, 48
Min peak height ratio 54
MSDS の入手 vii
```

Ρ

Panel インポート 19,41 エクスポート 65 解析の選択 12,34,55 採用 46 作成 19, 45 定義 7, 19, 41 リファレンス データの追加 23 Panel Manager 開始 19, 41 Plot Setting 17, 39, 60, 66 「Plot」ウィンドウ 「Genotypes Plot」ウィンドウを参照 「Samples Plot」ウィンドウを参照 Primer Focus Kit 3 Primer Focus Kit データ Kit への追加 43 説明 5, 31 Primer Focus サンプルの種類 34

R

Report Settings のエクスポート 66

S

「Samples Plot」 ウィンドウ 印刷 64 エクスポート 65 拡大表示 17,39 データの検証 18,40 表示 17,39 「Samples」タブ 34, 12, 52 Size Match Editor 15, 37 Size Quality (SQ) SQ を参照 Sizing テーブル 18, 40 SNaPshot Kit 2 SNaPshot Kit 解析 Marker 2 互換性のある装置 4 設定 49 定義 2 フローチャート 6 SQ 上書き 16,38 閲覧 14,36 原因調查 15,37 スコア 15,37

Т

Table Settings 12, 13, 34, 35, 52, 56, 66

Х

X軸 拡大表示 17, 39, 61 スケール 18, 40, 61

Y

```
Y軸
拡大表示 17, 39, 61
スケール 18, 40, 61
```

あ

アナリシス パラメータ 設定 12,34,52 定義 7,12,34,52 アレル コールの閲覧 60

い

印刷 「Genotypes Plot」ウィンドウ 64 「Samples Plot」ウィンドウ 64 結果 64 レポート 64 インポート Bin Set 19, 41 Panel 19, 41

え

エクスポート Analysis Method 66 Bin Set 65 「Genotypes Plot」ウィンドウ 65 Kit 65 Panel 65 Plot Settings 66 Report Settings 66 「Samples Plot」ウィンドウ 65 Size Standard 66 Table Settings 66 結果 65 プロジェクト 66 レポート 66

お

オンライン ヘルプへのアクセス vi, vii

か

解析 SNaPshot Kit 解析を参照 解析の設定 49 拡大表示 「Genotypes Plot」ウィンドウ 61 「Samples Plot」ウィンドウ 17, 39 X 軸 17, 39, 61 Y 軸 17, 39, 61 編小表示 17, 39, 61 関連資料 vi

け

蛍光色素ピークの非表示 18,40,61 蛍光色素ピークの表示と非表示 18,40,61 結果 印刷 64 エクスポート 65

2

ゴシック体での表記 v

さ

サイジングのみ 14,36 サイズコーリング曲線 16,38 サイズ スタンダード エクスポート 66 解析の選択 12,34,56 カスタム 16,38 検証 15,37 データ例 5 トラブルシューティング 16 ラベルの削除 16,38 ラベルの追加 16,38 ラベルの編集 16,38 サイズ スタンダードに関する トラブルシューティング 16 削除 Bin 30, 47 Marker 30, 48 サイズスタンダードラベル 16,38 作成 Analysis Method 12, 34, 52 Bin Set 21, 42 Bin (自動) 45 Bin (手動) 24 Kit 19, 41 Marker 24, 45 Panel 19, 45 プロジェクト 10,32 レポート 66 サンプル 拡大表示 17, 39, 61 ソート 15, 37, 59 サンプルの種類、Primer Focus 34 サンプルのソート 15,37,59 サンプル ファイル 拡大表示 17, 39, 61 情報の表示 16,38 ソート 15, 37, 59 場所 5 プロジェクトへの追加 10, 32, 50 プロット、表示 17, 39, 60 プロットの表示 17, 39, 60

サンプル プロット 拡大表示 17, 39, 61 表示 17, 39, 60

L

縮小表示 17, 39, 61

す

ステータスバー 14,36

そ

装置 SNaPshot Kit 解析と互換性 4 データ例 5 ソフトウェア 起動とログイン 7 用語の定義 7

つ

追加 Marker に Bin 29, 46 サイズ スタンダード ラベル 16, 38 プロジェクトへのサンプル ファイル 10, 32, 50

τ

データ
サンプルファイルを参照
データ例を参照
リファレンスデータを参照
データ例
Marker 情報 5, 24, 25, 26, 27, 45, 46
概要 5
サイズスタンダード 5
使用した装置 5
テクニカルサポートの連絡先 vii

と

トレーニングの情報 vii

ひ

ピーク検出アルゴリズム 54 表記法 v ゴシック体 v 重要! v 注意 v 太字 v マニュアル v メニュー コマンドの説明 v ユーザへの注意事項 v

ふ

ファイル サンプル ファイルを参照 太字での表記 v プロジェクト Table Settings の設定 52 アナリシスパラメータの設定 52 エクスポート 66 解析 58 解析(サイジングのみ) 14,36 作成 10,32 サンプルファイルの追加 10,32,50 保存 13, 35, 56 プロジェクトの解析 58 プロジェクトの解析(サイジングのみ) 14,36 プロット サンプル プロットを参照

説明 5,9

れ

```
レポート
印刷 64
エクスポート 66
作成 66
```

$\overline{}$

ヘルプへのアクセス vi, vii 編集 Bin 29, 47 Marker 30, 48 サイズ スタンダード ラベル 16, 38

ほ

保存 Analysis Method 55 プロジェクト 13, 35, 56

ま

マニュアルの使用前提条件 v

ଷ

メソッド Analysis Method を参照 メニュー コマンドの表記法 v

ゆ

ユーザへの注意事項 **v** ユーザ マニュアルに関する顧客フィードバック **vii**

5

ライセンス拒否 <mark>ii</mark> ライセンスの拒否 <mark>ii</mark>

り

リファレンス データ Kit への追加 21 Panel への追加 23