



マイクロサテライト解析

**Getting Started Guide** 



ご使用になる前に

1

2

マイクロサテライト解析の セットアップ

結果の解析と検証 3

4

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、構成物、試薬、キットの権利のいかなる権利(ライセンス)の取得も意味しません。

GeneMapper ソフトウェアは、HID(ヒト個人識別)アプリケーションについて特定の開発バリデーションを受けていません。単一ソースまたは親子サンプルを 解析するためのデータ解析に GeneMapper ソフトウェアを使用する HID(ヒト個人識別)研究施設や機関は、開発検認研究を独自に実行する必要があります。

#### 商標:

ABI PRISM, Applied Biosystems, GeneMapper, GeneScan, LIZ and SNaPshot are registered trademarks and AB(Design), Applera GeneMapper, GeneScan 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. All

rights reserved.

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

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.

# 目 次

	まえがき	v
	このガイドの使い方	v
	詳細情報の入手方法	vi
	サポートの入手方法	vii
第1章	ご使用になる前に	1
	マイクロサテライト解析について	2
	データ例について	3
	マイクロサテライト解析のワークフロー	4
	GeneMapper <sup>®</sup> ソフトウェア の用語	4
	ソフトウェアの起動とログイン	5
	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	5
	GeneMapper ソフトウェア v3.7 を使う場合のこのガイドの使用	5
	このガイドにおける手順のその他の方法	6
第2章	マイクロサテライト解析の設定	7
	概要	8
	Mit. Panel. Markerの作成	9

Kit、Panel、Markerの作成	9
新規プロジェクトの作成とサンプル ファイルの追加	12
プロジェクト用のアナリシス パラメータと Table Settings の設定	14
プロジェクトでのサイジングの実行	20
Auto Bin 機能を使用した Bin Set および Bin の作成	25

第3章	結果の解析と検証	33
	Analysis Method の編集	
	プロジェクトの解析	
	結果の検証	
第4章	結果の印刷とエクスポート	41

卓	結果の印刷とエクスホート	41
	結果の印刷	. 42
	結果のエクスポート	. 43

目 次

# まえがき

## このガイドの使い方

**このマニュアルの** GeneMapper<sup>®</sup> Software Version 4.0 マイクロサテライト解析 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 を使用してマイクロサテライト 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<sup>®</sup> Software Version 4.0 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<sup>®</sup> Winows<sup>®</sup> 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<sup>®</sup> 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<sup>®</sup> 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<sup>®</sup> 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<sup>®</sup> 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、メンテナンスする手順について記載 しています。
    - GeneMapper<sup>®</sup> 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<sup>®</sup> システム、 SNaPshot<sup>®</sup> キット、および SNPlex<sup>™</sup> 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<sup>®</sup> 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 ? をクリックする、という 3 つの方法があります。
    - GeneMapper<sup>®</sup> 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<sup>®</sup> 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 Software 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 からの情報の入手 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 ? 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





## マイクロサテライト解析について

- マイクロサテライト マーカは、別名ショート タンデム リピート (STR) とも呼ばれ、ヌクレ マーカ オチドの反復シーケンスで構成される多型 DNA 遺伝子座です。反復シーケンスの1単位は、 多くの場合2~7塩基対長です。また、反復単位の数はポピュレーション内で異なるため、1 つのマイクロサテライト遺伝子座に対して複数のアレルが作成されます。
- マイクロサテライト 解析 解析 単型的なマイクロサテライト解析では、マイクロサテライト遺伝子座(ローカス)は、蛍光標 識されたフォワードプライマーおよび蛍光標識されていないリバースプライマーを使用して PCRで増幅します。PCR産物は、電気泳動を使用して分子量(サイズ)ごとに分離されます。 次に蛍光検出によって、蛍光標識された産物が検出されます。これで、GeneMapper<sup>®</sup>ソフト ウェアを使用して、アレル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。
- **カスタム プライマー** Applied Biosystems では、マイクロサテライト マーカの PCR 増幅用カスタム プライマーを 提供しています。 詳細については、Applied Biosystems Web サイト (www.appliedbiosystems.co.jp) を参照 してください。
- 互換性のあるキットマイクロサテライト解析と互換性のある Applied Biosystems ケミストリ キットおよびシーケンサの詳細については、『GeneMapper® Software Version 4.0 Quick Reference Guide』<br/>(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** このガイドの演習を実行するには、コンピュータの次の場所に配置されている 3 つのサンプル の場所 ファイル (.fsa)を使用します。

</

**注:**上記の場所は、GeneMapper<sup>®</sup> 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 スタ サンプル ファイルは、PCR 増幅および蛍光標識された マイクロサテライトサンプルを ンダード GeneScan<sup>™</sup> 500 LIZ<sup>®</sup>サイズスタンダードとともにABI PRISM<sup>®</sup> 3100ジェネティックアナライ ザで泳動したものです。

マーカ情報

サンプルファイルには、次の17の蛍光標識されたマーカが含まれます。

マーカ	蛍光色素標識	アレルサイズレンジ (bp) 108 - 130 146 - 172 201 - 233 240 - 285 326 - 354 104 - 138 154 - 176 201 - 245 274 - 298 112 - 125 167 - 195 212 - 234 299 - 339 69 - 99 168 - 196			
D6S264	Blue	108 – 130			
D6S1574	Blue	146 – 172			
D6S276	Blue	201 – 233			
D5S408	D5S408 Blue 240 – 285				
D6S308	Blue	326 – 354			
D6S287	Green	104 – 138			
D6S292 Green		154 – 176			
D6S434 Green		201 – 245			
D5S426 Green		274 – 298			
D5S1981 Yellow		112 – 125			
D6S257 Yellow		167 – 195			
D6S446 Yellow		212 – 234			
D5S641	Yellow	299 – 339			
D5S433 Red		69 – 99			
D5S406	Red	168 – 196			
D5S400	Red	221 – 243			
D6S309 Red		307 – 333			

**注**: 上記の表の情報は、第2章「マイクロサテライト解析の設定」で Panel と Marker を作成する際に使用します。



## マイクロサテライト解析のワークフロー

次のフローチャートは、GeneMapper<sup>®</sup> ソフトウェア を使用してマイクロサテライト解析を実行する手順を要約したものです。

 マイクロサテライト解析の設定 (第2章)
 1. プロジェクト用の Kit、Panel、Marker を作成します。
 2. 新規プロジェクトを作成し、サンプルファイルを追加します。
 3. 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。
 4. サイジングを実行します。
 5. Bin Set を作成してから、Bin を作成します (Auto Bin 機能を使用)。

結果の解析と検証 (第3章)

 Analysis Method を編集して、Bin Set を指定します。
 プロジェクトのサンプルを解析します。
 結果を検証します。

結果の印刷とエクスポート (オプション)
(第4章)
・ 作用ますのいと フキレーナキ

- 結果をプリントアウトします。
- 結果をエクスポートします。

## GeneMapper<sup>®</sup> 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)。プロジェクトのすべてのサンプルの解析用に GeneMapper <sup>®</sup> ソ フトウェア で使用するサイジング 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。			
ビン Bin	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。			
ビン セット Bin Set	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 を作成します。			
マーカ Marker	マイクロサテライト マーカは、名前(ローカス名など)、フラグメント サ イズ レンジ(bp)、蛍光色素色、反復長によって定義されます。			
パネル Panel	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。			
キット Kit	Panel のグループ。			

## ソフトウェアの起動とログイン

#### GeneMapper<sup>®</sup> 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マイクロサテライト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<sup>®</sup> 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 解析する際のマイクロサテライト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<sup>®</sup> ソフトウェア を使用してマイクロサテライト 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。これを利用すればマイクロ サテライト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「マイクロ サテライト解析の設定」のほとんどが、マイクロサテライト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マイクロサテライトプロジェクトを設定する方 法の詳細については、『GeneMapper<sup>®</sup> ソフトウェア Version 4.0 Installation and Administration Guide』(PN 4363080)を参照してください。

コマンド ライン インタフェースを 使用した プロジェクトの設定 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、第 2章 「マ イクロサテライト解析の設定」で説明されている多くのタスクを自動化できるため、マイクロ サテライト プロジェクトを解析する際非常に便利なツールです。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 ソフトウェア の機能を自動化する方法の詳細 については、『GeneMapper<sup>®</sup> ソフトウェア Version 4.0 Installation and Administration Guide』 (PN 4363080) を参照してください。







## 概要

章の構成

この章では、次の方法について説明します。

- Kit、Panel、Marker の作成
- 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 Display Settings の設定
- プロジェクトでのサイジングの実行
- AutoBin 機能を使用した Bin Set および Bin の作成

詳細情報 この章では、基本的な手順を記載しています。GeneMapper ソフトウェアのすべての機能とパ ラメータについての詳細な説明はありません。この章に含まれる項目の詳細については、 GeneMapper<sup>®</sup> Software Online Help の次の項目を参照してください。

- Creating a New Kit (新規 Kit の作成)
- Creating a Custom Panel (カスタムパネルの作成)
- Creating Markers (Marker の作成)
- Creating a Project (プロジェクトの作成)
- Adding Samples (サンプルの追加)
- Applying Analysis Settings (Analysis Settings の適用)
- Starting Analysis (アナリシスの開始)
- Creating a New Bin Set (Bin Set の新規作成)
- Using the Auto Bin Function (Auto Bin 機能の使用)

「Help」メニューからオンライン ヘルプにアクセスするには、 🕜 をクリックするか、[F1] キーを押します。



## Kit、Panel、Marker の作成

概要 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)、蛍光色素色、反復長

**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。たとえば、Applied Biosystems が提供している Linkage Mapping Set キットの一部に対しては、GeneMapper<sup>®</sup> ソ フトウェア で Panel ファイルが利用できます。Panel のインポートに関する詳細については、 GeneMapper<sup>®</sup> Software Online Help を参照してください。

Kit、Panel、 Marker の作成

- Kit、Panel、Marker を作成するには、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」)。
- 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」)。

🖨 Panel Manager							
File Edit Bins View							
	📕 🛄 📕 Bin Set:						
Panel Manager							

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

New Kit	×
123	Missonstallija 1/4
Kit name	Microsaleine Ail
Kit type:	Microsatellite 💌
	OK Cancel





- 4. ナビゲーション ペインで Microsatellite 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 Microsatellite Panel と入力してから、「Enter」キーを押します。

L 1	Panel Name	Comment
1	Microsatellite Panel	none

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

🗑 Panel Manager					
File Edit Bins View					
🔐 🗙 📲 🔳 🔳	Bin Set:				
Anal Manager     Airosatellite Kit     Airosatellite Panel     K					

- 6. ナビゲーション ペインで Microsatellite Panel を選択し、 「File」 ▶ 「New Marker」)。
- 7. Panel Manager の右側ペインで、次の Marker 情報を入力します。

	Marker Name	Dye Color	Min Size	Max Size	Control Alleles	Marker Repeat
1	D6S264	blue	108.0	130.0		2



Marker	Dye Color	Minimum Size (bp)	Maximum Size (bp)	Marker Repeat
D6S1574	Blue	146	172	2
D6S276	Blue	201	233	2
D5S408	Blue	240	285	2
D6S308	Blue	326	354	2
D6S287	Green	104	138	2
D6S292	Green	154	176	2
D6S434	Green	201	245	2
D5S426	Green	274	298	2
D5S1981	Yellow	112	125	2
D6S257	Yellow	167	195	2
D6S446	Yellow	212	234	2
D5S641	Yellow	299	339	2
D5S433	Red	69	99	2
D5S406	Red	168	196	2
D5S400	Red	221	243	2
D6S309	Red	307	333	2

8. 手順 6~7 を繰り返し、次の Marker を作成します。

9. 変更を適用して、Panel Manager を閉じるには、「OK」をクリックします。

次の手順

12 ページ の説明に従って、新規プロジェクトを作成し、サンプル ファイル (.fsa) を追加します。



## 新規プロジェクトの作成とサンプル ファイルの追加

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。

Project の新規作成 とサンプル ファイ ルの追加 Project 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	
<ul> <li>Microsatellite</li> </ul>	
🔘 SNaPshot®	
🔘 OLA Analysis	
O SNPlex™	
O AFLP	

- 2. 「New Project」ダイアログ ボックスで、Microsatellite を選択し、「OK」をクリックします。
- 3. 「GeneMapper」ウィンドウで、 **し** 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

**注:**上記のパスは、GeneMapper<sup>®</sup> 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. Microsatellite 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、Microsatellite 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:
W Computer       Image: Computer (A)         Image: CD-RW Drive (E:)       Image: CD-RW Drive (E:)         Image: CD-RW Drive (E:)       Image: CD-RW Drive (E:)	<ul> <li>Microsatelline</li> <li>Co06_E019_Run_C006_2002-05-10_205_07.fsa</li> <li>C006_E039_Run_C006_2002-05-10_205_03.fsa</li> <li>C006_E109_Run_C006_2002-05-10_205_01.fsa</li> </ul>
Add To List >>  Options	Clear Add Add & Analyze Eance



Microsatellite 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 社製のシーケンサの 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。

\varTheta GeneMapper - *Un	titled -	gm Is	Logged In Data	base FRMINFO	DEVD01				
File Edit Analysis Viev	File Edit Analysis View Tools Help								
😂 🗀 📙 🗓 🚰 🔲 🔛 🖾 📖 🗱 📰 📖 🖉 🌐 👘 👘 🛛 Table Setting: 🛛 Microsatellite Default 🔤 📩 🔎 🌺 🖉 🥥									
😑 🚠 Project	□ 品Project Samples Genotypes								
🗄 🧰 Microsatellite		Status	Sample File	Sample Name	Comments	Sample Type	Analysis Method	Panel	Size Standard
	1	Т,	C006_E019_Run_	E019	None	Sample	None	None	None
	2	<b>N</b>	C006_E039_Run_	E039	None	Sample	None	None	None
	3	1	C006_E109_Run_	E109	None	Sample	None	None	None
								1	>
< · · · >									
Progress Status								0%	Stop

次の手順

14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<sup>®</sup> 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」タブからアナリシス メソッドを新規 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 Microsatellite を選 択し、「OK」をクリックします。

New Analysis Method			
Select analysis type: Microsatellite SNaPshot®	◯ OLA Analysis ◯ SNPlex™	O AFLP	
	OK Cancel		

- 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、アナリシスメソッドに関する参照情報が表示されます。 「Name」に「Microsatellite 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 - Microsatellite 🛛 🗙
General Allele Pe	ak Detector Peak Quality Quality Flags
Analysis Method De	escription
Name:	Microsatellite Analysis Method
Description:	
Instrument:	
Analysis Type:	Microsatellite
	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 None を選択し、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 Edi	tor - Mi	icrosate	ellite			X		
General Allele Peak Detector Peak Quality Quality Flags								
Bin Set: None					~			
Marker Repeat Type								
Use marker-specific stutter ratio if available								
Values for dinucleot	ide repea	nts are ca	alculated autor	natically.				
		Mono	Tri	Tetra	Penta	Hexa		
Cut-off value		0.25	0.2	0.25	0.25	0.25		
PlusA ratio		0.0	0.95	0.95	0.95	0.95		
PlusA distance		0.0	1.6	1.6	1.6	1.6		
Stutter ratio		0.0	0.95	0.15	0.15	0.15		
Stutter distance	From	0.0	0.0	0.0	0.0	0.0		
	To	0.0	3.5	4.5	5.5	6.5		
Range Fitter				Fac	tory Defaults			
	,							
					K Ca	incel		

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	is Method	Editor - Micr	osatellite			×
Genera	I Allele Per	ak Detector Pe	ak Quality Qua	ality Flags		
Peak D	etection Algo	rithm: Basic		*		
_ <sup>Minimu</sup>	m Peak Heigh	t				
	~					
	<ul> <li>Automat</li> </ul>	tic				
	User sp	ecified (rtu):				
	Blue	Green	Yellow	Red	Orange	
	50	50	50	50	50	
					Eactory Defaults	
					Con Cancer	

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。

Homozygous min peak height に 140.0 と入力します。

Heterozygous min peak height に 85.0 と入力します。

Min peak height ratio に 0.4 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 - Microsatellite	
General Allele Peak Detector Peak Quality	Quality Flags
Signal level	
Homozygous min peak height	140.0
Heterozygous min peak height	85.0
- Heterozygote balance	
Min peak height ratio	0.4
Peak morphology	
Max peak width (basepairs)	1.5
⊢Pull-up peak	
Pull-up ratio	0.1
Pull-up scan	1
Allele number	
Max expected alleles	2
Cross-talk peak	
Cross-talk ratio	0.05
	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 -	Microsal	tellite					X
General Allele Peak	Detector	Peak	Quality Qu	ality Flags			
Quality weights are be	tween 0	and 1.					
Quality Flag Settings							
Spectral Pull-up (SPU)	)	0.5		Control Concorda	nce (CC)	0.5	
Broad Peak (BD)		0.5		_ow Peak Height	(LPH)	0.5	
Single Peak Artifact (	SPA)	0.5		Off-scale (OS)		0.5	
Sharp Peak (SHP)		0.5		Peak Height Ratio	(PHR)	0.5	
Cross Talk (XTLK)		0.5		One Basepair Alle	ele (OBA)	0.5	
Out of Bin Allele (BIN)		0.8		Split Peak (SP)		0.5	
PQV Thresholds							
	Pa	ss Ranj	ge:	Low G	uality Ran	ge:	
Sizing Quality:	From	0.75	to 1.0	From 0.	D to 0.2	25	
Genotype Quality:	From	0.75	to 1.0	From 0.0	) to 0.2	25	
					Factory De	efaults	
							_
					<u>o</u> κ	Cancel	

すべての設定をデフォルト値のままにします。

Analysis Method パラメータの詳細については、GeneMapper<sup>®</sup> 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、 Microsatellite Kit を展開し、Microsatellite Panel をダブルクリックします(これは、9 ページ で作成したパネルです)。



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 GS500 (-250)LIZ を選択します。

**注**: GeneMapper ソフトウェアでは、GeneScan<sup>™</sup> 500 LIZ<sup>®</sup> サイズ スタンダードによる サンプル ラン用に次のサイズ スタンダードが使用できます。

- GS500LIZ 実際の GeneScan<sup>™</sup> 500 LIZ<sup>®</sup> サイズ スタンダードに存在するすべて のピークを包含
- GS500(-250)LIZ 250-bp ピークを除外
- GS500LIZ\_3730-35-、250-、340-bp ピークを除外
- GS500LIZ\_3130-35-、250-、340-bp ピークを除外

使用する装置、ポリマー種類、プライマーによっては、特定のピークを除外する他のサ イズスタンダードの1つを選択することが必要な場合もあります。具体的には、35-bp ピークが隣接するプライマー ピークによって隠れてしまう場合、または250-bp および 340-bp ピークがキャピラリシーケンサ上で異常に移動する場合などです。また、独自の カスタム サイズスタンダードを作成することもできます。カスタム サイズスタンダー ドの作成に関する詳細については、GeneMapper<sup>®</sup> Software Online Help を参照してくだ さい。

- 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 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
Microsatellite Analysis Method	Microsatellite Panel	GS500(-250)LIZ
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<br/>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<br/>Microsatellite Default を選択します。

Та	ble Setting:	Microsatellite Default	2
		New	hr
ts	Sample Type	AELD Defeut	þ
	Sample	AFLP Default Microsatellite Default	
	Sample	SNPlex_v3	þ
	Sample	SNaPshot Default	ի

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。Microsatellite 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<sup>®</sup> 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、Microsatellite Project と入力し、「OK」をクリックします。

Save Pr	oject	×
٢	Project name: Microsatellite Project	

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

次の手順 20ページの説明に従って、プロジェクトでのサイジングを実行します。



## プロジェクトでのサイジングの実行

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、データをサイジングします。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、ビン(アレル定義)を作成できるようになります。

**注**: プロジェクトにおけるサンプル ファイル用の Panel が選択されているため、 GeneMapper<sup>®</sup> ソフトウェア は、データをサイジングするだけでなく、データのジェノタイピ ングも行います。しかし、Analysis Method の Bin Set が指定されていないため、Genotype Quality (GQ) は失敗します。

最初の解析を実行するには、次の順序で行います。

- 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  $\triangleright$  をクリックします (「Analysis」  $\bullet$  「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 - Mici	rosatell	ite 4b Project	- gm Is Logg	ged In Database	FRMINFOD	EVD01							-	
File Edit Analysis View	v Tools	Help												
😂 😂 📗   🏂 🖻	P   M	Ш 📰 🛄		🗏 🕨 💕 🛛	Table Setting	: Micro	osatell	ite Defi	ault	~ [		P 8	2	🛯 🕐
😑 🚠 Project	Sample	s Genotypes												
🕀 🧀 Microsatellite		Matrix	Run Name	Instrument Type	Run Date &	Time	REF	SQI	SFNF	MNF	SNF	OS	SQ	UD1
	1		Microsatellite	ABI3100	2002-05-10	22:23:44								
	2		Microsatellite	ABI3100	2002-05-10	22:23:44								
	3		Microsatellite	ABI3100	2002-05-10	22:23:44								
		<											j	>
< >									_					
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。



**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、または解析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、21 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」)。



- 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。

ー方、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、22 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、22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 ん (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 <sup>®</sup> 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 <sup>®</sup> 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<sup>®</sup> 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。

	🛛 GeneMapper - Microsatellite	e 4b Project - gm Is Logged In Database FRMINFODEVD01	
	File Edit Analysis View Tools H	Help	
サンプル・フランル	😂 😂 📕   🍢 🛃   🌆	📰 📴 🛄 🛅 📰 📄 🕨 💕 Table Setting: 🛛 Microsatellite Default 🛛 🖬 🔎 🍇	, 🖪 🕜
リンフル ファイル	■ ♣Project	Info Raw Data EPT Data	
を選択し、	Microsatellite     COOS 5019 Run COOS	Sample Information	<u> </u>
サンブル情報と Raw Data を表示	C006_E039_Run_C006_ C006_E109_Run_C006_ C006_E109_Run_C006_	Sample File : C006_E019_Rum_C006_2002-05-10_205_07.fsa Sample Origin Path : C:\AppliedBiosystems\GeneMapper\Example Data\Microsatellite\C006_E019_Rum_C006_2002-05-10_205_07.fsa Status Message : Good File Source : Disk media Effor Message	
		Nessage : None Current Settings	<b>·</b>
	Analysis Completed.		Stop

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. Image Schule Schul

「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 Microsatellite 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。Microsatellite 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<sup>®</sup> 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 左右へ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。



第2章 マイクロサテライト解析の設定 プロジェクトでのサイジングの実行

「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

24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<sup>®</sup> 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

25 ページの説明に従い、Auto Bin 機能を使用して、Bin Set および Bin を作成します。



図 2-2 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



## Auto Bin 機能を使用した Bin Set および Bin の作成

概要

Panel Manager を使用して、Bin Set および Bin (アレル定義)を作成します。

Bin Set を作成する前に、Kit を選択する必要があります。Kit を選択後、その Kit 内の任意の Panel と Bin Set を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。リファレンス データ として選択した Panel に追加し、Bin 作成に利用できるのは、その Panel を使って解析された プロジェクト内のサンプル ファイルのみです。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: このガイドでは、リファレンス データ、および Auto Bin 機能を使用して Bin を作成す る方法について説明します。ただし、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。また、Bin を手動で作成することもできます。Bin Set のインポートと 手動による Bin 作成の詳細については、GeneMapper<sup>®</sup> Software Online Help を参照してくだ さい。

#### 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. 🔠 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、9ページで作成した Microsatellite Kit を選択します。
- 4. 「New Bin Set」ダイアログ ボックスで、Bin Set Name に Microsatellite Bin Set と入力 し、「OK」をクリックします。

New Bi	n Set	×
٩	Please enter name of new Bin Set. Microsatellite Bin Set	

Panel Manager 最上部の「Bin Set」ドロップダウン リストに、Microsatellite Bin Set が 追加されます。これで、Microsatellite Bin Set を Microsatellite Panel(または Microsatellite 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。



Panel および Bin Set への リファレンス データ の追加 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。

Microsatellite Panel と Microsatellite Bin Set 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。

「Bin Set」ドロップダウンリストで、Microsatellite Bin Set が選択されていることを確認します。

Bin Set:	Microsatellite Bin Set

**2.** 左側のナビゲーション ペインで、Microsatellite Kit を展開し、9 ページ で作成した Microsatellite Panel を選択します。

😁 Panel M	Manager	
File Edit	Bins View	
Щ 🗙	Bin Set:	
⊟ 🚠 Pane ⊟ 🧀 M	el Manager Microsatellite Kit Microsatellite Panel	

作成したすべての Marker が Panel Manager の右側のペインに表示されます (27 ページの図 2-3)。

3. 😼 をクリックします (「Bins」 ) 「Add Reference Data」)。

「Add Microsatellite Reference Data」ダイアログ ボックスが開き、左下のペインに、作 成した Microsatellite Project が表示されます。

注: 左下のペインには常に、選択した Panel を使用して解析されたすべてのプロジェクトが表示されます。

	🗢 Add Microsatellite Reference Data Edit	×
■ をクリック して、12ページ で作成した Microsatellite Project を展開	CM Database Search Criteria All Project Sample Analysis Genotypes SNP SNP Search Results Search Cancel	Samples To Add: → Microsatelite • C006_E019_Run_C006_2002-05-10_205_07.tsa • C006_E039_Run_C006_2002-05-10_205_03.tsa • C006_E109_Run_C006_2002-05-10_205_01.tsa
	Add To Lint >>	Clear Add Eance



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

Microsatellite Project 内のすべてのサンプル ファイルが、リファレンス サンプルとして Microsatellite Panel 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 (27 ページの図 2-3)。

🖶 Panel Manager										
File Edit Bins View										
		Bin Set: Microsate	illite Bin S	Set		~	🖬 🍒			
🖃 🚠 Panel Manager		Marker Name	Dye Color	Min Size	Max Size	Control Alleles	Marker Repeat	Marker Specific	Comments	Ladder Alleles
🗉 🚞 Microsatellite Kit	1	D6S264	blue	108.0	130.0		2	0.0	none	
🗄 🛄 Microsatellite Panel	2	D6S1574	blue	146.0	172.0		2	0.0	none	
	3	D6S276	blue	201.0	233.0		2	0.0	none	
	4	D5S408	blue	240.0	285.0		2	0.0	none	
	5	D6S308	blue	326.0	354.0		2	0.0	none	
	6	D6S287	GREEN	104.0	138.0		2	0.0	none	
	7	D6S292	GREEN	154.0	176.0		2	0.0	none	
	8	D6S434	GREEN	201.0	245.0		2	0.0	none	
A.T.	9	D5S426	GREEN	274.0	298.0		2	0.0	none	
😑 🚠 Reference Samples	10	D5S1981	YELLOW	112.0	125.0		2	0.0	none	
C006_E019_Run_C006_2002-05-	11	D6S257	YELLOW	167.0	195.0		2	0.0	none	
C006_E039_Run_C006_2002-05- C006_E109_Run_C006_2002-05-	12	D6S446	YELLOW	212.0	234.0		2	0.0	none	
	13	D5S641	YELLOW	299.0	339.0		2	0.0	none	
	14	D5S433	RED	69.0	99.0		2	0.0	none	
	15	D5S406	RED	168.0	196.0		2	0.0	none	
	16	D5S400	RED	221.0	243.0		2	0.0	none	
	17	D6S309	RED	307.0	333.0		2	0.0	none	
		<					Ш			>
<										
				ж 🗌	Cancel	Apply				

図 2-3 選択された Panel に対する Marker とリファレンス サンプル ファイルを表示する Panel Manager

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

- 1. ナビゲーション ペインで Microsatellite Panel が選択されており、「Bin Set」ドロップダウン リストで Microsatellite Bin Set が選択されていることを確認したら、 をクリックします (「Bins」 ▶ 「Auto Bin」)。
- 2. 「Auto Bin」ダイアログボックスで、次の手順を実行します。
  - Minimum quality value の設定を 0.1 のままにします。
  - Allele Naming Scheme の Number を選択します。
  - Auto Bin Options の Auto Bin (clear existing bins) を選択します。

Minimum quality value 0.1
Allele Naming Scheme
Number (1, 2, 3, 4,)
O Letter (A, B, C, D,)
Rounded basepair (100, 200, 300)
Auto Bin Options
Auto Bin (clear existing bins)
O Incremental Auto Bin (keep existing bins)

- 3. 「OK」をクリックします。
- 4. 「Autobinning completed」と表示されたら、「OK」をクリックします。
- 5. Panel Manager 下部にある「Apply」ボタンをクリックします。

Auto Bin を使用 した Bin の作成



Marker と Bin の 閲覧

#### リファレンス データから作成された Marker と Bin を閲覧するには、次の手順を実行します。

1. ■ をクリックして Microsatellite Panel を展開し、次に、ナビゲーション ペインの上半分で Marker を選択します。

プロット (図 2-4) に次の項目が表示されます。

- Marker (ピンクの線)
- その Marker の Bin (灰色の垂直のバー)
- 各 Bin のリファレンス アレル(赤い+印)
- 2. ナビゲーション ペインの上半分で Marker を選択し、ナビゲーション ペインの下半分で リファレンス サンプルを選択します。

プロットが更新され(図 2-5)、選択されたサンプルのエレクトロフェログラム ピークを示します。

**注**: Applied Biosystems では、各 Marker を選択して、Marker に対する Bin が作成されているか確認することをお勧めします。Bin が存在しない場合、原因を調査します。詳細については、『GeneMapper<sup>®</sup> Software Reference and Troubleshooting Guide』を参照してください。



Marker (bp レンジ)

図 2-4 Marker の選択



図 2-5 Marker およびリファレンス サンプルの選択

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 追加、編集、削除 (オプション)

このガイドの実験を完了するのに、Bin や Marker 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、Microsatellite Kit や Microsatellite 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!** Bin や Marker 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. 上部のナビゲーション ペインで Marker を選択し、 I をクリックします(「Bins」 ト 「Add Bin」)。
- 2. Bin を追加したい位置で、プロットをクリックします。
- **3.** 「Add Bin」ダイアログ ボックスで、Bin の Name、Location、Offsets を入力し、「OK」 をクリックします。

#### Bin を編集するには

- 1. Bin (灰色の垂直のバー)をクリックして、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 Bin」を選択します。
- **3**. 「Edit Bin」ダイアログ ボックスで、Bin の Name、Location、Offsets を編集し、「OK」 をクリックします。

#### Bin を視覚的に編集するには

- 1. Bin (灰色の垂直のバー)をクリックして、Bin を選択します (図 2-6)。
- 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。



図 2-6 Bin の視覚的編集

#### 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。

- Bin を選択し、 をクリックします (「Bins」 ▶ 「Delete Bin」)。 または
- Bin を選択して、Bin 上で右クリックし、「Delete Bin」を選択します。

#### Marker を編集するには

- 1. 上部のナビゲーション ペインで Marker を選択します (30 ページの図 2-7)。
- 2. 左右のハンドルをクリックしてドラッグし、Marker 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。



図 2-7 Marker の編集



#### Panel から Marker を削除するには

- 1. 上部のナビゲーション ペインで Marker を選択します。

次の手順 第3章の説明に従って、マイクロサテライトプロジェクトのデータを解析および検証します。









## Analysis Method の編集

#### 概要

14 ページ で、Microsatellite Analysis Method を作成したとき、最初の解析では、メソッド内 の Bin Set を選択しませんでした。Bin Set を作成したら次は、Analysis Method 内の Bin Set を選択して、データを解析します。Bin Set を使用すると、GeneMapper<sup>®</sup> ソフトウェア で、プ ロジェクトのサンプル ファイルを解析する際に、アレル コールできるようになります。

Analysis Method の編集

#### Microsatellite Analysis Method を編集するには、次の手順を実行します。

1. 
[i] をクリックして GeneMapper Manager を開きます(「Tools」 ▶ 「GeneMapper Manager」)。

Name         Last Saved         Owner         Instrument         Analysis Type         Description           AFLP Default         2004-07-16 19:21:1         gm         AFLP         AFLP         Factory Provided           Microsatellite Analysis Methy         2004-12-22 09:47:0         gm         Microsatellite         Microsatellite           Microsatellite Default         2003-07-25 21:09:0         gm         3730         Microsatellite           SNPlex_Model_3730         2004-09-21 14:38:4         gm         3730         SNPlex***         Default SNPlex set           SNPlex_Rules_3730         2004-09-21 14:39:0         gm         3730         SNPlex***         Default SNPlex set           SNAPshot Default         2003-07-25 21:10:0         gm         3730         SNPlex***         Default SNPlex set	ojects	Analysis Methods Table	Settings	Plot Settin	gs	Cluster Plot Settin	ngs Matrices	Size	Standards	SNP Sets	Report Settings
AFLP Default       2004-07-16 19:21:1       gm       AFLP       AFLP       Factory Provided         Microsatellite Analysis Meth       2004-12-22 09:47:0       gm       3730       Microsatellite         Microsatellite Default       2003-07-25 21:00:0       gm       3730       Microsatellite         SNPlex_Model_3730       2004-09-21 14:38:4       gm       3730       SNPlex.™       Default SNPlex set         SNPlex_Rules_3730       2004-09-21 14:39:0       gm       3730       SNPlex.™       Default SNPlex set         SNPlex_Rules_3730       2003-07-25 21:10:0       gm       3730       SNPlex.™       Default SNPlex set         SNAPshot Default       2003-07-25 21:10:0       gm       3730       SNaPshot®       SNaPshot®	N	lame	Last Save	ed	Ow	/ner	Instrument		Analysis Ty	pe D	escription
Microsatellite     Analysis Meth     2004-12-22 09:47:0     gm     Microsatellite       Microsatellite Default     2003-07-25 21:09:0     gm     3730     Microsatellite       SNPlex_Model_3730     2004-09-21 14:384     gm     3730     SNPlex™     Default SNPlex set       SNPlex_Rules_3730     2004-09-21 14:39:0     gm     3730     SNPlex™     Default SNPlex set       SNPlex_Rules_3730     2003-07-25 21:10:0     gm     3730     SNPlex™     Default SNPlex set       SNPshot Default     2003-07-25 21:10:0     gm     3730     SNaPshot®	A	\FLP Default	2004-07-	16 19:21:1	gm				AFLP	Fa	actory Provided
Microsatellite Default       2003-07-25 21:09:0       gm       3730       Microsatellite         SNPlex_Model_3730       2004-09-21 14:38:4       gm       3730       SNPlex™       Default SNPlex set         SNPlex_Rules_3730       2004-09-21 14:39:0       gm       3730       SNPlex™       Default SNPlex set         SNPlex_Rules_3730       2004-09-21 14:39:0       gm       3730       SNPlex™       Default SNPlex set         SNaPshot Default       2003-07-25 21:10:0       gm       3730       SNaPshot@	h	licrosatellite Analysis Meth	2004-12-	22 09:47:0	gm	s			Microsatellite		
SNPlex_Model_3730         2004-09-21 14:384         gm         3730         SNPlex™         Default SNPlex set           SNPlex_Rules_3730         2004-09-21 14:39:0         gm         3730         SNPlex™         Default SNPlex set           SNaPshot Default         2003-07-25 21:10:0         gm         3730         SNaPshot@	N	/licrosatellite Default	2003-07-	25 21:09:0	gm		3730		Microsatellite		
SNPlex_Rules_3730         2004-09-21 14:39.0         gm         3730         SNPlex™         Default SNPlex set           SNaPshot Default         2003-07-25 21:10:0         gm         3730         SNaPshot@	s	SNPlex_Model_3730	2004-09-	21 14:38:4	gm		3730		SNPlex™	D	efault SNPlex sett
SNaPshot Default 2003-07-25 21:10:0 gm 3730 SNaPshot®	s	SNPlex_Rules_3730	2004-09-	21 14:39:0	gm		3730		SNPlex™	D	efault SNPlex sett
	-		2003-07-	25 21-10-0	am		3730	-	SNaPshot®		
	2	Sharshol Dendul	2000-01-	20 21.10.0	gin	<u>.</u>					

2. 「Analysis Methods」タブを選択します。

3. Microsatellite Analysis Method を選択し、「Open」をクリックします。



4. Analysis Method Editor で、「Allele」タブを選択し、Bin Set に Microsatellite Bin Set を選択して、「OK」をクリックします。

Analysis Method Editor - M	icrosatel	lite			
General Allele Peak Detector	Peak Qual	ity Quality F	lags		
Bin Set: Microsatellite Bin S	et			*	
Marker Repeat Type					
Use marker-specific stu	tter ratioifa	rvailable			
Values for dinucleotide repe	ats are calc	ulated autom	atically.	<u> </u>	
	Mono	Tri	Tetra	Penta	Hexa
Cut-off value	0.25	0.2	0.25	0.25	0.25
PlusA ratio	0.0	0.95	0.95	0.95	0.95
PlusA distance	0.0	1.6	1.6	1.6	1.6
Stutter ratio	0.0	0.95	0.15	0.15	0.15
Stutter distance From	0.0	0.0	0.0	0.0	0.0
10	0.0	3.5	4.5	5.5	6.5
Range Filter			Fac	tory Defaults	
					_
			0	K Ca	ncel

5. 「Done」をクリックして GeneMapper Manager を閉じます。

**注**:「GeneMapper」ウィンドウの「Samples」タブにある「Analysis Method」カラムで任意 の列をダブルクリックすることによっても、Analysis Method Editor にアクセスできます(14 ページを参照してください)。

次の手順

36ページの説明に従って、プロジェクトを解析します。



## プロジェクトの解析

**概要** Analysis Method で Bin Set を指定後、サンプル ファイルを解析すると、GeneMapper<sup>®</sup> ソフ トウェア は、データのサイジングおよびジェノタイピングを実行します。

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

- 「Status」カラムに 
   で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。このアイコンは、Microsatellite Analysis Method に変更が加えら れたために表示されます。
- 「REF」カラムに ✓ アイコンが表示されていること。これは、そのサンプルが、Bin Set 作成用のリファレンス データとして使用されたことを示します。

**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 (20 ~ 24 ページ)
- GQ および関連する PQV の閲覧(37 ページ)
- 各サンプルに対するアレル コールの閲覧(38ページ)
- Genotype Plot の表示(38 ページ)
- 「Genotypes Plot」ウィンドウでのデータの検証(39ページ)
- プロジェクトアレルの表示(39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 右へスクロールします。

🕏 GeneMapper - Microsatellite	Project	gm Is Logge	ed In Databa	se FRMINFODE	/D01		(	
File Edit Analysis View Tools H	lelp							
😂 🗃 📕   🍒 🛃   🌆 🛙	Ш 🖪 🖩		1   🕨 💣	Table Setting:	Microsatellite Defa	ault 🔽 🗖	🔎 🍉	A 🛛
🖻 🚠 Panels	Samples Gen	otypes						
🗄 🦲 Microsatellite Panel	Data Po	iint 2 ADO AE	OS SHP	OBA SPA SP	BIN PHR LPH	SPU AN BE	XTLK GQ	UD1
1	2 3827							<u>^</u>
1	3 4375							
1	4 5389							
1	5 3028				NA 🔳			
1	6 3558				🔳 NA 🔳			
1	7 4184				NA 🔳			
1	8 4733				NA 🔳			
1	9 5625				NA 🔳			
2	20 4242				NA 🔳			
2	21 3511							
	<					ш		>
Analysis Completed.								Stop

「Genotypes」タブを選択

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、ほとんどのサンプルに対する GQ は (Pass) になります。GQ の算出に関連する Process Quality Values (PQV) (AN、BD、 BIN、LPH、OBA、OS、PHR、SHP、SP、SPA、SPU、XTLK) も () になります。

黄色の 📥 と赤色の 🔿 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 の作成に関連する PQV (AN、BD、BIN、CC、LPH、OBA、OS、PHR、SHP、SP、 SPA、SPU、XTLK) 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<sup>®</sup> Software Reference and Troubleshooting Guide』を参照し てください。



#### **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。

Samp	es Genotypes						1			
	Sample File	Sample Name	Panel	Marker	Dye	Allele 1	Allele 2	Size 1	Size 2	Height 1
12	C006_E019_Run_	E019	Microsatellite	D5S406	R	1	3	173.16	180.92	548
13	C006_E019_Run_	E019	Microsatellite	D5S400	R	1	2	224.96	227.02	403
14	C006_E019_Run_	E019	Microsatellite	D6S309	R	1	2	312.76	319.0	622
15	C006_E019_Run_	E019	Microsatellite	D6S264	в	2	2	113.8	113.8	1680
16	C006_E019_Run_	E019	Microsatellite	D6S1574	в	2	2	158.37	158.37	1481
17	C006_E019_Run_	E019	Microsatellite	D6S276	в	2	2	210.71	210.71	1041
18	C006_E019_Run_	E019	Microsatellite	D5S408	в	3	3	258.33	258.33	910
9	C006_E019_Run_	E019	Microsatellite	D6S308	в	1	1	341.13	341.13	665
20	C006 E019 Run	E019	Microsatellite	D6S446	Y	1	1	215.63	215.63	1804

#### 各 Marker に対するアレル コールを表示

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。

- 「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 Microsatellite 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。Microsatellite 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<sup>®</sup> 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。

「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<sup>®</sup> 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。

#### すべてのプロジェクト 各 Marker に対するサンプル データで検出されたすべてのアレルを表示するには、次の手順 アレルの表示 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.** ナビゲーション ペインで、Microsatellite Kit、Microsatellite Panel の順に展開し、左 側のナビゲーション ペインで Marker を選択します。
- 3. 「Bins」 ▶ 「Show Project Alleles」の順に選択します。

各ビンで、プロジェクト アレル (サンプル データで検出されたアレル) が青いアスタリ スク **\*** で表示されます。**\*** の Y 軸上での各位置は、その Marker とサンプルに対する GQ スコアを示します。





第3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<sup>®</sup> 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 よび「Genotypes」	「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。					
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。					
	<ul> <li>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」</li> <li>▶「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。</li> </ul>					
	<b>注:</b> 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 <sup>®</sup> 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。					
	<ul> <li>「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)。</li> </ul>					
Kitの	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。					
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. 「File」 ▶ 「Export All Kits」の順に選択します。					
Panel Ø	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。					



Project、メソッド、 Settings、 Size Standardの エクスポート Project、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. I 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 コンては、GeneMapper<sup>®</sup> 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 16 Analysis Method エクスポート 44 作成 14 ビンセットの選択 15,35 編集 34 保存 17 Analysis Method Editor 「Allele」タブ 15, 35 「General」タブ 15 「Peak Detector」タブ 16 「Peak Quality」タブ 16 「Quality Flags」タブ 17 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 Bin 27

## В

Bin Auto Bin を使用したビンの作成 27 Marker に追加 29 閲覧 28 削除 30 手動による作成 25 定義 4, 25 編集 29 Bin Set Analysis Method での選択 15, 35 インポート 25 エクスポート 43 採用 29 作成 25 定義 4, 25

## G

Genotype QualityGQ を参照 「Genotypes Plot」ウィンドウ 印刷 42 エクスポート 43 拡大表示 39 データの検証 39 表示 38 「Genotypes」タブ 37 Genotypes テーブル 24 GQ 閲覧 37 原因調査 37

### Η

Heterozygous min peak height16Homozygous min peak height16

### Κ

Kit エクスポート 43 作成 9 種類 9 定義 4,9 マイクロサテライト解析との互換性 2

### Μ

Marker Bin の追加 29 Panel からの削除 31 閲覧 28 作成 9 定義 2, 4, 9 データ例 3, 11 プロットの表示 38 編集 30 Min peak height ratio 16 MSDS の入手 vii

### Ρ

```
Panel
インポート 9
エクスポート 43
解析の選択 18
作成 9
定義 4,9
リファレンス データの追加 26
Panel Manager
開始 9
Plot Setting 23, 38, 44
「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

### R

Report 印刷 42 エクスポート 44 作成 44 Report Settings のエクスポート 44

### S

「Samples Plot」ウィンドウ 印刷 42 エクスポート 43 拡大表示 23 データの検証 24 表示 23 「Samples」タブ 14 Size Match Editor 21 Size Standard エクスポート 44 解析の選択 18 カスタム 22 検証 21 データ例 3 トラブルシューティング 22 ラベルの削除 22 ラベルの追加 22 ラベルの編集 22 Size Quality (SQ) SQ を参照 Sizing テーブル 24 SQ 上書き 22 閲覧 20 原因調查 21 スコア 21

### Т

Table Settings 14, 19, 44

### Х

X 軸 拡大表示 23,39 スケール 24,39

### Y

Y軸 拡大表示 23,39 スケール 24,39

### あ

アレルコールの閲覧 38

アレルのプロジェクト 39

### い

印刷 「Genotypes Plot」ウィンドウ 42 Report 42 「Samples Plot」ウィンドウ 42 結果 42 インポート Bin Set 25 Panel 9

## え

エクスポート Analysis Method 44 Bin Set 43 「Genotypes Plot」ウィンドウ 43 Kit 43 Panel 43 Plot Settings 44 Project 44 Report Settings 44 「Samples Plot」ウィンドウ 43 Size Standard 44 Table Settings 44 結果 43

### お

オンライン ヘルプへのアクセス vi

### か

```
解析
マイクロサテライト解析を参照
解析の設定 7
解析パラメータ
設定 14
定義 4,14
拡大表示 23,39
「Genotypes Plot」ウィンドウ 39
「Samples Plot」ウィンドウ 23
X 軸 23,39
Y 軸 23,39
関連資料 vi
```

### け

蛍光色素ピークの表示と非表示 24,39 結果 印刷 42 エクスポート 43

### 2

ゴシック体での表記 v

## さ

サイズ コーリング曲線 22 削除 Bin 30 Marker 31 サイズ スタンダード ラベル 22 作成 Analysis Method 14 Bin 27 Bin Set 25 Kit 9 Marker 9 Panel 9 Project 12 Report 44 サンプル 拡大表示 23,39 ソート 21,37 サンプル ファイル 拡大表示 23,39 情報の表示 22 ソート 21,37 場所 3 プロジェクトへの追加 12 プロットの表示 23,38 サンプル プロット 拡大表示 23,39 表示 23,38

## L

縮小表示 23,39

### す

ステータスバー 20

## そ

装置
データ例 3
マイクロサテライト解析との互換性 2
ソフトウェア
起動とログイン 5
用語の定義 4

### つ

追加 Marker に Bin 29 サイズ スタンダード ラベル 22 プロジェクトへのサンプル ファイル 12

## 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
Size Standard 3
概要 3
使用した装置 3
マーカ情報 3, 11
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 16 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 ビン 編集 30

### ふ

ファイルサンプル ファイルを参照 フィルダウン 18 太字での表記 v プライマー、カスタム 2 プロジェクト Table Settings の設定 14 エクスポート 44 解析 20,36 解析パラメータの設定 14 作成 12 サンプルファイルの追加 12 保存 19 プロジェクト アレル 表示 39 プロジェクトの解析 20,36 プロットサンプル プロットを参照

### ~

ヘルプへのアクセス vi 編集

```
Analysis Method 34
Bin 29, 30
Marker 30
サイズスタンダードラベル 22
```

#### ほ

```
保存
Analysis Method 17
プロジェクト 19
```

### ま

```
マイクロサテライト解析
互換性のあるキット 2
互換性のある装置 2
設定 7
定義 2
フローチャート 4
マーカ 2
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

ライセンスの拒否 ii

### り

リファレンス データ Panel への追加 26 アイコン 36