1. eArrayへログイン後画面の右上Application Typeが、ダウンロードしたいアレイデザインのアプリ ケーションになっているか確認します。

Home	Microarray	Probe Group	Probes	My Account			Application Type:	Expression	Switch Appli	cation Typ
Search © Microan	ray y Name:	O Probe Group	p	O Simple Prot	pe	Design Wizards © Create a Microarray Design by Uploa © Create a Microarray Design from Exi © Create a Microarray Design from Tar	ading Probes I <u>nfo</u> sting ProbeGroup(s) rget Transcripts I <u>nfo</u>) <u>Info</u>	<u>Refresh</u>	<u>View Al</u>
Species: Design Nu	ımber:			ielect and Add Ipload		Search Resulte: 0 matching results fo	und		Ne	жt >>

2. 異なるアプリケーションの場合は、Switch Application Typeをクリックし、目的のアプリケーションに 変更します。



 Expression、CGH、ChIP、microRNA、TargetEnrichmentの いずれかを選び、Changeをクリック

・Set As Default View Typeにチェックを入れ、Saveをクリックするとロ グイン時のアプリケーションとして選択されます。





3. HomeタブでMicroarrayにチェックを入れ、Design ID欄に<u>デザイン番号の接頭に0をつ</u> <u>けて</u>入力します。

Home	Microarray	Probe Group	Probes	My Functions	My Account		
Search							
Microarra	ay	C Probe G	roup	0	Probe		
Microarray	Name:						
Species:			Sele	ect and Add			
Design ID:	014	693	Upl	oad			
Search Reset							

*Design ID(AMADID)

各アレイのデザインに固有の番号です。デ ザイン番号はマイクロアレイのスライドにつ いている12桁のバーコード番号から調べる 事が出来ます。

バーコード番号

例:Human Genome CGH 244K 25<u>14693</u>XXXXX 25の次の5桁の数字がそのアレイのデザイ ンIDです。

<u>この場合は14693の接頭に0をつけ、</u> <u>Design IDに**014693**と入力</u>



デザインファイルのダウンロード

4. Searchをクリックする。

Design IDあるいはバーコード番号が不明な場合

HomeタブでMicroarrayにチェックを入れ、Microarray Name欄や、Species欄に検索条件を半角英字で入力し、Searchをクリックします(入力欄に何も入力しなくても検索することができます)。

Microarray	Probe Group	Probes	My Functions	My Account				
•	Probe Group	O F	Probe O	HD Probe				
Microarray Name: Human								
		Sele	ct and Add					
		Uplo	iad					
Search	Reset							
	Name: Hurr	Microarray Probe Group C Probe Group Name: Human Search Reset	O Probe Group O F Name: Human Sele Uplo Search Reset	Microarray Probe Group Probes My Functions O Probe Group O Probe O I Name: Human Select and Add Upload Search Reset				

検索結果中、目的のデザインのActions欄内のDownloadをクリックします。

Share Move								
	Microarray Name 🔺	Microarrayset name	Folder Name	<u>Status</u>	<u>Design ID</u>	<u>GPL</u> Identifier	Created Date	Actions
□ Human C 105A	Genome CGH Microarray		AgilentCatalog	Submitted	014698		20-Aug-2007	Copy <u>View</u> <u>Quote</u> <u>Download</u> <u>Attachment</u>
E Human C	Genome CGH Microarray		AgilentCatalog	Submitted	014693		20-Aug-2007	<u>Copy View Quote Download Attachment</u>
E Human C 244A-Su	Genome CGH Microarray upplemental		AgilentCatalog	Submitted	016266		20-Aug-2007	<u>Copy View Quote Download Attachment</u>
Human C 44A	Genome CGH Microarray		AgilentCatalog	Submitted	012750		08-Jul-2005	<u>View Download Attachment</u>
Human C 44B	Genome CGH Microarray		AgilentCatalog	Submitted	013282		08-Jul-2005	View Download Attachment
Human C 4×44K	Genome CGH Microarray		AgilentCatalog	Submitted	014950		20-Aug-2007	<u>Copy View Quote Download Attachment</u>





5.検索結果中、目的のデザインのActions欄内のDownloadをクリックします。







7. 目的のファイルにチェックを入れ"Download"をクリックします。ダウンロードができない場合は、<u>SaveあるいはOpenをクリックするまで</u>Ctrlキーを押しながらダウンロードしてください。

注意:インターネットの設定でPopupBplockを解除しておく必要があります。

Download	If you have difficulty downloading the desired file, hold down the <ctrl> key until a File Dowr appears. This bypasses pop-up blocking software.</ctrl>	load dialog box	
	Category	File Type	
ReadMe		Attachment	
LayoutAgilent		Attachment	
LayoutThirdParty		Attachment	Feature Extractiony田には
ALLANNOTATIONS		All Annotations	GFMI 1.0をダウン
BED		BED	ロードしてください。
EXTERNALFULLGEML		<u>GEML 1.0</u>	
EXTERNALFULLGEML2		<u>GEML 2.0</u>	
FASTA		FASTA	
GAL		GAL	

EXTERNALFULLGEML: Feature Extraction、Agilent Genomic Workbench等アジレントソフト ウェアにインポートできるデザインファイル

GAL: GenePixで使用できる形式のデザインファイル

SEQUENCELIST: プローブ配列のリスト

BED: UCSCのGenome BrowserでカスタムトラックとしてインポートできるBedファイル

TDT:アレイ上のスポット位置情報を含んだ、タブ区切りのテキストファイル



デザインファイルのダウンロード

eArrayからダウンロードされたデザインファイルファイルは2重にZIP化されています 解凍ソフトウェアの種類によってはファイルに障害を生じるため(FeatureExtraction で認識されない)、下記のいずれかで解凍してください

- ・Windows XPの解凍機能(右クリック>Open with(プログラムから開く)> Compressed(zipped)Folders
- Lhaz version 1.33
- WinZip

【Compressed(zipped)Foldersの解凍手順】

①ダウンロードされたzipファイルを右クリック>Open with(プログラムから

開く) >Compressed(zipped)Foldersを選択します

②右のボックスが表示されるのでExternal Full GEML フォルダをダブルクリックします。

③zipファイルが表示されるので ドラッグでデスク トップにzipファイルを移動させます





014850 D ...

💐 014850_1217211910343.zip

④移動したzipファイルを 右クリック>Open with (プログラムから開く)
 >Compressed(zipped)Foldersを選択します

xmlファイルが表示されるので ドラッグでデスクトップにzip





【FeatureExtractionへのImport方法】

①zip形式から解凍したファイルは、下記のようにxmlファイルになります Cドライブの中のProgram File フォルダの中(分かりやすい任意の場所)に入れてください (フォルダ名に全角文字は含まないようにしてください)。

🔮 012097_D_20050419.xml XML ドキュメント

② FeatureExtractionをたちあげてください

③ 下図のように一番上段にあります"Tools">"Grid Template"

>"Add"を選択します。

④で保存したxmlファイルを選択します



Importが完了すると、Grid Template Browserに グリッドが表示されます

