Thermal Cycler Dice Real Time Systemシリーズ 食品環境検査用ソフトウェアQuick Manual -CycleavePCRキットによる定性解析用-

——— 目次 ———

1	・起動と終了					
	(1)Thermal Cycler Dice Real Time System IIIの場合	•	•	•	•	2
	(2)Thermal Cycler Dice Real Time System <i>II / Lite</i> の場合			•	•	4
2.	初期画面と解析タイプの選択		•	•	•	5
3.	実験ファイルの画面構成	•	•	•	•	5
4.	反応条件設定					
	(1) PCR条件の設定	•	•	•	•	6
	(2) ランの開始と進行状況の確認	•	•	•	•	8
5.	サンプル設定					
	(1) サンプル情報の入力	•	•	•	•	9
	(2) 補足	•	•	•	•]	10
	(3) 設定例	•	•	•	•]	10
6.	結果/解析					
	(1) 基本操作	•	•	•	•	11
	(2) データの種類と解析法	•	•	•	•	11
7.	結果の出力	•	•	•	•]	16
8.	トラブルシューティング	•	•	•	•]	18
9.	リアルタイムPCR装置関連製品	•	•	•	•]	19

Thermal Cycler Dice Real Time System $\mathcal{V} \mathcal{V} - \mathcal{X}$

1. 起動と終了

(1) Thermal Cycler Dice Real Time System IIIの場合

■ 起動

① 本体とコンピューターがLANケーブルで接続されていることを確認します。

(注意)本体とコンピューターの接続はLANクロスケーブルをご使用ください。

- ② 本体背面の主電源をONにします。
- ③本体前面の電源ボタンを押します。
- ④ 本体が完全に起動後、LCDにホーム画面が表示されます。
- (補足) 本体の電源を入れるとリッドヒーターが約102℃を超えるまではウォーミングアップの状態となります。 ウォーミングアップ中は、LCDの"スタンバイ"状態表示が点滅します。ウォーミングアップが完了し 使用可能 な状態になると点灯に変わります。



- ⑤ コンピューターの電源をONにします。
- ⑥ デスクトップ上のアイコンをダブルクリックして、ソフトウェアを起動します。



⑦ 画面右下の 機器 と カメラ が"接続"になっていることを確認します。

	\sim	\sim	
リッド:閉温度:107.7℃	機器 接続	カメラ 接続	at
リッド: 温度:	機器:未接続	カメラ : 未接続	

※ 本体とコンピューターが接続されていないときは、機器とカメラに"未接続"と表示されます。

タカラバイオ

■ 終了

- ① ソフトウェアを終了します。
- ② コンピューターをシャットダウンします。
- 本体LCDのホーム画面に表示されている シャットダウンボタンをタップし、本体を待機 状態にします。

④本体背面の主電源をOFFにします。

(2) Thermal Cycler Dice Real Time System II / Liteの場合

∎ 起動

 本体の電源をONにします。
 本体の電源を入れると、ランプやリッドヒーターのウォーミングアップを行います。ウォー ミングアップ中は、"STANDBY"ランプが点滅し、ウォーミングアップが完了すると "STANDBY"ランプが点灯に変わります。



- ② コンピューターの電源をONにします。
- ③ デスクトップ上のアイコンをダブルクリックして、ソフトウェアを起動します。



■ 終了

- ① ソフトウェアを終了します。
- ② コンピューターの電源をOFFにします。
- ③ 本体の電源をOFFにします。

2. 初期画面と解析タイプの選択

■ 初期画面

ソフトウェアを立ち上げると右図のよう な初期画面となります。新しい実験を行う ときは実験ファイルを新規作成し、以前に 行った実験データの解析をするときは既 存の実験ファイルを開きます。

実験ファイルの新規作成

"新規"アイコンをクリック、またはメニュ ーバーから[ファイル]→[新規]を選択。 既存の実験ファイルを開く "開く"アイコンをクリック、またはメニュ ーバーから[ファイル]→[開く]を選択。

第三級連邦軍用ソフトウエア					and the second		HCH (O) = C
2943(7) 第6(1) ユーザー(1) へ62(4) 296 日 谷 市 市 市 中 市 日 日 今 🎙							
M							
	4						
s a 🛤 🚳 🛓		3	27 OK 92798040	1 12 416 1979	■ 主版:107410 総務 図 A 約 世 ● ● ●	aa 0 ⊒ î - ⊷ p	5/9 484 30 1 3011/02/73
🖉 食品環境検	査用ソフ	7トウ:	エア				
ファイル(F)	機器(I)	고-	-ザ-	(U)	へレプ(H)	
🗅 🗃 🔳	<i>6</i> P		ĸ		€ Q	٥	?
新規	開く						

■ 解析タイプの選択

実験ファイルを新規作成すると、"新規測定"ウィ ンドウが表示されます。解析タイプの中から、 +/-判定(CycleavePCR Kit)を選択します。

新規測定		_
解析タイプ 測定者名	+/-判定 (CycleavePCR Ki ▼	

3. 実験ファイルの画面構成

実験ファイルは、3つの画面で構成されており、画面左側のボタンで表示を切り換えます。



タカラバイオ

Thermal Cycler Dice Real Time System シリーズ

4. 反応条件設定

反応条件設定画面では、測定に使用する蛍光フィルターの選択とPCR反応条件の設定を行いま す。ラン開始もこの画面で行い、ラン実行中はランの進行状態を確認できます。

(1) PCR条件の設定

■ 蛍光検出フィルターの選択

+/-判定(CycleavePCR Kit)では、デフォルトで FAMとROXの両方のフィルターが選択されており、 変更の必要はありません。

検出フィルター √ FAM √ ROX	
パターン	Hold
セグメント	1
100-	

■ PCR条件の設定

デフォルトでは、以下の温度条件が表示されます。サイクル数などを必要に応じて変更します。

Hold (初期変性) 95℃、10 sec. 3 Step PCR: 45 cycles 95℃、5 sec. 55℃、10 sec. 72℃、20 sec. (データ取得)

 注: "Speed"は、Thermal Cycler Dice Real Time System III with PC (製品コード TP970)の場合は Normal に、Thermal Cycler Dice Real Time System II (製品コード TP900) やLite (製品コー ド TP700) ではFast に設定します。
 食品環境検査用ソフトウェアをご使用の場合、Speed の設定は各装置のデフォルト条件のままな ので、設定の変更は不要です。

【温度、時間、サイクル数の変更】

変更したい数字をダブルクリックして、数値を入力します。

サイクル数	1		*		
温度(℃)	95.0	95.0	55.0	72.0	
時間(分、秒)	00:10	00:05	00:10	00:20	
データ取得					

タカラバイオ

Thermal Cycler Dice Real Time System シリーズ

■ 既存ファイルからの設定読込み

以前と同じ条件でランを行う場合には、他のファイルから反応条件設定や後述のサンプル設定を 読み込むことができます。画面右上のウェル情報"読み込み"または、"反応条件読み込み"ボタン をクリックすると、ファイルを選択するブラウザが開きますので、目的のファイルを選択して、 "開く"をクリックします。

	後出フィルター IFAM □ROX Speed Fast ♥ Dissociation time 4.0 ♣ sec	表示 ・ Normal 〇 Extend 反応条件読込み
≥	、反応条件を読み込むと、PCR条件の他に蛍光フィルターの選択なども読み	、込まれます。

(検出フィルター)	表示切替	ウェル情報	\frown
インターナルコントロール <none> 💌</none>	● マーク ○ 名称 FAM	しんか 補助 🚺	読込み
			\smile

※ウェル情報を読み込むと、サンプル情報が読み込まれます。

(2) ランの開始と進行状況の確認

■ ランの開始

PCR反応液を分注したチューブ(またはプレート)を装置にセットし、画面右下の"反応開始"ボ タンをクリックします。ファイル名と保存場所を指定し、保存ボタンをクリックします。

パターン · · · · · · · · · · · · · · · · · · ·	反応開始
ランブ:On ランブ使用時間:66時間 リッド:閉 温度:88.9	9℃ 機器:接続 カメラ:接続

■ ラン進行状況の確認

ランを開始すると、画面左側にランの進行状況が表示されます。デフォルトでは、ランの残り時間と温度が表示されますが、"経過"に切り換えると、現在実行中のパターン、セグメント、サイクル数を確認できます。



■ ラン実行中の制御

"装置制御"には、ラン実行中に操作可能な制御ボタンがあります。

- ・ サイクル追加(サイクル数の追加を行います)
- ・ 一時停止 (ランを一時停止します)
- ・ 再開(一時停止中のランを再開します)
- ・ スキップ(実行中のパターンを途中で終了して、次のパターンに移ります)
- ・ ストップ (ランを強制的に終了します)

■ ラン終了後のランプ自動消灯

"自動ランプ Off"をチェックしておくと、ラン終了後にランプが自動的に消灯します。次にラン を行う予定のないときは、ここをチェックしておくとランプ寿命の節約になります。

※ 光源としてハロゲンランプを使用しているTP900シリーズのみの機能です。LEDランプを使用しているTP700/TP970シリーズには、この機能はありません。

タカラバイオ

5. サンプル設定

サンプル設定画面では、サンプル情報の入力を行います。サンプル設定画面での設定は、ラ ン開始前・ラン実行中・ラン終了後のいずれの時点でも行うことができ、設定内容を変更し て再解析することもできます。

(1)サンプル情報の入力

- (インターナルコントロール検出用フィルター)
 +/-判定(CycleavePCR Kit)では、予めROXに設定されており、変更は不要です。
- 画面右上のウェル情報の"入力"ボタンをクリッ クすると"ウェル情報設定"が表示されます。"ウェ ル情報設定"の上の項目から順番に設定していく と入力作業がスムーズに行えます。
- ② サンプルタイプの設定 該当するウェルを選択し、サンプルタイプをプルダウンメニ ューから選択します。
 UNKN (Unknown): 測定対象である未知サンプル PC (Positive Control): ポジティブコントロール
 NC (Negative Control): ネガティブコントロール
- (インターナルコントロールの設定)

+/-判定(CycleavePCR Kit)では、IC(+)に固定されており、設定は不要です。

- ③ レプリケートの設定
 同じ検体サンプルでの反応を複数ウェルで行う場合、同一サンプルを測定するウェルを選択して、レプリケートマーク 1, 2, 3, 4・・・を指定します。
 連続設定機能により、連続入力が可能です。
 - (1) 最初のレプリケートのウェルを選択し、"マー ク"プルダウンメニューから番号を選択。
 - (2) "連続設定"ボタンをクリック。
 - (3) 次のレプリケートのウェルを選択すると、次の 数字が繰り上げ入力される。
 - (4) 設定を解除するには、再度"連続設定"ボタンを クリックする。

タカラバイオ

Thermal Cycler Dice Real Time System シリーズ







ウェル情報設定	×
サンプル名	
サンプルタイプ	UNKN -
インターナルコントロ・	<blank> UNKN</blank>
ターゲット設定	PC NC
マーク <none th="" ▼<=""><th>連続設定</th></none>	連続設定
レブリケート設定	連続設定
🔲 Omit	

④ Omitの設定

反応に使用しないウェルは、Omitにチェックを入れることで、 解析から除外することができます。

ウェル情報設定
サンプル名
サンプルタイプ UNKN ・
インターナルコントロール IC(+) マ
ターゲット設定 📃 複数
マーク <none td="" マ="" 連続設定<=""></none>
レブリケート設定
マーク 1 🥣 連続設定
🔲 Omit

(2) 補足

■ ウェル情報 補助による名称設定

画面右上のウェル情報"補助"ボタンをクリックすると "補助設定"が表示されます。レプリケートのタブをク リックすると、サンプル設定画面で設定したレプリケ ートの名称を入力できます。

補助設	淀				x
g-	ゲット	ノブリケート]		
	<u>–л</u>	之称		 伯	
-	1	A 体			
	2	本 年			
	3	検体3			
	4	検体4			
	5	検体5			
			ОК	キャンセル	

(3) 設定例

9検体サンプル、陰性コントロール、陽性コントロールを各2連で行う反応について、以上の項 目を設定すると以下のようになります。



タカラバイオ

Thermal Cycler Dice Real Time System シリーズ

6. 結果/解析

結果/解析画面では、結果の表示や解析パラメータの設定を行います。画面は、上下二画面に分 かれており、それぞれに任意の図を表示させることができます。

(1) 基本操作

- ① 検出フィルターボタンをクリックするとそのフィルターのグラフが表示されます。
- 2 データ解析のプルダウンメニューからデータの種類を選択します。
- ③ 表示セレクトで表示/解析するウェルを選択します。表示セレクトは、表示セレクトタブの クリックで表示/非表示を切り替えます。

🔟 ファイル(E) 編集(E) セ	かっいの 解析(A) 機器(D ユーザー(U) <u>いって 日</u> (C) <u>の</u> ?	データ解析 増幅曲線	表示セレクト
+ /- 判定 (CycleavePCR Kit)	検出フィルター FAM ROX	クラフ表示 Primary Curve ▼ 解析設定	 ウェル クターゲット / タイプ 1 2 3 4 5 6 7 8 9 10 11 12 A N N U
サンプル設定	70 - 60 -	フィルター FAM ▼ 2回用 ペースライン 瞬値	
反応条件設定 結果/解析	(a) 50 20 20 20 40 E E 20 20 20 20 20 20 20 20 20 20	● Auto 個別設定 ● Manual Start 8 End 45	
①検出フ	ィルターの選択	②データの種類の選択	③表示するウェルの選択

(2) データの種類と解析法

■ 増幅曲線

増幅曲線には、グラフ表示の種類としてPrimary CurveやRawがあります。

Primary Curve: 通常の一次曲線です。

Crossing Point法(CP法)によるCt値の算出に用います。

Raw: 蛍光値の生データです。

バックグラウンドなど、測定結果の確認が必要なときに参照します。

【解析パラメータの設定変更の仕方】

増幅曲線を表示させると、グラフの右側には解析パラメータの設定項目が表示されます。閾値の タブをクリックするとグラフに閾値が表示されます。デフォルトのAuto設定が適切でない場合 は、Manualを選択し、解析パラメータを変更してから"適用"ボタンをクリックしてください。

データ解析	增幅曲線 🔽	データ解析 は	曾幅曲線	
グラフ表示	Primary Curve 🔽	グラフ表示 F	Primary Curve	-
解析設定		解析設定		
フィルター FAI	M 🖌 適用	フィルター FAM	▶ 適用	
ベースライン	閾値	ベースライン 隆	國值	
💿 Auto	Fluorescence Cvcle	 Auto 	個別設定	1
🔘 Manual	7.028 🗘 💈	OManual	Start 8 End 45	
				1

タカラバイオ

Thermal Cycler Dice Real Time System $\checkmark \mathbb{Y} - \vec{x}$

【解析の手順】

① NCの反応結果確認

表示セレクトで《N》のウェルを選択し、検出フィルター《FAM》においてベースラインの 蛍光シグナルに変化がなく閾値を超えないこと、および検出フィルター《ROX》で増幅曲 線が描かれ閾値を超えていることを確認します。《FAM》においてベースラインが安定せ ず、閾値を超えた場合は、トラブルシューティングを参照ください。



② PCの反応結果確認

表示セレクトで《P》のウェルを選択し、検出フィルター《FAM》で増幅曲線が描かれ閾値 を超えていること、および検出フィルター《ROX》で増幅曲線が描かれ閾値を超えている ことを確認します。



タカラバイオ

Thermal Cycler Dice Real Time System $\mathfrak{>} \mathbb{V} - \mathfrak{X}$

③ 検体サンプルの反応結果確認

表示セレクトで反応を行ったウェルをすべて選択し、検出フィルター《FAM》で増幅曲線 やベースラインが、正常に描かれていることを確認します。



■ 測定値分布

データ解析から"測定値分布"を選択し、+/-判定基準の閾値を確認します。

初期設定では45となっています。この設定値は45サイクルのリアルタイムPCRにおいてCt 値が算出されたときに+判定となることを意味します。



■ 判定結果

データ解析から"判定結果"を選択し、判定結果を確認します。

検出フィルターの"総合判定"ボタンが選択された状態では、インターナルコントロールの結果を 踏まえた総合判定結果が表示されます。検出フィルター《FAM》または《ROX》を選択すると、 各々の検出フィルターでの検出結果が+、-で表示されます。

+/-判定	検	出フィルター	F/	AM (R0X		<u>**</u> -	判定										2画面 全画面
(CycleavePCR Kit) サンブル設定 反応条件設定	A	1		2		3		4	5	6 OK OK	7 OK OK	8 Posi. Posi	9	10	11	12	データ解析 判定方法	判定結果 ♥ Cf値(CP) ♥
結果/解析	C D	-								Posi. Posi.	Posi. Posi.	Posi. Posi.					 ・ウェル 1 2 3 	表示セレクト 図 クターゲット / タイプ 4 5 6 7 8 9 10 11 12
	E									Posi.	Posi.	Nega.					AB	
	F									Posi.	Posi.	Nega.					C	
	G									Posi.	Posi.						F	
	н									Posi.	Posi.						GH	

コントロール反応の判定結果

- **OK**: コントロール反応:正常
- **OUT**: コントロール反応:異常

検体サンプルの判定結果

Posi.: 陽性

- Nega.: 陰性(検出限界以下)
- **ND**: 偽陰性、判定不能

(インターナルコントロール、ターゲット遺伝子ともに検出されていない場合) ERROR: エラー(同一レプリケート内で判定結果が異なる場合)

判定結果についての注意

- ・ 陰性コントロール反応で、FAMフィルターにおいて増幅が認められた場合
 → コンタミネーションの疑いがあるため、反応液の調製場所および使用する機器を除染す
 る。
- ・ 陽性コントロール反応で、FAM、ROX両方のフィルターで増幅が認められなかった場合
 → 何らかの原因でPCR反応またはサイクリングプローブ検出が正常に行われていない。
 → サンプル中に反応阻害物質が含まれている可能性もあるので、サンプルを希釈して再反応を行う。または検体の再調製を行い、再反応を行う。
- 陽性コントロール反応で、ROXフィルターでは増幅が認められ、FAMフィルターでは増幅 が認められなかった場合
 - → Primer/Probe Mixに問題があるか、陽性コントロールが分解している可能性がある。

タカラバイオ

(補足) グラフ表示形式の変更

グラフ上でダブルクリックすると、"プロパティ"ウィンドウが表示されます。軸目盛りの変更や Linear、Log スケールの切り換え、ラインやシンボルのデザインの変更ができます。なお、これら のグラフ表示の変更は、右クリックのショートカットでも選択できます。変更内容がひとつだけの 場合には、この方法が便利です。



7. 結果の出力

■ 総合判定結果の出力

総合判定の結果は、数値データあるいはレポートとして出力できます。

【データ出力】

右クリックでショートカットを表示させ、[データ出力]→[CSV]または[Excel]を選択してくだ さい。

【レポート出力】

右クリックでショートカットを表示させ、[レポート出力]→[Word]または[Power Point]を選択して ください。

検	出フィルター 🗌	FAM	ROX 🖁	给判定								
	1	2	3	4	5	6	7	8	9	10	11	12
1	4					ОК	OK	Posi.				
E	3					ок	ОК	Posi.				
						Posi.	Posi.	Posi.	画像出力 データ出力	CSV		
0	<mark>></mark>					Posi.	Posi.	Posi.	レポート出た 印刷	ל וּ Exce		
E	E					Posi.	Posi.	Nega.		_		
F	-					Posi.	Posi.	Nega.				
0	à					Posi.	Posi.					
ŀ	4					Posi.	Posi.					

■ テキストレポートの出力

テキストレポートの内容は、CSV形式またはExcel形式のファイルとして出力できます。 テキストレポートを表示させ、表示項目では"CP法データ"に、詳細項目では必要な項目に☑を 入れます。右クリックでショートカットを表示させ、[データ出力]→[CSV]または[Excel]を選択 してください。

					データ解析	テキストレポート	*
					表示形式	ウェル	~
サンブル 名	サンブルタイプ	レプリケートマ・ 📥	検出フィルター	Ct值(CP)		2746 m //	
滅菌水	NC	1	FAM		表示項目	解析条件	
滅菌水	NC	1	FAM			✓ UP法テータ	
コントロール DNA	PC	- 	L FAM		詳細項目		
コントロール DNA	PC —		- FAM				
サンプル1	UNKN	ウェルの並べ方	FAM				
サンプル1	UNKN	列を元の並びに戻す	FAM			(-)	
サンプル2	UNKN	データ出力	CSV		▼ レプリケート	, ン 、マーカ	
サンブル2	UNKN -	4	Excel		レブリケート	, ·名	
サンプル3	UNKN	5	FAM		▶ 検出フィル	_ &-	
サンブル3	UNKN	5	FAM		Image: Ct(值(CP)		
サンプル4	UNKN	6	FAM		✓ 結果(CP)		
サンプル4	UNKN	6	FAM		✓ 判定(CP)		
サンプルち	UNKN	7	FAM				

※ 他のグラフ等も上記と同様な操作で出力できます。出力したいグラフ等の上で右クリックでショート カットを表示させ、出力形式を選択してください。

タカラバイオ

■ レポート作成機能

いくつかの図をまとめてレポートを作成することもできます。

【レポート作成(Word, Power Point)】

ファイルメニューからフルレポート作成を選 択すると、"フルレポート設定"画面が表示さ れます。必要な図とファイル形式を選択して、 "OK"ボタンをクリックするとレポートが作 成されます。

フルレポート設定	
出力データ選択	
☑ サンブル設定	
📝 反応条件設定	
☑ 結果/解析	
グラフ表示 Primary	Curve *
☑ 増幅曲線	関値 💿 表示 💿 非表示
☑ 測定値分布	関値 💿 表示 💿 非表示
📝 判定結果	🗌 最終蛍光値 📝 Ct值
出力形式 Word	 OK キャンセル

【レポート印刷(PDF ファイル)】

レポートをPDFファイルとして保存することもできます。ファイルメニューから印刷を選択すると、 "フルレポート"画面が表示されます。必要な図とファイル形式を選択して、"OK"ボタンをクリック すると"Print Preview"画面が表示され、ここのFileメニューからsaveを選択するとPDFファイルと して保存されます。

8. トラブルシューティング

- ◆ 陰性コントロール反応においてベースラインが閾値を超えた場合
- 1 データ解析から"増幅曲線"を選択し、グラフ表示から"Raw"を選択する。表示セレクトで ≪N≫のウェルを選択し、検出フィルター≪FAM≫を選択する。
- 2 増幅曲線(Raw)の形状を確認する。
 - 2.1 ベースラインが不安定で閾値を超えたと判断される場合 閾値をManual設定に変更し、閾値をベースラインを超えない位置に設定して"適用"ボ タンを押す。



2.2 PCR増幅によるシグナル増加と判断される場合

コンタミネーションの疑いがあるため、反応液の調製場所および使用機器を除染する。

45000	
40000 -	
2 ³⁵⁰⁰⁰	
변 30000 -	
帮 25000 -	
10 20000	
15000	
	42 44 46
サイクル 数	

9. リアルタイムPCR装置関連製品

■消耗品

製品名	用途	製品コード	容量	価格
0.1 ml 8-strip tube, individual Flat Caps	独立型キャップ付き 8 連チューブ	NJ902	120 strips	¥19,000
0.1 ml 8-strip -neo- tube & cap Set	8 連チューブ	NJ907	120 strips	¥18,500
0.2 ml 8-strip tube, individual Flat Caps	独立型キャップ付き 8 連チューブ	NJ600	120 strips	¥21,000
0.2 ml Hi-8-Tube	8 連チューブ	NJ300	125 strips	¥17,500
0.2 ml Hi-8-Flat Cap	8 連チューブキャップ	NJ302	125 strips	¥4,500
FrameStar® 0.1ml 96 well qPCR plate	96well プレート	NJ904	10 枚	¥9,000
Sealing Film for Real Time	96 well プレート用のシー ル(圧着タイプ)	NJ500	100 枚	¥33,500
Plate Sealing Pads	圧着用パッド	9090	5 個	¥9,500
Sealing Film for Real Time (Adhesive) Ver.2	96 well プレート用のシー ル	NJ502	100 枚	¥27,500
48 well snap plate	48wel プレート	NJ700	20plates	¥9,500
Flat cap for snap plate	48wel プレート用のフラッ トキャップ	NJ720	120 strips	¥7,000

表示価格はすべて税別です。

・最新のライセンス情報に関しては弊社ウェブサイトにてご確認下さい。
 ・本冊子の内容の一部または全部を無断で転載あるいは複製することはご遠慮ください。
 ・本冊子に記載されている会社名および商品名などは、各社の商号、または登録済みもしくは未登録の商標であり、これらは各

所有者に帰属します。

●本冊子記載の価格は2021年4月7日現在の希望小売価格です。価格に消費税は含まれておりません。

タカラバイオ株式会社

テクニカルサポートライン

	TEL. 077-565-6999	FAX. 077-565-6995
首都圈支店	TEL. 03-3271-8553	FAX. 03-3271-7282
関西支店	TEL. 077-565-6969	FAX. 077-565-6995

202104

タカラバイオ