Agilent 1200 シリーズ インスタントパイロットクイックガイド

1. Setup Wizard 画面



インスタントパイロットが接続されている装置の電源を入れると、Setup Wizard が表示されます。 ここではデータと分析時間、装置名、LAN インターフェースの設定/変更が行えます。 Cont.キー、Next キーでそれぞれの設定画面に移動します。最後に Welcome 画面に切り替わります。

2. Welcome 画面



この画面から Method 設定画面等に移動できます。

接続されている全てのモジュールの製品番号/シリアル番号が表示されています。

各モジュールのステータス状態がカラーコードされています。

黄色:Not ready

グレー:Ready

緑:分析中

赤:エラー

Control:分析関連以外の装置に関する機能にアクセス。

下記が Control キーでアクセスできる機能です。

	1 System 2 System 3 System 4 Autosa 5 VWD S 6 DAD SI	n : On / Off n : Get Ready n : Clear Errors mpler : Needle Cleaning L : Balance : Balance
1 System : Set Defau 2 Autosampler : Rele 3 Autosampler : Hom 4 Autosampler : Park 5 Autosampler : Res 6 VWD SL : Take Sca 7 DAD SL : Take Sca	ults ase Vial Gripper Gripper et Sampler in	9

3. Method を設定します。

Welcome 画面で Method キーを押します。

	Me	thod - UNNAM	ED	
				-
Setting	Valu	10		
	S	ystem		A Edit
Stoptime	20.0	0 min		
Posttime	OFF			
Quate	ernary Pu	ump : DE239231	124	Control
Stoptime	20.0	0 min		Control
Posttime	OFF			
Flow	0.00	0 ml/min		Tarrela
%B,%C,%D	OFF	, OFF , OFF		Toggle
Min. Pressure	OFF			
Max. Pressure	400	bar		
Minimum Stroke	AUT	0		
Compressibility	100	* 10E-6/bar		
Max. Flow Gradien	t 100.	0 ml/min²		
Primary Channel	Auto			🗉 Exit
Specifies a time limit !	or your a	naluele		13:45
specifies a ciffe mine	or your a	rialysis.		1 10,40
Filter Co	mpare	Timetable	Properties	File

各モジュールのメソッドを設定します。スクロールすると接続しているモジュールの設定項目が表示されますので、 分析条件を設定してください。Exit キーで Welcome 画面に戻ります。

Welcome 画面の Control キーから System On を選択します。システムが On になり、ポンプ送液が開始され、検出器のランプが点灯します。



4. 分析を開始します。

	Welcome			00	1	○ 🛉 🖲
				Control	\square	€.
Quat	Pump	G1311	A - DE23923124			1
Autosa	mpler	G1329	A - DE91603245	Details	\square	
Col	Comp) G1316	A - DE14923865		\bigcirc	789
	DAD	G1315	B - DE03010634			(4) (5) (8
12/	13 Val.	G1160	A - DE14900157		\square	1 2 3
7	ディス	プレイ		89	\bigcirc	• • •
Select a buttor	to continue.			09:58		
Method	Sequence	Status	Logbook	More 🛆		Start Sto

キーボードの Start ボタンを押します。下記の画面が表示されます。



Sample Range に分析するサンプルのオートサンプラ ローケーションを入力します。

From(どこから)to(どこまで)で設定します。サンプルが1検体の場合は、FormとToに同じロケーションを入力します。

#Injに注入回数/サンプルを入力します。Cont.キーで分析が開始します。

Status を確認し、クロマトグラムを表示します。
Welcome 画面で、Status キーを押します。



Plot キーを押します。



クロマトグラムが表示されます。

Select:表示するシグナルを設定/変更ができます。

Rescale:表示するクロマトグラムのスケールが変えられます。