Agilent 7890 GC 用 診断・モニタリングソフト Lab Advisor の操作方法

1. ソフトの起動

 ①はじめに、各クライアント PC から右の アイコンをクリックします。



②Lab Advisor ソフトが立ち上がり、以下の画面が出ます。

※立ち上げ時は、機器と通信していないため(モニタされていません)と表示されます。

Agilent Lab Advisor	۵	My Agilent Lab Advisor	ユーザー:	AGILENT\hianazaw [管	理者]	
 		 	選択されている機器: 109090621 (モニタされて)	() 名称未設定Agill () 名称未設定Agill () 名称未設定Agill	ent 7890 [CN10909062] 1個の機器のづち1個を表	·示
 後篇 ジステム情報	*	Agilent 7890 / CN10909062 Hidetaka Anazawa				
 ヘルブ ② 全体表示 ③ 設定 ③ ソフトウェアについて 	۲					
タスク ② 全体表示をカスタマイズする ③ 機器の追加または変更 ④ 機器のモニタの開始	۲					
使利なリンク	\$					

③機器表示エリアにポインタを移動させ右クリックするとメニューが出ます。

メニューから「モニタ開始」をクリックします。



④機器の通信が完了すると以下のような画面になります。

Agilent Lab Advisor Agilent	۲	Agilent Lab Advisor バージョン B.1.03.10053.0521 全体表示	ユーザー: 選択されている機器:	AGILENT\hianazaw [管 ³ []]: 名称末設定Agile	理者] nt 7890 [CN10909062] 📲
 マテームウェア更新 計算ツール 	•	 	09062]		1個の機器のうち1個を表示
 ● システム情報 ○ システム情報 ○ カレンダー △ クロマトグラフの詳細 ○ テスト ① テスト ① EMF ○ ステータスレポート ○ ログと結果 	×	Agilent 7890 / CN10909062			ステータス: アイドル 準備状況: ノットレディ
 ヘルブ ② 全体表示 ③ 設定 ③ ソフトウェアについて 	۲				
タスク ② 全体表示をカスタマイズする ③ 機器の追加または変更 ④ 機器のモニタの開始	۲				
便利なリンク	*				

2. メンテナンス予告機能(EMF)の操作

モジュールの各稼働状況を確認できます。「警告」「限界」のリミット値設定により、メンテナンス時期をアナウンスする機能です。

①「EMF 機能」の起動

画面右のメニューから「EMF」をクリックすると、以下の画面になります。

Agilent Lab Advisor	🍌 Agilent	Lab Advisor			ユーザー:		AGILENT\hianazaw [管理	里者]		
🏦 全体表示	🧥 バージョン 🛚	.1.03.10053.0521			選択されて	いる機器:	() 名称未設定Agile	nt 7890 [CN10	909062]	~
🔒 コンフィグレーション	EMF									
 ドキュメント パー・・・・・・・・・・・・・・・・・・・・・・・・・・・・・・・・・・・・	多称未設定∆cilent	7890 [CN10909062]								
ファーム・フェア 足利 計算ツール	リリソース		単位	値	警告	限界	経過			
	機器の分析実行	行力ウンタ	注入	0					0%	Π
	カラム1の長さ		m	30					0%	
	カラム1注入サイ	つん	注入	0	1000	1000			0%	=
	70ンド注入口:	ライナー	注入	0	500	500			0%	
// ジロマドジジンの#単和曲 「34 テフト	70ンド注入口:	ライナーの使用時間	時間	00:09:48	1.00:00:00	1.00:00:00			1%	
	70ンド注入口・	セプタム	注入	0	100	100			0%	
	70ント注入口:	カラムアダプタ/シルバーシール	注入	0	100	100			0%	
 ログと結果 	70ンド注入口:	スプリットベントトラップ 	注入	0	1000	1000			0%	
	バック注入ロライ		注入	0	500	500			0%	-
∧ <i>µ</i> ,7	バック/主人ロライ	(ナーの使用時間	時間	00:09:47	1.00:00:00	1.00:00:00			1%	\mathbf{v}
O EMF										
emfのしくみ						□ ► レ =	えきれていないインジケータ	ヌを非表示	削除	
😢 注意動作	詳細 新しいEMF	インジケータの追加								
タスク 🗇	モジュール	GC7890.CN10909062								
🕜 リソースの限界値を設定、変更、または消	名前	機器の分析実行カウンタ						カウン	トの停止	
② UV−2 値のUセット	≣兑 ^日 月									
	値	0 ^{(±} λ						URY	トのクリア	
指示切新規追加	10			28 88			24-7	//`tht	なのリセット	
🕐 注意 Eメールの送信				P氏环			i±Λ			
😢 注意テキストメッセージの送信	注意の通知	□ ノットレディに設定 □ 9	Service Due	オン	メール	□ テキスト;	いセージ	変更	を保存	
便利なリンク 😵										

②画面の紹介

単位	値	警告	限界	径過	
m	30				0%
注入	0	1000	1000		0%
注入	0	500	500		0%
時間	00:11:18	1.00:00:00	1.00:00:00		1%
注入	0	100	100		0%
注入	0	100	100		0%
注入	0	1000	1000		0%
注入	0	500	500		0%
時間	00:11:17	1.00:00:00	1.00:00:00		1%
注入	0	1000	1000		0%
時間	00:11:16	1.00:00:00	1.00:00:00		1%
注入	0	500	500		0%
時間	00:11:16	1.00:00:00	1.00:00:00		1%
注入	0	1000	1000		0%
注入	0	100	100		0%
	単位 注入 m 注入 時間 注入	単位 値 注入 0 m 30 注入 0 注入 0 時間 00:11:18 注入 0 注入 0 注入 0 時間 00:11:17 注入 0 時間 00:11:17 注入 0 時間 00:11:16 注入 0 時間 00:11:16 注入 0 時間 00:11:16 注入 0	単位 値 警告 注入 0 第二、 1000 注入 0 1000 注入 0 500 時間 00:11:18 1.00:00:00 注入 0 100 注入 0 100 注入 0 100 注入 0 1000 注入 0 1000 注入 0 500 時間 00:11:17 1.00:00:00 注入 0 500 時間 00:11:16 1.00:00:00 注入 0 500 時間 00:11:16 1.00:00:00 注入 0 500 時間 00:11:16 1.00:00:00 注入 0 1000 注入 0 1000 注入 0 1000	単位値<警告限界注入0m30注入0注入0注入0500500時間00:11:181.00:00:00注入0100100注入0100100注入0100100注入0100100注入0100100注入0時間00:11:171.00:00:001.00:00:00注入0500500時間00:11:161.00:00:00500注入0注入0注入01001000注入01001000	単位 値 警告 限界 経過 注入 0

稼働状況が数値表示されます。 現在の値とリミット値(「警告」と「限界」) の値を表示します。

稼働状況がグラフで表示されます。

予め設定された「警告」「限界」のリミットに対して、 警告リミット以下 =緑表示 警告リミットを超える=黄色表示

限界リミットを超える=赤表示となります。

③各種設定変更

設定変更したいリソースをクリックします。

リソース	単位	値	警告	限界	経過	
機器の分析実行カウンタ	注入	0				0%
カラム1の長さ	m	30				0%
カラム1注入サイクル	注入	0	1000	1000		0%
フロント注入ロライナー	注入	0	500	500		0%
フロント注入ロライナーの使用時間	時間	00:15:48	1.00:00:00	1.00:00:00		1%
フロンド注入ロセプタム	注入					0%
フロント注入ロカラムアダプタ/シルバーシール	注入	0	100	100		0%
フロント注入ロスプリットベントトラップ	注入	0	1000	1000		0%
バック注入ロライナー	注入	0	500	500		0%
バック注入ロライナーの使用時間	時間	00:15:47	1.00:00:00	1.00:00:00		1%
バック注入ロゴールドシール	注入	0	1000	1000		0%
バック注入ロゴールドシールの使用時間	時間	00:15:46	1.00:00:00	1.00:00:00		1%
バック注入口ロリング注入回数	注入	0	500	500		0%
バック注入口ロリングの使用時間	時間	00:15:46	1.00:00:00	1.00:00:00		1%
バック注入ロスプリットベントトラップ	注入	0	1000	1000		0%
バック注入ロセプタム	注入	0	100	100		0%

画面下が選択したリソースに切り替わるので、このエリアで設定変更します。

「詳細 新しいEM	Fインジケータの追加	
モジュール	GC7890.CN10909062	
名前	フロンド注入口セプタム	カウントの停止
記日月		
値	0 注入 B	リミットのクリア
警告	100 注入 限界 100 注入	A カウンタのリセット
注意の通知	□ ノットレディに設定 Service Due オン □ メール □ テキストメッセージ	変更を保存

A. カウンタのリセット

消耗部品交換などメンテナンスを実行した後、カウンタをゼロにしたい場合は「カウンタのリセット」 ボタンを押します。

B. リミット値の変更

リミット値を変更したい場合は「警告」「限界」に数値を入力し、「変更を保存」ボタンを押します。

3. テスト機能の操作

各モジュールの診断テストを実施し、良否の判定ができる機能です。

①テスト画面の起動

•

画面右のメニューから「テスト」をクリックすると以下のようなテスト画面になります。

「テスト」をクリック

Agilent Instrument Utilities ② 計 機器の表示 ③ コンフィグレーション ③ コンフィグレーション ③ ドキュメント ③ ファームウェア更新	Agilent Instrument Utilities バージョン B.1.03.10012.0334 テスト モジュール表示
 計算ツール 機器 ※ 	注入ロ:スプリット/スプリットレス 位置:バック
 ミステム情報 テスト ステータスレポート 	検出器: FID 位置: フロント

②テスト実行例

②-1:テストしたいモジュール表示(この場合「注入口」)をクリックします
 ②-2:テストしたい項目(この場合「リーク検査」)をクリックします
 ②-3:「テストを今すぐ実行」ボタンを押します



②-4:画面の指示に従い、操作を実行します。

テスト名 概算時間 ステータ	司 ス	バックのスプリット/スプリットレス注入口 7890 リ・ 5 分 分析中	説明	- 177	スプリット/スプリットレスリーク検査を行うと、メンテナンスを行った 多、または、グロスリークの疑いがある場合に、注入口の漏れを素 早く検査できます。結果はすぐ出ます。分解は不要です。信頼 生は圧減衰テストほど高くありません。
「 「テスト手ル」	ļā				
1. 2. 3. 4. 5. 6.	GCコンフ 選択した 覚択した プレランの システム(オリジナル	ィグレーションを検査します。 G C のコントロールを確立します。 G C (Jリーク検査メソッドをダウンロードします。 ステータスを入力します。 tニューマティック条件を確立しようとします。 、メソッドを元に戻し、G C から切断します。		אידעב א	びレーションの確認
カテゴリ	ソース		時間		メッセージ
	バックのス: バックのス:	グリット/スプリットレス注入口 7890 リーク検査 グリット/スプリットレス注入口 7890 リーク検査	2010/03/	28 12:36:26 28 12:36:26	バックのスプリット/スプリットレス注入口 7890 リーク検査 が Administrator (より 名称未設定Agilent 7890 [CN10924073] 開始されました GCのコンフィグレーションを調べ、選択したテストの要件を満たして いることを確認します。

②-6:メニューから「結果の印刷」をクリックすると、レポート出力できます。

診断結果	۲
 ■ テストの停止 ● 結果の印刷 	
ヘルプ	۲
🥑 FIDジェット制限テストの実行	
テスト結果を改善するには	۲
🥑 FIDジェットを変更するには	
🕑 FIDジェットを選択するには	

③実施可能なテスト項目

7890A GC 検出器テスト

●検出器チェックアウト

検出器およびその他の GC 部品が古くなると、検出器の性能に影響を及ぼすことがあります。Lab Advisor は、GC と検出器の性能が工場出荷時の状態に保たれていることを確認するために、FID、TCD、および uECD の検出器チェックアウトに関連する自動作業を実行します。

●ジェット制限テスト

FID の点火に関連した問題の最も一般的な原因は、ジェットが完全に詰まっているか、部分的に詰まって いることです。 ジェットが完全に詰まっていてもフレームが点火する場合は、ピークリテンションタイ ムが長くなるという二次的な症状が現れます。 ジェットの詰まりは、膜厚カラム、高ブリードカラム、 またはパックドカラムを高温の用途で使用した場合に起こりやすくなります。

FID ジェット制限テストでは、水素流量を増やすと同時にメークアップガス流量を遮断することによって、 完全または部分的に詰まったジェットがないかどうかがチェックされます。

7890A GC 注入ロテスト

●注入口の圧減衰テスト

注入口本体およびフローモジュールベントを通るGCフローモジュールからの注入ロシステムに漏れがな いかを調べるために圧減衰テストが実行されます。 注入ロカラムフィッティングや検出器は、このテス トの対象外です。

各圧減衰テストでは、十分なガスの供給があること、およびテストを実行する前に注入口が冷却されて いることを前提にしています。 テストの実行には約30分かかります。

●注入口スプリットベント制限テスト

注入ロスプリットベントは時間の経過と共に制限され、圧力設定値に達しないエラーや併行精度に関連 した問題を招くことがあります。 また、ライナーのタイプによっては背圧が生じ、低流量/低圧メソッ ドがいつまでも設定値に達しなくなる場合もあります。 スプリットベント制限検査を行うと、注入ロス プリットベントが制限されていないかどうか、およびライナーが現在の用途に対応しているかを判定で きます。 ●注入口プレランリーク検査

プレランリーク検査は、フィッティング締め付け時に漏れがないかどうかを確認したり、メンテナンス 実行時の事前/事後テストとして実行します。

※ヘルプの「テスト」より、手順、各注入口ごとの詳しい説明を閲覧することができます。



4. システム情報の確認

「システム情報」をクリックすることによりモニターしている GC の詳細情報(注入口の種類や、カラム、GC のコントロールパネルなど)を確認することができます。

Agilent Lab Advisor 🏾 💲	Agilent Lab Advisor	ユーザー:	AGILENT\hianazaw [管理者]
🏦 全体表示	レンジョン B.1.03.10053.0521	選択されている機器:	[] 名称未設定Agilent 7890 [CN10909062]
 □ コンフィグレーション ◇ ドキュメント ◇ アームウェア更新 □ 計算ツール 	システム情報 GC7890 様況の詳細	機器の)表示:
機器			
 システム情報 ゴカレンダー ヘ、クロマトグラフの詳細 マスト 	名称未設定Agilent 7890 [CN10909062]		COLUMN 1 (He) Dimensions 60.0 m 250 u Pressure 17.686 17.686< Flow 1.200
 MF ステータスレポート ログと結果 	モデル: Agilent 7890 シリアル: CN10909062 ファームウェアバ ネットワークアドレス: 146.208.29.129 ス	担当者: <u>Hidetaka Anazawa</u> ベージョン: A.01.10.3 、テータス: アイドル - 準備状i	<u>。</u> 兄: ノットレディ
ヘルブ ※	説明: Agilent 7890 [CN10909062] が 146.20	8.29.129 で検出されました	
🕖 システム情報			
タスク 余	コンフィグレーション メッセージ		
 ジッセージの解決 (ジ) 根器コンフィグレーションを表示する 	またつ: PTV 位置: 7ロント 注入ロライナー: 不明		
 	注入ロ:スプリット/スプリットレス 位置:パック		=
② 全体表示	1 1 注入ロライナー: 不明		
 砂 機器コンフィグレーション 砂 EMFインジケータ 	カラム 1: J&W 122-5562 長さ(現在): 60.0 m 内径: 250µm 長さ(オリジナル): 0.0 m 膜厚: 0.25µm	注入口接続: フロント注入 出口接続: 真空	