Agilent University ePass

ユーザーガイド





Contents

1.	Agilent University ePass	3
2.	<u>ePass利用手続き</u>	4
3.	Agilent University アカウントにログインする	8
4.	<u>コース情報を見る・登録する</u>	9
5.	<u>問い合わせ先</u>	15



Agilent University ePass

Agilent University ePass は、24時間365日すべての Agilent University e-Learningコンテンツに何度でも アクセスいただけます。期間は3か月間、または 12か月間から選択いただけます。

Agilent University ePassをご利用いただくと、ブラウ ザ上でのソフトウェアやデータ解析を実践できるクラ ウドラボコースをはじめ、3か月間または12か月間の 希望する期間中、プロダクトや言語に関わらず全ての オンラインコンテンツを視聴可能です。

(600時間以上に及ぶ、500コース以上のコンテンツ、 日本語版コンテンツは一部のみに限定されます)

同じコースを何度でも繰り返しご覧いただけます。



Information

Product Number	Description	Price
R5002A	3か月間 ePass	600 トレーニングクレジット
R5003A	12か月間 ePass	1200 トレーニングクレジット

ePassを開始する前に

日本国内で、初めてePassサービスをご利用いただく場合(まだePassを購入していない場合) 利用開始のために、あらかじめePassを購入していただきます。 購入希望の場合は<u>こちらから</u>お問い合わせをお願いします。 ※オンライン購入はできません



1. ePassの購入が完了すると、登録されたメールアドレス宛にクレジットキーが記載された 購入確認メールが配信されます。

III クレジットキーが不明な場合、トレーニングクレジットが不足している場合は、<u>お問い合わせ</u>ください。

- 2. Agilent Universityのアカウントにログインします。ログイン方法は<u>こちら</u>をご参照くださ い。画面の表示言語が英語になっている場合は、画面右上部より言語を日本語に変更して ください。
- 3. <u>こちら</u>をクリックし、スケジュールと受講料をクリックします。

Agilent Tusted Answers Revealed I and Answers
Agilent University TOP 選択中のコース トレーニングクレジット残高確認 マイページ ログイン/ログアウト
希 All Courses Classroom ▼ e-Learning ▼ ラーニングパス ▼ Live e-Learning ▼ ヘルプ ▼
コース概要
戻る
Agilent University ePass - 何度でもアクセスできる、自己学習型オンラインコンテンツ e-Learning ePass
ePassは、600時間以上あるオンラインコンテンツに3か月または12か月の希望する期間中、何度でもアクセスでき、必須の ラボスキルを向上することができます。 クロマトグラフィー、質望分析、元素分析、前処理など、アジレントのエキスパートがアジレントの機器に合わせて構成し た分析や操作方法を学ぶことができます。 トラブルシューティング、メンテナンス、高度な分析手法のコースを受講し、知識を高め、即戦力となるスキルを向上させ ることができます。
 ePassは、3か月間または12か月間のサブスクリブションサービス*です。 3か月間または12か月間の希望する期間中、プロダクトや言語に開わらず全ての オンラインコンテンツを視聴可能です。 目己ペースで学習できるオンラインコースは、主に英語で提供されます(注:日本語版コンテンツは一部のみに限定されます)。 600時間以上に及ぶ、500コース以上のコンテンツがご利用いただけます。 コースは、単独、または20以上のラーニングパズ(学習ガイド)に基づいて受講することができます。ラーニングパスは、特定装置の分析方法や操作を体系的に習得できるよう設計されています。 クラウドラボコースでは実践的な操作を注意を発見いただけます。このオンライントレーニングでは、ブラウザ上で装置のソフトウェアを実践するラボ(資源が含まれており、学びを深めることができます。
*ePassサブスクリプションサービスは、前払いの一括払いになり、自動更新はありません。
日本国内で、初めてePassサービスをご利用いただく場合(まだePassを購入していない場合) ご利用開始のために、あらかじめePassを購入していただきます。 購入希望の場合は Japan_training@agilent.com までお問い合わせをお願いします。 ※オンライン購入はできません
日本国内で、すでにePassサービスをご購入いただいたお客様 1.弊社より、メールにてご案内したクレジットキーを準備してください。 2.ごちらのユーザーガイドに沿ってお客様目身でアカウントを作成してください。 3.ユーザーガイドに沿ってePassの利用手続きを行ってください。 4. ePassの利用手続き完了後、弊社にてePassのアグティベーションを行います(1~3営業日以内)。 5. ePassのアグティベーションが完了しましたら、お客様にメールで通知が届きます。 6. 通知が届いたらアカウントにログインしてください。3か月間または12か月間の希望する期間中、全てのオンラインコン テンツの視聴が可能となります。
概要 スケジュールと受講料
'ジレントトレーニング - アジレント装置を最大限活用できるトレーニング 抑なトレーニングは不具合を減らし、効率と生産性を向上させ、装置のダウンタイムを低減します。また継続的なトレーニングは、自信と生産性のあるス ?ッフを育成し、離職率を減らすことができる研究結果もあります。
'ジ レントの自己ペース型オンライントレーニング・出張コストを抑えて学習できます 'ンライントレーニングは、いつでも、どこでも自己ペースで学習することができます。アジレントの自己ペース型オンラインコースは、幅広い内容と専門 職レベルを網羅していますので、スキルを向上できるコースが必ず見つかります。



4. 購入した ePass のプラン(3か月または12か月)を選択し、コースの追加をクリックします。



5. 登録内容を確認し、次へをクリックします。

Agilent	Trusted Anowers						表示言語:日本語 -	
Agilent Univer	rsity			TOP 選択中のC	1-2 NU-III	ングクレジット残蔵網路	לעליםערילים פֿיזעידי⊳	
All Courses C	lassroom +	e-Learning +	ラーニングバス・	Live e-Learning +	~いしづ ・			
			fi	ックアウト				
Agilent Universityトレー	ニングのお申込	込みにあたり以下の)ものが必須になります。					
クレジットキー: 個別の 装置シリアル番号	のクレジットキ	_						
 クレジットキーは、 装置付属の受講権をお クレジットキーをおお クレジットキーをおお 	トレーニングク お持ちの場合に 持ちの場合は、 持ちでない場合	レジットを購入す も、クレジットキー 最後の入力画面で は、最後の入力画面	ると発行される、英数字 ーが発行されます。 お持ちのクレジットキー 面で『JPBUND17』をご	で組み合わされたキー をご入力ください。 入力ください。	:例、ABC300	0000000_1000_xxxx	xです。	
※eLearningまたは、Li ※eLearningまたは、Li ※償別のクレジットキー ※シリアル番号は装置作 ※「トレーニングクレジ	※eLearningまたは、Live e-Learningをご希望のお客様は、クレジットキー『JPBUND17』 は使用できません。 ※eLearningまたは、Live e-Learningをご希望のお客様は、弊社からメールでお伝えする個別のクレジットキーが必要です。 ※個別のクレジットキーをお持ちでない場合は、担当営業またはトレーニング受付窓口までお問合せください。 ※シリアル番号は装置付属の受請権利、および該当コースの確認のため必ず入力ください。 ※「トレーニングクレジットの購入」に関して、現在日本では取扱いがございません。							
		7-74			トレーニン	・グクレジット(受講		
						料)		
Agilent University ePase コース詳細	s - Unlimited ad	cess to self-paced	online courses			600	MIR:	
			合計トレー	ニングクレジットの購	Л	600.00		
and the fail and a								
**石田(満子) ふりがた (せい・めい)			24					
**姓 (漢字)			友理恵					
**個人メールアドレス			xxxxxxxxx	agilent.com				
*電話番号								
*会社名			アジレント	・テクノロジー株式会	社			
部署								
役職								
*住所(都道府県)								
*住所(市町村・番地・)	建物)							
州(未選択のままでお適	(みください)			~				
*鄭使蕃号								
7 12			民	20	¥			



6. お支払い方法で、トレーニングクレジット **(**受講料**)** を選択し、クレジットキーを入力します。

!!! クレジットキーが不明な場合、トレーニングクレジットが不足している場合は、<u>お問い合わせ</u>ください。

Agilent Trusted Answers	裏示言語:日本語・
Agilent University	TOP 選択中のコース トレーニングクレジット残高確認 マイページ ログイン/ログアウト
# All Courses Classroom • e-Learning • ラーニングパス	・ Live e-Learning ・ ヘルプ・
雪友理恵	
「お支払方法」から「トレーニングクレジット」を選択後、クレジットキ・	チェックアウト -を入力し、
「キーを適用」を必ずクリックしてください。	
 クレジットキーは、トレーニングクレジットを購入すると発行される、 装置付属の受講権をお持ちの場合にも、クレジットキーが発行されます クレジットキーをお持ちの場合は、最後の入力画面でお持ちのクレジッ クレジットキーをお持ちでない場合は、最後の入力画面で『JPBUND17 	英数字で組み合わされたキー:例、ABC30000000_1000_xxxxxです。 ・ トキーをご入力ください。 』をご入力ください。
※e-Learningまたは、Live e-Learningをご希望のお客様は、クレジットキー ※e-Learningまたは、Live e-Learningをご希望のお客様は、弊社からメール ※個別のクレジットキーをお持ちでない場合は、担当営業またはトレーニ ※シリアル番号は装置付属の受講権利、および該当コースの確認のため必 ※「トレーニングクレジットの購入」に関して、現在日本では取扱いがご	『JPBUND17』 は使用できません。 でお伝えする個別のクレジットキーが必要です。 ング受付窓口までお問合せください。 ず入力ください。 ざいません。
コース名	トレーニングクレジット(受講料)
Agilent University ePass - Unlimited access to self-paced online courses コース詳細	600
合計	トレーニングクレジットの購入 600.00
お支払方法 トレーニングクレジット(受講科) ✔	
クレジットキー	

7. クレジットキーを入力し、申し込みの完了をクリックします。

‡ –	利用可能	ニューバランス	アクション				
ABC123 456789	600	0	取り去る				
				戻る	クリア	申し込みの完了	



8. ePassの登録が完了しました。

🖶 All Courses	Classroom •	Self-paced 🔻	Virtual Instructor-led 🔹	Keeping You Safe					
Inhi Kim									
			Registration	Completed					
Confirmation	Number		Registratio	n Details				Training Credits	Status
549-18857-166251 410863	-221- Agiler cours	nt University ePa es	ss - Unlimited access to s	elf-paced online	Launch	Course Details		600	Confirmed
							Total	600	
You will receive a First Name: Last Name: Email: Preferred Phone Company: Street 1: City: Zip:	an email notificat Kelly Kim xxxxxxxxx @agi 2: xxx-xxx-xxxx Agilent 11F 369 Gang Seoul 137858	ion with your regi lent.com nam-daero, Seocl	stration information. 10-gu						
Country:	South Korea								
Payment Metho	d: Training Credit	ts							

- 9. ご登録いただいたメールアドレスに、Agilent University から登録確認メールが配信されま す。
- 10. 弊社にてePass のアクティベーション手続きをします。手続き完了後、1~3営業日以内に アクティベーション完了通知がメールで配信されます。
- **11.** 完了通知配信後よりePassの利用が可能です。Agilent Universityに<u>ログイン</u>し視聴してくだ さい。アカウントにログインすると、すべてのe-Learningコースに登録ボタンが表示されま す。個別のコースを検索する方法は<u>こちら</u>をご覧ください。

■ アカウントログイン前、ePass アクティベーション前





Agilent Universityアカウントにログインする

1. 言語を日本語にします。



2. Agilent University Webサイトにログインします。<u>Agilent Technologies (viewcentral.com)</u>. ログインまたは、 初めてご利用される場合は新規にアカウント作成します。

Agilent Trusted Answers	表示言語:日本語・
Agilent University	
All Courses Classroom • e-L	earning ・ ラーニングパス ・ Live e-Learning ・ ヘルプ ・
ユーザー名の確認 パスワードリセット アカウント作成	ユーザー名 パスワード クリア ズヘ
🖨 All Courses Classroom 🔹	e-Learning ▼ ラーニングパス ▼ Live e-Learning ▼ ヘルプ ▼
	*名前(漢字)



 以下のリンクをクリックいただくと、各装置のラーニングパスが表示されます。アジレント 推奨のコンテンツに沿って受講いただくと、効率的にスキルアップいただくことができます。

備考:各種ラーニングパスで紹介するコンテンツの多くは英語で提供されます。 一部、日本語字幕付きで視聴が可能なコンテンツも用意しております。

Instrument	Links
	<u>HPLC ラーニングパス</u>
	<u>LC/MS ラーニングパス</u>
	<u>GC ラーニングパス</u>
	<u>GC/MS ラーニングパス</u>
	<u>ICP-MS ラーニングパス</u>
	<u>元素分析 ラーニングパス</u>
	<u>データマネジメントラーニングパス</u>





個々のコースを見るには、メニューのトレーニングとイベントから、トレーニングを クリックし、オンライントレーニング項目内の、お申し込みはこちらをクリックします。







2. トレーニング形態からe-Learningを選択し、必要に応じて言語から日本語を選択し、 コースを検索します。



検索キーワードや、以下の条件で検索いただけます 言語: コースの配信言語を選択します 機種: 装置の種類を選択します 学習内容:学びたい学習分野を選択します ソフトウェアプラットフォーム: ソフトウェアを選択します

上記を選択後、適用をダブルクリックします



 コース概要をクリックすると、トレーニングコースの詳細をご確認いただけます。 スケジュールと受講料からコースを登録します。

Agilent Trusted Ans				表示言語:日本語 -
Agilent University			TOP 選択中のコース トレーニングクレジット残高確認 マイベ	ージ ログイン/ログアウト
All Courses Classroom	• • e-Learning •	ラーニングパス・	Live e-Learning ・ ヘルプ・	
		トレーニン	ッグコースカタログ	
フィルター				
フィルター: <mark>通用</mark> すべてクリア (2)	0			
キーワード へ	(e) X-L		▶レーニングが感 ④	日数
B ~	A IS	Agilent 5977B MassHunter Software e-	e-Learning	 ③ コース概要 圖 スケジュールと受
会場 く		Introduction (Acq 10.1; Qual 10.0 / Quant 10.2) GCMS-MH-1100zsV4	L	= 講料

4. コース情報を確認します



Agilent OpenLab CDS ChemStation – Sequences (C.01.09 and C.01.10) e-Learning HPLC-OL-1180rV2

This interactive module describes how single runs can be grouped into sequences of runs. The runs are acquired, evaluated and stored together.

- * This course may be accessed over a 90-day period.
- ****** This is a certificate course.

★ 一部有効期限が設定されているコースもあります。
★★ 完了すると、完了メールが届きます。



5. スケジュールと受講料を選択し、登録をクリックします。

概要 スケジュールと受講料		
言語: 日本語 日数: 0.75 時間		91 9
ePass が利用できない場合 * :	概要 スケジュールと受講科	
登録ボタンが表示されません	重要: 日本縣 日勤: 0.75時間	0.13
ePassが利用できる場合・	概要 スケジュールと受達料	
登録ボタンが表示されます	重語:日本語 日歌: 0.75 時間	20
*登録ボタンが表示されない場合は、	アカウントにログインができているか、	
ePass の利用手続きが全て完了して	いるかをご確認ください。	

6. コース登録完了

	Re	gistration Completed	
Confirmatio	n Number	Registration Details	Status
549-35707-1662	251-3-403099 HPLC-OLII-1140eV2 - Agilent Infinity LC systems with Open	ab CDS - Integration and viewing results (Ver. 2.6)	Course Details Confirmed
You will receive a	n email notification with your registration information.		
First Name:	Kelly		
Last Name:	Kim		
Email:	xxxxxxx@agilent.com		
Preferred Phone	XXX-XXX-XXXX		
Company:	Agilent		
Street 1:	11F 369 Gangnam-daero, Seocho-gu		
City:	Seoul		
Zip:	137858		
Country:	South Korea		





7. 右上のマイページをクリックします。

Agilent	Trusted Answers			表示言	語:日本語 -
Agilent Uni	versity			TOP 選択中のコース トレーニングクレジット残高確認 マイページ コグイ	ン/ログアウト
All Courses	Classroom •	e-Learning 🕇	ラーニングパス・	Live e-Learning ・ ヘルプ・	

8. マイページに新しいe-Learningが追加されます。 開始するをクリックするとコースが開始します。

		アカウント情報		
申込済みコース 受講履歴	アカウント情報			
- <i>∠</i> €	会場④	ステータス・	日程④	
Agilent 1260 LC シ ステム - 装置の握 礎 549-2758-233774-188- 50555	eLearning - Japanese	Confirmed		① □ス概要③ 開始する



問い合わせ先

E-mail japan_training@agilent.com

Website https://www.chem-agilent.com/contents.php?id=1005849

THAILAND

Agilent Technologies (Thailand) Ltd. U Chu Liang Building 22/F Unit A.D, 968 Rama IV Rd, Silom, Bangrak, Bangkok 10500 (66) 2 6376363 Option 3 (Sales & Marketing) Option 1 (Instrument & Service)

SINGAPORE

Agilent Technologies Singapore (Sales) Pte Ltd 1 Yishun Avenue 7 Singapore 768923 1800 276 2622 (Select Option 3)

SOUTH KOREA

한국애질런트테크놀로지스(주) 대한민국 서울특별시 서초구 강남대로 369 에이플러스 에셋타워 9층 06621 • 080 004 5090 (옵션 3, 영업 및 마케팅)

JAPAN

アジレント・テクノロジー株式会社 東京都八王子市高倉町9-1 ⑤ 0120-477-111

MALAYSIA

Agilent Technologies Sales (Malaysia) Sdn Bhd Unit 201, Level 2, Uptown 2, 2 Jalan SS21/37 Damansara Uptown 47400 Petaling Jaya, Selangor 1800 88 0805 Option 3 (Sales & Marketing) Option 1 (Instrument & Service)

AUSTRALIA and New Zealand Agilent Technologies Australia 679 Springvale Road, Mulgrave, Victoria 3170 Australia : 1800 802 402 New Zealand : 0508555344 Option 2 (Sales) Option 3 (Service)

Agilent

記載内容は事前通知なく変更される可能性があります

DE29778367 © Agilent Technologies, Inc. 2024 Published in the Japan JJuly 2024