

OncoMine™ Dx Express Test

Pub. No. MAN0025897 Rev. D.0

User guides for this product and updates to Genexus™ Dx Software and Assay Definition Files (ADFs) for the OncoMine™ Dx Express Test are provided in a Thermo Fisher™ Connect Platform resource library.

Download user documentation

- Go to apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, then sign in to thermofisher.com.
Note: If you do not have a thermofisher.com account, click **Create Account** in the sign in screen, then follow the instructions.
- In the resource library, locate the OncoMine™ Dx Express Test user guides in the language of interest, then download and extract the ZIP folder containing the user guides.
- Open the user guides with a PDF viewer such as Adobe Reader™ or equivalent.
Note: Adobe Reader™ can be downloaded at [Adobe Reader](#).

Download Genexus™ Dx Software and ADF updates


Link a Thermo Fisher™ Connect Platform account with a QR code (manager/administrator)

To access software updates, you must first link an instrument to a Connect Platform account. Linking an instrument is a one-time action.


- In Genexus™ Dx Software, click  **(Settings)** ▶ **Thermo Fisher Account**.
- Click **+ Link Account to Device**.
The **Link Account to Device** dialog box opens and a QR code image and QR code text are shown.
- Link the Connect Platform account with the software using one of two following options.


Options for linking a Connect Platform account

Option	Description
Link with the Instrument Connect mobile app	Note: To use this option, you must have the InstrumentConnect mobile app installed on a mobile device. <ol style="list-style-type: none"> Open the InstrumentConnect app, then sign in. In the app, click +. Click QR Code, then scan the QR code image shown in the Link Account to Device dialog box in Genexus™ Dx Software.

Option	Description
Link through the Connect Platform website	<ol style="list-style-type: none"> Copy the QR code text given in the dialog box, then click Thermo Fisher Connect to open the Connect Platform website. Note: Sign in is required. If you do not have a Thermo Fisher™ Connect Platform account, click Create Account in the sign in screen, then follow the instructions. In the Connect Platform website, click  InstrumentConnect in the left sidebar. Click Link instrument in the upper right corner of the screen. Enter the QR code text shown in the dialog box in the Link Instrument screen. Click Link.

Download and install software and ADF updates

When the Thermo Fisher™ Connect Platform account is linked, available software and ADF updates are visible in the Genexus™ Dx Software  **(Settings)** ▶ **Software Updates** screen. Manager- and administrator-level users can download and install the updates.

- In Genexus™ Dx Software, click ▶  **(Settings)** ▶ **Software Updates**.
- In the **Installed Software** screen, click **Software Updates** to view updates to software and ADFs that are available at the Connect Platform. If not selected, click the **APP STORE** button.
- In the **Actions** column for the software or ADF package of interest, click **Download**.
- After the download is complete, click **Install**.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

Oncomine™ Dx Express Test

№ на публикацията MAN0025897 Ред. D.0

Ръководства за потребителя на този продукт и актуализации за Genexus™ Dx Software, и файлове с дефиниции на анализи в (ADF) за Oncomine™ Dx Express Test са налични в библиотеката с ресурси на Thermo Fisher™ Connect Platform.

Изтегляне на документация за потребителя

1. Отворете apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, след което влезте в thermofisher.com.

Забележка: Ако нямате акаунт в thermofisher.com, щракнете върху **Create Account (Създаване на акаунт)** на екрана за вписване, след което следвайте инструкциите.


2. В библиотеката с ресурси намерете ръководствата за потребителя на Oncomine™ Dx Express Test на желанния език, след което изтеглетте и разархивирайте ZIP папката, съдържаща ръководствата за потребителя.
3. Отворете ръководствата за потребителя с PDF визуализатор като Adobe Reader™ или еквивалентен.

Забележка: Adobe Reader™ може да се изтегли от [Adobe Reader](#).

Изтеглетте Genexus™ Dx Software и актуализации на ADF

Свържете акаунт в Thermo Fisher™ Connect Platform с QR код (ръководител/администратор)


За да получите достъп до актуализации на софтуера, трябва да свържете инструмент с акаунт в Connect Platform. Свързването на инструмент е еднократно действие.

1. В Genexus™ Dx Software щракнете върху  **(Settings) (Настройки) ▶ Thermo Fisher Account (Акаунт в Thermo Fisher)**.
2. Щракнете върху **+ Link Account to Device (Свързване на акаунта с устройството)**.


Отваря се диалоговият прозорец **Link Account to Device (Свързване на акаунта с устройството)** и се показва изображение на QR код и текст на QR код.


3. Свържете акаунта в Connect Platform със софтуера с помощта на една от следните две опции.

Опции за свързване на акаунт в Connect Platform

Опция	Описание
Връзка с мобилно приложение InstrumentConnect	<p>Забележка: За да използвате тази опция, трябва да имате мобилното приложение InstrumentConnect инсталирано на мобилно устройство.</p> <ol style="list-style-type: none"> 1. Отворете приложението InstrumentConnect и се впишете. 2. В приложението щракнете върху +. 3. Щракнете върху QR Code (QR код) и сканирайте изображението на QR кода, показано в диалоговия прозорец Link Account to Device (Свързване на акаунта с устройството) в Genexus™ Dx Software.
Връзка чрез уебсайта Connect Platform	<ol style="list-style-type: none"> 1. Копирайте текста от QR кода, посочен в диалоговия прозорец, след което щракнете върху Thermo Fisher Connect, за да отворите уебсайта на Connect Platform. <p>Забележка: Изисква се вписване. Ако нямате акаунт в Thermo Fisher™ Connect Platform, щракнете върху Create Account (Създаване на акаунт) на екрана за вписване, след което следвайте инструкциите.</p> <ol style="list-style-type: none"> 2. В уебсайта на Connect Platform щракнете върху  InstrumentConnect в лявата странична лента. 3. Щракнете върху Link instrument (Свързване на инструмента) в горния десен ъгъл на екрана. 4. Въведете текста на QR кода, показан в диалоговия прозорец на екрана Link Instrument (Свързване на инструмента). 5. Щракнете върху Link (Свързване).

Изтегляне и инсталиране на актуализации на софтуера и ADF

Когато е свързан акаунт в Thermo Fisher™ Connect Platform, наличните актуализации на софтуера и ADF се виждат на екрана Genexus™ Dx Software  **(Settings) (Настройки) ▶ Software Updates (Актуализации на софтуера)**. Потребителите на ниво ръководител и администратор могат да изтеглят и инсталират актуализациите.

1. В Genexus™ Dx Software щракнете върху **▶  (Settings) (Настройки) ▶ Software Updates (Актуализации на софтуера)**.
2. На екрана **Installed Software (Инсталиран софтуер)** щракнете върху **Software Updates (Актуализации на софтуера)**, за да прегледате актуализациите на софтуера и ADF, налични от Connect Platform. Ако не е избран, щракнете върху бутона **APP STORE (МАГАЗИН ЗА ПРИЛОЖЕНИЯ)**.
3. В колоната **Actions (Действия)** на желанния софтуерен или ADF пакет, щракнете върху **Download (Изтегляне)**.
4. Когато изтеглянето завърши, щракнете върху **Install (Инсталиране)**.

Ограничена гаранция на продукта

Life Technologies Corporation и/или нейните свързани дружества гарантират своите продукти, както е посочено в Общите условия за продажба на Life Technologies, намиращи се на www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Ако имате въпроси, моля, свържете се с Life Technologies на адрес www.thermofisher.com/support.

Oncomine™ Dx Express Test

Pub. č. MAN0025897 Rev. D.0

Uživatelské příručky k tomuto produktu a aktualizace softwaru Genexus™ Dx Software a souborů s definicemi analýz (Assay Definition File, ADF) pro test Oncomine™ Dx Express Test jsou k dispozici v knihovně zdrojů Thermo Fisher™ Connect Platform.

Stažení uživatelské dokumentace

1. Přejděte na stránku apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, poté se přihlaste na webu thermofisher.com.

Poznámka: Pokud účet na webu thermofisher.com nemáte, klikněte na přihlašovací obrazovce na možnost **Create Account (Vytvořit účet)** a postupujte podle pokynů.


2. V knihovně zdrojů vyhledejte uživatelské příručky pro test Oncomine™ Dx Express Test v jazyce, který vás zajímá, a poté si stáhněte a rozbalte složku ZIP obsahující uživatelské příručky.
3. Otevřete uživatelské příručky v prohlížeči PDF, například Adobe Reader™ nebo jiném obdobném programu.

Poznámka: Program Adobe Reader™ si můžete stáhnout zde: [Adobe Reader](#).

Stažení aktualizací softwaru Genexus™ Dx Software a souboru ADF

Propojení účtu Thermo Fisher™ Connect Platform s QR kódem (manažer/správce)

Chcete-li získat přístup k aktualizacím softwaru, musíte nejprve propojit přístroj s účtem Connect Platform. Propojení přístroje je jednorázová záležitost.


1. V softwaru Genexus™ Dx Software klikněte na možnost  **(Settings) (Nastavení) ▶ Thermo Fisher Account (Účet Thermo Fisher)**.
2. Klikněte na možnost **+ Link Account to Device (Propojit účet se zařízením)**.

Otevře se dialogové okno **Link Account to Device (Propojit účet se zařízením)** a zobrazí se obrázek a text QR kódu.

3. Propojte účet Connect Platform se softwarem pomocí jedné ze dvou následujících možností.

Možnosti propojení účtu Connect Platform

Možnost	Popis
Propojení s mobilní aplikací InstrumentConnect	<p>Poznámka: Chcete-li tuto možnost použít, musíte mít v mobilním zařízení nainstalovanou mobilní aplikaci InstrumentConnect.</p> <ol style="list-style-type: none"> 1. Otevřete aplikaci InstrumentConnect a přihlaste se. 2. V aplikaci klikněte na možnost +. 3. Klikněte na možnost QR Code (QR kód) a poté naskenujte snímek QR kódu zobrazený v dialogovém okně Link Account to Device (Propojit účet se zařízením) v softwaru Genexus™ Dx Software.

Možnost	Popis
Propojení přes webovou stránku Connect Platform	<ol style="list-style-type: none"> 1. Zkopírujte text QR kódu uvedený v dialogovém okně a poté kliknutím na Thermo Fisher Connect otevřete webovou stránku Connect Platform. <p>Poznámka: Je nutné se přihlásit. Pokud účet Thermo Fisher™ Connect Platform nemáte, klikněte na přihlašovací obrazovce na možnost Create Account (Vytvořit účet) a postupujte podle pokynů.</p> <ol style="list-style-type: none"> 2. Na webové stránce Connect Platform klikněte v levém postranním panelu na možnost  InstrumentConnect. 3. V pravém horním rohu obrazovky klikněte na možnost Link instrument (Připojit přístroj). 4. Zadejte text QR kódu zobrazený v dialogovém okně na obrazovce Link Instrument (Propojit přístroj). 5. Klikněte na možnost Link (Propojit).

Stažení a instalace aktualizací softwaru a souborů ADF

Po propojení účtu Thermo Fisher™ Connect Platform jsou na obrazovce Genexus™ Dx Software  **(Settings) (Nastavení) ▶ Software Updates (Aktualizace softwaru)** viditelné dostupné aktualizace softwaru a souborů ADF. Uživatelé na úrovni manažera a správce mohou aktualizace stahovat a instalovat.

1. V softwaru Genexus™ Dx Software klikněte na možnost  **(Settings) (Nastavení) ▶ Software Updates (Aktualizace softwaru)**.
2. Na obrazovce **Installed Software (Nainstalovaný software)** klikněte na položku **Software Updates (Aktualizace softwaru)** a zobrazte aktualizace softwaru a souborů ADF, které jsou k dispozici na platformě Connect Platform. Pokud není vybrána, klikněte na tlačítko **APP STORE**.
3. Ve sloupci **Actions (Akce)** u softwarového nebo ADF balíčku, který vás zajímá, klikněte na tlačítko **Download (Stáhnout)**.
4. Po dokončení stažení klikněte na možnost **Install (Instalovat)**.

Omezená záruka na výrobek

Společnost Life Technologies Corporation a/nebo její pobočky poskytují záruku na své výrobky podle ustanovení ve smluvních podmínkách pro prodej společnosti Life Technologies na adrese www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Máte-li jakékoli dotazy, kontaktujte prosím společnost Life Technologies na adrese www.thermofisher.com/support.

Oncomine™ Dx Express Test

Publ. nr. MAN0025897 Rev. D.0

Brugervejledninger til dette produkt og opdateringer til Genexus™ Dx Software og analysedefinitionsfil (ADF'er) til Oncomine™ Dx Express Test findes i et Thermo Fisher™ Connect Platform-ressourcebibliotek.

Download brugerdokumentation

1. Gå til apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, og log derefter ind på thermofisher.com.

Bemærk: Hvis du ikke har en konto på thermofisher.com account, skal du klikke på **Create Account (Opret konto)** på logonskærmen og derefter følge instruktionerne.


2. Find brugervejledningerne til Oncomine™ Dx Express Test på det ønskede sprog i ressourcebiblioteket, og download og udtræk derefter ZIP-mappen med brugervejledningerne.
3. Åbn brugervejledningerne med en PDF-fremviser som Adobe Reader™ eller lignende.

Bemærk: Adobe Reader™ kan downloades på [Adobe Reader](#).

Download Genexus™ Dx Software- og ADF-opdateringer

Tilknyt en Thermo Fisher™ Connect Platform-konto til en QR-kode (chef/administrator)


For at få adgang til softwareopdateringer skal du først knytte et instrument til en Connect Platform-konto. Tilknytning af et instrument skal kun gøres én gang.

1. I Genexus™ Dx Software skal du klikke på  (**Settings (Indstillinger)**) ▶ **Thermo Fisher Account (Thermo Fisher-konto)**.
2. Klik på **+ Link Account to Device (Tilknyt konto til enhed)**. Dette åbner dialogboksen **Link Account to Device (Tilknyt konto til enhed)**, som indeholder en QR-kode i billed- og tekstformat.
3. Knyt Connect Platform-kontoen til softwaren ved hjælp af en af de følgende to muligheder.

Valgmuligheder for tilknytning af en Connect Platform-konto

Valgmulighed	Beskrivelse
Tilknytning med mobilappen InstrumentConnect	<p>Bemærk: For at bruge denne valgmulighed skal du have en mobilenhed med mobilappen InstrumentConnect installeret.</p> <ol style="list-style-type: none"> 1. Åbn appen InstrumentConnect, og log ind. 2. Klik på + i appen. 3. Klik på QR Code (QR-kode), og scan derefter det QR-kodebillede, der vises i dialogboksen Link Account to Device (Knyt konto til enhed) i Genexus™ Dx Software.

Valgmulighed	Beskrivelse
Tilknytning via webstedet Connect Platform	<ol style="list-style-type: none"> 1. Kopiér den QR-kodetekst, der er angivet i dialogboksen, og klik derefter på Thermo Fisher Connect for at åbne webstedet Connect Platform. Bemærk: Login er påkrævet. Hvis du ikke har en konto på Thermo Fisher™ Connect Platform, skal du klikke på Create Account (Opret konto) i på login-skærmbilledet og derefter følge anvisningerne. 2. På webstedet Connect Platform skal du klikke på  InstrumentConnect i venstre sidepanel. 3. Klik på Link instrument (Tilknyt instrument) i øverste højre hjørne af skærmbilledet. 4. Indtast QR-kodeteksten, der vises i dialogboksen på skærmbilledet Link Instrument (Tilknyt instrument). 5. Klik på Link (Tilknyt).

Download og installation af software- og ADF-opdateringer
Når Thermo Fisher™ Connect Platform-kontoen er tilknyttet, vises de tilgængelige software- og ADF-opdateringer på skærmbilledet Genexus™ Dx Software  (**Settings (Indstillinger)**) ▶ **Software Updates (Softwareopdateringer)**. Disse opdateringer kan downloades og installeres af brugere med chef- og administratorrettigheder.

1. I Genexus™ Dx Software skal du klikke på ▶  (**Settings (Indstillinger)**) ▶ **Software Updates (Softwareopdateringer)**.
2. På skærmbilledet **Installed Software (Installeret software)** skal du klikke på **Software Updates (Softwareopdateringer)** for at få vist opdateringer til software og ADF'er, som er tilgængelig i Connect Platform. Klik på knappen **APP STORE**, hvis det ikke er valgt.
3. Klik på **Download** for den relevante software- eller ADF-pakke i kolonnen **Actions (Handlinger)**.
4. Klik på **Install (Installer)**, når overførslen er færdig.

Begrænset produktgaranti

Life Technologies Corporation og/eller deres datterselskab(er) yder garanti på deres produkter i henhold til Life Technologies' generelle vilkår og betingelser for salg (General Terms and Conditions of Sale), som kan ses på adressen www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Kontakt Life Technologies på adressen www.thermofisher.com/support i tilfælde af spørgsmål.

Oncomine™ Dx Express Test

Pub.-Nr. MAN0025897 Vers. D.0

Gebrauchsanweisungen für dieses Produkt, Aktualisierungen für die Genexus™ Dx Software und Assaydefinitionsdateien (ADF) für den Oncomine™ Dx Express Test werden in einer Thermo Fisher™ Connect Platform Ressourcenbibliothek bereitgestellt.


Benutzerdokumentation herunterladen

- Gehen Sie zu apps.thermofisher.com/apps/spa/#/publiclib/genexusdx und melden Sie sich bei thermofisher.com an.
Hinweis: Wenn Sie kein thermofisher.com Konto haben, klicken Sie im Anmeldebildschirm auf **Create Account (Konto anlegen)**, und folgen Sie dann den Anweisungen.
- Suchen Sie in der Ressourcenbibliothek nach den Oncomine™ Dx Express Test Benutzerhandbüchern in der gewünschten Sprache. Laden Sie dann das ZIP-Verzeichnis mit den Benutzerhandbüchern herunter und entpacken Sie es.
- Öffnen Sie die Benutzerhandbücher mit einem PDF-Viewer wie Adobe Reader™ (oder einem gleichwertigen Programm).
Hinweis: Adobe Reader™ kann unter [Adobe Reader](#) heruntergeladen werden.


Genexus™ Dx Software und ADF-Aktualisierungen herunterladen

Ein Thermo Fisher™ Connect Platform Konto mit einem QR-Code verknüpfen (Manager/Administrator)

Um auf Software-Updates zuzugreifen, müssen Sie zuerst ein Gerät mit einem Connect Platform Konto verknüpfen. Geräte müssen nur einmal verknüpft werden.


- Klicken Sie in der Genexus™ Dx Software auf  (**Settings (Einstellungen)**) ▶ **Thermo Fisher Account (Thermo Fisher Konto)**.
- Klicken Sie auf **+ Link Account to Device (Konto mit Gerät verbinden)**.
Das Dialogfeld **Link Account to Device (Konto mit Gerät verbinden)** wird geöffnet und ein QR-Code-Bild und QR-Code-Text werden angezeigt.
- Verknüpfen Sie das Connect Platform Konto mit einer der beiden folgenden Optionen mit der Software.

Optionen zum Verknüpfen eines Connect Platform Kontos

Option	Beschreibung
Mit der Instrument-Connect Mobile App verknüpfen	<p>Hinweis: Um diese Option zu nutzen, müssen Sie die InstrumentConnect Mobile App auf einem Mobilgerät installiert haben.</p> <ol style="list-style-type: none"> Öffnen Sie die InstrumentConnect App und melden Sie sich an. Klicken Sie in der App auf +. Klicken Sie auf QR Code (QR-Code) und scannen Sie dann die QR-Code-Grafik, die im Dialogfeld Link Account to Device (Konto mit Gerät verbinden) in der Genexus™ Dx Software angezeigt wird.
Über die Connect Platform Website verknüpfen	<ol style="list-style-type: none"> Kopieren Sie den im Dialogfeld angezeigten QR-Code-Text, und klicken Sie dann auf Thermo Fisher Connect, um die Connect Platform Website zu öffnen. Hinweis: Eine Anmeldung ist erforderlich. Wenn Sie kein Thermo Fisher™ Connect Platform Konto haben, klicken Sie im Anmeldebildschirm auf Create Account (Konto anlegen), und folgen Sie dann den Anweisungen. Klicken Sie in der Connect Platform Website auf  InstrumentConnect in der linken Randleiste. Klicken Sie in der oberen rechten Ecke des Bildschirms auf Link instrument (Gerät verknüpfen). Geben Sie den QR-Code-Text ein, der in dem Dialogfeld im Bildschirm Link Instrument (Gerät verknüpfen) angezeigt wird. Klicken Sie auf Link (Verknüpfen).

Software- und ADF-Aktualisierungen herunterladen und installieren

Wenn das Thermo Fisher™ Connect Platform Konto verknüpft ist, werden verfügbare Software- und ADF-Aktualisierungen in der Genexus™ Dx Software im Bildschirm  (**Settings (Einstellungen)**) ▶ **Software Updates (Software-Updates)** angezeigt. Benutzer mit Manager- und Administratorrechten können die Aktualisierungen herunterladen und installieren.

- Klicken Sie in der Genexus™ Dx Software auf ▶  (**Settings (Einstellungen)**) ▶ **Software Updates (Software-Updates)**.
- Klicken Sie im Bildschirm **Installed Software (Installierte Software)** auf **Software Updates (Software-Updates)**, um Aktualisierungen für Software und ADFs anzuzeigen, die in der Connect Platform verfügbar sind. Falls nicht ausgewählt, klicken Sie auf die Schaltfläche **APP STORE (App-Store)**.
- Klicken Sie in der Spalte **Actions (Aktionen)** des gewünschten Software- oder ADF-Pakets auf **Download (Herunterladen)**.
- Nachdem der Download abgeschlossen ist, klicken Sie auf **Install (Installieren)**.

Eingeschränkte Produktgarantie

Life Technologies Corporation und/oder seine Tochtergesellschaft(en) bieten für ihre Produkte eine Gewährleistung gemäß den allgemeinen Verkaufs- und Lieferbedingungen von Life Technologies, die unter www.thermofisher.com/us/en/home/global/terms-and-conditions.html eingesehen werden können. Bitte wenden Sie sich bei Fragen unter www.thermofisher.com/support an Life Technologies.

Oncomine™ Dx Express Test

Αρ. δημ. MAN0025897 Αναθ. D.0

Οι οδηγοί χρήστη για αυτό το προϊόν και οι ενημερώσεις για το Genexus™ Dx Software και τα αρχεία ορισμού δοκιμασίας (ADF) για το Oncomine™ Dx Express Test παρέχονται στη βιβλιοθήκη πόρων Thermo Fisher™ Connect Platform.

Λήψη τεκμηρίωσης χρήστη

1. Μεταβείτε στη διεύθυνση apps.thermofisher.com/apps/spa/#/publiclib/genexusdx και, στη συνέχεια, συνδεθείτε στον ιστότοπο thermofisher.com.

Σημείωση: Εάν δεν έχετε λογαριασμό στον ιστότοπο thermofisher.com, κάντε κλικ στο **Create Account (Δημιουργία λογαριασμού)** στην οθόνη σύνδεσης και, στη συνέχεια, ακολουθήστε τις οδηγίες.


2. Στη βιβλιοθήκη πόρων, εντοπίστε τους οδηγούς χρήστη του Oncomine™ Dx Express Test στη γλώσσα που σας ενδιαφέρει και, στη συνέχεια, προβείτε σε λήψη και εξαγωγή του φακέλου ZIP που περιέχει τους οδηγούς χρήστη.
3. Ανοίξτε τους οδηγούς χρήστη με ένα πρόγραμμα προβολής PDF, όπως Adobe Reader™ ή άλλο αντίστοιχο.

Σημείωση: Μπορείτε να πραγματοποιήσετε λήψη του Adobe Reader™ από τον ιστότοπο [Adobe Reader](https://www.adobe.com/reader).

Λήψη Genexus™ Dx Software και ενημερώσεων ADF

Σύνδεση ενός λογαριασμού Thermo Fisher™ Connect Platform με κωδικό QR (υπεύθυνος/διαχειριστής)


Για να αποκτήσετε πρόσβαση σε ενημερώσεις λογισμικού, πρέπει πρώτα να συνδέσετε ένα όργανο με έναν λογαριασμό Connect Platform. Η σύνδεση ενός οργάνου πραγματοποιείται μία φορά.

1. Στο Genexus™ Dx Software, κάντε κλικ στο  **[Settings (Ρυθμίσεις)]** ▶ **Thermo Fisher Account (Λογαριασμός Thermo Fisher)**.
2. Κάντε κλικ στο **+ Link Account to Device (Σύνδεση λογαριασμού με συσκευή)**.

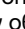
Το παράθυρο διαλόγου **Link Account to Device (Σύνδεση λογαριασμού με συσκευή)** ανοίγει και εμφανίζονται μια εικόνα κωδικού QR και κείμενο κωδικού QR.

3. Συνδέστε τον λογαριασμό Connect Platform με το λογισμικό χρησιμοποιώντας έναν από τους δύο ακόλουθους τρόπους.

Επιλογές για τη σύνδεση ενός λογαριασμού Connect Platform

Επιλογή	Περιγραφή
Σύνδεση με την εφαρμογή για κινητά Instrument Connect	<p>Σημείωση: Για να χρησιμοποιήσετε αυτήν την επιλογή, πρέπει να διαθέτετε την εφαρμογή για κινητά InstrumentConnect εγκατεστημένη σε μια κινητή συσκευή.</p> <ol style="list-style-type: none"> 1. Ανοίξτε την εφαρμογή InstrumentConnect και, στη συνέχεια, συνδεθείτε. 2. Στην εφαρμογή, κάντε κλικ στο +. 3. Κάντε κλικ στην επιλογή QR Code (Κωδικός QR) και, στη συνέχεια, σαρώστε την εικόνα κωδικού QR που εμφανίζεται στο παράθυρο διαλόγου Link Account to Device (Σύνδεση λογαριασμού με συσκευή) στο Genexus™ Dx Software.
Σύνδεση μέσω του ιστότοπου Connect Platform	<ol style="list-style-type: none"> 1. Αντιγράψτε το κείμενο κωδικού QR που παρέχεται στο παράθυρο διαλόγου και, στη συνέχεια, κάντε κλικ στο Thermo Fisher Connect για να ανοίξετε τον ιστότοπο Connect Platform. Σημείωση: Απαιτείται σύνδεση. Εάν δεν έχετε λογαριασμό Thermo Fisher™ Connect Platform, κάντε κλικ στο Create Account (Δημιουργία λογαριασμού) στην οθόνη σύνδεσης και, στη συνέχεια, ακολουθήστε τις οδηγίες. 2. Στον ιστότοπο Connect Platform, κάντε κλικ στην επιλογή  InstrumentConnect στην αριστερή πλευρική μπάρα. 3. Κάντε κλικ στην επιλογή Link instrument (Σύνδεση οργάνου) στην επάνω δεξιά γωνία της οθόνης. 4. Καταχωρίστε το κείμενο κωδικού QR που εμφανίζεται στο παράθυρο διαλόγου της οθόνης Link instrument (Σύνδεση οργάνου). 5. Κάντε κλικ στην επιλογή Link (Σύνδεση).

Λήψη και εγκατάσταση ενημερώσεων λογισμικού και ενημερώσεων ADF

Όταν υπάρχει συνδεδεμένος λογαριασμός Thermo Fisher™ Connect Platform, οι διαθέσιμες ενημερώσεις λογισμικού και ADF εμφανίζονται στην οθόνη Genexus™ Dx Software  **[Settings (Ρυθμίσεις)]** ▶ **Software Updates (Ενημερώσεις λογισμικού)**. Οι χρήστες επιπέδου υπεύθυνου και διαχειριστή μπορούν να πραγματοποιήσουν λήψη και εγκατάσταση των ενημερώσεων.

1. Στο Genexus™ Dx Software, κάντε κλικ στην επιλογή  **[Settings (Ρυθμίσεις)]** ▶ **Software Updates (Ενημερώσεις λογισμικού)**.
2. Στην οθόνη **Installed Software (Εγκατεστημένο λογισμικό)**, κάντε κλικ στην επιλογή **Software Updates (Ενημερώσεις λογισμικού)** για να δείτε τις ενημερώσεις λογισμικού και ADF που είναι διαθέσιμες στο Connect Platform. Εάν δεν έχει επιλεγεί, κάντε κλικ στο κουμπί **APP STORE (ΚΑΤΑΣΤΗΜΑ ΕΦΑΡΜΟΓΩΝ)**.
3. Στη στήλη **Actions (Ενέργειες)** για το πακέτο λογισμικού ή ADF που σας ενδιαφέρει, κάντε κλικ στην επιλογή **Download (Λήψη)**.
4. Όταν η λήψη ολοκληρωθεί, κάντε κλικ στην επιλογή **Install (Εγκατάσταση)**.

Περιορισμένη εγγύηση προϊόντος

Η Life Technologies Corporation ή/και οι συνδεδεμένες με αυτήν εταιρείες παρέχουν εγγύηση για τα προϊόντα τους, όπως ορίζουν οι Γενικοί Όροι και Προϋποθέσεις Πώλησης της Life Technologies στον ιστότοπο www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Εάν έχετε ερωτήσεις, επικοινωνήστε με τη Life Technologies στον ιστότοπο www.thermofisher.com/support.

Oncomine™ Dx Express Test

N.º de pub. MAN0025897 Rev. D.0


Tanto las guías de usuario de este producto y las actualizaciones para el Genexus™ Dx Software como los archivos de definiciones de los ensayos (Assay Definition Files, ADF) para el Oncomine™ Dx Express Test se proporcionan en una biblioteca de recursos en Thermo Fisher™ Connect Platform.

Descargar la documentación del usuario

- Vaya a apps.thermofisher.com/apps/spa/#/publiclib/genexusdx e inicie sesión en thermofisher.com.
Nota: Si no tiene una cuenta de thermofisher.com, haga clic en **Create Account (Crear cuenta)** en la pantalla de inicio de sesión y siga las instrucciones.
- En la biblioteca de recursos, busque las guías de usuario del Oncomine™ Dx Express Test en el idioma que desee y, a continuación, descargue y extraiga la carpeta ZIP que contiene las guías de usuario.
- Abra las guías de usuario con un programa adecuado para archivos PDF, como Adobe Reader™ o equivalente.
Nota: Puede descargar Adobe Reader™ en [Adobe Reader](#).


Descargar el Genexus™ Dx Software y actualizaciones de ADF


Vincular una cuenta de Thermo Fisher™ Connect Platform con un código QR (director/administrador)
Para acceder a las actualizaciones de software, primero debe vincular un instrumento a una cuenta de Connect Platform. Solo es necesario vincular un instrumento una vez.


- En Genexus™ Dx Software, haga clic en  **(Settings [Ajustes]) ▶ Thermo Fisher Account (Cuenta de Thermo Fisher)**.
- Haga clic en **+ Link Account to Device (Vincular cuenta al dispositivo)**.
Se abre el cuadro de diálogo **Link Account to Device (Vincular cuenta al dispositivo)** y se muestra tanto la imagen como el texto del código QR.
- Use una de las dos siguientes opciones para vincular la cuenta de Connect Platform con el software.

Opciones para vincular una cuenta de Connect Platform

Opción	Descripción
Vincular con la aplicación móvil Instrument Connect	Nota: Para usar esta opción, debe tener la aplicación móvil InstrumentConnect instalada en un dispositivo móvil. <ol style="list-style-type: none"> Abra la aplicación InstrumentConnect e inicie sesión. En la aplicación, haga clic en +. Haga clic en QR Code (Código QR) y escanee la imagen del código QR que se muestra en el cuadro de diálogo Link Account to Device (Vincular cuenta al dispositivo) en Genexus™ Dx Software.

Opción	Descripción
Vincular a través del sitio web de Connect Platform	<ol style="list-style-type: none"> Copie el texto del código QR que se proporciona en el cuadro de diálogo y haga clic en Thermo Fisher Connect para abrir el sitio web de Connect Platform. Nota: Es necesario iniciar sesión. Si no tiene una cuenta de Thermo Fisher™ Connect Platform, haga clic en Create Account (Crear cuenta) en la pantalla de inicio de sesión y siga las instrucciones. En el sitio web de Connect Platform, haga clic en  InstrumentConnect en la barra lateral izquierda. Haga clic en Link instrument (Vincular instrumento) en la esquina superior derecha de la pantalla. Introduzca el texto del código QR que se muestra en el cuadro de diálogo de la pantalla Link Instrument (Vincular instrumento). Haga clic en Link (Vincular).

Descargar e instalar actualizaciones de ADF y software
Cuando la cuenta de Thermo Fisher™ Connect Platform está vinculada, las actualizaciones de ADF y software son visibles en la pantalla Genexus™ Dx Software  **(Settings [Ajustes]) ▶ Software Updates (Actualizaciones de software)**. Los usuarios de nivel director y administrador pueden descargar e instalar las actualizaciones.

- En Genexus™ Dx Software, haga clic en **▶  (Settings [Ajustes]) ▶ Software Updates (Actualizaciones de software)**.
- En la pantalla **Installed Software (Software instalado)**, haga clic en **Software Updates (Actualizaciones de software)** para ver las actualizaciones de software y ADF que están disponibles en Connect Platform. Si no está seleccionado, haga clic en el botón **APP STORE**.
- En la columna **Actions (Acciones)** del paquete de ADF o software de interés, haga clic en **Download (Descargar)**.
- Cuando termine la descarga, haga clic en **Install (Instalar)**.

Garantía limitada del producto

Life Technologies Corporation y/o sus filiales garantizan sus productos tal y como se establece en los términos y condiciones de venta de Life Technologies en www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Si tiene cualquier duda, póngase en contacto con Life Technologies en www.thermofisher.com/support.

Oncomine™ Dx Express Test

Väljaande nr MAN0025897 Red D.0

Selle toote kasutusjuhendid ja tarkvara Genexus™ Dx Software ning analüüside definitsiooni failide (ADF-id) värskendused analüüsi Oncomine™ Dx Express Test kohta leiate Thermo Fisher™ Connect Platformi ressursside pangast.

Kasutajadokumentatsiooni allalaadimine

1. Avage veebileht apps.thermofisher.com/apps/spa/#/publiclib/genexusdx ja seejärel logige lehel thermofisher.com sisse.

Märkus: Kui teil lehel thermofisher.com kontot ei ole, siis klõpsake sisselogimiskuvale nuppu **Create Account (Loo konto)** ja järgige juhiseid.


2. Leidke ressursside pangast üles sekveerija Oncomine™ Dx Express Test soovitud keeles kasutusjuhendid, laadige neid sisaldav ZIP-kaust alla ja seejärel pakkige lahti.
3. Avage kasutusjuhendid PDF-vaaturiga nagu Adobe Reader™ või muu sarnase vaaturiga.

Märkus: Vaatur Adobe Reader™ saab alla laadida veebilehelt [Adobe Reader](#).

Tarkvara Genexus™ Dx Software ja ADF-i värskenduste allalaadimine

Thermo Fisher™ Connect Platformi konto sidumine QR-koodiga (haldur/administraator)

Tarkvaravärskenduste installimiseks peate esmalt siduma instrumendi Connect Platformi kontoga. Instrumendi sidumine on ühekordne toiming.

1. Klõpsake tarkvaras Genexus™ Dx Software nuppu  **(Settings) (Sätted)** ▶ **Thermo Fisher Account (Thermo Fisheri konto)**.
2. Klõpsake nuppu **+ Link Account to Device (Konto sidumine seadmega)**.

Avaneb dialoogiaken **Link Account to Device (Konto sidumine seadmega)** ja kuvatakse QR-koodi pilt või QR-koodi tekst.


3. Siduge Connect Platformi konto tarkvaraga ühel järgmisest kahest viisist.


Konto Connect Platform sidumise võimalused

Suvand	Kirjeldus
Sidumine mobiilrakendusega InstrumentConnect	<p>Märkus: Selle võimaluse kasutamiseks peab teil olema mobiilseadmesse installitud mobiilirakendus InstrumentConnect.</p> <ol style="list-style-type: none"> 1. Avage rakendus InstrumentConnect ja logige sisse. 2. Klõpsake rakenduses nuppu +. 3. Klõpsake nuppu QR Code (QR-kood) ja seejärel skannige tarkvaras Genexus™ Dx Software dialoogiaknas Link Account to Device (Konto sidumine seadmega) kuvatav QR-kood.

Suvand	Kirjeldus
Sidumine veebilehe Connect Platform kaudu	<ol style="list-style-type: none"> 1. Kopeerige dialoogiaknas kuvatava QR-koodi tekst ja seejärel klõpsake nuppu Thermo Fisher Connect, et avada veebileht Connect Platform. Märkus: Vajalik on sisselogimine. Kui teil Thermo Fisher™ Connect Platformi kontot ei ole, siis klõpsake sisselogimiskuvale nuppu Create Account (Loo konto) ja järgige juhiseid. 2. Klõpsake Connect Platformi veebilehel vasakul külgribal kuvatavat nuppu InstrumentConnect. 3. Klõpsake kuva ülemises paremas nurgas nuppu Link instrument (Instrumendi sidumine). 4. Sisestage kuvale Link Instrument (Instrumendi sidumine) dialoogiaknas näidatud QR-koodi tekst. 5. Klõpsake nuppu Link (Seo).

Tarkvara ja ADF-i värskenduste allalaadimine ning installimine

Kui Thermo Fisher™ Connect Platformi konto on seotud, on saadaval tarkvaravärskendused ja ADF-i värskendused nähtavad tarkvara Genexus™ Dx Software kuvale  **(Settings) (Sätted)** ▶ **Software Updates (Tarkvaravärskendused)**. Halduri ja administraatori taseme kasutajad saavad värskendusi alla laadida ning installida.

1. Klõpsake tarkvaras Genexus™ Dx Software nuppu ▶  **(Settings) (Sätted)** ▶ **Software Updates (Tarkvaravärskendused)**.
2. Klõpsake kuvale **Installed Software (Installitud tarkvara)** nuppu **Software Updates (Tarkvaravärskendused)**, et näha tarkvara ja ADF-i värskendusi, mis on veebilehelt Connect Platform kättesaadavad. Kui see pole veel valitud, klõpsake nuppu **APP STORE (RAKENDUSTE POOD)**.
3. Klõpsake huvipakkuva tarkvara või ADF-i värskenduste paketi veerus **Actions (Toimingud)** nuppu **Download (Laadi alla)**.
4. Pärast allalaadimise lõppu klõpsake nuppu **Install (Installi)**.

Piiratud tootegarantii

Life Technologies Corporation ja/või selle tütarettevõtte/tütarettevõtte annavad oma toodetele garantii, nagu on kirjeldatud ettevõtte Life Technologies üldistes müügisätetes ja -tingimustes veebilehel www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Palun võtke küsimuste korral ühendust ettevõttega Life Technologies järgmisel saidil: www.thermofisher.com/support.

Oncomine™ Dx Express Test

Pub. n° MAN0025897 Rév. D.0

Les guides de l'utilisateur pour ce produit et les mises à jour du Genexus™ Dx Software et des fichiers de définition de dosage (ADF) pour l'Oncomine™ Dx Express Test sont fournis dans une bibliothèque de ressources Thermo Fisher™ Connect Platform.

Télécharger la documentation destinée à l'utilisateur

1. Rendez-vous sur apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, puis connectez-vous sur thermofisher.com.

Remarque : Si vous ne disposez pas d'un compte sur thermofisher.com, cliquez sur **Create Account (Créer un compte)** sur l'écran de connexion, puis suivez les instructions.


2. Dans la bibliothèque des ressources, recherchez les guides de l'utilisateur de l'Oncomine™ Dx Express Test dans la langue qui vous intéresse, puis téléchargez et extrayez le dossier ZIP contenant les guides de l'utilisateur.
3. Ouvrez les guides de l'utilisateur avec un lecteur de fichiers PDF tel qu'Adobe Reader™ ou un équivalent.

Remarque : Adobe Reader™ peut être téléchargé sur [Adobe Reader](https://www.adobe.com/fr/reader).

Télécharger le Genexus™ Dx Software et les mises à jour (ADF)

Lier un compte Thermo Fisher™ Connect Platform avec un code QR (responsable / administrateur)


Pour accéder aux mises à jour du logiciel, vous devez d'abord lier un instrument à un compte Connect Platform. La liaison d'un instrument est une action ponctuelle.

1. Dans Genexus™ Dx Software, cliquez sur  **(Settings) (Paramètres) ▶ Thermo Fisher Account (Compte Thermo Fisher)**.
2. Cliquez sur **+ Link Account to Device (Lier un compte à un appareil)**.

La boîte de dialogue **Link Account to Device (Lier un compte à un appareil)** s'ouvre et une image de code QR et du texte de code QR s'affichent.


3. Liez le compte Connect Platform avec le logiciel à l'aide d'une des deux options suivantes.

Options pour lier un compte Connect Platform

Option	Description
Liaison avec l'application mobile Instrument Connect	<p>Remarque : Pour utiliser cette option, l'application mobile InstrumentConnect doit être installée sur un de vos appareils mobiles.</p> <ol style="list-style-type: none"> 1. Ouvrez l'application InstrumentConnect, puis connectez-vous. 2. Dans l'application, cliquez sur +. 3. Cliquez sur QR Code (Code QR), puis scannez le code QR indiqué dans la boîte de dialogue Link Account to Device (Lier un compte à un appareil) dans Genexus™ Dx Software.
Liaison par le biais du site Web de Connect Platform	<ol style="list-style-type: none"> 1. Copiez le texte du code QR dans la boîte de dialogue, puis cliquez sur Thermo Fisher Connect pour ouvrir le site Web de Connect Platform. <p>Remarque : L'authentification est requise. Si vous ne disposez pas d'un compte Thermo Fisher™ Connect Platform, cliquez sur Create Account (Créer un compte) sur l'écran de connexion, puis suivez les instructions.</p> 2. Sur le site Web Connect Platform, cliquez sur  InstrumentConnect dans la barre latérale de gauche. 3. Cliquez sur Link instrument (Lier un instrument) dans le coin supérieur droit de l'écran. 4. Saisissez le texte du code QR affiché dans la boîte de dialogue sur l'écran Link Instrument (Lier un instrument). 5. Cliquez sur Link (Lier).

Télécharger et installer les mises à jour du logiciel et des ADF

Lorsque le compte Thermo Fisher™ Connect Platform est lié, les mises à jour du logiciel et des ADF disponibles sont visibles sur l'écran Genexus™ Dx Software  **(Settings) (Paramètres) ▶ Software Updates (Mises à jour logicielles)**. Les utilisateurs de niveau responsable et administrateur peuvent télécharger et installer les mises à jour.

1. Dans Genexus™ Dx Software, cliquez sur **▶  (Settings) (Paramètres) ▶ Software Update (Mise à jour du logiciel)**.
2. Sur l'écran **Installed Software (Logiciel installé)**, cliquez sur **Software Updates (Mises à jour du logiciel)** pour voir les mises à jour du logiciel et des ADF qui sont disponibles dans la Connect Platform. S'il n'est pas sélectionné, cliquez sur le bouton **APP STORE**.
3. Dans la colonne **Actions** pour le pack logiciel ou ADF d'intérêt, cliquez sur **Download (Télécharger)**.
4. Lorsque le téléchargement est terminé, cliquez sur **Install (Installer)**.

Garantie limitée du produit

Life Technologies Corporation et ses filiales garantissent leurs produits selon les termes et conditions générales de ventes disponibles sur le site www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Si vous avez des questions, vous pouvez prendre contact avec Life Technologies à l'adresse web suivante : www.thermofisher.com/support.

Oncomine™ Dx Express Test

Br. pub. MAN0025897 Rev. D.0


Korisnički priručnici za ovaj proizvod i ažuriranja softvera Genexus™ Dx Software te datoteke s definicijama testova (ADF-ovi) za Oncomine™ Dx Express Test nalaze se u biblioteci resursa platforme Thermo Fisher™ Connect Platform.

Preuzimanje korisničke dokumentacije

- Idite na apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, a zatim se prijavite na thermofisher.com.
Napomena: Ako nemate račun na stranici thermofisher.com, kliknite **Create Account (Izradi račun)** na zaslonu za prijavu i slijedite upute.
- U biblioteci resursa pronađite korisničke priručnike za Oncomine™ Dx Express Test na željenom jeziku, a zatim preuzmite i izdvojite datoteke ZIP mape u kojoj se nalaze korisnički priručnici.
- Otvorite korisničke priručnike u PDF pregledniku kao što je Adobe Reader™ ili sličan preglednik.
Napomena: Adobe Reader™ možete preuzeti na [Adobe Reader](#).


Preuzimanje softvera Genexus™ Dx Software i ažuriranja ADF-ova


Povezivanje računa za platformu Thermo Fisher™ Connect Platform s pomoću QR koda (upravitelj/administrator)
Da biste pristupili ažuriranjima softvera, najprije morate povezati instrument s računom za platformu Connect Platform. Povezivanje instrumenta obavlja se samo jednom.


- U softveru Genexus™ Dx Software kliknite  **(Settings) (Postavke) ▶ Thermo Fisher Account (Račun Thermo Fisher)**.
- Kliknite **+ Link Account to Device (Poveži račun s uređajem)**.
Otvora se dijaloški okvir **Link Account to Device (Poveži račun s uređajem)** te se prikazuju slika QR koda i tekst QR koda.
- Povežite račun za platformu Connect Platform sa softverom koristeći se jednom od dviju opcija navedenih u nastavku.

Opcije za povezivanje računa za platformu Connect Platform

Opcija	Opis
Povezivanje s mobilnom aplikacijom InstrumentConnect	Napomena: Da biste mogli upotrijebiti ovu opciju, morate imati instaliranu mobilnu aplikaciju InstrumentConnect na mobilnom uređaju. <ol style="list-style-type: none"> Otvorite aplikaciju InstrumentConnect, a zatim se prijavite. U aplikaciji kliknite +. Kliknite QR Code (QR kod), a zatim skenirajte sliku QR koda koji se prikazuje u dijaloškom okviru Link Account to Device (Poveži račun s uređajem) u softveru Genexus™ Dx Software.

Opcija	Opis
Povezivanje putem web-mjesta Connect Platform	<ol style="list-style-type: none"> Kopirajte tekst QR koda naveden u dijaloškom okviru, a zatim kliknite Thermo Fisher Connect da biste otvorili web-mjesto Connect Platform. Napomena: Potrebna je prijava. Ako nemate račun za platformu Thermo Fisher™ Connect Platform, kliknite Create Account (Izradi račun) na zaslonu za prijavu, a zatim slijedite upute. Na web-mjestu Connect Platform kliknite  InstrumentConnect u lijevoj bočnoj traci. Kliknite Link instrument (Poveži instrument) u gornjem desnom kutu zaslona. Unesite tekst QR koda koji se prikazuje u dijaloškom okviru na zaslonu Link Instrument (Poveži instrument). Kliknite Link (Poveži).

Preuzimanje i instaliranje ažuriranja softvera i ADF-ova
Kada povežete račun za Thermo Fisher™ Connect Platform, dostupna ažuriranja softvera i ADF-ova vidljiva su na zaslonu softvera Genexus™ Dx Software  **(Settings) (Postavke) ▶ Software Updates (Ažuriranja softvera)**. Korisnici na razini upravitelja i administratora mogu preuzimati i instalirati ažuriranja.

- U softveru Genexus™ Dx Software kliknite **▶  (Settings) (Postavke) ▶ Software Updates (Ažuriranja softvera)**.
- Na zaslonu **Installed Software (Instalirani softver)** kliknite **Software Updates (Ažuriranja softvera)** da biste prikazali ažuriranja softvera i ADF-ova koja su dostupna na platformi Connect Platform. Ako opcija nije odabrana, kliknite gumb **APP STORE (Trgovina aplikacijama)**.
- U stupcu **Actions (Radnje)** za željeni softverski paket ili paket ADF-ova kliknite **Download (Preuzmi)**.
- Kada preuzimanje završi, kliknite **Install (Instaliraj)**.

Ograničeno jamstvo proizvoda

Društvo Life Technologies Corporation i/ili njegova pridružena društva jamče za svoje proizvode kako je navedeno u Općim uvjetima prodaje društva Life Technologies na stranici www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Ako imate bilo kakvih pitanja, obratite se društvu Life Technologies na www.thermofisher.com/support.

Oncomine™ Dx Express Test

Kiadv.szám MAN0025897 D.0 átdolg.

A termék felhasználói útmutatói, valamint a Genexus™ Dx Software és az Oncomine™ Dx Express Test assay definíciós fájlok (ADF-ek) a Thermo Fisher™ Connect Platform egyik forráskönyvtárában található.

A felhasználói dokumentumok letöltése

- Lépjen a apps.thermofisher.com/apps/spa/#/publiclib/genexusdx weboldalra, majd jelentkezzen be a thermofisher.com fiókba.

Megjegyzés: Ha nem rendelkezik thermofisher.com fiókkal, kattintson a bejelentkező képernyőn a **Create Account (Fiók létrehozása)** lehetőségre, és kövesse az utasításokat.


- A forráskönyvtárban keresse meg a Oncomine™ Dx Express Test felhasználói útmutatóját a megfelelő nyelven, majd tölts le és csomagolja ki a felhasználói útmutatókat tartalmazó ZIP mappát.
- Nyissa meg a felhasználói útmutatókat egy PDF olvasóval, például az Adobe Reader™ vagy más, ennek megfelelő programmal.

Megjegyzés: Az Adobe Reader™ letölthető az alábbi helyről: [Adobe Reader](#).

A Genexus™ Dx Software és az ADF frissítéseinek letöltése

A Thermo Fisher™ Connect Platform fiók összekapcsolása QR-kód segítségével (menedzser/rendszergazda)

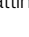
A szoftverfrissítések eléréséhez először egy készüléket kell hozzákapcsolnia a Connect Platform fiókhoz. A készülék összekapcsolását csak egyszer kell elvégezni.

- A Genexus™ Dx Software szoftverben kattintson a  **(Settings [Beállítások]) ▶ Thermo Fisher Account (Thermo Fisher felhasználói fiók)** lehetőségre.
- Kattintson a **+ Link Account to Device (Fiók összekapcsolása eszközzel)** lehetőségre.


Megnyílik a **Link Account to Device (Fiók összekapcsolása eszközzel)** párbeszédablak, amelyen megjelenik egy QR-kód képe és a QR-kód szövege.

- Az alábbi két lehetőség egyikével kapcsolja össze a Connect Platform fiókot a szoftverrel.

A Connect Platform fiók összekapcsolásának lehetőségei

Opció	Leírás
Összekapcsolás az Instrument-Connect mobil alkalmazással	<p>Megjegyzés: Ezen opció használatához telepítenie kell az InstrumentConnect mobil alkalmazást egy mobil eszközre.</p> <ol style="list-style-type: none"> Nyissa meg az InstrumentConnect alkalmazást, és jelentkezzen be. Az alkalmazásban kattintson a + jelre. Kattintson a QR Code (QR-kód) lehetőségre, majd olvassa be a Genexus™ Dx Software Link Account to Device (Fiók összekapcsolása eszközzel) párbeszédablakában látható QR-kód képét.
Összekapcsolás a Connect Platform weboldalon keresztül	<ol style="list-style-type: none"> Másolja a QR-kód megadott szövegét a párbeszédablakba, majd kattintson a Thermo Fisher Connect (Thermo Fisher kapcsolódás) lehetőségre a Connect Platform weboldalon megnyitásához. <p>Megjegyzés: Bejelentkezés szükséges. Ha nem rendelkezik Thermo Fisher™ Connect Platform fiókkal, kattintson a bejelentkező képernyőn a Create Account (Fiók létrehozása) lehetőségre, és kövesse az utasításokat.</p> <ol style="list-style-type: none"> A Connect Platform weboldalon kattintson a bal oldalsáv  InstrumentConnect elemére. Kattintson a képernyő jobb felső sarkában található Link instrument (Készülék összekapcsolása) gombra. Írja be a QR-kód szövegét a Link Instrument (Készülék összekapcsolása) képernyőn megjelenő párbeszédablakba. Kattintson a Link (Összekapcsolás) gombra.

A szoftverek és az ADF frissítéseinek letöltése és telepítése

Amikor megtörténik a Thermo Fisher™ Connect Platform fiókjának összekapcsolása, láthatóvá válnak a szoftverek és az ADF frissítései a Genexus™ Dx Software  **(Settings [Beállítások]) ▶ Software Updates (Szoftverfrissítések)** képernyőn. A menedzser és rendszergazda szintű felhasználók letölthetik és telepíthetik a frissítéseket.

- A Genexus™ Dx Software szoftverben kattintson a **▶  (Settings [Beállítások]) ▶ Software Updates (Szoftverfrissítések)** gombra.
- Az **Installed Software (Telepített szoftver)** képernyőn kattintson a **Software Updates (Szoftverfrissítések)** lehetőségre, hogy megtekintse a szoftver és az ADF-ek Connect Platform platformon elérhető frissítéseit. Ha nincs kiválasztva, kattintson az **APP STORE (Alkalmazásbolt)** gombra.
- Az adott szoftver- vagy ADF-csomaghoz tartozó **Actions (Tevékenységek)** oszlopban kattintson a **Download (Letöltés)** gombra.
- A letöltés befejeződésekor kattintson az **Install (Telepítés)** lehetőségre.

Korlátozott termékszavatosság

A Life Technologies Corporation és/vagy leányvállalata(i) a www.thermofisher.com/us/en/home/global/

[terms-and-conditions.html](#) weboldalon található, „A Life Technologies általános értékesítési feltételei” című dokumentumban meghatározott módon vállalnak garanciát termékeikre. Ha bármilyen kérdése van, vegye fel a kapcsolatot a Life Technologies vállalattal a www.thermofisher.com/support webcímen.

Oncomine™ Dx Express Test

No. Pub. MAN0025897 Rev. D.0

Panduan pengguna untuk produk ini dan pembaruan untuk Genexus™ Dx Software dan File Definisi Uji Kadar (ADF) untuk Oncomine™ Dx Express Test diberikan dalam pustaka sumber daya Thermo Fisher™ Connect Platform.


Mengunduh dokumentasi pengguna

- Buka apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, lalu masuk ke thermofisher.com.
Catatan: Jika Anda tidak memiliki akun thermofisher.com, klik **Create Account (Buat Akun)** di layar masuk, lalu ikuti petunjuknya.
- Di pustaka sumber daya, buka panduan pengguna Oncomine™ Dx Express Test dalam bahasa yang diinginkan, kemudian unduh dan ekstrak folder ZIP yang berisi panduan pengguna.
- Buka panduan pengguna dengan penampil PDF seperti Adobe Reader™ atau yang setara.
Catatan: Adobe Reader™ dapat diunduh di [Adobe Reader](#).

Mengunduh Genexus™ Dx Software dan pembaruan ADF

Menautkan akun Thermo Fisher™ Connect Platform dengan kode QR (manajer/administrator)

Untuk mengakses pembaruan perangkat lunak, Anda harus terlebih dahulu menautkan instrumen ke akun Connect Platform. Menautkan instrumen adalah tindakan satu kali.

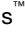
- Di Genexus™ Dx Software, klik  **(Settings) (Pengaturan) ▶ Thermo Fisher Account (Akun Thermo Fisher)**.
- Klik **+ Link Account to Device (Tautkan Akun ke Perangkat)**.
Kotak dialog **Link Account to Device (Tautkan Akun ke Perangkat)** terbuka dan gambar kode QR dan teks kode QR ditampilkan.
- Tautkan akun Connect Platform dengan perangkat lunak menggunakan salah satu dari dua opsi berikut.


Opsi untuk menautkan akun Connect Platform

Opsi	Deskripsi
Tautkan dengan aplikasi seluler InstrumentConnect	Catatan: Untuk menggunakan opsi ini, Anda harus menginstal aplikasi seluler InstrumentConnect di perangkat seluler. <ol style="list-style-type: none"> Buka aplikasi InstrumentConnect, lalu masuk. Di aplikasi, klik +. Klik QR Code (Kode QR), lalu pindai gambar kode QR yang ditampilkan di kotak dialog Link Account to Device (Tautkan Akun ke Perangkat) di Genexus™ Dx Software.

Opsi	Deskripsi
Tautkan melalui situs web Connect Platform	<ol style="list-style-type: none"> Salin teks kode QR yang diberikan di kotak dialog, lalu klik Thermo Fisher Connect (Koneksi Thermo Fisher) untuk membuka situs web Connect Platform. Catatan: Anda perlu melakukan sign-in. Jika Anda tidak memiliki akun Thermo Fisher™ Connect Platform, klik Create Account (Buat Akun) di layar sign in, lalu ikuti panduannya. Di situs web Connect Platform, klik  InstrumentConnect di bilah sisi kiri. Klik Link instrument (Tautkan Instrumen) di sudut kanan atas layar. Masukkan teks kode QR yang ditampilkan di kotak dialog di layar Link Instrument (Tautkan Instrumen). Klik Link (Tautkan).

Mengunduh dan memasang pembaruan perangkat lunak dan ADF

Saat akun Thermo Fisher™ Connect Platform ditautkan, perangkat lunak yang tersedia dan pembaruan ADF dapat dilihat di layar Genexus™ Dx Software  **(Settings) (Pengaturan) ▶ Software Updates (Pembaruan Perangkat Lunak)**. Pengguna tingkat manajer dan administrator dapat mengunduh dan menginstal pembaruan.

- Di Genexus™ Dx Software, klik **▶  (Settings) (Pengaturan) ▶ Software Updates (Pembaruan Perangkat Lunak)**.
- Di layar **Installed Software (Perangkat Lunak yang Dipasang)**, klik **Software Updates (Pembaruan Perangkat Lunak)** untuk melihat pembaruan perangkat lunak dan ADF yang tersedia di Connect Platform. Jika tidak dipilih, klik tombol **APP STORE**.
- Di kolom **Actions (Tindakan)** untuk perangkat lunak atau paket ADF yang diinginkan, klik **Download (Unduh)**.
- Setelah unduhan selesai, klik **Install (Instal)**.

Garansi produk terbatas

Life Technologies Corporation dan/atau afliasinya menjamin produk mereka sebagaimana diatur dalam Persyaratan dan Ketentuan Umum Penjualan Life Technologies di www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Jika ada pertanyaan, silakan hubungi Life Technologies di www.thermofisher.com/support.

Oncomine™ Dx Express Test

N. di pub. MAN0025897 Rev. D.0

Le guide dell'utente per questo prodotto e gli aggiornamenti del Genexus™ Dx Software e dei file di definizione del saggio (ADF) per Oncomine™ Dx Express Test sono disponibili in una libreria risorse su Thermo Fisher™ Connect Platform.



Scaricare la documentazione dell'utente

1. Accedere a apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, quindi eseguire l'accesso a thermofisher.com.
Nota: Se non si possiede un account thermofisher.com, nella schermata di accesso fare clic su **Create Account (Crea account)** e seguire le istruzioni.
2. Nella libreria risorse, individuare le guide dell'utente Oncomine™ Dx Express Test nella lingua desiderata, quindi scaricare ed estrarre la cartella ZIP contenente le guide dell'utente.
3. Aprire le guide dell'utente con un visualizzatore di PDF come Adobe Reader™ o equivalente.
Nota: Adobe Reader™ può essere scaricato da [Adobe Reader](#).



Download del Genexus™ Dx Software e aggiornamenti ADF

Collegare un account Thermo Fisher™ Connect Platform con un codice QR (manager/amministratore)

Per poter accedere agli aggiornamenti software, occorre prima collegare uno strumento a un account Connect Platform. Il collegamento di uno strumento è un'azione una tantum.

1. Nel Genexus™ Dx Software, fare clic su  (**Settings [Impostazioni]**) ▶ **Thermo Fisher Account**.
2. Fare clic su  **Link Account to Device (Collega account al dispositivo)**.
La finestra di dialogo **Link Account to Device (Collega account al dispositivo)** si apre e vengono mostrati un'immagine del codice QR e un testo del codice QR.
3. Collegare l'account Connect Platform al software utilizzando una delle due opzioni che seguono.


Opzioni per il collegamento a un account Connect Platform

Opzione	Descrizione
Collegamento tramite l'app mobile InstrumentConnect	<p>Nota: Per utilizzare questa opzione è necessario aver installato l'app mobile InstrumentConnect su un dispositivo mobile.</p> <ol style="list-style-type: none"> 1. Aprire l'applicazione InstrumentConnect ed eseguire l'accesso. 2. Nell'applicazione, fare clic su . 3. Fare clic su QR Code (Codice QR), quindi scansionare il codice QR mostrato nella finestra di dialogo Link Account to Device (Collega account al dispositivo) in Genexus™ Dx Software.
Collegamento tramite il sito Web Connect Platform	<ol style="list-style-type: none"> 1. Copiare il testo del codice QR visualizzato nella finestra di dialogo, quindi fare clic su Thermo Fisher Connect per aprire il sito Web Connect Platform. Nota: È richiesta la registrazione. Se non si possiede un account Thermo Fisher™ Connect Platform, fare clic su Create Account (Crea account) nella schermata di registrazione e seguire le istruzioni. 2. Nel sito web Connect Platform, fare clic su  InstrumentConnect nella barra laterale sinistra. 3. Nell'angolo in alto a destra della schermata, fare clic su Link instrument (Collega strumento). 4. Inserire il testo del codice QR mostrato nella finestra di dialogo Link Instrument (Collega strumento). 5. Fare clic su Link (Collega).

Download e installazione del software e degli aggiornamenti ADF

Una volta collegato l'account Thermo Fisher™ Connect Platform, gli aggiornamenti di software e ADF sono visualizzati nella schermata Genexus™ Dx Software  (**Settings (Impostazioni)**) ▶ **Software Updates (Aggiornamenti software)**.

Gli utenti con privilegi manager e amministratore possono scaricare e installare gli aggiornamenti.

1. In Genexus™ Dx Software, fare clic su  (**Settings (Impostazioni)**) ▶ **Software Updates (Aggiornamenti software)**.
2. Nella schermata **Installed Software (Software installato)**, fare clic su **Software Updates (Aggiornamenti software)** per visualizzare gli aggiornamenti di software e ADF disponibili nella Connect Platform. Se non è selezionato, fare clic sul pulsante **APP STORE**.
3. Nella colonna **Actions (Azioni)** per il pacchetto software o ADF desiderato, quindi fare clic su **Download (Scarica)**.
4. Una volta completato il download, fare clic su **Install (Installa)**.

Garanzia limitata del prodotto

Life Technologies Corporation e/o le sue affiliate garantiscono i loro prodotti come indicato nel documento Termini generali e condizioni di vendita di Life Technologies disponibile all'indirizzo www.thermofisher.com/us/en/home/global/terms-and-conditions.html. In caso di domande, contattare Life Technologies all'indirizzo www.thermofisher.com/support.

Oncomine™ Dx Express Test

出版物番号 MAN0025897 改訂版 D.0

本製品のユーザーガイドおよび Genexus™ Dx Software の更新と Oncomine™ Dx Express Test のアッセイ定義ファイル (ADF) は Thermo Fisher™ Connect Platform リソースライブラリで提供されています。


ユーザー文書のダウンロード

1. apps.thermofisher.com/apps/spa/#/publiclib/genexusdx に移動し、thermofisher.com にサインインします。
注記: thermofisher.com アカウントがない場合には、サインイン画面で **Create Account** (アカウントの作成) をクリックし、指示に従ってください。
2. リソースライブラリで、対象とする言語の Oncomine™ Dx Express Test ユーザーガイドを探し、そのユーザーガイドを含む ZIP フォルダをダウンロードして解凍します。
3. Adobe Reader™などの PDF ビューアーでユーザーガイドを開きます。
注記: Adobe Reader™は、[Adobe Reader](#) でダウンロードできます。

Genexus™ Dx Software および ADF アップデートのダウンロード


QR コードを使った Thermo Fisher™ Connect Platform アカウントのリンク (マネージャー/管理者)

ソフトウェア更新にアクセスするには、まず装置を Connect Platform アカウントにリンクする必要があります。装置へのリンクは 1 回だけの操作です。

1. Genexus™ Dx Software で、 (Settings) (設定) ▶ Thermo Fisher Account (Thermo Fisher アカウント) をクリックします。
2. **+** Link Account to Device (アカウントをデバイスにリンク) をクリックします。
Link Account to Device (アカウントをデバイスにリンク) ダイアログボックスが開き、QR コードの画像と QR コードテキストが表示されます。
3. 以下の 2 つのオプションのいずれかを使用して、Connect Platform アカウントをソフトウェアとリンクします。

Connect Platform アカウントをリンクするオプション

オプション	説明
Instrument Connect モバイルアプリとリンクする	注記: このオプションを使用するには、モバイルデバイスに InstrumentConnect モバイルアプリがインストールされている必要があります。 1. InstrumentConnect アプリを開いてサインインします。 2. アプリで、 + をクリックします。 3. QR Code (QR コード) をクリックし、Link Account to Device (デバイスにアカウントをリンク) ダイアログボックス (Genexus™ Dx Software 内) に表示されている QR コードの画像をスキャンします。

オプション	説明
Connect Platform ウェブサイトからリンクする	1. ダイアログボックスに表示される QR コードテキストをコピーし、Thermo Fisher Connect (Thermo Fisher Connect) をクリックして Connect Platform ウェブサイトを開きます。 注記: サインインが必要です。Thermo Fisher™ Connect Platform アカウントがない場合には、サインイン画面で Create Account (アカウントの作成) をクリックし、指示に従ってください。 2. Connect Platform の Web サイトで、左側のサイドバーにある  InstrumentConnect (InstrumentConnect) をクリックします。 3. 画面の右上隅にある Link instrument (機器のリンク) をクリックします。 4. Link Instrument (装置のリンク) 画面のダイアログボックスに QR コードのテキストを入力します。 5. Link (リンク) をクリックします。

ソフトウェアおよび ADF 更新のダウンロードとインストール

Thermo Fisher™ Connect Platform アカウントがリンクされると、利用可能なソフトウェアと ADF 更新が Genexus™ Dx Software  (Settings) (設定) ▶ Software Updates (ソフトウェアの更新) 画面に表示されます。マネージャーレベルおよび管理者レベルのユーザーは、更新をダウンロードしてインストールできます。

1. Genexus™ Dx Software で  (Settings) (設定) ▶ Software Updates (ソフトウェア更新) をクリックします。
2. Installed Software (インストールされるソフトウェア) 画面で、Software Updates (ソフトウェア更新) をクリックして Connect Platform で利用可能なソフトウェアと ADF の更新を表示します。選択されていない場合には、APP STORE ボタンをクリックします。
3. ダウンロードしたいソフトウェア/ADF パッケージの Actions (アクション) 列で **Download** (ダウンロード) をクリックします。
4. ダウンロードが完了したら、**Install** (インストール) をクリックします。

限定製品保証

Life Technologies Corporation またはその関連会社は、ウェブサイト (www.thermofisher.com/us/en/home/global/terms-and-conditions.html) にある Life Technologies の General Terms and Conditions of Sale (販売に関する一般条件) の定めにしたがって製品を保証します。ご質問のある方は、www.thermofisher.com/support で Life Technologies までお問い合わせください。



体外診断用。

ThermoFisher
SCIENTIFIC

Oncomine™ Dx Express Test

공개 번호 MAN0025897 개정 D.0

이 제품에 대한 사용자 가이드 및 Genexus™ Dx Software 업데이트와 Oncomine™ Dx Express Test에 대한 분석 정의 파일(ADF)은 Thermo Fisher™ Connect Platform 리소스 라이브러리에 제공됩니다.

사용자 문서 다운로드

1. apps.thermofisher.com/apps/spa/#/publiclib/genexusdx 로 이동한 후 thermofisher.com에 로그인하십시오.
참고: thermofisher.com 계정이 없는 경우, 로그인 화면에서 **Create Account(계정 생성)**를 클릭한 후 지침에 따르십시오.
2. 리소스 라이브러리에서 원하는 언어의 Oncomine™ Dx Express Test 사용 설명서를 찾은 후 사용 설명서를 포함하는 ZIP 폴더를 다운로드 및 추출하십시오.
3. Adobe Reader™ 또는 동등한 뷰어 등 PDF 뷰어로 사용 설명서를 여십시오.
참고: Adobe Reader™는 [Adobe Reader](#)에서 다운로드할 수 있습니다.

Genexus™ Dx Software 및 ADF 업데이트 다운로드

QR 코드로 Thermo Fisher™ Connect Platform 계정 연결 (매니저/관리자)

소프트웨어 업데이트에 액세스하려면 먼저 Connect Platform 계정에 기기를 연결해야 합니다. 한 번의 작업으로 기기를 연결할 수 있습니다.

1. Genexus™ Dx Software에서 **(Settings)((설정)) ▶ Thermo Fisher Account(Thermo Fisher 계정)**를 클릭합니다.
2. **+ Link Account to Device(장치에 계정 연결)**를 클릭합니다.
Link Account to Device(장치에 계정 연결) 대화 상자가 열리고 QR 코드 이미지와 QR 코드 텍스트가 나타납니다.
3. 다음 두 가지 옵션 중 하나를 사용하여 소프트웨어로 Connect Platform 계정을 연결하십시오.

Connect Platform 계정 연결 옵션

옵션	설명
Instrument Connect 모바일 앱 연결	참고: 이 옵션을 사용하려면 모바일 장치에 InstrumentConnect 모바일 앱이 설치되어 있어야 합니다. 1. InstrumentConnect 앱을 연 후 로그인합니다. 2. 앱에서 + 를 클릭합니다. 3. QR Code(QR 코드) 를 클릭한 다음 Genexus™ Dx Software의 Link Account to Device(장치에 계정 연결) 대화 상자에 표시되는 QR 코드 이미지를 스캔합니다.

옵션	설명
Connect Platform 웹사이트를 통해 연결	1. 대화 상자에 제공된 QR 코드 텍스트를 복사한 다음 Thermo Fisher Connect 를 클릭하여 Connect Platform 웹사이트를 엽니다. 참고: 로그인이 필요합니다. Thermo Fisher™ Connect Platform 계정이 없는 경우 로그인 화면에서 Create Account(계정 생성) 를 클릭한 다음 지침을 따르십시오. 2. Connect Platform 웹사이트에서 왼쪽 사이드바의 InstrumentConnect 를 클릭합니다. 3. 화면의 오른쪽 상단 모서리에 있는 Link instrument(기기 연결) 를 클릭합니다. 4. Link Instrument(기기 연결) 화면의 대화 상자에 표시되는 QR 코드 텍스트를 입력합니다. 5. Link(연결) 를 클릭합니다.

소프트웨어 및 ADF 업데이트 다운로드 및 설치

Thermo Fisher™ Connect Platform 계정이 연결되면 이용 가능한 소프트웨어 및 ADF 업데이트가 Genexus™ Dx Software **(Settings)((설정)) ▶ Software Updates(소프트웨어 업데이트)** 화면에 표시됩니다. 매니저 및 관리자 수준 사용자는 업데이트를 다운로드하여 설치할 수 있습니다.

1. Genexus™ Dx Software에서 **(Settings)((설정)) ▶ Software Updates(소프트웨어 업데이트)**를 클릭합니다.
2. **Installed Software(설치된 소프트웨어)** 화면에서 **Software Updates(소프트웨어 업데이트)**를 클릭하여 Connect Platform에서 이용 가능한 소프트웨어 및 ADF 업데이트를 보십시오. 선택하지 않은 경우, **APP STORE(앱 스토어)** 버튼을 클릭합니다.
3. 원하는 소프트웨어 또는 ADF 패키지에 대한 **Actions(작업)** 열에서 **Download(다운로드)**를 클릭합니다.
4. 다운로드가 완료된 후 **Install(설치)**를 클릭합니다.

제한적인 제품 보증

Life Technologies Corporation 및/또는 그 계열사는 www.thermofisher.com/us/en/home/global/terms-and-conditions.html의 Life Technologies의 일반 판매 약관에 명시된 대로 제품을 보증합니다. 질문 사항이 있을 경우, www.thermofisher.com/support를 통해 Life Technologies에 연락해 주십시오.



체외 진단용.

Oncomine™ Dx Express Test

Leid. nr. MAN0025897 D.O perž.

Šio gaminio naudotojo vadovai ir Genexus™ Dx Software programinės įrangos naujiniai bei Oncomine™ Dx Express Test „Assay Definition Files“ (Analizės apibrėžties failai, ADF) pateikiami Thermo Fisher™ Connect Platform išteklių bibliotekoje.

Atsisiųskite naudotojo dokumentaciją

- Eikite į apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, tada prisijunkite prie thermofisher.com.
Pastaba: Jei neturite thermofisher.com paskyros, prisijungimo ekrane spustelėkite **Create Account (Sukurti paskyrą)** ir vykdykite nurodymus.
- Išteklių bibliotekoje raskite Oncomine™ Dx Express Test naudotojo vadovus dominančia kalba, tada atsisiųskite ir išskleiskite ZIP aplanką su naudotojo vadovais.
- Atidarykite naudotojo vadovus naudodami PDF naršyklę, pavyzdžiui, Adobe Reader™ arba lygiavertę programą.
Pastaba: Adobe Reader™ galite atsisiųsti adresu [Adobe Reader](#).

Genexus™ Dx Software ir ADF naujinių atsiuntimas

Susiekite Thermo Fisher™ Connect Platform paskyrą su QR kodu (vadovas / administratorius)

Norėdami naudotis programinės įrangos atnaujinimais, pirmiausia turite susieti prietaisą su Connect Platform paskyra. Prietaiso susiejimas yra vienkartinis veiksmas.

- Genexus™ Dx Software programinėje įrangoje spustelėkite **(Settings) (Nustatymai) ▶ Thermo Fisher Account („Thermo Fisher“ paskyra)**.
- Spustelėkite **+ Link Account to Device (Susieti paskyrą su prietaisu)**.
Atveriamas dialogo langas **Link Account to Device (Susieti paskyrą su prietaisu)** ir rodomas QR kodo paveikslėlis bei QR kodo tekstas.
- Susiekite Connect Platform paskyrą su programine įranga naudodami vieną iš dviejų toliau nurodytų parinkčių.

Connect Platform paskyros susiejimo parinktys

Parinktis	Aprašymas
Sąsaja su „Instrument Connect“ mobiliąja programėle	Pastaba: Kad galėtumėte naudotis šia parinktimi, mobiliajame įrenginyje turi būti įdiegta „InstrumentConnect“ mobilioji programėlė. <ol style="list-style-type: none"> Atidarykite „InstrumentConnect“ programėlę, tada prisijunkite. Programėlėje spustelėkite +. Spustelėkite QR Code (QR kodas), tada nuskaitykite QR kodo vaizdą, rodomą Genexus™ Dx Software dialogo lange Link Account to Device (Susieti paskyrą su prietaisu).

Parinktis	Aprašymas
Sąsaja per Connect Platform svetainę	<ol style="list-style-type: none"> Nukopijuokite dialogo lange pateiktą QR kodo tekstą, tada spustelėkite Thermo Fisher Connect, kad atidarytumėte Connect Platform svetainę. Pastaba: Būtina prisijungti. Jei neturite Thermo Fisher™ Connect Platform paskyros, prisijungimo ekrane spustelėkite Create Account (Sukurti paskyrą) ir vykdykite nurodymus. Svetainėje Connect Platform kairėje šoninėje juostoje spustelėkite InstrumentConnect. Viršutiniame dešiniajame ekrano kampe spustelėkite Link instrument (Susieti prietaisą). Dialogo lange įveskite QR kodo tekstą, rodomą Link instrument (Susieti prietaisą) ekrane. Spustelėkite Link (Susieti).

Atsisiųskite ir įdiekite programinės įrangos ir ADF naujinius Kai Thermo Fisher™ Connect Platform paskyra susieta, galimi programinės įrangos ir ADF atnaujinimai matomi Genexus™ Dx Software programinės įrangos **(Settings) (Nustatymai) ▶ Software Updates (Programinės įrangos atnaujinimai)** (Nustatymai / Programinės įrangos atnaujinimai) ekrane. Vadovo ir administratoriaus lygmens naudotojai gali atsisiųsti ir įdiegti naujinius.

- Genexus™ Dx Software programinėje įrangoje spustelėkite **(Settings) (Nustatymai) ▶ Software Updates (Programinės įrangos atnaujinimai)**.
- Ekrane **Installed Software (Įdiegta programinė įranga)** spustelėkite **Software Updates (Programinės įrangos naujiniai)**, kad peržiūrėtumėte Connect Platform platformoje prieinamus programinės įrangos ir ADF naujinius. Jei nepasirinkta, spustelėkite mygtuką **APP STORE**.
- Norimos programinės įrangos arba ADF paketo **Actions (Veiksmų)** stulpelyje spustelėkite **Download (Atsisiųsti)**.
- Pasibaigus atsisiuntimui, spustelėkite **Install (Įdiegti)**.

Ribotoji produkto garantija

Bendrovė „Life Technologies Corporation“ ir (arba) jos patronuojamoji (-osios) įmonė (-ės) suteikia produktams garantiją, kaip aprašyta bendrovės „Life Technologies“ bendrosiose pardavimo sąlygose: www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Kilus klausimų, susisiekite su bendrove „Life Technologies“: www.thermofisher.com/support.

Oncomine™ Dx Express Test

Pub. nr. MAN0025897 Red. D.0

Šī izstrādājuma lietotāja rokasgrāmatas un Genexus™ Dx Software un Oncomine™ Dx Express Test analīzes definīcijas failu (Assay Definition Files, ADF) atjauninājumi ir sniegti Thermo Fisher™ Connect Platform resursu bibliotēkā.


Lietotāja dokumentācijas lejupielāde

1. Dodieties uz apps.thermofisher.com/apps/spa/#/publiclib/genexusdx un tad pierakstieties vietnē thermofisher.com.
Piezīme: Ja jums nav thermofisher.com konta, pierakstīšanās ekrānā noklikšķiniet uz **Create Account (Izveidot kontu)** un tad izpildiet norādījumus.
2. Resursu bibliotēkā sameklējiet Oncomine™ Dx Express Test lietotāja rokasgrāmatas vēlamajā valodā un pēc tam lejupielādējiet un izvelciet ZIP mapes, kas satur lietotāja rokasgrāmatas.
3. Atveriet lietotāja rokasgrāmatas ar PDF skatītāju, piemēram, Adobe Reader™ vai līdzīgu.
Piezīme: Adobe Reader™ var lejupielādēt vietnē [Adobe Reader](#).

Genexus™ Dx Software un ADF atjauninājumu lejupielāde


Thermo Fisher™ Connect Platform konta saistīšana ar QR kodu (vadītājs/administrators)

Lai piekļūtu programmatūras atjauninājumiem, jums vispirms ir jāsaista iekārta ar Connect Platform kontu. Iekārtas saistīšana veicama vienu reizi.

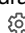
1. Programmatūrā Genexus™ Dx Software noklikšķiniet uz  **(Settings) (Iestatījumi) ▶ Thermo Fisher Account (Thermo Fisher konts)**.
2. Noklikšķiniet uz **+ Link Account to Device (Saistīt kontu ierīcei)**.
Tiek atvērts dialoglodziņš **Link Account to Device (Saistīt kontu ierīcei)** un tiek parādīts QR koda attēls ar QR koda tekstu.
3. Saistiet Connect Platform kontu ar programmatūru, izmantojot vienu no divām tālāk norādītajām opcijām.


Connect Platform konta saistīšanas opcijas

Opcija	Apraksts
Saistīt ar Instrument Connect mobilās ierīces lietotni	Piezīme: Lai izmantotu šo opciju, jums mobilajā ierīcē jābūt instalētai InstrumentConnect lietotnei. <ol style="list-style-type: none"> 1. Atveriet InstrumentConnect lietotni un piesakieties profilā. 2. Lietotnē noklikšķiniet uz +. 3. Noklikšķiniet uz QR Code (QR kods) un tad noskenējiet QR koda attēlu, kas tiek parādīts dialoglodziņā Link Account to Device (Saistīt kontu ierīcei) programmatūrā Genexus™ Dx Software.

Opcija	Apraksts
Saistīt, izmantojot Connect Platform tīmekļa vietni	<ol style="list-style-type: none"> 1. Iekopējiet parādīto QR kodu dialoglodziņā un tad noklikšķiniet uz Thermo Fisher Connect, lai atvērtu Connect Platform tīmekļa vietni. Piezīme: Nepieciešams pierakstīties. Ja jums nav Thermo Fisher™ Connect Platform konta, pierakstīšanās ekrānā noklikšķiniet uz Create Account (Izveidot kontu) un tad izpildiet norādījumus. 2. Connect Platform tīmekļa vietnē noklikšķiniet uz  InstrumentConnect kreisajā sāņņoslā. 3. Noklikšķiniet uz Link instrument (Saistīt iekārtu) ekrāna augšējā labajā stūrī. 4. Ievadiet ekrāna Link Instrument (Saistīt iekārtu) dialoglodziņā parādīto QR kodu. 5. Noklikšķiniet uz Link (Saistīt).

Programmatūras un ADF atjauninājumu lejupielāde un instalēšana

Kad Thermo Fisher™ Connect Platform konts ir saistīts, pieejamie programmatūras un ADF atjauninājumi ir redzami ekrānā Genexus™ Dx Software  **(Settings) (Iestatījumi) ▶ Software Updates (Programmatūras atjauninājumi)**. Vadītāja un administratora līmeņa lietotāji var lejupielādēt un instalēt atjauninājumus.

1. Programmatūrā Genexus™ Dx Software noklikšķiniet uz  **(Settings) (Iestatījumi) ▶ Software Updates (Programmatūras atjauninājumi)**.
2. Ekrānā **Installed Software (Instalētā programmatūra)** noklikšķiniet uz **Software Updates (Programmatūras atjauninājumi)**, lai skatītu programmatūras un ADF atjauninājumus, kas ir pieejami Connect Platform. Ja nav atlasīta, noklikšķiniet uz pogas **APP STORE**.
3. Interesējošās programmatūras vai ADF pakotnes kolonnā **Actions (Darbības)** noklikšķiniet uz **Download (Lejupielādēt)**.
4. Kad instalēšana ir pabeigta, noklikšķiniet uz **Install (Instalēt)**.

Ierobežota produkta garantija

Life Technologies Corporation un/vai tās filiāle(-es) sniedz savu produktu garantiju, kā noteikts Life Technologies Vispārējos pārdošanas noteikumos un nosacījumos, kas atrodami vietnē www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Jautājumu gadījumā lūdzam sazināties ar Life Technologies vietnē www.thermofisher.com/support.

Oncomine™ Dx Express Test

Pub.nr. MAN0025897 Rev. D.0

Brukerveiledninger for dette produktet og oppdateringer for Genexus™ Dx Software og analysedefinisjonsfiler (ADF-er) for Oncomine™ Dx Express Test finnes i et Thermo Fisher™ Connect Platform-ressursbibliotek.


Laste ned brukerdokumentasjon

- Gå til apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, og logg deretter på thermofisher.com.
Merk: Hvis du ikke har en konto på thermofisher.com, klikker du på **Create Account (Opprett konto)** i påloggings skjermbildet og følger instruksjonene.
- I ressursbiblioteket finner du brukerveiledningene for Oncomine™ Dx Express Test på ønsket språk. Last deretter ned og pakk ut ZIP-mappen som inneholder brukerveiledningene.
- Åpne brukerveiledningene i en PDF-fremviser, for eksempel Adobe Reader™ eller tilsvarende.
Merk: Adobe Reader™ kan lastes ned på [Adobe Reader](#).

Last ned oppdateringer for Genexus™ Dx Software og ADF


Koble en Thermo Fisher™ Connect Platform-konto med en QR-kode (leder/administrator)

For å få tilgang til programvareoppdateringer må du først koble et instrument til en Connect Platform-konto. Kobling av instrument gjøres bare én gang.

- I Genexus™ Dx Software klikker du på  **(Settings (Innstillinger))** ▶ **Thermo Fisher Account (Thermo Fisher-konto)**.
- Klikk på **+ Link Account to Device (Koble konto til enhet)**.
Dialogboksen **Link Account to Device (Koble konto til enhet)** åpnes og et QR-kodebilde og QR-kodetekst vises.
- Koble Connect Platform-kontoen med programvaren ved hjelp av ett av disse to alternativene.

Alternativer for å koble en Connect Platform-konto

Alternativ	Beskrivelse
Koble med mobilappen InstrumentConnect	<p>Merk: For å bruke dette alternativet må du ha mobilappen InstrumentConnect installert på en mobilenhet.</p> <ol style="list-style-type: none"> Åpne InstrumentConnect-appen og logg på. I appen klikker du på +. Klikk på QR Code (QR-kode), og skann deretter QR-kodebildet som er vist i dialogboksen Link Account to Device (Koble konto til enhet) i Genexus™ Dx Software.

Alternativ	Beskrivelse
Koble gjennom Connect Platform-nettstedet	<ol style="list-style-type: none"> Kopier QR-kodeteksten i dialogboksen, og klikk deretter på Thermo Fisher Connect for å åpne nettstedet Connect Platform. Merk: Pålogging er påkrevd. Hvis du ikke har en Thermo Fisher™ Connect Platform-konto, klikker du på Create Account (Opprett konto) i påloggings skjermbildet og følger deretter instruksjonene. På Connect Platform-nettstedet klikker du på  InstrumentConnect i venstre sidelinje. Klikk på Link instrument (Koble instrument) i øvre høyre hjørne av skjermbildet. Skriv inn QR-kodeteksten som er vist i dialogboksen i skjermbildet Link Instrument (Koble instrument). Klikk på Link (Koble).

Laste ned og installere programvare- og ADF-oppdateringer

Når Thermo Fisher™ Connect Platform-kontoen er koblet, vises tilgjengelige programvare- og ADF-oppdateringer i skjermbildet Genexus™ Dx Software  **(Settings (Innstillinger))** ▶ **Software Updates (Programvareoppdateringer)**. Brukere på leder- og administratornivå kan laste ned og installere oppdateringene.

- I Genexus™ Dx Software klikker du på ▶  **(Settings (Innstillinger))** ▶ **Software Updates (Programvareoppdateringer)**.
- I skjermbildet **Installed Software (Installert programvare)** klikker du på **Software Updates (Programvareoppdateringer)** for å vise oppdateringer for programvaren og ADF-er som er tilgjengelig på Connect Platform. Hvis ikke det er valgt, klikker du på knappen **APP STORE**.
- I kolonnen **Actions (Handlinger)** for ønsket programvare- eller ADF-pakke klikker du på **Download (Last ned)**.
- Når nedlastingen er fullført, klikker du på **Install (Installer)**.

Begrenset produktgaranti

Life Technologies Corporation og/eller dets tilknyttede selskap(er) garanterer sine produkter i henhold til Life Technologies' General Terms and Conditions of Sale (generelle vilkår og betingelser for salg) www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Hvis du har spørsmål, kan du ta kontakt med Life Technologies på www.thermofisher.com/support.

OncoMine™ Dx Express Test

Pub.nr. MAN0025897 Rev. D.0

Gebruikshandleidingen en updates voor de Genexus™ Dx Software en Assay Definition Files (Assaydefinitiebestanden, ADF's) voor de OncoMine™ Dx Express Test worden geleverd in een Thermo Fisher™ Connect Platform bronnenbibliotheek.

Gebruikersdocumentatie downloaden

- Ga naar apps.thermofisher.com/apps/spa/#/publiclib/genexusdx en meld u aan bij thermofisher.com.
Opmerking: Als u geen thermofisher.co-account hebt, klik dan op **Create Account (Account aanmaken)** op het aanmeldscherm en volg de instructies.
- Zoek in de bronnenbibliotheek de OncoMine™ Dx Express Test-gebruikershandleidingen in de gewenste taal, en download de ZIP-map met de gebruikershandleidingen en pak deze uit.
- Open de gebruikershandleidingen met een pdf-viewer zoals Adobe Reader™ of een gelijkwaardig programma.
Opmerking: Adobe Reader™ kan worden gedownload op [Adobe Reader](#).

Download Genexus™ Dx Software en ADF-updates

Een Thermo Fisher™ Connect Platform-account koppelen met een QR-code (manager/beheerder)

Om software-updates te kunnen gebruiken, moet u eerst een instrument aan een Connect Platform-account koppelen. Het koppelen van een instrument is een eenmalige actie.


- In Genexus™ Dx Software, klik op  **(Settings) (Instellingen)** ▶ **Thermo Fisher Account**.
- Klik op **+ Link Account to Device (Account koppelen aan toestel)**.

Het dialoogvenster **Link Account to Device (Account koppelen aan toestel)** verschijnt en een QR-codebeeld en QR-codetext worden getoond.

- Koppel de Connect Platform-account aan de software met behulp van een van de twee volgende opties.

Opties voor het koppelen van een Connect Platformaccount

Optie	Beschrijving
Koppelen met de mobiele app InstrumentConnect	<p>Opmerking: Als u deze optie wilt gebruiken, moet de mobiele app InstrumentConnect op een mobiel apparaat zijn geïnstalleerd.</p> <ol style="list-style-type: none"> Open de InstrumentConnect-app en meld u aan. In de app klikt u op +. Klik op QR Code (QR-code) en scan de QR-code die wordt weergegeven in het dialoogvenster Link Account to Device (Account koppelen aan toestel) in Genexus™ Dx Software.

Optie	Beschrijving
Koppelen via de Connect Platform-website	<ol style="list-style-type: none"> Kopieer de tekst van de QR-code in het dialoogvenster en klik vervolgens op Thermo Fisher Connect om de Connect Platform-website te openen. Opmerking: Aanmelden is vereist. Als u geen Thermo Fisher™ Connect Platform-account hebt, klik dan op Create Account (Account aanmaken) op het aanmeldscherm en volg de instructies. Op de Connect Platform-website klikt u op  InstrumentConnect in de linker zijbalk. Klik op Link instrument (Instrument koppelen) in de rechterbovenhoek van het scherm. Voer de QR-codetext in die wordt weergegeven in het dialoogvenster op het scherm Link Instrument (Instrument koppelen). Klik op Link (Koppelen).

Software- en ADF-updates downloaden en installeren

Wanneer de Thermo Fisher™ Connect Platform-account gekoppeld is, worden beschikbare software- en ADF-updates weergegeven op het Genexus™ Dx Software  **(Settings) (Instellingen)** ▶ **Software Updates (Software-updates)**-scherm. Gebruikers op het niveau van manager en beheerder kunnen de updates downloaden en installeren.

- In Genexus™ Dx Software, klik op ▶  **(Settings) (Instellingen)** ▶ **Software Updates (Software-updates)**.
- Klik in het scherm **Installed Software (Geïnstalleerde software)** op **Software Updates (Software-updates)** om de updates voor software en ADF weer te geven die beschikbaar zijn op het Connect Platform. Als dit niet is geselecteerd, klik dan op de knop **APP STORE**.
- Selecteer in de kolom **Actions (Acties)** het gewenste software- of ADF-pakket en klik vervolgens op **Download (Downloaden)**.
- Klik op **installeren (Voltooien)** als de download beëindigd is.

Beperkte productgarantie

Life Technologies Corporation en/of zijn filialen geven garantie op hun producten volgens de Algemene verkoopvoorwaarden van Life Technologies op www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Indien u nog vragen hebt, neem dan contact op met Life Technologies op www.thermofisher.com/support.

Oncomine™ Dx Express Test

Nr publ. MAN0025897 Wer. D.0

Podręczniki użytkownika niniejszego produktu, aktualizacje oprogramowania Genexus™ Dx Software oraz pliki definicji oznaczeń (Assay Definition File, ADF) dla testu Oncomine™ Dx Express Test są dostępne w bibliotece zasobów platformy Thermo Fisher™ Connect Platform.

Pobieranie dokumentacji użytkownika

1. Przejdź do strony apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, a następnie zalogować się do serwisu thermofisher.com.

Uwaga: W przypadku nieposiadania konta w serwisie thermofisher.com należy na ekranie logowania kliknąć opcję **Create Account (Utwórz konto)**, a następnie postępować zgodnie z instrukcjami.

2. W bibliotece zasobów odszukać podręczniki użytkownika sekwencjonatora Oncomine™ Dx Express Test w wybranym języku, a następnie pobrać i wypakować folder ZIP zawierający dane podręczniki.
3. Otworzyć podręczniki w przeglądarce plików PDF, np. Adobe Reader™ lub jej odpowiedniku.

Uwaga: Program Adobe Reader™ można pobrać ze strony [Adobe Reader](https://www.adobe.com/reader).

Pobieranie aktualizacji oprogramowania Genexus™ Dx Software i plików ADF

Łączenie konta w platformie Thermo Fisher™ Connect Platform przy użyciu kodu QR (menedżer/administrator)

Aby uzyskać dostęp do aktualizacji oprogramowania, należy najpierw połączyć urządzenie z kontem w platformie Connect Platform.

Łączenie urządzenia to działanie jednorazowe.

1. W oprogramowaniu Genexus™ Dx Software kliknąć kolejno opcję **(Settings (Ustawienia)) ▶ Thermo Fisher Account (Konto w serwisie Thermo Fisher)**.
2. Kliknąć opcję **+ Link Account to Device (Połącz konto z urządzeniem)**.

Otwarte zostanie okno dialogowe **Link Account to Device (Połącz konto z urządzeniem)** wraz z obrazem przedstawiającym kod QR oraz tekstem kodu QR.

3. Połączyć konto w platformie Connect Platform z oprogramowaniem, korzystając z jednej z poniższych dwóch opcji.

Opcje połączenia konta w platformie Connect Platform

Opcja	Opis
Połączenie z aplikacją mobilną Instrument Connect	<p>Uwaga: Aby skorzystać z tej opcji, należy zainstalować aplikację mobilną InstrumentConnect na urządzeniu mobilnym.</p> <ol style="list-style-type: none"> 1. Otworzyć aplikację InstrumentConnect, a następnie zalogować się. 2. W aplikacji kliknąć ikonę +. 3. Kliknąć opcję QR Code (Kod QR), następnie zeskanować kod QR wyświetlany w oknie dialogowym Link Account to Device (Połącz konto z urządzeniem) w oprogramowaniu Genexus™ Dx Software.
Połączenie przez stronę internetową platformy Connect Platform	<ol style="list-style-type: none"> 1. Skopiować tekst kodu QR widoczny w oknie dialogowym, a następnie kliknąć opcję Thermo Fisher Connect, aby otworzyć stronę internetową platformy Connect Platform. Uwaga: Wymagane jest zalogowanie. W przypadku nieposiadania konta w platformie Thermo Fisher™ Connect Platform należy na ekranie logowania kliknąć opcję Create Account (Utwórz konto), a następnie postępować zgodnie z instrukcjami. 2. Na stronie internetowej platformy Connect Platform kliknąć opcję InstrumentConnect na lewym pasku bocznym. 3. Kliknąć opcję Link instrument (Połącz urządzenie) w prawym górnym rogu ekranu. 4. Wprowadzić tekst kodu QR widoczny w oknie dialogowym Link Instrument (Połącz urządzenie). 5. Kliknąć opcję Link (Połącz).

Pobieranie i instalowanie aktualizacji oprogramowania oraz plików ADF

Jeśli konto w platformie Thermo Fisher™ Connect Platform jest połączone, dostępne aktualizacje oprogramowania i plików ADF są widoczne na ekranie **(Settings (Ustawienia)) ▶ Software Updates (Aktualizacje oprogramowania)** oprogramowania Genexus™ Dx Software. Użytkownicy o roli menedżera i administratora mogą pobierać i instalować aktualizacje.

1. W oprogramowaniu Genexus™ Dx Software kliknąć kolejno opcje **▶ (Settings (Ustawienia)) ▶ Software Updates (Aktualizacje oprogramowania)**.
2. Na ekranie **Installed Software (Zainstalowane oprogramowanie)** kliknąć opcję **Software Updates (Aktualizacje oprogramowania)**, aby wyświetlić aktualizacje oprogramowania i plików ADF dostępne w platformie Connect Platform. Jeśli nie został wybrany, kliknąć przycisk **APP STORE (Sklep z aplikacjami)**.
3. W kolumnie **Actions (Działania)** żadanego pakietu oprogramowania lub plików ADF kliknąć opcję **Download (Pobierz)**.
4. Po zakończeniu pobierania kliknąć opcję **Install (Instaluj)**.

Ograniczona gwarancja na produkt

Firma Life Technologies Corporation i/lub jej podmioty stowarzyszone udzielają gwarancji na produkty, jak opisano w dokumencie Life Technologies' General Terms and Conditions of Sale, dostępnym w witrynie internetowej www.thermofisher.com/us/en/home/global/terms-and-conditions.html. W razie pytań należy skontaktować się z firmą Life Technologies pod adresem www.thermofisher.com/support.

Oncomine™ Dx Express Test

N.º de pub. MAN0025897 Rev. D.0

Os guias do utilizador deste produto e as atualizações do Genexus™ Dx Software e dos ficheiros de definição do ensaio (ADF) do Oncomine™ Dx Express Test são fornecidos numa biblioteca de recursos da Thermo Fisher™ Connect Platform.

Transferir documentação do utilizador

- Vá para apps.thermofisher.com/apps/spa/#/publiclib/genexusdx e depois inicie sessão em thermofisher.com.
Nota: Se não tiver uma conta thermofisher.com, clique em **Create Account (Criar conta)** no ecrã de início de sessão e depois siga as instruções.
- Na biblioteca de recursos, localize os guias de utilizador do Oncomine™ Dx Express Test no idioma de interesse e depois transfira e extraia a pasta ZIP que contém os guias do utilizador.
- Abra os guias do utilizador com um visualizador de PDF como o Adobe Reader™ ou equivalente.
Nota: O Adobe Reader™ pode ser transferido em [Adobe Reader](#).

Transferir atualizações do Genexus™ Dx Software e do ADF

Associar uma conta Thermo Fisher™ Connect Platform a um código QR (gestor/administrador)

Para aceder a atualizações de software, primeiro tem de associar um instrumento a uma conta Connect Platform. Associar um instrumento é uma ação que só se faz uma vez.

- No Genexus™ Dx Software, clique em **(Settings (Definições))** ▶ **Thermo Fisher Account (Conta Thermo Fisher)**.
- Clique em **Link Account to Device (Associar conta a dispositivo)**.

A caixa de diálogo **Link Account to Device (Associar conta a dispositivo)** abre-se e apresenta uma imagem com o código QR e o texto do código QR.

- Associe a conta Connect Platform ao software utilizando uma das seguintes duas opções.

Opções de associação de uma conta Connect Platform

Opção	Descrição
Associar à aplicação móvel Instrument Connect	<p>Nota: Para utilizar esta opção, a aplicação móvel InstrumentConnect tem de estar instalada num dispositivo móvel.</p> <ol style="list-style-type: none"> Abra a aplicação InstrumentConnect e depois inicie sessão. Na aplicação, clique em . Clique no QR Code (Código QR) e depois leia a imagem do código QR apresentada na caixa de diálogo Link Account to Device (Associar conta a dispositivo) no Genexus™ Dx Software.

Opção	Descrição
Associar através do site da Connect Platform	<ol style="list-style-type: none"> Copie o texto do código QR que é apresentado na caixa de diálogo e depois clique em Thermo Fisher Connect para abrir o site Connect Platform. Nota: É necessário iniciar sessão. Se não tiver uma conta Thermo Fisher™ Connect Platform, clique em Create Account (Criar conta) no ecrã de início de sessão e depois siga as instruções. No site Connect Platform, clique em InstrumentConnect na barra lateral esquerda. Clique em Link instrument (Associar instrumento) no canto superior direito do ecrã. Introduza o texto do código QR que é apresentado na caixa de diálogo no ecrã Link Instrument (Associar instrumento). Clique em Link (Associar).

Transferir e instalar atualizações de software e de ADF
Quando a conta Thermo Fisher™ Connect Platform está associada, as atualizações de software e de ADF disponíveis ficam visíveis no ecrã do Genexus™ Dx Software **(Settings (Definições))** ▶ **Software Updates (Atualizações de software)**. Os utilizadores com nível de gestor e de administrador podem transferir e instalar as atualizações.

- Em Genexus™ Dx Software, clique em ▶ **(Settings (Definições))** ▶ **Software Updates (Atualizações de software)**.
- No ecrã **Installed Software (Software instalado)**, clique em **Software Updates (Atualizações de software)** para ver as atualizações do software e de ADF que estão disponíveis na Connect Platform. Se não estiver selecionado, clique no botão **APP STORE (LOJA DE APLICAÇÕES)**.
- Na coluna **Actions (Ações)** do pacote de software ou ADF de interesse, clique em **Download (Transferir)**.
- Depois de concluída a transferência, clique em **Install (Instalar)**.

Garantia limitada do produto

A Life Technologies Corporation e/ou a(s) sua(s) filial(ais) concede(m) uma garantia relativamente aos seus produtos conforme estabelecido nos Termos e condições de venda gerais da Life Technologies em www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Em caso de dúvidas, contacte a Life Technologies em www.thermofisher.com/support.

Oncomine™ Dx Express Test

Nr. pub. MAN0025897 Rev. D.0

Ghidurile de utilizare pentru acest produs și actualizările pentru Genexus™ Dx Software, precum și fișierele de definire a testului (ADF) pentru Oncomine™ Dx Express Test sunt furnizate într-o bibliotecă de resurse Thermo Fisher™ Connect Platform.


Descărcarea documentației de utilizare

- Accesați apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, apoi conectați-vă pe thermofisher.com.
Notă: Dacă nu aveți cont pe thermofisher.com, faceți clic pe **Create Account (Creare cont)** în ecranul de conectare, apoi urmați instrucțiunile.
- În biblioteca de resurse, localizați ghidurile de utilizare Oncomine™ Dx Express Test în limba de interes, apoi descărcați și extrageți folderul ZIP care conține ghidurile de utilizare.
- Deschideți ghidurile de utilizare cu un vizualizator PDF, precum Adobe Reader™ sau echivalent.
Notă: Adobe Reader™ se poate descărca de pe [Adobe Reader](#).

Descărcarea actualizărilor pentru Genexus™ Dx Software și ADF

Asociați un cont Thermo Fisher™ Connect Platform cu un cod QR (manager/administrator)

Pentru a accesa actualizări de software, mai întâi trebuie să asociați un instrument la un cont Connect Platform. Asocierea unui instrument este o acțiune pe care trebuie să o efectuați o singură dată.

- În Genexus™ Dx Software, faceți clic pe  **(Settings) (Setări)** ▶ **Thermo Fisher Account (Cont Thermo Fisher)**.
- Faceți clic pe **+** **Link Account to Device (Asociere cont la dispozitiv)**.


Se deschide caseta de dialog **Link Account to Device (Asociere cont la dispozitiv)** și sunt afișate o imagine cu codul QR și textul codului QR.


- Asociați contul Connect Platform cu software-ul utilizând una dintre următoarele două opțiuni.

Opțiuni pentru asocierea unui cont Connect Platform

Opțiune	Descriere
Asocierea cu aplicația mobilă Instrument Connect	<p>Notă: Pentru a utiliza această opțiune, trebuie să aveți instalată aplicația mobilă InstrumentConnect pe un dispozitiv mobil.</p> <ol style="list-style-type: none"> Deschideți aplicația InstrumentConnect, apoi conectați-vă. În aplicație, faceți clic pe +. Faceți clic pe QR Code (Cod QR), apoi scanați imaginea cu codul QR afișată în caseta de dialog Link Account to Device (Asociere cont la dispozitiv) din Genexus™ Dx Software.

Opțiune	Descriere
Asocierea prin intermediul unui site web Connect Platform	<ol style="list-style-type: none"> Copiați textul cu codul QR indicat în caseta de dialog, apoi faceți clic pe Thermo Fisher Connect pentru a deschide site-ul web Connect Platform. Notă: Conectarea este obligatorie. Dacă nu aveți cont Thermo Fisher™ Connect Platform, faceți clic pe Create Account (Creare cont) în ecranul de conectare, apoi urmați instrucțiunile. Pe site-ul web Connect Platform, faceți clic pe  InstrumentConnect (Conectare instrument) din bara laterală din stânga. Faceți clic pe Link instrument (Asociere instrument) în colțul din dreapta sus al ecranului. Introduceți textul codului QR afișat în caseta de dialog în ecranul Link Instrument (Asociere instrument). Faceți clic pe Link (Asociere).

Descărcarea și instalarea actualizărilor de software și ADF
Atunci când este asociat contul Thermo Fisher™ Connect Platform, actualizările de software și ADF disponibile sunt vizibile în ecranul Genexus™ Dx Software  **(Settings) (Setări)** ▶ **Software Updates (Actualizări de software)**. Utilizatorii cu nivel de manager și administrator pot descărca și instala actualizările.

- În Genexus™ Dx Software, faceți clic pe ▶  **(Settings) (Setări)** ▶ **Software Updates (Actualizări software)**.
- În ecranul **Installed Software (Software instalat)**, faceți clic pe **Software Updates (Actualizări de software)** pentru a vizualiza actualizările de software și ADF care sunt disponibile la Connect Platform. Dacă nu este selectat, faceți clic pe butonul **APP STORE**.
- În coloana **Actions (Acțiuni)** pentru pachetul software sau ADF de interes, faceți clic pe **Download (Descărcare)**.
- După ce instalarea este încheiată, faceți clic pe **Install (Instalare)**.

Garanție limitată a produsului

Life Technologies Corporation și/sau filialele sale garantează produsele sale conform termenilor și condițiilor generale de vânzare ale Life Technologies disponibile la www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Dacă aveți întrebări, contactați Life Technologies la www.thermofisher.com/support.

Oncomine™ Dx Express Test

№ публ. MAN0025897 Ред. D.0

Руководства пользователя для этого продукта и обновления программного обеспечения Genexus™ Dx Software и файлов определения тест-системы (ADF) для Oncomine™ Dx Express Test предоставляются через библиотеку ресурсов Thermo Fisher™ Connect Platform.


Загрузка пользовательской документации

1. Перейдите к apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, затем выполните вход в thermofisher.com.
Примечание: При отсутствии учетной записи thermofisher.com, нажмите **Create Account (Создать учетную запись)** на экране входа в систему и следуйте инструкциям.
2. Найдите руководства пользователя Oncomine™ Dx Express Test на нужном языке в библиотеке ресурсов, затем загрузите и распакуйте папку ZIP, содержащую руководства пользователя.
3. Откройте руководства пользователя в средстве просмотра PDF, например в Adobe Reader™ или эквивалентной программе.
Примечание: Adobe Reader™ можно загрузить из [Adobe Reader](#).


Загрузка обновлений программного обеспечения Genexus™ Dx Software и ADF

Связывание учетной записи Thermo Fisher™ Connect Platform с помощью QR-кода (менеджер/администратор)

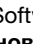
Для получения доступа к обновлениям программного обеспечения вначале нужно связать прибор с учетной записью Connect Platform. Связывание прибора это действие, выполняемое однократно.

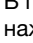
1. В программном обеспечении Genexus™ Dx Software нажмите  **(Settings) (Настройки) ▶ Thermo Fisher Account (Учетная запись Thermo Fisher)**.
2. Нажмите **+ Link Account to Device (Связать учетную запись с устройством)**.
Откроется диалоговое окно **Link Account to Device (Связать учетную запись с устройством)**, будут отображены изображение и текст QR-кода.
3. Свяжите учетную запись Connect Platform с программным обеспечением одним из двух следующих способов.

Возможности связывания с учетной записью Connect Platform

Функция	Описание
Связывание с мобильным приложением Instrument Connect	<p>Примечание: Чтобы использовать эту возможность, необходимо установить на мобильное устройство мобильное приложение Instrument Connect.</p> <ol style="list-style-type: none"> 1. Откройте приложение InstrumentConnect и выполните вход в систему. 2. В приложении нажмите +. 3. Нажмите на QR Code (QR-код), затем отсканируйте изображение QR-кода в диалоговом окне Link Account to Device (Связать учетную запись с устройством) в программном обеспечении Genexus™ Dx Software.
Связывание через веб-страницу Connect Platform	<ol style="list-style-type: none"> 1. Скопируйте текст QR-кода, приведенный в диалоговом окне, затем нажмите Thermo Fisher Connect, чтобы открыть веб-страницу Connect Platform. Примечание: Потребуется войти в систему. При отсутствии учетной записи Thermo Fisher™ Connect Platform, нажмите Create Account (Создать учетную запись) на экране входа в систему и следуйте инструкциям. 2. На веб-странице Connect Platform нажмите  InstrumentConnect на левой боковой панели. 3. Нажмите Link instrument (Связать прибор) в правом верхнем углу экрана. 4. Введите текст QR-кода, отображаемый в диалоговом окне на экране Link Instrument (Связать прибор). 5. Нажмите Link (Связать).

Загрузка и установка обновлений программного обеспечения и ADF

При установлении связи с учетной записью Thermo Fisher™ Connect Platform доступные обновления программного обеспечения и ADF становятся видны на экране Genexus™ Dx Software  **(Settings) (Настройки) ▶ Software Updates (Обновление программного обеспечения)**. Пользователи уровня администратора и менеджера смогут загружать и устанавливать обновления.

1. В программном обеспечении Genexus™ Dx Software нажмите **▶  (Settings) (Настройки) ▶ Software Updates (Обновления программного обеспечения)**.
2. На экране **Installed Software (Установленное программное обеспечение)** нажмите **Software Updates (Обновления программного обеспечения)**, чтобы просмотреть обновления программного обеспечения и ADF, доступные в Connect Platform. Нажмите на кнопку **APP STORE**, если она не выбрана.
3. В столбце **Actions (Действия)** нужного пакета для программного обеспечения или ADF нажмите **Download (Загрузить)**.
4. Когда загрузка будет закончена, нажмите **Install (Установить)**.

Ограниченная гарантия на продукт

Life Technologies Corporation и/или ее филиал(ы) предоставляют гарантию на свою продукцию в порядке, изложенном в документе «General Terms and Conditions of Sale» (Общие положения и условия продажи) компании Life Technologies, который опубликован на веб-сайте по адресу: www.thermofisher.com/us/en/home/global/terms-and-conditions.html. С любыми вопросами обращайтесь в компанию Life Technologies на веб-сайте по адресу: www.thermofisher.com/support.

Oncomine™ Dx Express Test

Č. pub. MAN0025897 Rev. D.0

Používateľské príručky pre tento produkt a aktualizácie Genexus™ Dx Software a súbory definície analýz (ADF) pre Oncomine™ Dx Express Test sú uvedené v knižnici zdrojov Thermo Fisher™ Connect Platform.



Stiahnutie používateľskej dokumentácie

1. Prejdite na apps.thermofisher.com/apps/spa/#/publiclib/genexusdx a prihláste sa na stránke thermofisher.com.
Poznámka: Ak nemáte účet thermofisher.com, na prihlasovacej obrazovke kliknite na možnosť **Create Account (Vytvoriť účet)** a potom postupujte podľa pokynov.
2. V knižnici zdrojov vyhľadajte používateľské príručky Oncomine™ Dx Express Test v požadovanom jazyku a potom si stiahnite a rozbaľte priečinok ZIP obsahujúci používateľské príručky.
3. Otvorte používateľské príručky pomocou prehliadača PDF, ako je napríklad Adobe Reader™ alebo ekvivalentného.
Poznámka: Adobe Reader™ si môžete stiahnuť na adrese [Adobe Reader](https://www.adobe.com/reader).


Stiahnite si Genexus™ Dx Software a aktualizácie ADF


Prepojenie účtu Thermo Fisher™ Connect Platform pomocou QR kódu (manažér/správca)

Ak chcete získať prístup k aktualizáciám softvéru, musíte najprv prepojiť prístroj s účtom Connect Platform. Prepojenie prístroja je jednorazová akcia.


1. V Genexus™ Dx Software kliknite na možnosť  **(Settings) (Nastavenia) ▶ Account (Účet Thermo Fisher)**.
2. Kliknite na možnosť  **Link Account to Device (Prepojiť účet so zariadením)**.
Otvorí sa dialógové okno **Link Account to Device (Prepojiť účet so zariadením)** a zobrazí sa obrázok QR kódu a text QR kódu.
3. Prepojte účet Connect Platform so softvérom pomocou jednej z dvoch nasledujúcich možností.


Možnosti prepojenia účtu Connect Platform

Možnosť	Popis
Prepojenie s mobilnou aplikáciou Instrument Connect	<p>Poznámka: Ak chcete použiť túto možnosť, musíte mať v mobilnom zariadení nainštalovanú mobilnú aplikáciu Instrument Connect.</p> <ol style="list-style-type: none"> 1. Otvorte aplikáciu InstrumentConnect a potom sa prihláste. 2. V aplikácii kliknite na . 3. Kliknite na možnosť QR Code (Kód QR), potom naskenujte obrázok s kódom QR zobrazený v dialógovom okne Link Account to Device (Prepojiť účet so zariadením) v Genexus™ Dx Software.

Možnosť	Popis
Prepojenie prostredníctvom webovej stránky Connect Platform	<ol style="list-style-type: none"> 1. Skopírujte text QR kódu uvedený v dialógovom okne a potom kliknutím na Thermo Fisher Connect otvorte webovú stránku Connect Platform. Poznámka: Vyžaduje sa prihlásenie. Ak nemáte účet Thermo Fisher™ Connect Platform, na prihlasovacej obrazovke kliknite na možnosť Create Account (Vytvoriť účet) a potom postupujte podľa pokynov. 2. Na webovej lokalite Connect Platform kliknite na  InstrumentConnect na ľavom bočnom paneli. 3. Kliknite na položku Link instrument (Prepojiť prístroj) v pravom hornom rohu obrazovky. 4. Zadajte text QR kódu zobrazený v dialógovom okne na obrazovke Link Instrument (Prepojiť prístroj). 5. Kliknite na možnosť Link (Prepojiť).

Stiahnutie a inštalácia aktualizácií softvéru a ADF

Keď je prepojený účet Thermo Fisher™ Connect Platform, dostupné aktualizácie softvéru a ADF sú viditeľné na obrazovke Genexus™ Dx Software  **(Settings) (Nastavenia) ▶ Software Updates (Aktualizácie softvéru)**. Stiahnuť a nainštalovať aktualizácie môžu používatelia na úrovni manažéra a správcu.

1. V Genexus™ Dx Software kliknite na možnosť  **(Settings) (Nastavenia) ▶ Software Updates (Aktualizácie softvéru)**.
2. Na obrazovke **Installed Software (Nainštalovaný softvér)** kliknite na tlačidlo **Software Updates (Aktualizácie softvéru)**, aby ste zobrazili aktualizácie softvéru a ADF, ktoré sú dostupné na Connect Platform. Ak táto možnosť nie je vybraná, kliknite na tlačidlo **APP STORE**.
3. V stĺpci **Actions (Akcie)** pre požadovaný softvér alebo balík ADF kliknite na možnosť **Download (Stiahnuť)**.
4. Po dokončení sťahovania kliknite na možnosť **Install (Inštalovať)**.

Obmedzená záruka na výrobok

Spoločnosť Life Technologies Corporation a/alebo jej pridružené spoločnosti poskytujú záruku na svoje výrobky tak, ako je uvedené vo všeobecných predajných zmluvných podmienkach spoločnosti Life Technologies na webovej lokalite www.thermofisher.com/us/en/home/global/terms-and-conditions.html. V prípade akýchkoľvek otázok kontaktujte spoločnosť Life Technologies na adrese www.thermofisher.com/support.

Oncomine™ Dx Express Test

Št. izd. MAN0025897 Rev. D.0

Uporabniški priročniki za ta izdelek in posodobitve za Genexus™ Dx Software ter datoteke z definicijami testa (ADF) za Oncomine™ Dx Express Test so na voljo v knjižnici virov platforme Thermo Fisher™ Connect Platform.

Prenos uporabniške dokumentacije

1. Pomaknite se na apps.thermofisher.com/apps/spa/#/publiclib/genexusdx in se nato prijavite v thermofisher.com.

Opomba: Če nimate računa thermofisher.com, kliknite **Create Account (Ustvari račun)** na zaslону za prijavo in upoštevajte navodila.

2. V knjižnici virov poiščite uporabniške priročnike za Oncomine™ Dx Express Test v jeziku, ki vas zanima, nato pa prenesite in ekstrahirajte mapo ZIP, ki vsebuje uporabniške priročnike.
3. Odprite uporabniške priročnike s pregledovalnikom PDF, kot je Adobe Reader™ ali enakovrednim pregledovalnikom.

Opomba: Adobe Reader™ je mogoče prenesti na [Adobe Reader](https://www.adobe.com/reader).

Prenos posodobitev za Genexus™ Dx Software in ADF

Povezovanje računa Thermo Fisher™ Connect Platform s kodo QR (upravljavce/skrbnik)

Za dostop do posodobitev programske opreme morate instrument najprej povezati z računom Connect Platform. Povezovanje instrumenta je enkratno dejanje.


1. V Genexus™ Dx Software kliknite  **(Settings (Nastavitve)) ▶ Thermo Fisher Account (Račun Thermo Fisher Account)**.
2. Kliknite **+ Link Account to Device (Poveži račun z napravo)**.

Odpre se pogovorno okno **Link Account to Device (Poveži račun z napravo)** ter prikazeta se slika kode QR in besedilo kode QR.

3. Račun Connect Platform povežite s programsko opremo z eno izmed naslednjih dveh možnosti.

Možnosti za povezovanje računa Connect Platform

Možnost	Opis
Povezava z mobilno aplikacijo Instrument Connect	<p>Opomba: Za uporabo te možnosti morate imeti v mobilni napravi nameščeno mobilno aplikacijo InstrumentConnect.</p> <ol style="list-style-type: none"> 1. Odprite aplikacijo InstrumentConnect in se prijavite. 2. V aplikaciji kliknite +. 3. Kliknite QR Code (Koda QR), nato skenirajte sliko kode QR, prikazano v pogovornem oknu Link Account to Device (Poveži račun z napravo) v Genexus™ Dx Software.

Možnost	Opis
Povezava prek spletnega mesta Connect Platform	<ol style="list-style-type: none"> 1. Kopirajte besedilo kode QR, navedeno v pogovornem oknu, nato kliknite Thermo Fisher Connect, da odprete spletno mesto Connect Platform. Opomba: Prijava je obvezna. Če nimate računa Thermo Fisher™ Connect Platform, kliknite Create Account (Ustvari račun) na zaslону za prijavo in upoštevajte navodila. 2. Na spletnem mestu Connect Platform kliknite  InstrumentConnect v levi stranski vrstici. 3. Kliknite Link instrument (Poveži instrument) v zgornjem desnem kotu zaslona. 4. Vnesite kodo QR, prikazano v pogovornem oknu na zaslону Link Instrument (Poveži instrument). 5. Kliknite Link (Poveži).

Prenos in namestitev posodobitev programske opreme in ADF

Ko je račun Thermo Fisher™ Connect Platform povezan, so na zaslону Genexus™ Dx Software  **(Settings (Nastavitve)) ▶ Software Updates (Posodobitve programske opreme)** vidne razpoložljive posodobitve za programsko opremo in ADF. Uporabniki na ravni upravljavca in skrbnika lahko prenesejo in namestijo posodobitve.

1. V Genexus™ Dx Software kliknite **▶  (Settings (Nastavitve)) ▶ Software Updates (Posodobitve programske opreme)**.
2. Na zaslону **Installed Software (Nameščena programska oprema)** kliknite **Software Updates (Posodobitve programske opreme)**, da si ogledate posodobitve za programsko opremo in ADF, ki so na voljo na Connect Platform. Če gumb **APP STORE** ni izbran, ga kliknite.
3. V stolpcu **Actions (Dejanja)** kliknite **Download (Prenesi)** za paket programske opreme ali ADF, ki vas zanima.
4. Ko je prenos dokončan, kliknite **Install (Namesti)**.

Omejena garancija za izdelek

Družba Life Technologies Corporation in/ali njene podružnice zagotavljajo garancijo za izdelke, kot je določeno v splošnih prodajnih pogojih (Life Technologies' General Terms and Conditions of Sale) na spletni strani www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Če imate kakršna koli vprašanja, se obrnite na družbo Life Technologies prek spletne strani www.thermofisher.com/support.

Oncomine™ Dx Express Test

Utg.nr MAN0025897 Rev. D.0

Användarguider för denna produkt och uppdateringar till Genexus™ Dx Software och Assay Definition Files (ADFs, analysdefinitionsfiler) för Oncomine™ Dx Express Test finns i ett Thermo Fisher™ Connect Platform-resursbibliotek.


Ladda ned användardokumentation

- Gå till apps.thermofisher.com/apps/spa/#/publiclib/genexusdx och logga in på thermofisher.com.
Obs: Om du inte har ett thermofisher.com-konto, klicka på **Create Account (Skapa konto)** på inloggningsskärmen, följ sedan anvisningarna.
- I resursbiblioteket, lokalisera Oncomine™ Dx Express Test-användarhandböckerna för det språk du vill använda, ladda sedan ner och extrahera på ZIP-mappen som innehåller användarhandböckerna.
- Öppna användarhandböckerna med en PDF-läsare såsom Adobe Reader™ eller motsvarande.
Obs: Adobe Reader™ kan laddas ned på [Adobe Reader](#).

Ladda ned Genexus™ Dx Software och ADF-uppdateringar


Länka ett Thermo Fisher™ Connect Platform-konto med en QR-kod (chef/administratör)

För att komma åt programuppdateringar måste du först länka ett instrument till ett Connect Platform-konto. Länkning av ett instrument är en engångsåtgärd.

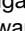
- På Genexus™ Dx Software, klicka på  (**Settings**) (**Inställningar**) ▶ **Thermo Fisher Account (Thermo Fisher-konto)**.
- Klicka på **+ Link Account to Device (Länka konto till enhet)**. **Link Account to Device (Länka konto till enhet)**-dialogrutan öppnas och en QR-kodbild och en QR-kodtext visas.
- Länka Connect Platform-kontot till programvaran med ett av två följande alternativ.

Alternativ för att länka ett Connect Platform-konto

Alternativ	Beskrivning
Länka med Instrument Connect-mobilappen	<p>Obs: För att använda det här alternativet måste du ha installerat InstrumentConnect-mobilappen på en mobil enhet.</p> <ol style="list-style-type: none"> Öppna appen InstrumentConnect, och logga sedan in. Klicka på + i appen. Klicka på QR Code (QR-kod) och skanna sedan QR-kod-bilden som visas i Link Account to Device (Länka konto till enhet)-dialogrutan i Genexus™ Dx Software.

Alternativ	Beskrivning
Länka via Connect Platform-webbplatsen	<ol style="list-style-type: none"> Kopiera QR-koden i dialogrutan och klicka sedan på Thermo Fisher Connect för att öppna Connect Platform-webbplatsen. Obs: Inloggning krävs. Om du inte har ett Thermo Fisher™ Connect Platform-konto, klicka på Create Account (Skapa konto) på inloggningsskärmen, följ sedan anvisningarna. Klicka på  InstrumentConnect på den vänstra sidopanelen på Connect Platform-webbplatsen. Klicka på Link instrument (Länka instrument) i skärmbildens övre högra hörn. Ange QR-koden som visas i dialogrutan på skärmbilden Link instrument (Länka instrument). Klicka på Link (Länka).

Ladda ned och installera ADF-uppdateringar

När Thermo Fisher™ Connect Platform-kontot är länkat kan tillgängliga programvaru- och ADF-uppdateringar ses på Genexus™ Dx Software  (**Settings**) (**Inställningar**) ▶ **Software Updates (Programvaruuppdateringar)**-skärmen. Användare på chefs- och administratörsnivå kan ladda ned och installera uppdateringarna.

- I Genexus™ Dx Software, klicka på ▶  (**Settings**) (**Inställningar**) ▶ **programvaruuppdateringar**.
- På skärmen **Installed Software (Installerad programvara)**, klicka på **Software Updates (Programvaruuppdateringar)** för att visa uppdateringar till programvara och ADF:er som är tillgängliga på Connect Platform. Om ej markerad, klicka på knappen **APP STORE (APP STORE)**.
- I kolumnen **Actions (Åtgärder)** för programvaru- eller ADF-paket av intresse, klicka på **Download (Ladda ner)**.
- När nedladdning är slutförd, klicka på **Install (Installera)**.

Begränsad produktgaranti

Life Technologies Corporation och/eller dess dotterbolag garanterar sina produkter i enlighet med Life Technologies allmänna villkor och försäljningsvillkor på www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Vid frågor, var god kontakta Life Technologies på www.thermofisher.com/support.

Oncomine™ Dx Express Test

Yay. No. MAN0025897 Rev. D.0

Bu ürüne yönelik kullanım kılavuzları ve Genexus™ Dx Software ve Oncomine™ Dx Express Test Tahlil Tanımlama Dosyaları (ADF'ler) güncellemeleri bir Thermo Fisher™ Connect Platform kaynak kütüphanesinde sağlanır.


Kullanıcı belgelerini indirme

1. apps.thermofisher.com/apps/spa/#/publiclib/genexusdx adresine gidin ve ardından thermofisher.com üzerinde oturum açın.
Not: thermofisher.com hesabınız yoksa oturum açma ekranında **Create Account (Hesap Oluştur)** ögesine tıklayın ve ardından talimatları izleyin.
2. Kaynak kütüphanesinde, ilgilenilen dildeki Oncomine™ Dx Express Test kullanım kılavuzlarını bulun, ardından kullanım kılavuzlarını içeren ZIP klasörünü indirin ve çıkartın.
3. Kullanım kılavuzlarını Adobe Reader™ veya eşdeğer bir PDF görüntüleyici ile açın.
Not: Adobe Reader™, [Adobe Reader](#) adresinden indirilebilir.

Genexus™ Dx Software ve ADF güncellemelerini indirme

QR kodu ile bir Thermo Fisher™ Connect Platform hesabını bağlama (yönetici/sistem yöneticisi)

Yazılım güncellemelerine erişmek için öncelikle bir cihazı bir Connect Platform hesabına bağlamanız gerekir. Cihaz bağlama tek seferlik bir eylemdir.


1. Genexus™ Dx Software yazılımında,  **(Settings (Ayarlar))** ▶ **Thermo Fisher Account (Thermo Fisher Hesabı)** ögesine tıklayın.
2. **+ Link Account to Device (Hesabı Cihaza Bağla)** ögesine tıklayın.
Link Account to Device (Hesabı Cihaza Bağla) iletişim kutusu açılır ve bir QR kodu görüntüsü ve QR kodu metni gösterilir.
3. Aşağıdaki iki seçeneğin birini kullanarak Connect Platform hesabını yazılıma bağlayın.


Bir Connect Platform hesabını bağlama seçenekleri

Seçenek	Açıklama
Instrument Connect mobil uygulaması ile bağlama	Not: Bu seçeneği kullanmak için InstrumentConnect mobil uygulamasının bir mobil cihazda kurulu olması gerekir. 1. InstrumentConnect uygulamasını açın ve oturum açın. 2. Uygulamada + ögesine tıklayın. 3. QR Code (QR Kodu) ögesine tıklayın ve ardından, Link Account to Device (Hesabı Cihaza Bağla) iletişim kutusunda gösterilen QR kodu resmini Genexus™ Dx Software yazılımında tarayın.

Seçenek	Açıklama
Connect Platform web sitesi aracılığıyla bağlama	1. İletişim kutusunda verilen QR kodu metnini kopyalayın, ardından Connect Platform web sitesini açmak için Thermo Fisher Connect ögesine tıklayın. Not: Oturum açma gereklidir. Thermo Fisher™ Connect Platform hesabınız yoksa oturum açma ekranında Create Account (Hesap Oluştur) ögesine tıklayın ve ardından talimatları izleyin. 2. Connect Platform web sitesinde, sol kenar çubuğundaki InstrumentConnect ögesine tıklayın. 3. Ekranın sağ üst köşesindeki Link instrument (Cihazı bağla) ögesine tıklayın. 4. Link Instrument (Cihazı Bağla) ekranındaki iletişim kutusunda gösterilen QR kodu metnini girin. 5. Link (Bağla) ögesine tıklayın.

Yazılım ve ADF güncellemelerini indirme ve kurma

Thermo Fisher™ Connect Platform hesabı bağlandığında, mevcut yazılım ve ADF güncellemeleri Genexus™ Dx Software  **(Settings (Ayarlar))** ▶ **Software Updates (Yazılım Güncellemeleri)** ekranında görünür. Yönetici ve sistem yöneticisi seviyesindeki kullanıcılar, güncellemeleri indirebilir ve kurabilirler.

1. Genexus™ Dx Software yazılımında, ▶  **(Settings (Ayarlar))** ▶ **Software Updates (Yazılım Güncellemeleri)** ögesine tıklayın.
2. **Installed Software (Kurulu Yazılım)** ekranında, Connect Platform içinde bulunan yazılım ve ADF güncellemelerini görüntülemek için **Software Updates (Yazılım Güncellemeleri)** ögesine tıklayın. Seçili değilse **APP STORE** düğmesine tıklayın.
3. İlgilenilen yazılım veya ADF paketinin **Actions (Eylemler)** sütununda **Download (İndir)** ögesine tıklayın.
4. İndirme tamamlandıktan sonra **Install (Kur)** ögesine tıklayın.

Sınırlı ürün garantisi

Life Technologies Corporation ve/veya bağlı kuruluşları, ürünlerini www.thermofisher.com/us/en/home/global/terms-and-conditions.html adresindeki Life Technologies Genel Satış Şartları ve Koşulları içinde ifade edildiği şekilde garantilemektedir. Sorularınız varsa lütfen www.thermofisher.com/support adresinden Life Technologies ile iletişime geçin.

Oncomine™ Dx Express Test

Số Ấn phẩm MAN0025897 Bản sửa đổi D.0

Hướng dẫn sử dụng cho sản phẩm này, các bản cập nhật Genexus™ Dx Software và các Tập Định nghĩa Xét nghiệm (ADF) cho Oncomine™ Dx Express Test được cung cấp trong thư viện tài nguyên Thermo Fisher™ Connect Platform.


Tải xuống tài liệu người dùng

- Đi đến apps.thermofisher.com/apps/spa/#/publiclib/genexusdx, sau đó đăng nhập vào thermofisher.com.
Lưu ý: Nếu bạn không có tài khoản thermofisher.com, hãy nhấp vào **Create Account (Tạo Tài khoản)** trong màn hình đăng nhập, sau đó làm theo các hướng dẫn.
- Trong thư viện tài nguyên, hãy xác định vị trí của các hướng dẫn sử dụng Oncomine™ Dx Express Test bằng ngôn ngữ được quan tâm, sau đó tải xuống và giải nén thư mục ZIP chứa các hướng dẫn sử dụng.
- Mở hướng dẫn sử dụng bằng trình xem PDF chẳng hạn như Adobe Reader™ hoặc tương đương.
Lưu ý: Adobe Reader™ có thể được tải xuống tại [Adobe Reader](#).

Tải xuống Genexus™ Dx Software và các bản cập nhật ADF


Liên kết tài khoản Thermo Fisher™ Connect Platform với mã QR (người quản lý/quản trị viên)

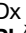
Để truy cập các bản cập nhật phần mềm, trước tiên bạn phải liên kết một công cụ với tài khoản Connect Platform. Liên kết một thiết bị là hành động một lần.


- Trong Genexus™ Dx Software, nhấp vào  **(Settings) (Cài đặt) ▶ Thermo Fisher Account (Tài khoản Thermo Fisher)**.
- Nhấp vào **+ Link Account to Device (Liên kết Tài khoản với Thiết bị)**.
Hộp thoại **Link Account to Device (Liên kết Tài khoản với Thiết bị)** mở ra và hình ảnh mã QR và văn bản mã QR được hiển thị.
- Liên kết tài khoản Connect Platform với phần mềm bằng cách sử dụng một trong hai tùy chọn sau.

Các tùy chọn để kết nối một tài khoản Connect Platform

Tùy chọn	Mô tả
Liên kết với ứng dụng di động InstrumentConnect	<p>Lưu ý: Để sử dụng tùy chọn này, bạn phải cài đặt ứng dụng di động InstrumentConnect trên thiết bị di động.</p> <ol style="list-style-type: none"> Mở ứng dụng InstrumentConnect, sau đó đăng nhập. Trong ứng dụng, nhấp vào +. Nhấp vào QR Code (Mã QR), sau đó quét hình ảnh mã QR hiển thị trong hộp thoại Link Account to Device (Liên kết Tài khoản với Thiết bị) trong Genexus™ Dx Software.

Tùy chọn	Mô tả
Liên kết qua trang web Connect Platform	<ol style="list-style-type: none"> Sao chép văn bản mã QR được cung cấp trong hộp thoại, sau đó nhấp vào Thermo Fisher Connect để mở trang web Connect Platform. Lưu ý: Bạn cần phải đăng nhập. Nếu bạn không có tài khoản Thermo Fisher™ Connect Platform, hãy nhấp vào Create Account (Tạo Tài khoản) trong màn hình đăng nhập, sau đó làm theo các hướng dẫn. Trong trang web Connect Platform, hãy nhấp vào  InstrumentConnect ở thanh bên trái. Nhấp vào Link instrument (Liên kết thiết bị) ở góc trên bên phải màn hình. Nhập văn bản mã QR được hiển thị trong hộp thoại trong màn hình Link Instrument (Liên kết Thiết bị). Nhấp vào Link (Liên kết).

Tải xuống và cài đặt các bản cập nhật phần mềm và ADF Sau khi liên kết tài khoản Thermo Fisher™ Connect Platform, phần mềm và các bản cập nhật ADF có sẵn sẽ được hiển thị trong màn hình Genexus™ Dx Software  **(Settings) (Cài đặt) ▶ Software Updates (Cập nhật Phần mềm)**. Người dùng cấp quản lý và cấp quản trị viên có thể tải xuống và cài đặt các bản cập nhật.

- Trong Genexus™ Dx Software, nhấp vào  **(Settings) (Cài đặt) ▶ Software Updates (Cập nhật Phần mềm)**.
- Trong màn hình **Installed Software (Phần mềm đã Cài đặt)**, nhấp vào **Software Updates (Cập nhật Phần mềm)** để xem các bản cập nhật cho phần mềm và ADF có sẵn tại Connect Platform. Nếu không được chọn, hãy nhấp vào nút **APP STORE (CỬA HÀNG ỨNG DỤNG)**.
- Trong cột **Actions (Tác vụ)** của phần mềm hoặc gói ADF được quan tâm, hãy nhấp vào **Download (Tải xuống)**.
- Sau khi quá trình tải xuống hoàn tất, hãy nhấp vào **Install (Cài đặt)**.

Bảo hành sản phẩm giới hạn

Life Technologies Corporation và/hoặc (các) chi nhánh của họ bảo hành sản phẩm như quy định trong các Điều khoản và Điều kiện Bán hàng Chung của Life Technologies tại www.thermofisher.com/us/en/home/global/terms-and-conditions.html. Nếu bạn có thắc mắc, vui lòng liên hệ Life Technologies theo www.thermofisher.com/support.

Oncomine™ Dx Express Test

出版编号 MAN0025897 修订版 D.0

本产品的用户指南、Genexus™ Dx Software 更新和 Oncomine™ Dx Express Test 的检测定义文件 (ADF) 见 Thermo Fisher™ Connect Platform 资源库。


下载用户文档

1. 打开 apps.thermofisher.com/apps/spa/#/publiclib/genexusdx，然后登录 thermofisher.com。
注释: 如果您没有 thermofisher.com 帐户，请在登录屏幕单击 **Create Account (创建帐户)**，然后按照说明操作。
2. 在资源库中，找到相关语言的 Oncomine™ Dx Express Test 用户指南，然后下载并解压包含用户指南的 ZIP 文件夹。
3. 使用 PDF 查看器 (如 Adobe Reader™ 或同等软件) 打开用户指南。
注释: Adobe Reader™ 可在 [Adobe Reader](#) 下载。

下载 Genexus™ Dx Software 软件和 ADF 更新

使用 QR 代码关联 Thermo Fisher™ Connect Platform 帐户 (经理/管理员)

要访问软件更新，必须首先将仪器关联到 Connect Platform 帐户。关联仪器是一次性操作。

1. 在 Genexus™ Dx Software 中，单击  (设置) ▶ **Thermo Fisher Account (Thermo Fisher 帐户)**。
2. 单击 **+ Link Account to Device (关联帐户与设备)**。
系统会打开 **Link Account to Device (关联帐户与设备)** 对话框，并会显示二维码图像和二维码文本。
3. 使用以下两种选项之一将 Connect Platform 帐户与软件关联。

关联 Connect Platform 帐户的选项

选项	说明
关联 Instrument Connect 移动应用	<p>注释: 若要使用此选项，必须在移动设备上安装 InstrumentConnect 移动应用。</p> <ol style="list-style-type: none"> 1. 打开 InstrumentConnect 应用，然后登录。 2. 在该应用中，单击 +。 3. 单击 QR Code (二维码)，然后扫描 Genexus™ Dx Software 中 Link Account to Device (关联帐户与设备) 对话框中显示的二维码图像。
通过 Connect Platform 网站进行关联	<ol style="list-style-type: none"> 1. 复制对话框中的二维码文本，然后单击 Thermo Fisher Connect 打开 Connect Platform 网站。 注释: 需要登录。如果您没有 Thermo Fisher™ Connect Platform 帐户，请在登录屏幕单击 Create Account (创建帐户)，然后按照说明操作。 2. 在 Connect Platform 网站，单击左侧边栏中的  InstrumentConnect。 3. 单击屏幕右上角的 Link instrument (关联仪器)。 4. 输入 Link Instrument (关联仪器) 屏幕中对话框上显示的二维码文本。 5. 单击 Link (关联)。

下载并安装软件和 ADF 更新

关联 Thermo Fisher™ Connect Platform 帐户后，可用的软件和 ADF 更新会显示在 Genexus™ Dx Software  (**Settings (设置)**) ▶ **Software Updates (软件更新)** 屏幕中。经理和管理员级别用户可以下载并安装这些更新。

1. 在 Genexus™ Dx Software 中，单击 ▶  (**Settings (设置)**) ▶ **Software Updates (软件更新)**。
2. 在 **Installed Software (安装的软件)** 屏幕中，单击 **Software Updates (软件更新)** 以查看 Connect Platform 上可用的软件更新和 ADF。如果没有选中，请单击 **APP STORE (APP 商店)** 按钮。
3. 在相关软件或 ADF 软件包的 **Actions (操作)** 列中，单击 **Download (下载)**。
4. 下载完成后，单击 **Install (安装)**。

有限产品保修

Life Technologies Corporation 和/或其附属公司根据 www.thermofisher.com/us/en/home/global/terms-and-conditions.html 上提供的 Life Technologies 一般销售条款和条件为其产品提供质保。如果有任何问题，请联系 Life Technologies，网址：www.thermofisher.com/support。



Life Technologies Europe B.V.
Kwartsweg 2, 2665 NN Bleiswijk
The Netherlands

UKRP Life Technologies Ltd
7 Kingsland Grange
Woolston
Warrington
WA1 4SR
UK



Manufacturer:

Life Technologies Corporation | 7335
Executive Way | Frederick, Maryland 21704
USA

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0025897

Revision	Date	Description
D.0	19 January 2023	More translations were added to the document.
C.0	12 July 2022	More translations were added to the document.
B.0	20 June 2022	Translations were added to the document.
A.0	11 March 2022	New document for the OncoPrint™ Dx Express Test

©2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Adobe Reader is a trademark of Adobe Systems, Inc.