



SSL Certificate Update Instructions

As part of our commitment to data security, updated SSL certificates are available for your MSD[®] instrument computer. The Methodical Mind[®] software on your MSD instrument computer uses SSL certificates to securely connect to [methodicalmind.com](https://www.methodicalmind.com). To ensure the software continues to operate without interruption, follow these instructions to install updated certificates. You need administrator privileges on the computer to install the certificates. For assistance, contact Scientific Support: 1-240-314-2798 or scientificsupport@mesoscale.com.

Update Certificates

1. Download the following five root SSL certificates:

<https://www.amazontrust.com/repository/SFSRootCAG2.cer>

<https://www.amazontrust.com/repository/AmazonRootCA1.cer>

<https://www.amazontrust.com/repository/AmazonRootCA2.cer>

<https://www.amazontrust.com/repository/AmazonRootCA3.cer>

<https://www.amazontrust.com/repository/AmazonRootCA4.cer>

If prompted, select the appropriate option to save each file.

2. Go to your *Downloads* folder to confirm you have downloaded five CER files.
3. Install each certificate:
 - a. Right click on the certificate file and select **Install Certificate**. If prompted, click **Open**.
 - b. Select **Local Machine** and click **Next**. If prompted, click **Yes**.
 - c. Select **Automatically select the certificate store based on the type of certificate** and click **Next**.
 - d. Click **Finish**.
 - e. Click **OK**.
4. Repeat this process on each instrument computer.

Confirm Certificate Installation

1. Open the Windows program **Manage Computer Certificates**.
2. Go to the folder *Trusted Root Certificates Authority > Certificates*.
3. Confirm the following certificates are installed:

Starfield Services Root Certificate Authority – G2 with an Expiration Date of 12/31/2037

Amazon Root CA 1 with an Expiration Date of 1/16/2038

Amazon Root CA 2 with an Expiration Date of 5/25/2040

Amazon Root CA 3 with an Expiration Date of 5/25/2040

Amazon Root CA 4 with an Expiration Date of 5/25/2040

Meso Scale Discovery

A division of Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA
www.mesoscale.com

Customer Service
1-240-314-2795
CustomerService@mesoscale.com

Scientific Support
1-240-314-2798
ScientificSupport@mesoscale.com

MSD, Methodical Mind, and MSD (design) are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners.

© 2025 Meso Scale Diagnostics, LLC. All rights reserved.