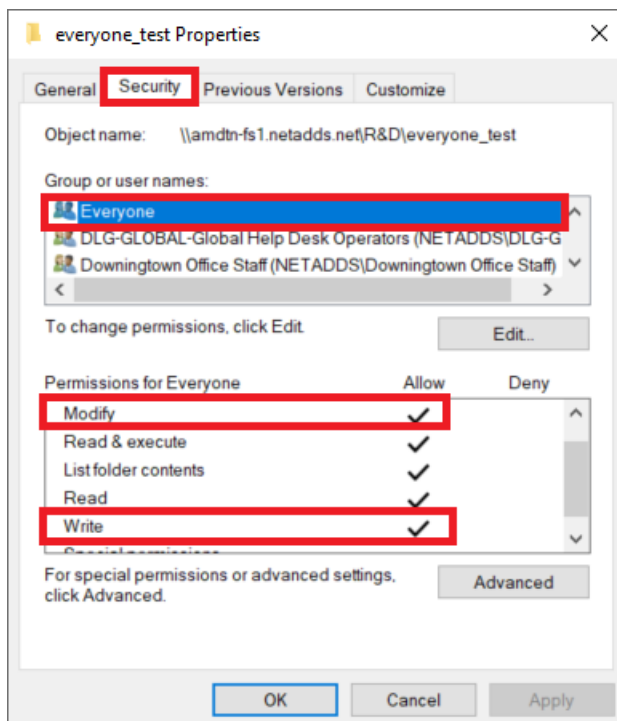


Using UNC Paths with CellReporterXpress Software Version 2.9.3

This document describes how to use UNC paths with CellReporterXpress Software Version 2.9.3. To map a UNC path as a network storage location in the software, the following must be true:

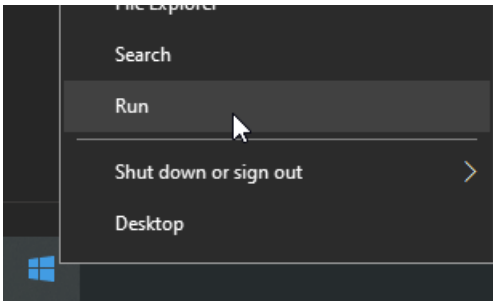
- The CellReporterXpress host computer must be running all software services, including the Analysis Service and Location Service. No software services should be running on any remote computer.
- The full UNC path (not a mapped drive letter) must be accessible on the host computer. You may need to follow the Windows prompts to log in manually before mapping the UNC path as a folder for storage in the CellReporterXpress software.
- The UNC path properties must include Modify and Write permissions for the Everyone group.



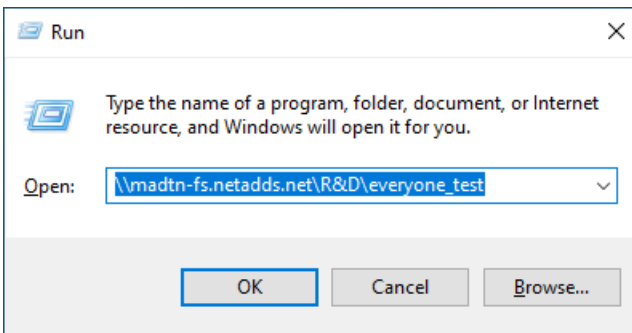
- Network access and throughput must be consistent and reliable. We recommend using a wired Ethernet connection (instead of a Wi-Fi connection). Note that network latency, which can result from heavy network traffic, can cause issues during acquisition, analysis, and exporting.

Mapping a UNC Path as a Storage Location

1. In Windows, right-click  **Start**, and select **Run**.

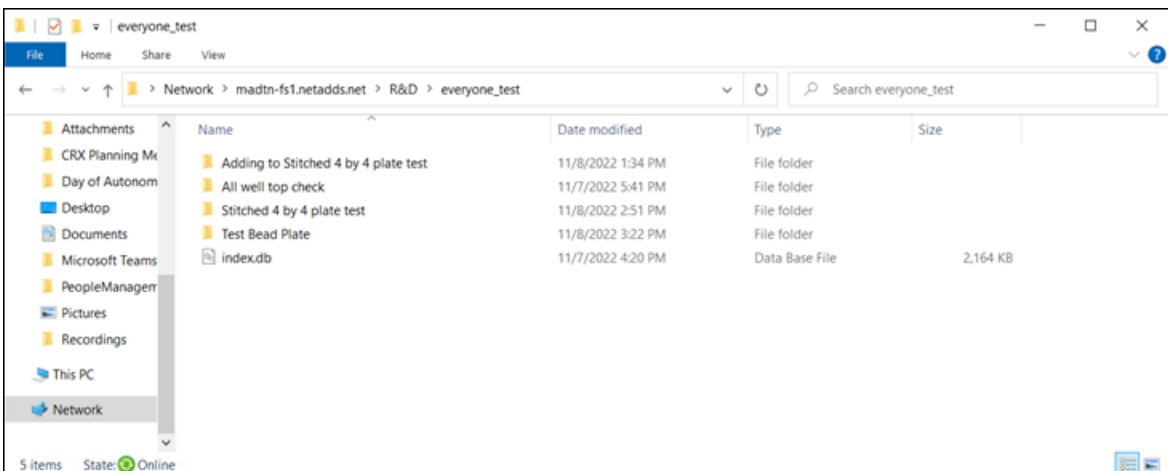



2. In the **Run** window, enter the UNC Path starting with two backslashes (\\), as shown below, and click **OK**.



You may need to enter login credentials if prompted by Windows.

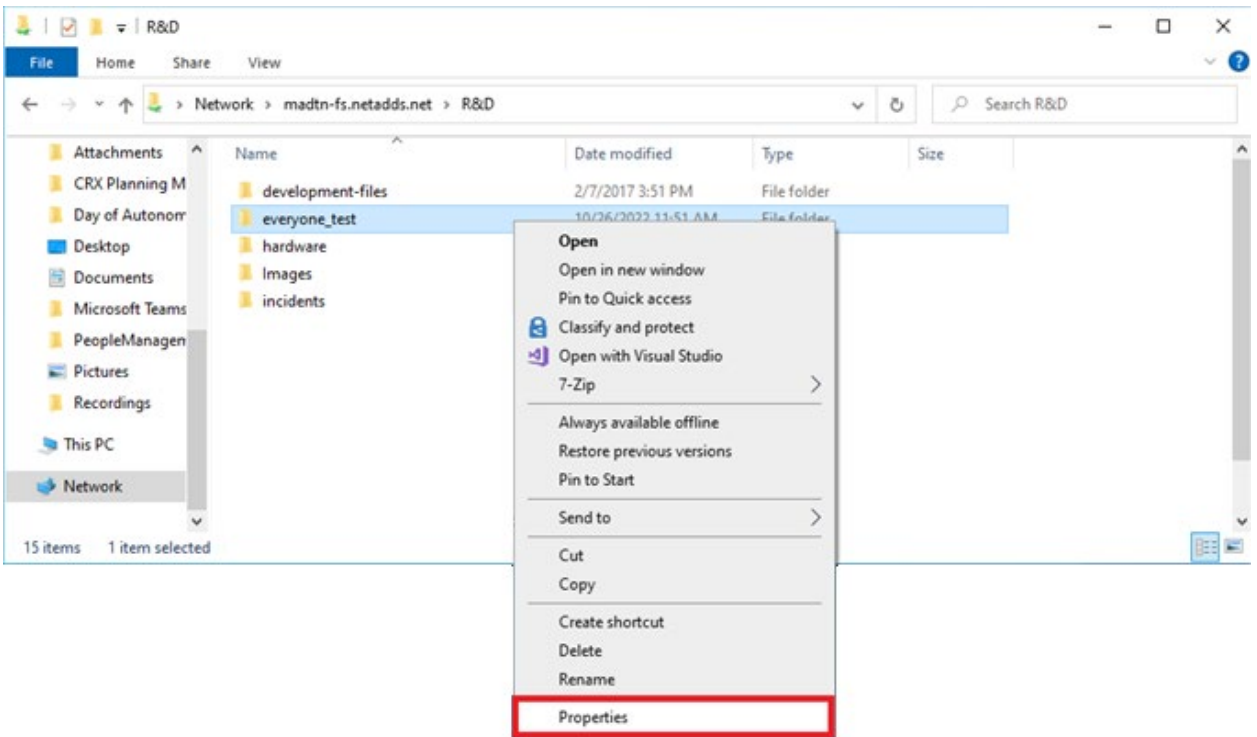
When you are logged in, Windows File Explorer opens and shows the UNC path location, which indicates that you can access the network location.



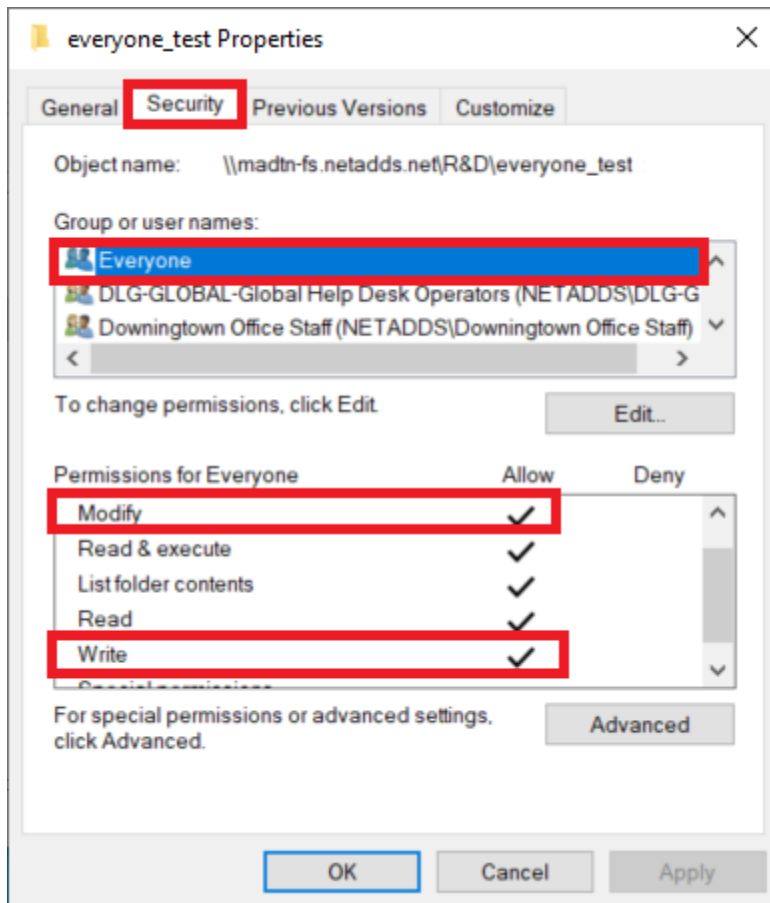
3. Click  **Up** to move up one level in the directory structure.



4. Right-click on the folder to be used as a storage location, and select **Properties**.

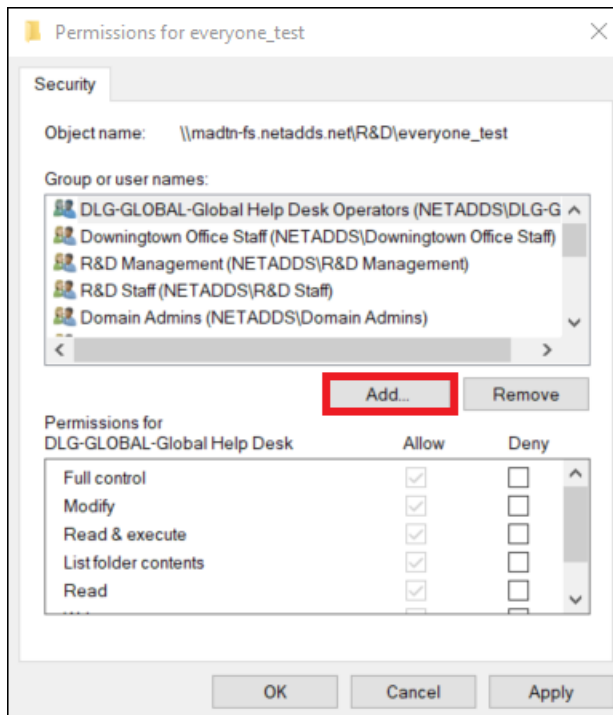


5. Select the **Security** tab.

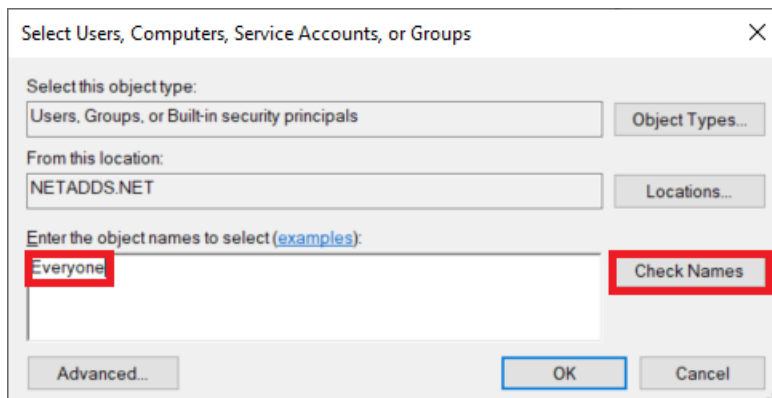


6. Confirm that **Everyone** is listed in the **Groups or user names** list.
If it is not listed, continue to the next step to add it. If it is listed, continue to step 13 of this procedure.
7. If **Everyone** is not listed, click **Edit**.

8. In the **Permissions** dialog, click **Add**.

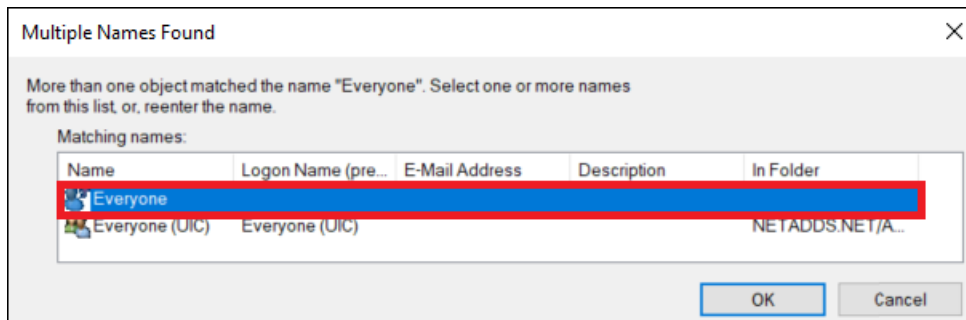


9. In the **Select Users or Groups** dialog, in the **Enter the object names to select** field, enter **Everyone**.



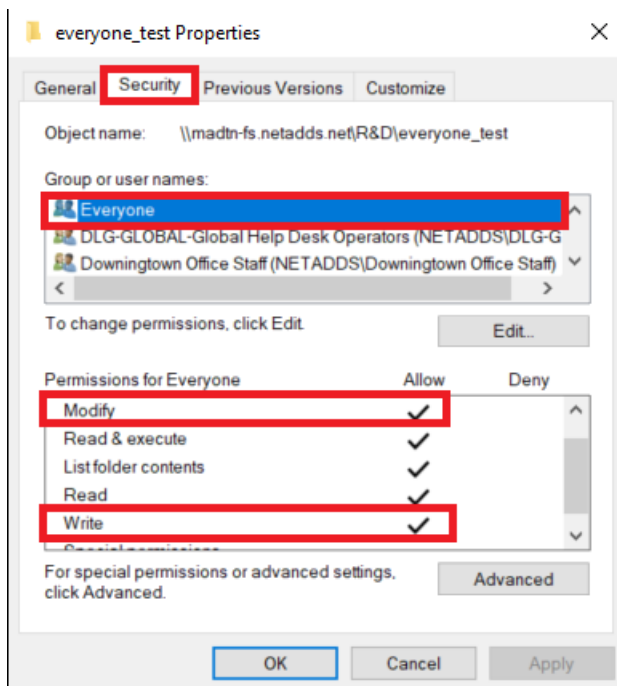
10. Click **Check Names**.

11. If multiple names are found, select **Everyone**, and click **OK**.



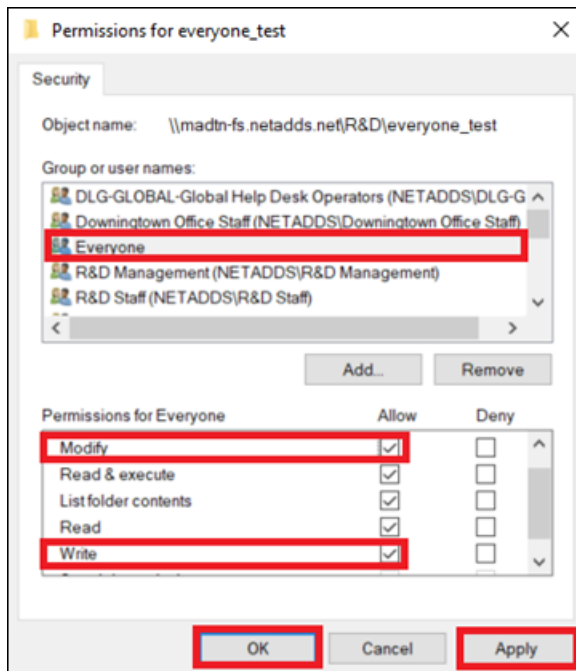
In the **Enter the object names to select** field, the word “Everyone” becomes underlined.

12. Click **OK**.
13. In the **Permissions for Everyone** section, confirm that the **Modify** and **Write** permissions both have checkmarks under **Allow**.



14. If both permissions are allowed, you have completed this procedure. Otherwise, continue to the next step.

15. If one or both permissions are not allowed, click **Edit**.



16. In the **Permissions for Everyone** section, select the **Modify** and **Write** check boxes under **Allow**.

17. Click **OK**.

18. Click **OK**.

Next, you need to map the UNC path for storage in software. See the procedure titled “Mapping a Folder for Storage” in the CellReporterXpress Help under **Configuration Mode > Data Storage**.