

Remote Tools for Simple Western

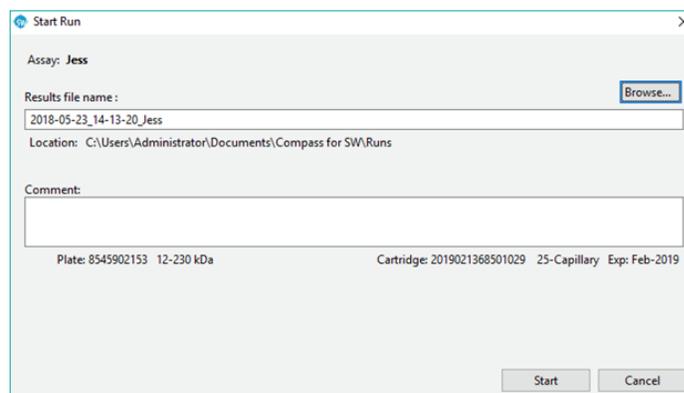
Introduction

The instructions below will enable users of Simple Western™ instruments to direct Compass for Simple Western software to automatically save data in a shared folder on a network drive to be accessed remotely. This will allow users to remotely monitor, access and analyze data after starting the run without having to return to the lab. These instructions are applicable to all supported Simple Western platforms (Jess™, Wes™, Sally Sue™, Peggy Sue™, NanoPro™ 1000) provided the following conditions are met:

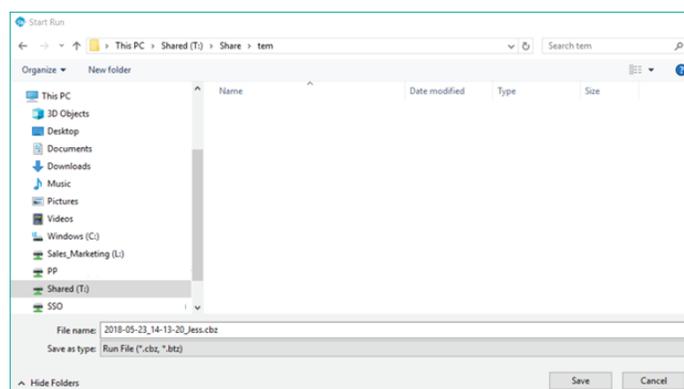
- The local computer for the Simple Western instrument is connected to the institution or company network and has a stable network connection.
- Users have writing permission to save data to the specified shared folder on the network drive.
- Users have access to the shared folder on the network drive when working from home. Users have permission to copy data from the network drive to a computer at home for data analysis, or have permission to perform data analysis directly on the network drive.
- Compass for Simple Western remembers the last location run data is saved. Data will be saved in the same folder on the network drive provided other users do not manually change the file save location from a previous run.
- Data saved on the network drive may be corrupted if network connectivity is interrupted during a run. However, a copy of the original run data is securely saved on the instrument. Users can retrieve the run data from the instrument via the Compass for SW>Runs folder on the local computer in the lab should the network be interrupted during a run.

INSTRUCTIONS FOR SAVING DATA IN A NETWORK DRIVE FOR REMOTE ACCESS

1. When setting up a run, change the directory where the data file will be stored by clicking **Browse...** in the **Start Run** window.



2. Navigate to the desired folder on the network drive, enter your desired **File name** and click **Save**.



3. Click **Start**. When the run is complete, data will be automatically saved to the specified folder location on the network drive.

INSTRUCTIONS FOR ENABLING TWITTER ACCESS

Users may link a Twitter account to any supported Simple Western platform to tweet the status of the run. Tweets may indicate that the run has started, the run has completed, or the run has an error. We recommend setting up a separate Twitter account for Jess, Wes, Sally Sue, Peggy Sue or NanoPro 1000. This lets multiple people in the lab follow run progress. It also lets you send tweets directly from the instrument to all users, for example to notify others when the instrument is available or when an error has been reset, etc.

1. Select **Edit** in the **Main Menu**.
2. Select **Preferences**.
3. Select **Twitter**.
4. Fill in the appropriate Twitter settings and select **OK**.
5. Refer to pages 438–443 of the [Compass for Simple Western User Guide](#) for more detailed instructions on configuring the Twitter settings.

