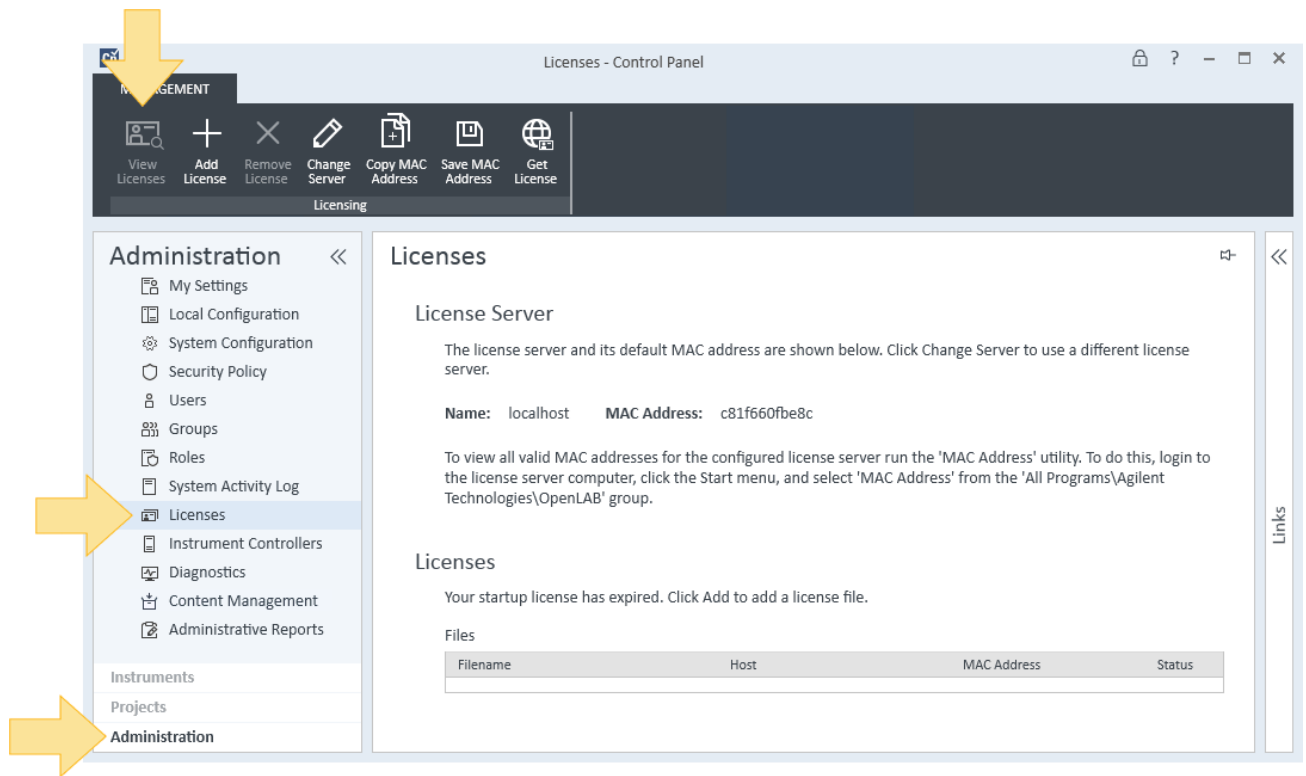




# View licenses

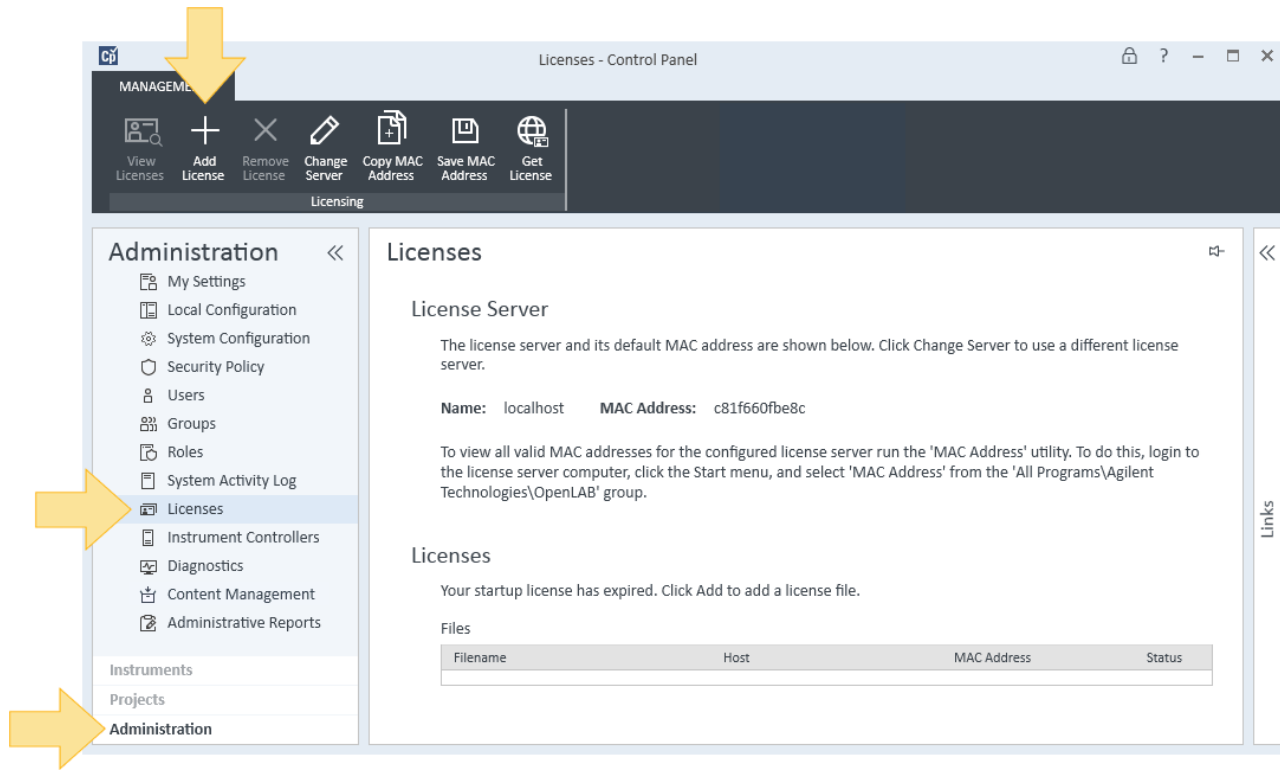
Click **Administration > Licenses > View Licenses**.



The information will display in an internet window.

# Add a license

1. Click **Administration > Licenses > Add Licenses**.

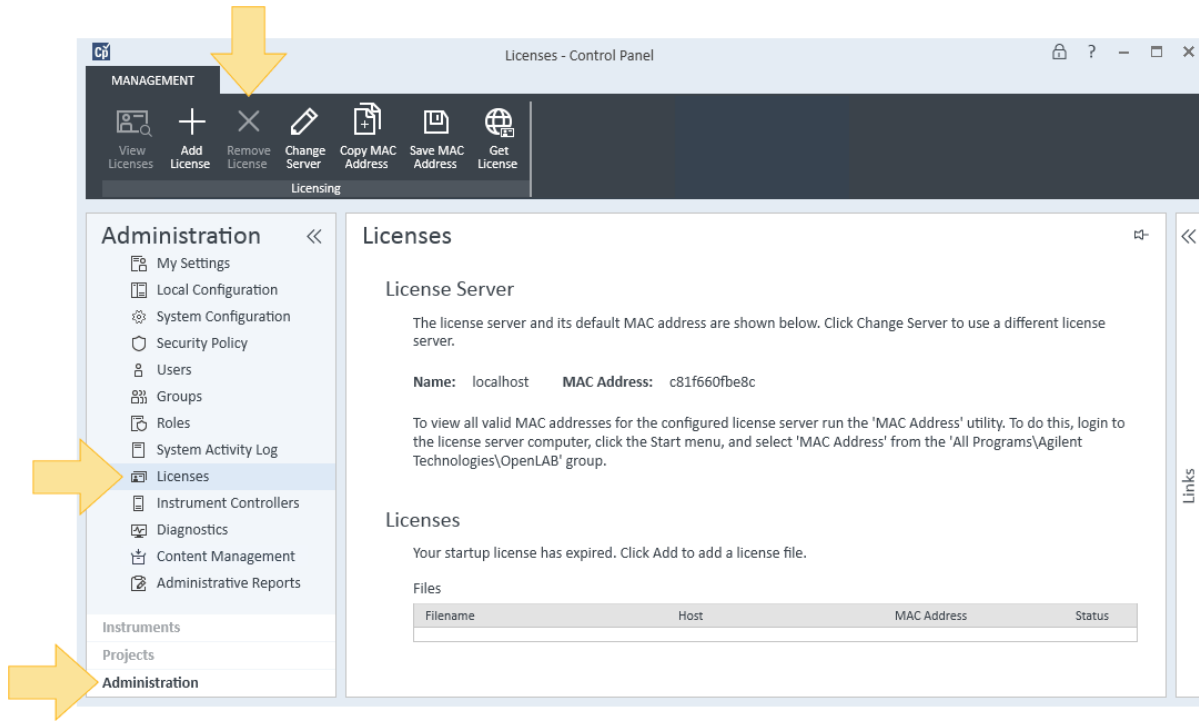


2. You can:

- **Select a license file**
  - a. Click **Browse**.
  - b. Navigate to and select the file.
  - c. Click **Open**.
- **Paste the contents of a license file**
  - a. Enter the license text.
  - b. Click **OK**.

## Remove a license

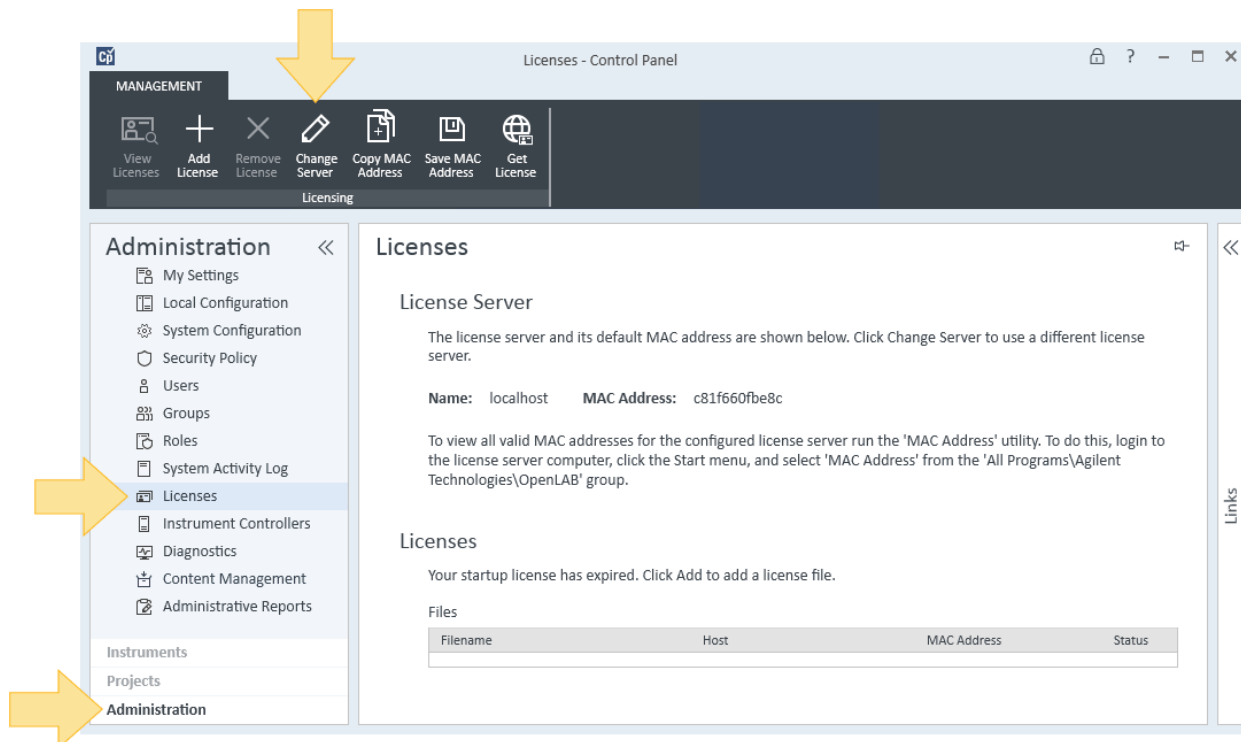
1. Click **Administration > Licenses**.
2. Select a license to remove and click **Remove License**.



3. Click **Yes**.

## Change the license server

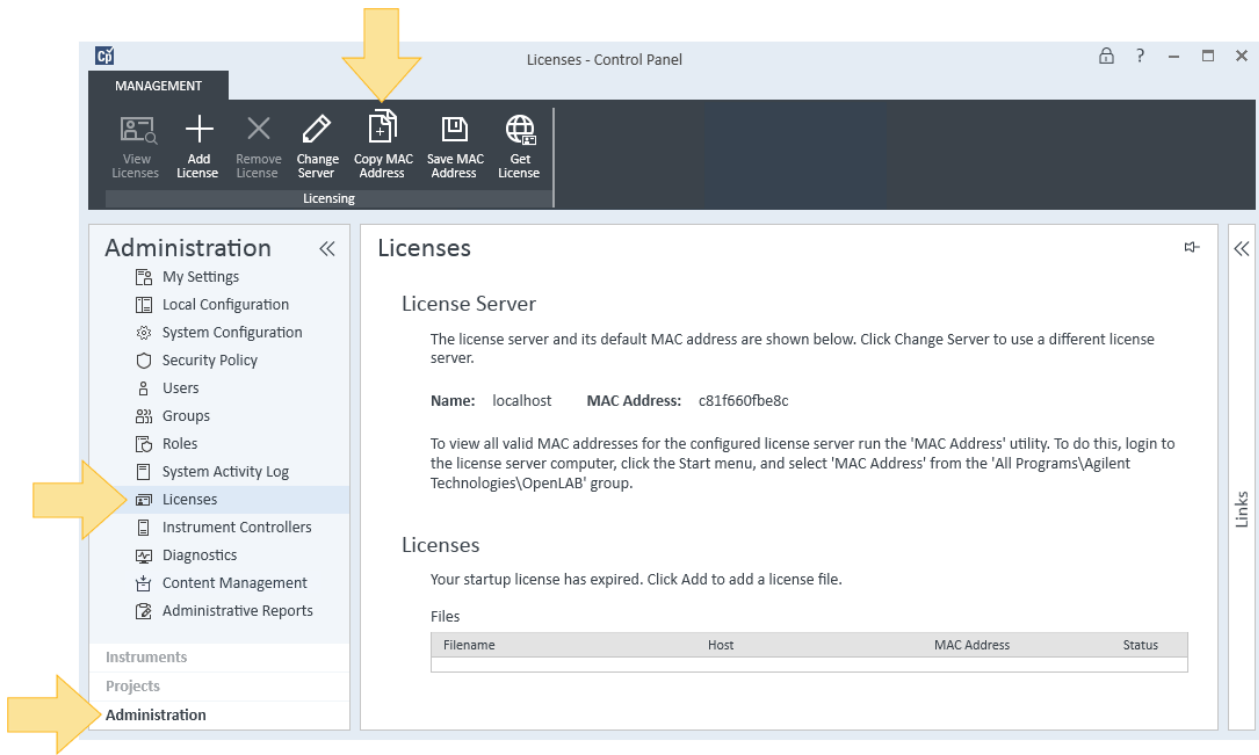
1. Click **Administration > Licenses > Change Server**.



2. Enter the computer name for your **License server**.
3. Click **Ping** to verify the connection.
4. Click **OK**.

## Copy the MAC address

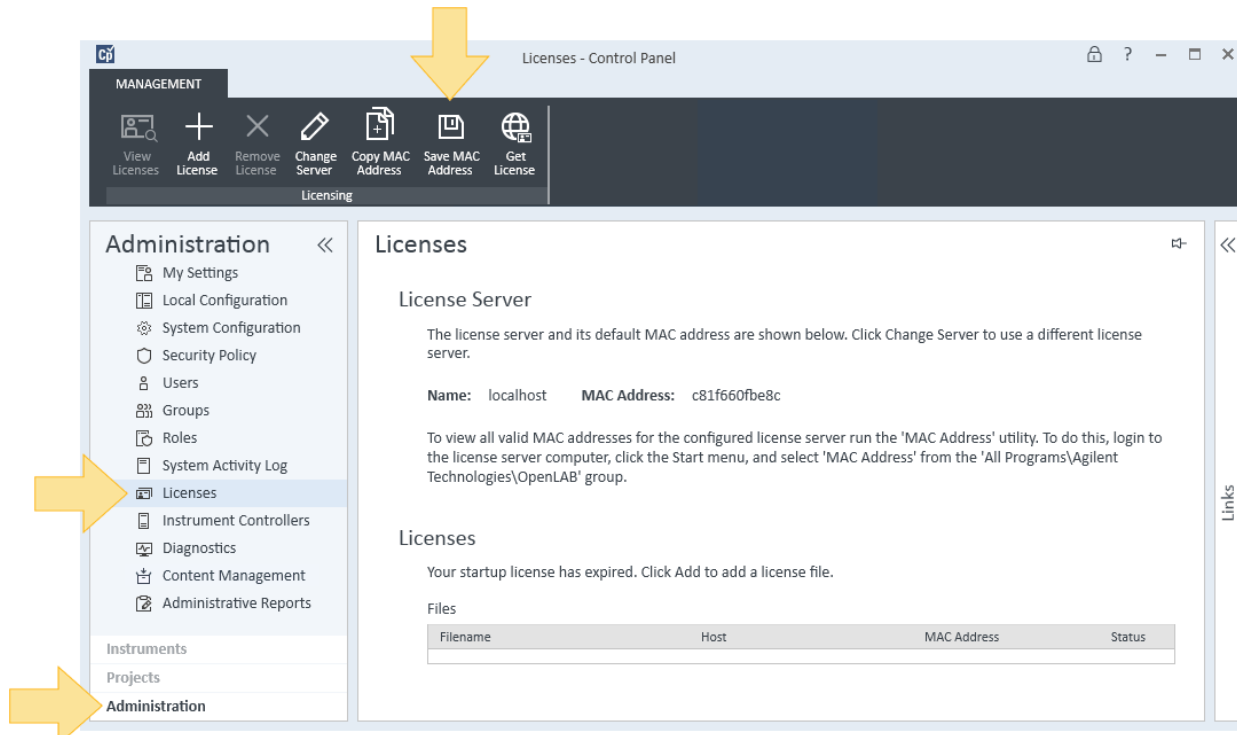
1. Click **Administration > Licenses > Copy MAC Address**.



2. Click **Okay**.

## Save the MAC address

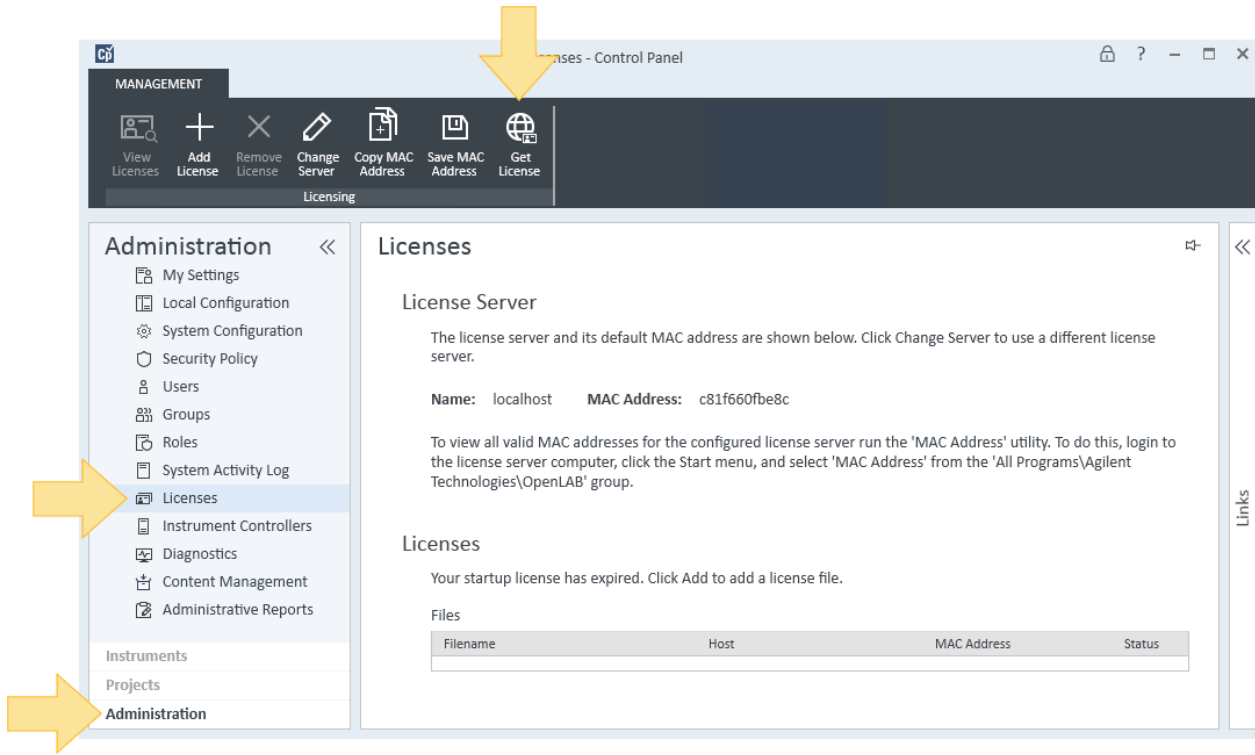
1. Click **Administration > Licenses > Save MAC Address**.



2. Select a location to save the License Server MAC Address as a text file.
3. Click **Save**.

# Get a license

1. Click **Administration > Licenses > Get License**.



Agilent.subscribenet.com opens in your browser.

2. Log on to Electronic Software and License Delivery



[www.agilent.com/chem/sample-scheduler-for-openlab](http://www.agilent.com/chem/sample-scheduler-for-openlab)

DE44343.3230208333

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021  
Published in the USA, June 1, 2021  
5994-3647EN

