

Compass Authorization Server for iCE Installation Guide

Before You Start

This document assumes that you are installing **Compass Authorization Server (CAS) for iCE 3.0.0** on a PC that has an older version of CAS for iCE already installed. If this is a new installation, all you need to do is run the installer and follow the prompts.

Upgrading the Compass Authorization Server

1. Why do I need to upgrade to Compass Authorization Server for iCE 3.0.0?

Compass Authorization Server for iCE 3.0.0 provides enhanced security features. Previous version of CAS were built on Python-Version 2.7. As of January 1, 2020, Python.org announced that Python 2 versions are no longer supported. CAS for iCE 3.0.0 is built on Python 3.x, and thus avoids any security vulnerabilities reported in older Python versions.

CAS for iCE 3.0.0 also supports HTTPS, which protects the communication between your Compass for iCE software and the server from being intercepted and tampered with.

2. Do I also need to update Compass for iCE software?

Upgrade to Compass for iCE version 3.0.0 prior to upgrading the Compass Authorization Server. You can download the latest Compass for iCE installer from the Bio-Techne website or contact our Technical Support team.

- 3. How do I upgrade to Compass Authorization Server for iCE 3.0.0?
 - a. Close Compass for iCE software if it is already running. Log out of the Compass Authorization Server and close the browser window.

- b. If you are using LDAP, Compass Authorization Server for iCE requires a **secure** connection.
 - If you are using LDAP but it is not secure, do not upgrade to CAS for iCE 3.0.0 as it will not support insecure connections.
 - If you have a secure LDAP connection, go to your LDAP settings and change the port number to 636 prior to upgrading or installing CAS for iCE 3.0.0.

Note: You may use CAS for iCE 3.0.0 without LDAP.

- Download the Compass Authorization Server for iCE 3.0.0.exe installer from the Bio-Techne website or contact our Technical Support team.
- d. Double click on the installer and follow the installation prompts.
- e. After a successful installation, the following window will open in your browser:



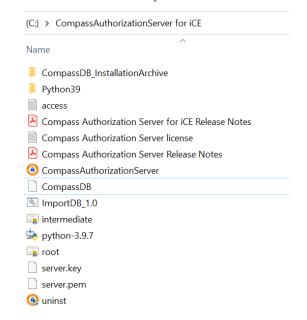
Note: Ignore any browser warnings – we recommend using Google Chrome.

f. Log in with the administrator account.

Note: default server administrator: Username: admin Password: admin.

4. What happens during the installation?

- a. Any older versions of Python installed on the PC will be removed.
- Python 3.9, along with CAS for iCE 3.0.0, will be installed.
- c. The existing CompassDB database file will be migrated automatically to the new folder named 'CompassAuthorizationServer for iCE'.
- d. A copy of the CompassDB database file will also be stored in the installation archive in a separate sub-folder named 'CompassDB_InstallationArchive'.



Frequently Asked Questions

1. Can two different Python versions be installed on the same PC?

No. During the installation of CAS for iCE 3.0.0, the installer checks for any previous Python versions and will remove them automatically.

2. Can CAS 2.2.x and CAS for iCE 3.0.0 be installed on the same computer?

No. When CAS 3.0.0 is installed, CAS 2.2.x will be overwritten.

3. Can I copy and paste the database file between different versions of Compass for iCE?

- No. The database file for CAS 2.2.x will not be recognized by CAS for iCE 3.0.0. When CAS 3.0.0 is installed over an existing database, an updated database file will be created
- The contents of older databases created on CAS versions prior to 3.0.0 cannot be accessed by CAS for iCE 3.0.0. To access this data, we recommend that you install the version of CAS that the database was last modified with on a separate PC (installation of an older version of CAS on the same PC would require uninstalling CAS for iCE 3.0.0).

bio-techne[®]

