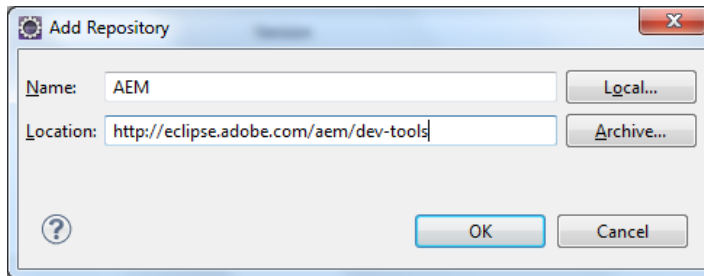


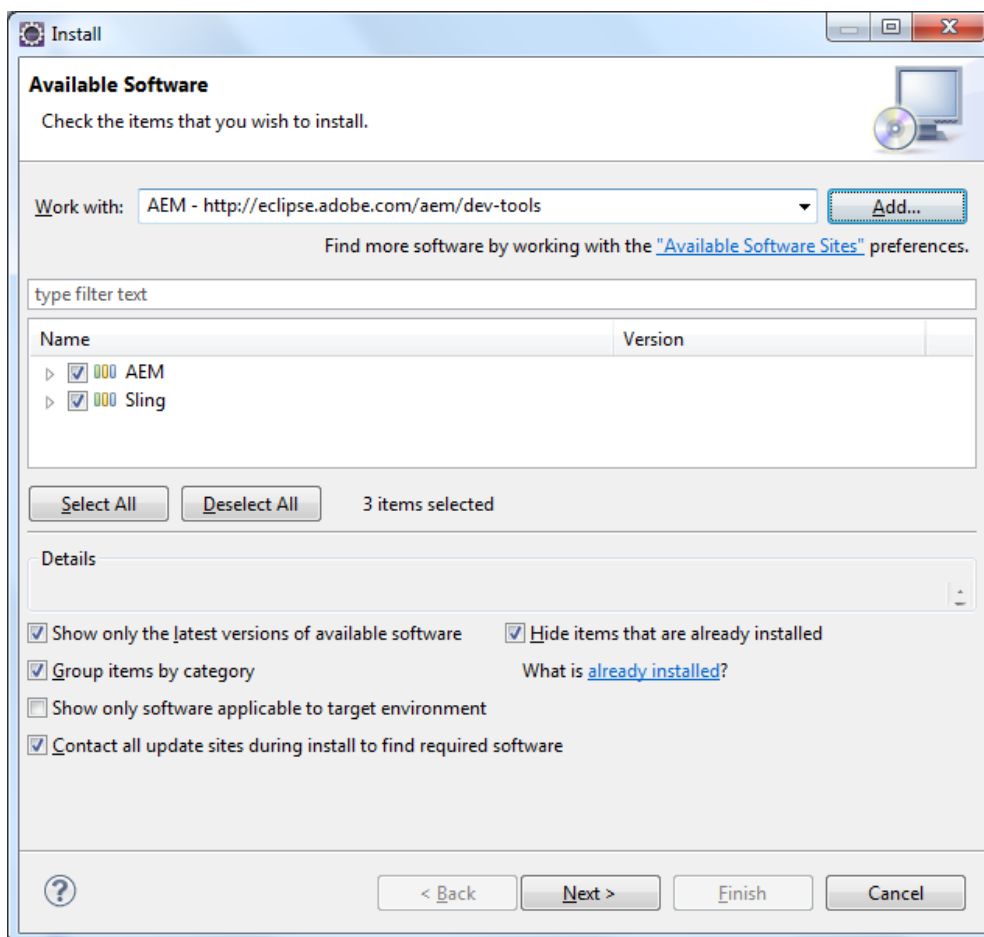
AEM plugin for eclipse

To install the plugin follow the below step.

- In Eclipse, open the Help menu.
- Click Install New Software then click on **Add**.

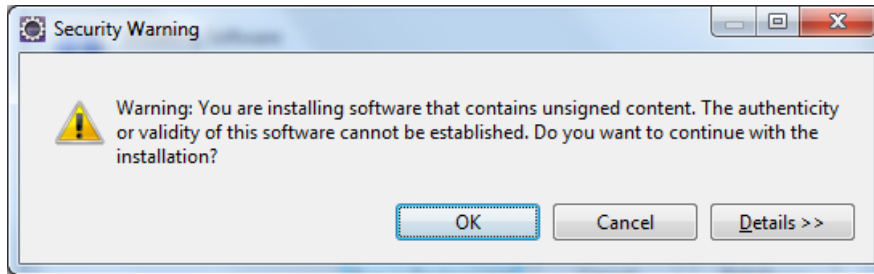


- In Name type AEM
- In Location copy the installation URL: **http://eclipse.adobe.com/aem/dev-tools**
- Click **OK**.
- Check both AEM and Sling plugins.



- Click **Next**.
- Click **Next**.

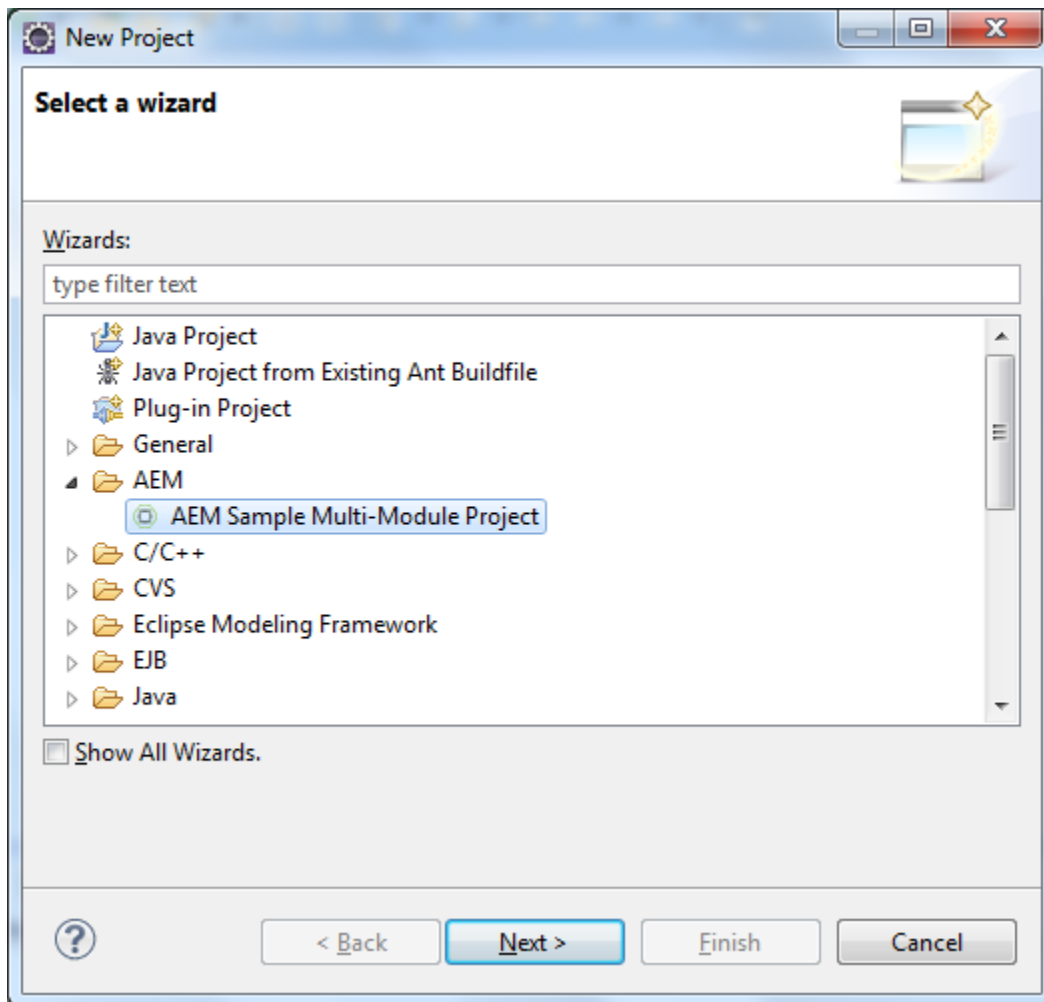
- ✚ Accept the license agreements and click **Finish**.
- ✚ If any security warning popup, Click **OK**.



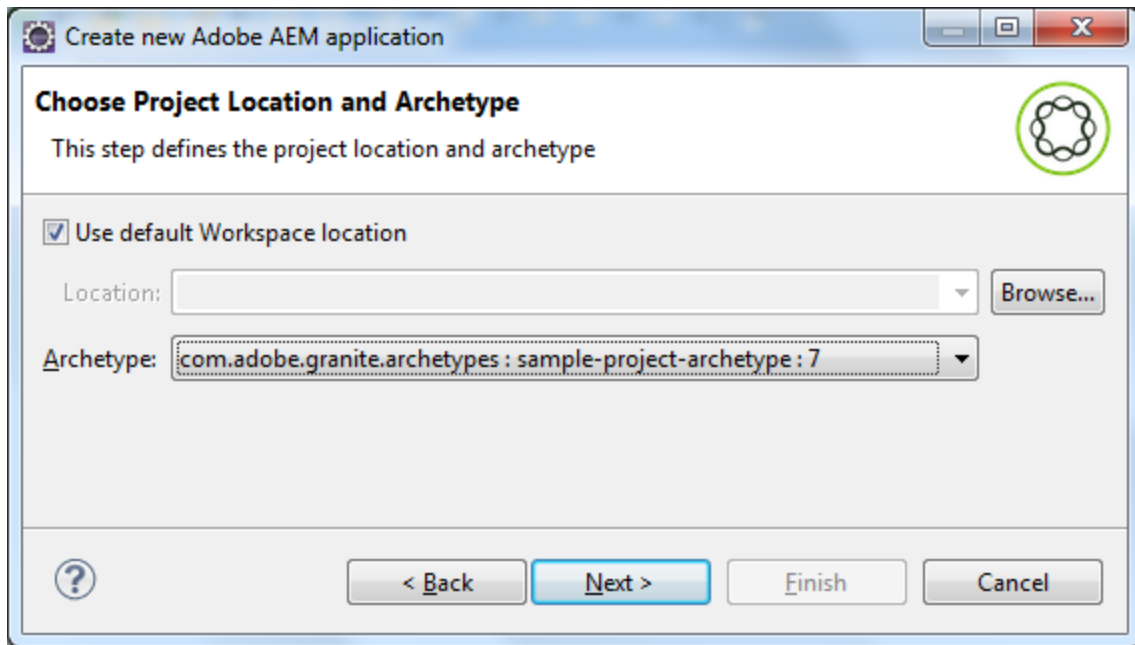
- ✚ Click **Yes** in order to restart Eclipse.

To create project follow the below steps.

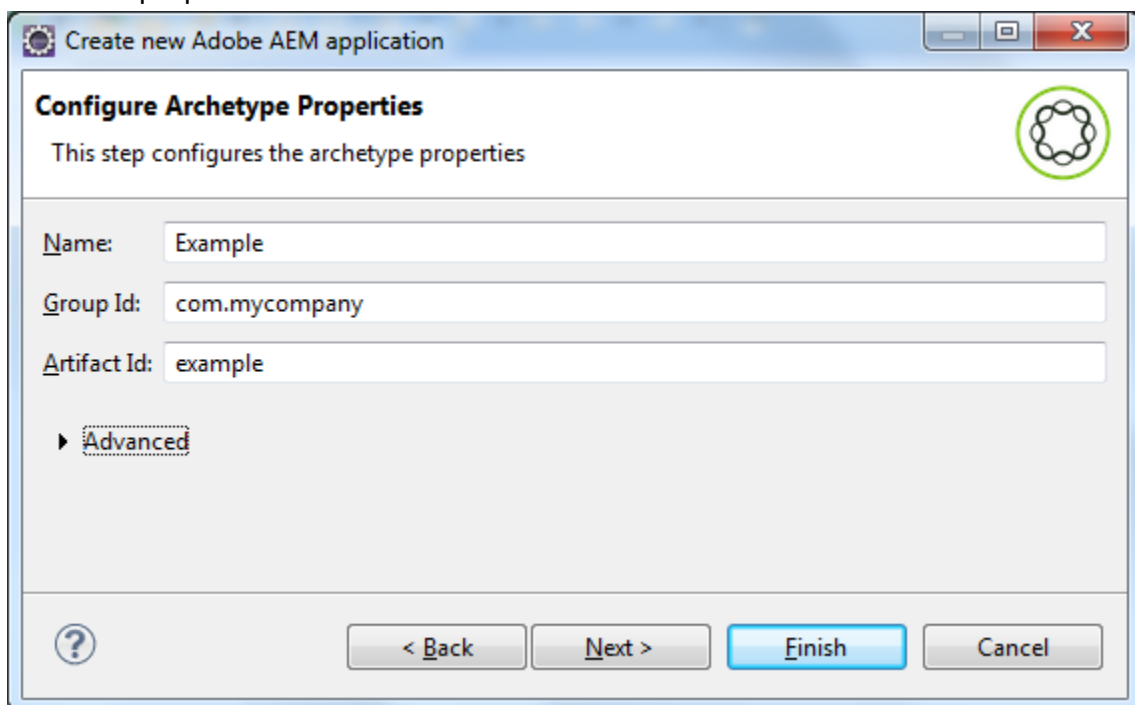
- In the File > New > Project menu, browse to the AEM section and select AEM Sample Multi-Module Project.



- Click **Next**.
- Choose **com.adobe.granite.archetypes : sample-project-archetype : 7** from the menu, then click **Next**.



- Click **Next**.
- Fill in a **Name**, **Group id** and an **Artifact id** for the project. You can also opt to set some advanced properties.



- You should then configure an AEM server to which Eclipse will connect.

Create new Adobe AEM application

Select or Create Server

This step defines which server to use with the new project.

☐ Add to existing server

Location:

☒ Start server after project creation (if server not yet started).

☒ Setup new server

Server name:

Host name:


Port:

Username:

Password:

Debug Port:

☒ Check/Install org.apache.sling.tooling.support.install bundle



Click **Finish**. The project structure is created.