

SkaniIt™ Software Installation for Windows™ 10 Operating System

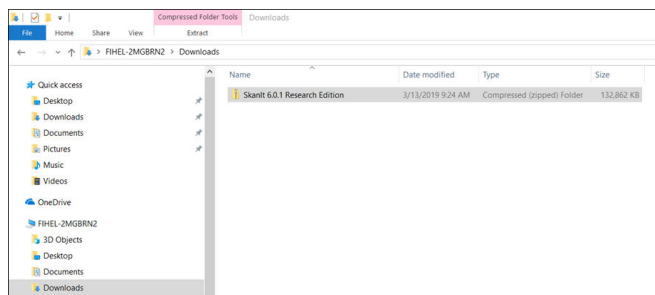
Pub. No. MAN0018651 Rev. A.0

When this installation package is downloaded onto your computer, the .zip file is stored in the Downloads folder. When Thermo Scientific™ SkaniIt™ Software is being installed from the .zip file using the Windows™ 10 operating system:

- Contents of the .zip folder **must** be extracted before running the installation. See **step-by-step** instructions below.
- Running the software installation from the .zip file without extraction may lead failure of the SkaniIt™ Software installation process.

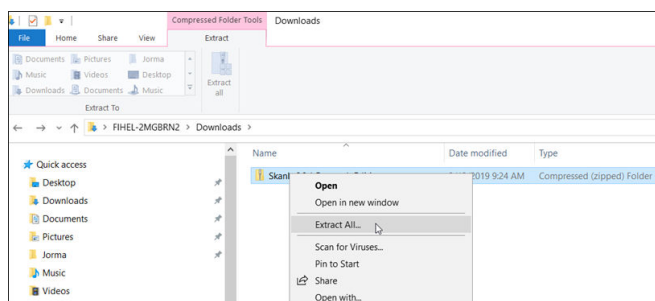
Install the software

1. Download the SkaniIt™ Software package. It will be stored in the **Downloads** folder.

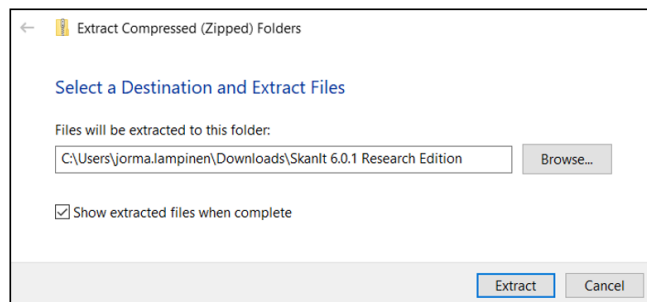


As shown, the folder type is compressed, so files need to be extracted before running the installation.

2. Select the compressed folder, right-click using the mouse and select **Extract All**.

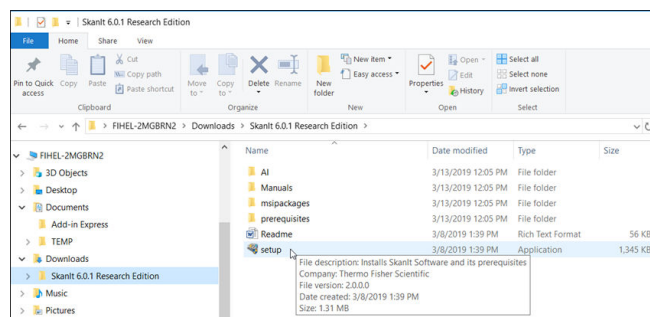


3. Select the destination folder where compressed folders and files are extracted and click **Extract**.



Files and folders are extracted and shown in the destination folder.

4. Select setup from this folder to start the SkaniIt™ Software installation.



Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

thermofisher.com/support | thermofisher.com/askaquestion



Thermo Fisher Scientific Oy | Ratatie 2 | FI-01620 Vantaa | Finland

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For Research Use Only. Not for use in diagnostic procedures.