

Smart Deep™ Basecaller License

Activate the Smart Deep™ Basecaller license in Sequencing Analysis Software 8

Catalog Numbers A53212, A53213, A53215, A53216

Publication Number MAN0025944 Revision A.0

■ About Smart Deep™ Basecaller	1
■ About the Smart Deep™ Basecaller license	2
■ Activate the Smart Deep™ Basecaller license	3

This Quick Reference provides brief information on the Smart Deep™ Basecaller license. For more details about Sequencing Analysis Software 8 and the Smart Deep™ Basecaller license, see the *Sequencing Analysis Software 8 User Guide* (Pub. No. MAN0019595).

About Smart Deep™ Basecaller

IMPORTANT! A Smart Deep™ Basecaller license is required to use Smart Deep™ Basecaller in Sequencing Analysis Software 8. See “About the Smart Deep™ Basecaller license”.

Smart Deep™ Basecaller is compatible with AB1 (.ab1) files generated on the 3500/3500xL Genetic Analyzer, 3730/3730xL DNA Analyzer, SeqStudio™ Genetic Analyzer, and SeqStudio™ Flex Series Genetic Analyzer. Smart Deep™ Basecaller is not compatible with 3130, 3100, or 310 instrument data.

Smart Deep™ Basecaller outperforms KB™ Basecaller in the following ways.

- Longer read length.
- More accurate pure and mixed base basecalls.
- Improved basecalling accuracy through heterozygous insertion deletion variants.
- More robust basecalling results through dye blob, mobility shift, and n-1 peak artifacts.
- Optional view for plasmid sequences with cleaned up baseline and improved 3' resolution of the trace.

About the Smart Deep™ Basecaller license

Smart Deep™ Basecaller demonstration license

A 90-day Smart Deep™ Basecaller demonstration license is included with your purchase of Sequencing Analysis Software 8.

IMPORTANT! A license key will be sent to your email. The license key is used to register the Smart Deep™ Basecaller license on your computer.

The demonstration license expires 90 days after the license is activated. Each Smart Deep™ Basecaller license is tied to the computer ID.

You can purchase the Smart Deep™ Basecaller license through the Thermo Fisher Scientific website (<http://thermofisher.com>) or your local sales office. For more information, see “Purchase options for the Smart Deep™ Basecaller license”. If you have questions or need help accessing your license key, email Support at:

seqa.support@thermofisher.com

Purchase options for the Smart Deep™ Basecaller license

To continue using Smart Deep™ Basecaller after the demonstration license expires, purchase the Smart Deep™ Basecaller license through the Thermo Fisher Scientific website (<http://thermofisher.com>) or your local sales office.

You have the option to purchase a 1-year or a 3-year license. After you purchase a license, a license key will be sent to your email. The license key is used to activate the Smart Deep™ Basecaller on your computer. If you have questions or need help accessing your license key, email Support at:

seqa.support@thermofisher.com

Table 1 Smart Deep™ Basecaller license options

Cat. No.	Description
A53212	New Smart Deep™ Basecaller license (1-year)
A53213	New Smart Deep™ Basecaller license (3-year)
A53215	Renew Smart Deep™ Basecaller license (1-year)
A53216	Renew Smart Deep™ Basecaller license (3-year)

Note: The license is valid until the expiration date. If you need to uninstall, then reinstall the software on the same computer, use the same license until the expiration date. Each Smart Deep™ Basecaller license is tied to the computer ID.

Activate the Smart Deep™ Basecaller license

Note: For detailed information, see the *Sequencing Analysis Software 8 User Guide* (Pub. No. MAN0019595).

To activate the Smart Deep™ Basecaller license, you must complete the following.

1. Install and register Sequencing Analysis Software 8.
2. Add .ab1 files, select **SDB.bcp** from the **BaseCaller** dropdown menu, then start sample analysis.
The **Warning: Smart Deep Basecaller** window is displayed.
3. In the **Warning: Smart Deep Basecaller** window, click **Yes**.
The license activation window appears, automatically displaying the Computer ID, which cannot be edited.
4. In the activation window, input the license key you received by email.
If you have questions or need help accessing your license key, email Support at:
seqa.support@thermofisher.com
5. Complete the activation of the Smart Deep™ Basecaller license based on your Internet connection.

Internet connection	Steps to activate the license
(Recommended) The computer is connected to the Internet	<ol style="list-style-type: none">1. In the activation window, click Yes. Connected. The Success: Smart Deep Basecaller window appears, confirming successful application of the Smart Deep™ Basecaller license.2. In the Success: Smart Deep Basecaller window, click OK.
The computer is not connected to the Internet	<ol style="list-style-type: none">1. In the activation window, click No. Not connected.2. When prompted, save the SampleRequest.bin file.3. Email the saved SampleRequest.bin file to Support at: seqa.support@thermofisher.com A CapabilityResponse.bin file will be emailed to you. Download this file.4. In the activation window, upload the CapabilityResponse.bin file, then select Install and Validate License.

The Smart Deep™ Basecaller license is activated. The Smart Deep™ Basecaller is ready to be used for analysis.

Note: The license is valid until its expiration date. If you need to uninstall, then reinstall the software on the same computer, use the same license until its expiration date.



Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256

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

Revision history: Pub. No. MAN0025944

Revision	Date	Description
A.0	14 November 2022	New document for the Smart Deep™ Basecaller license.

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.

NOTICE TO PURCHASER: DISCLAIMER OF LICENSE: Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under patent claims owned or otherwise controlled by Thermo Fisher Scientific, either expressly, or by estoppel.

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2022 Thermo Fisher Scientific Inc. All rights reserved.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

14 November 2022

ThermoFisher
SCIENTIFIC