

Security Bulletin

SUMMARY

CVE: Unknown File Modification Risk for Applied Biosystems™ Human Identification Software

PUBLISHED DATE:	CVE NUMBER:	CVSS V4.0 SCORE:	SEVERITY:
JULY 31, 2026	CVE-2026-17583	8.2	High

DESCRIPTION

Thermo Fisher Scientific has been made aware of a risk for nearly undetectable modification of .fsa/.hid file types generated by select Applied Biosystems™ Human Identification software. The risk is that nearly undetectable modifications to .fsa/.hid file outputs could occur before the file is loaded into analysis software if laboratory controls are circumvented.

SOLUTION

Thermo Fisher has developed security updates to address the risk for the software identified below (see Table 1). The security updates implement the use of digital signatures to add an extra layer of protection that, moving forward, will help customers verify that data files have not been modified.

For customers who are unable to implement the updates or use another third-party analysis platform, Thermo Fisher Scientific continues to recommend, consistent with best practices, evaluating the following measures alongside other established laboratory procedures, in alignment with your organization's security measures and policies:

- Maintain a secure chain of custody for files generated by the HID instrumentation throughout the analysis workflow.
- Store generated files on encrypted, password-protected storage media (for example, encrypted USB drives or encrypted hard drives).

- Restrict access to generated files to authorized personnel in accordance with your laboratory's access control policies.
- Apply the principle of least privilege by limiting user permissions on systems operating the HID instrumentation or hosting associated data analysis and secondary analysis software.
- Leverage firewall rules and network access control lists (NACLs) to restrict internet connectivity to only trusted sources.

Please contact your local Thermo Fisher Scientific support representative for any additional questions or contact product.security@thermofisher.com for any questions related to the issue.

Thermo Fisher Scientific is committed to maintaining the security, integrity and reliability of our products. We appreciate your partnership and recommend implementing the applicable updates as soon as practical.

ACKNOWLEDGEMENTS

Thermo Fisher Scientific would like to thank Nathan Adams, Kevin Dyer and Laura Gaydosh Combs as well as the Cybersecurity and Infrastructure Security Agency (CISA) for their identification and coordinated disclosure of the issue covered in this security bulletin.

Table 1. Software List & Updated Software

Software	Updated Software Versions
Applied Biosystems™ 3500/3500xL Series Data Collection Software version 4.0.2 and earlier.	Download the Applied Biosystems™ 3500/3500xL Series Data Collection Software version 4.0.3 update. Please refer to the specific software release notes for more information.
Applied Biosystems™ 3730/3730xL Series Data Collection Software version 5.0.2 and earlier.	Download the Applied Biosystems™ 3730/3730xL Series Data Collection Software version 5.0.3 update. Please refer to the specific software release notes for more information.
Applied Biosystems™ SeqStudio™ Genetic Analyzer Data Collection Software version 1.2.5 and earlier.	Download the Applied Biosystems™ SeqStudio™ Genetic Analyzer Data Collection Software version 1.2.6 update. Please refer to the specific software release notes for more information.
Applied Biosystems™ SeqStudio™ Flex Series Instrument Software version 1.2.0 and earlier.	Download the Applied Biosystems™ SeqStudio™ Flex Series Instrument Software version 1.2.1 update. For customers using the SeqStudio™ Flex system with SAE enabled, the user must first install the latest SAE profile on the SAE Admin Console before installing the update. Download the latest SAE profile using the link here. Please refer to the specific software release notes for more information.
Applied Biosystems™ GeneMapper™ ID-X Software version v1.7.3 and earlier	Download the Applied Biosystems™ GeneMapper™ ID-X Software version v1.7.4 update. Please refer to the specific software release notes for more information.
Applied Biosystems™ 3130 Series Data Collection Software version 4.1 and earlier.	As this software has reached its end-of-life and is no longer supported by Thermo Fisher Scientific, no update will be provided.
ABI PRISM® 3100/3100-Avant™ Data Collection Software version 2.0 and earlier.	As this software has reached its end-of-life and is no longer supported by Thermo Fisher Scientific, no update will be provided.
ABI PRISM® 310 Data Collection Software version 3.1 and earlier.	As this software has reached its end-of-life and is no longer supported by Thermo Fisher Scientific, no update will be provided.