

# Agilent Overview Management Server: Generating Reports

## Introduction

Agilent Overview management server facilitates monitoring and administration of a fleet of Agilent Insight200M liquid explosive detection systems (LEDs). Real-time information for system checks, maintenance, and overall alarm breakdown is readily available on the dashboard. Further analysis can be provided by scheduling and generating regular reports from the report wizard. The data can be filtered by machine, group, or location. These reports can be used to demonstrate European Civil Aviation Conference (ECAC) compliance, provide data on false alarm occurrences, and offer information on the general Insight health across airports.

## Easy management of data

The report wizard allows easy management of data and report generation. Generated reports are visible in the interface and are readily available to print to PDF and CSV. For example, the ECAC audit report is quickly populated with the relevant information in the dashboard, and can be printed to PDF for distribution, or downloaded as a CSV sheet for internal logging (Figure 1).

## Investigate false alarm rate

False alarm rates provide an indication of how many items are alarming incorrectly. These alarms still require an escalation process, which can cost time and money, and can cause passenger dissatisfaction. The Agilent Insight200M has the lowest false alarm rate on the market, and offers 100% threat detection. Reviewing the false alarm rate of an Insight can also provide information in case refresher training of personnel is required, or to inform the user if maintenance of the system is needed. False alarm rate reports can be obtained through the Overview report wizard using a prepopulated template (Figure 2) or by using the custom reporting option (Figure 3).

In both instances, the report can be scheduled for creation on certain dates, shared with others in Overview, and exported via PDF or CSV.

From the Dashboard...

**Report - Oct21 Feb22 ECAC**  
Generated at 2023-07-07 12:04:00

Serial	LastSync	SoftVerSystem	HardVerDetection	DetectedModelVer	ModelNum
INS200_82062	2025-12-06 09:41:23	3.2.0.150	LAGS Type B Insight200M	--	
INS200_83007	2025-12-04 13:50:16	3.2.0.150	LAGS Type B Insight200M	--	
INS200_82103	--	3.0.2.275	LAGS Type B Insight200M	--	

Download as an Excel spreadsheet

	A	B	C	D	E	F
1	Serial	LastSync	SoftVerSystem	HardVerDetection	DetectedModelVer	ModelNum
2	INS200_82062	2025-12-06 09:41:23	3.2.0.150	LAGS Type B Insight200M	--	
3	INS200_83007	2025-12-04 13:50:16	3.2.0.150	LAGS Type B Insight200M	--	
4	INS200_82103	--	3.0.2.275	LAGS Type B Insight200M	--	

Figure 1. (ECAC) report, printed to PDF and downloaded as a CSV.

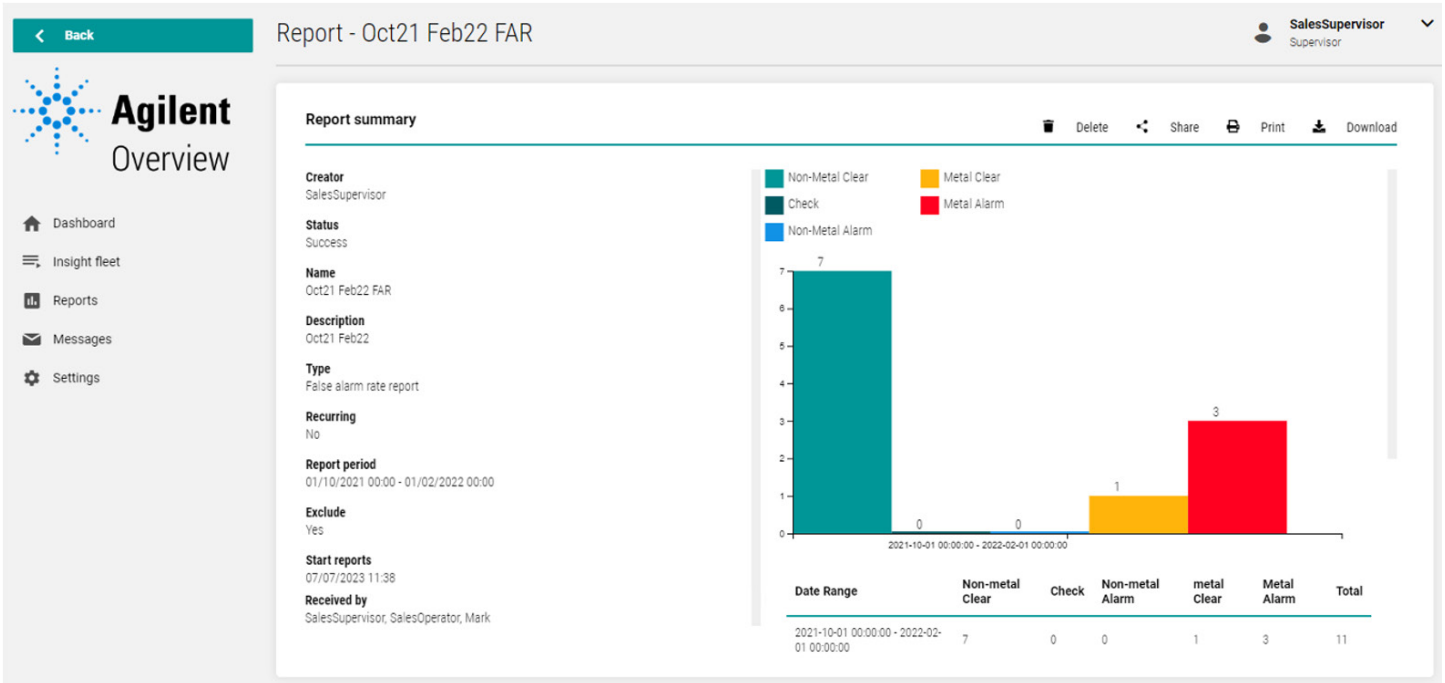


Figure 2. Prepopulated template false alarm report (FAR).

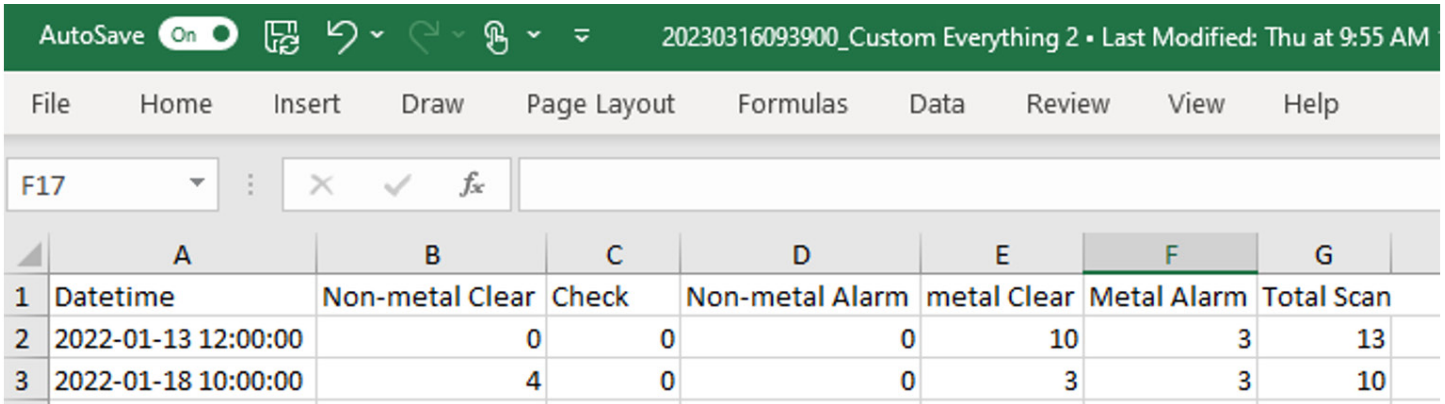


Figure 3. Example of a custom report.

## Engineering tracking

Maintenance of Insights is important to ensure both optimal usage and the prevention of downtime of a unit on a lane. Overview allows in-depth engineer assessment of the following:

- Location of Insights
- Software versions
- Configuration management
- ECAC auditing
- Disk space

- User setup and password management (if using Microsoft Windows 10 on the Insight200M and cybersecurity mode)
- System checks
- System health and diagnostics
- When system checks are due to be performed

Overview is also designed with preventative maintenance in mind (alerting engineers before maintenance is required), which provides opportunities for both optimal scheduling and cost-savings.

## Summary of reports

Agilent Overview helps to generate a variety of useful reports. These include reports for compliance with ECAC audits, plus reports for false alarm rate, and monitoring the health of the Insight fleet overall. The reporting wizard can offer both prepopulated and custom reports. All reports come readily populated, and can be printed to PDF for easy distribution, or downloaded to CSV for customized data analysis. Table 1 details the custom reports available in Overview.

**Table 1.** Summary table of all prepopulated reports.

Report Type	Information Provided in Report					Function
	Insight Location	Information by Specific Insight	Insight Information	Breakdown of Alarms	Operator Response Information	
ECAC Audit Report	No	Yes	Yes	No	No	Easily check software, hardware, and model version compliance against ECAC standard, per unit
False Alarm Report (FAR)	No	Yes	No	Yes	No	Compare metals and nonmetals as cause of false alarms
Custom Report	No	Yes	No	Yes	No	FAR report can be further broken down by timeframe
Configuration of Liquids, Aerosols, and Gels (LAG) Unit Report	Yes	Yes	Yes	Yes	Yes	Compare health of Insight with the number of false alarm occurrences and location within airport
Postponed Test Piece Report	No	Yes	No	No	Yes	Review operator training

## Conclusion

The Agilent Overview management server provides key capabilities for data tracking, false alarm rate investigation, monitoring of user training, and system health. Overview can be installed easily and securely, following the highest cybersecurity standards. It offers management of other cybersecurity features such as user login and password management (Windows 10 Insight200M only). Overview is a key tool for many engineering departments for managing and tracking the health of the fleet, both simply and cost-effectively.

[www.agilent.com](http://www.agilent.com)

DE17960523

This information is subject to change without notice.

© Agilent Technologies, Inc. 2023  
Printed in the USA, August 25, 2023  
5994-6703EN