

Environment Management Policy Statement



Approved by: Tom Mulder

Owner: Nichola Harrison

Date: 01/05/2025

Environmental Policy Statement

Malvern Panalytical recognises our responsibility to minimise the negative effect the Company's Business activities have on the environment. The Company has developed an Environmental Management System (EMS) to meet the requirements of ISO14001 and reinforce the Company's commitment to the prevention of pollution and the continual improvement of our environmental performance.

As a Spectris company, we share the same drive towards Net Zero. Malvern Panalytical commits to the following:

- Monitor and reduce the use of energy and raw materials by improved process efficiencies and by reviewing the opportunities of using recycled materials
- Reduce the production of waste and the release of harmful substances into the environment through elimination or substitution of materials where possible
- Managing, monitoring and measuring environmental performance and establishing objectives and targets to minimise these impacts.
- Comply with relevant environmental local and international regulations and legislations
- Communicating this policy to all employees, contractors and other stakeholders, as well as making this policy available to the public.
- Reporting on the company's environmental performance in both internal and external communications, where relevant.
- The EcoVadis rating system with a target to become a global sustainable organization and improve the level of sustainability.

Malvern Panalytical ensures that all people that work for or on behalf of the company are aware of and trained in their environmental responsibilities.

A handwritten signature in black ink, appearing to read 'Tom Mulder', positioned above a horizontal blue line.

Tom Mulder (May 1, 2026 13:20:21 GMT+2)

Tom Mulder
EVP Corporate Services

May 1, 2026

Review date 01/01/2027






Global Environment Policy Statement

Final Audit Report

2026-05-01

Created:	2026-05-01
By:	Nichola Harrison (nichola.harrison@malvernpanalytical.com)
Status:	Signed
Transaction ID:	CBJCHBCAABAA3IPM08EH1R-DDZbM-A3AfXlrkE4YOQej

"Global Environment Policy Statement" History

-  Document created by Nichola Harrison (nichola.harrison@malvernpanalytical.com)
2026-05-01 - 10:34:50 AM GMT
-  Document emailed to Tom Mulder (tom.mulder@malvernpanalytical.com) for signature
2026-05-01 - 10:35:43 AM GMT
-  Email viewed by Tom Mulder (tom.mulder@malvernpanalytical.com)
2026-05-01 - 11:19:32 AM GMT
-  Document e-signed by Tom Mulder (tom.mulder@malvernpanalytical.com)
Signature Date: 2026-05-01 - 11:20:21 AM GMT - Time Source: server
-  Agreement completed.
2026-05-01 - 11:20:21 AM GMT