

# Environment Management System Policy Statement



Author: Paul Prosser

File Name: Environmental Management System Policy Manual

Date: 24.10.2022

Page: 1 of 1

## Environmental Policy Statement

Malvern Panalytical recognises the responsibility it has to reduce the negative effect the Company's business activities have on the environment. The company has developed and integrated an Environmental Management System (EMS) to meet the requirements of BS EN ISO14001 and to reinforce the company's commitment to the **prevention of pollution** and the continual improvement in environmental performance. The EMS is used as a vehicle for the setting, measuring and monitoring of environmental objectives and operates in parallel with the Business Management System (BMS) introduced to ensure compliance with the demands of BS EN ISO 9001.

Malvern Panalytical complies with all relevant environmental legislation and regulations and other requirements to which the organisation subscribes related to its environmental aspects. The company is committed to the consideration of environmental issues being an integral part of our business strategy, and to ensuring that all activities which carrying a risk of pollution, in any form, are identified and eliminated, or controlled effectively within statutory limits.

Wherever possible the company monitors and **reduces** the use of energy and **raw materials** by improved process efficiencies and by reviewing the opportunities of using recycled materials. Through continual improvement objectives, the company improves processes to **reduce the production of waste**, and the release of harmful substances into the environment.

Malvern Panalytical meets the requirements of the WEEE legislation and ensures that all **waste is recycled** or incinerated to save waste going into land fill.

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

Environmental considerations form a fundamental part of the Company's future investment decisions.

The company supports the adoption of similar standards and commitment by its customers and suppliers.

Copies of this policy statement are freely available to the public and other interested bodies.

DocuSigned by:  
  
49C6BAD9C0104AB...

**Tom Mulder**  
Vice President Business Enablement Partners