

S0501SPUK

## Global Health & Safety Policy Statement

Approved by:

Terry Kelly

Owner:

Nichola Harrison

Date:

01/01/2025

**We at Malvern Panalytical are committed to ensuring our staff are not exposed to unnecessary risk which may impact their health or safety. We are also committed to ensuring the health and safety of those who may be impacted by our activities and operations.**

### Our Commitments

Our business purpose is to enable our customers to create a better world using our technologies. We want to do that in a way which is safe and healthy for our staff and those who may be impacted by our activities around the globe. By looking after our people, we can support the growth targets of our organisation.

Our values underpin our corporate strategy and create a framework in which we do business. Our values equally support us to look after our people.

#### Be true

- We are empowered to speak up when something isn't as safe or healthy as it could be. If you think there is an immediate risk to your safety or health or to that of someone else, stop work and speak to your line manager.

#### Own it

- We are all responsible for creating a healthy and safe work environment, whether we are at home, at a customer site or in our premises. We are accountable for our actions.

#### Aim high

- By continually improving, our culture moves forward and we take care of ourselves and each other. We share better ways of working and find solutions to hazards that could be controlled better.

- As a minimum we will meet mandatory local health and safety legislation and other relevant regulatory requirements. We aim high by looking to go beyond these and implement a global standard where possible.
- We will ensure health and safety is a key part of our decision-making process. This will ensure we continual look for ways to improve our control measures of hazards.
- We empower everyone to be true and speak up if we believe an unsafe or unhealthy situation exists. We will stop work if we believe there is an imminent and material risk to our staff and others.
- We will own it by following processes and procedures and wearing the necessary personal protective equipment. If we don't know or understand, we will ask someone who does.
- Our global health and safety team are our subject matter experts. They will work with the business to provide advice and guidance on adhering to legislation and internal requirements. We will actively seek their opinions.
- We operate a no blame culture – when things go wrong, we will use that to openly look for ways to improve.

*Terry Kelly*

Terry Kelly (Jan 9, 2025 13:44 EST)

**Terry Kelly**

**President**

*'We share responsibility for creating an environment that is healthy and safe and we take ownership of our actions.'*

Review Date: 01/01/2026


# MP Global Health & Safety Policy English 2025

Final Audit Report

2025-01-09

Created:	2025-01-06
By:	Nichola Harrison (nichola.harrison@malvernpanalytical.com)
Status:	Signed
Transaction ID:	CBJCHBCAABAANKnPoKoEO6BXQKE-ahI2NPXM3UEFXfbWG


## "MP Global Health & Safety Policy English 2025" History

 Document created by Nichola Harrison (nichola.harrison@malvernpanalytical.com)

2025-01-06 - 2:28:01 PM GMT

 Document emailed to terry.kelly@malvernpanalytical.com for signature


2025-01-06 - 2:29:35 PM GMT

 Email viewed by terry.kelly@malvernpanalytical.com

2025-01-09 - 6:40:31 PM GMT

 Signer terry.kelly@malvernpanalytical.com entered name at signing as Terry Kelly

2025-01-09 - 6:44:19 PM GMT

 Document e-signed by Terry Kelly (terry.kelly@malvernpanalytical.com)

Signature Date: 2025-01-09 - 6:44:21 PM GMT - Time Source: server

 Agreement completed.

2025-01-09 - 6:44:21 PM GMT