

Carbon Reduction Plan

Supplier name	Promega UK Limited
Publication date	15th July 2025
Commitment	Promega UK Ltd is committed to achieving Net Zero by 2045

Baseline emissions footprint

Baseline year	2022
Additional details relating to the baseline emissions calculations	<p>Aligned with policy requirements, baseline emissions are estimated, representing the greenhouse gases produced before implementing any emission reduction strategies. For Promega UK, the baseline year is set as 2022, serving as a reference point against which future reduction efforts will be measured. However, it is essential to acknowledge that Promega UK had already implemented several measures to reduce emissions, ranging from the utilisation of Renewable & Nuclear energy to recycling waste generation. This core calculation adheres to the Greenhouse Gas Protocol corporate standard and serves as the baseline and current emissions reporting year for our progress assessment towards the stated emissions reduction targets. Initially, we include only the required subset of emissions, but we have plans to expand the scope of our calculations for a full carbon footprint assessment in accordance with the Greenhouse Gas Protocol. Throughout the calculations, direct activity data is primarily used, with spend data applied specifically for business travel purposes.</p>

Emissions Category	CO ₂ e
Scope 1	126.94 t
Scope 2	0 t
Scope 3	<p>Purchased goods and services 30.76t</p> <p>Downstream transportation & distribution 19t</p> <p>Business travel 12.49t</p> <p>Employee Commuting 3.79t</p> <p>Waste generated in operations 220.69kg</p> <p>Total Scope 3 = 66.26t</p>
Total Emissions	193.2 t

Current Emissions Reporting

Reporting Year	2024
Emissions Category	CO2e
Scope 1	79.9 t
Scope 2	Market Based 3.2 t Location based 75.0 t
Scope 3	Purchased goods and services 60.8 t Upstream Distribution 377.8 t Business travel 57.4 t Waste generated in operations 0.1 t Employee Commuting 17.80 t Downstream transportation & distribution 43.9 t Total Scope 3 = 557.7 t
Total Emissions	Market Based 640.8 t Location Based: 712.6 t

Emissions reduction targets

In order to continue our progress to achieving Net Zero, we have adopted the following carbon reduction targets.

- Scope 1 - 100% reduction by 2030
- Scope 2 - maintain the 100% reduction achieved in 2022
- Scope 3 - 100% reduction by 2045

Carbon Reduction Projects

At our premises, measures have been implemented to reduce our carbon footprint:

- Installed solar panels in the staff car park to generate renewable energy.
- Installed EV chargers in the staff car park to facilitate EV uptake.
- Replaced bathroom paper towels for hand towels to reduce waste and electricity consumption.
- Certificate our environmental management system to ISO 14001
- Utilised dry ice suppliers with carbon neutral feedstock
- Increased the proportion of fully electric and hybrid electric vehicles in the company vehicle fleet

In the future we will review, and if feasible implement, a number of the following initiatives:

1. Scope 1
 - a. Reduce the intensity of dry ice use by 15% per £m revenue by 2030
 - b. Identifier carbon neutral suppliers of dry ice
 - c. Transition to 95% of company cars to EVs by 2029
 - d. Implement EV adoption policy for employees
 - e. Transition to the use of biofuel for the onsite backup generator
2. Scope 2
 - a. purchase carbon offsets to neutralise emissions generated to produce grid mix electricity when using EV charging facilities not under our control
3. Scope 3
 - a. Identify highest impact suppliers by 2025
 - b. Prepare supplier communication strategy by 2025
 - c. Engage Tier 1 impact suppliers by end 2025

Declaration Sign Off

This Carbon Reduction Plan has been completed in accordance with PPN 06/21 and associated guidance and reporting standards.

Emissions have been reported and recorded in accordance with the published reporting standard for Carbon Reduction Plans and the GHG Reporting Protocol corporate standard and uses the appropriate Government emission conversion factors for greenhouse gas company reporting.

Scope 1 and Scope 2 emissions have been reported in accordance with SECR requirements, and the required subset of Scope 3 emissions have been reported in accordance with the published reporting standard for Carbon Reduction Plans and the Corporate Value Chain (Scope 3) Standard.

This Carbon Reduction Plan has been reviewed and signed off by the board of directors (or equivalent management body).


For and on behalf of Promega UK Limited

Signed:

Name:

Title:

Date:


DEREK McCALL (MD)
28/07/2025

