

Carbon Reduction Plan

Supplier name	Promega UK Limited
Publication date	30th April 2026
Commitment	Promega UK Ltd is committed to achieving Net Zero by 2045

Baseline emissions footprint

Baseline year	2023
Additional details relating to the baseline emissions calculations	<p>In 2026 the decision was taken to recalculate the baseline using the 2023 data set (as published in the 2024 CRP). This recalculation reflects material improvements in the accuracy of both emission factors and activity data applied across the inventory, which would otherwise compromise the consistency and comparability of year-on-year reporting. The recalculation is consistent with the GHG Protocol Corporate Standard, which requires base year emissions to be retroactively recalculated where changes in calculation methodology, or improvements in the accuracy of emission factors or activity data, result in a significant impact on the base year emissions data.</p> <p>It should be acknowledged that Promega UK had already implemented a number of emissions reduction measures prior to the baseline year, including the procurement of electricity from renewable and nuclear sources and established waste segregation and recycling practices. The baseline therefore reflects an operational position in which several efficiencies were already in place.</p> <p>The underlying calculation adheres to the GHG Protocol Corporate Standard and serves as both the baseline and current emissions reporting year for assessing progress against our stated emissions reduction targets. At this stage, reporting is limited to the required subset of emissions defined under PPN 06/21, with plans to extend the scope of our calculations toward a full organisational carbon footprint in line with the GHG Protocol. Direct activity data is used throughout, with spend-based data applied specifically to business travel.</p>

Emissions Category	CO2e
Scope 1	100.6t
Scope 2	Market Based 0.8t Location based 74.6t Total Scope 2 = 75.4 t
Scope 3	Purchased goods and services 29.0t Upstream Distribution 417.4t Business travel 61.0t Waste generated in operations 0.21t Employee Commuting 2.8t Downstream transportation & distribution 98.2t Total Scope 3 = 608.6t
Total Emissions	784.6 t

Current Emissions Reporting

Reporting Year	2025
Emissions Category	CO2e
Scope 1	79.9 t
Scope 2	Market Based 3.1 t Location based 61.3 t Total Scope 2 = 64.4 t
Scope 3	Purchased goods and services 56.4 t Upstream Distribution 319.6 t Business travel 54.2 t Waste generated in operations 0.2 t Employee Commuting 18.61 t Downstream transportation & distribution 36 t Total Scope 3 = 485.1 t
Total Emissions	629.5 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 - 100% reduction by 2030
- 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:

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 2026
 - b. Prepare supplier communication strategy by 2026
 - c. Engage Tier 1 impact suppliers by end 2026

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 McALL
MJ
14/05/2026

