



This certificate acknowledges that

Promega UK Ltd

has contributed to global Net Zero goals by funding the mitigation of

10 tonnes of carbon dioxide (CO₂)

through the following verified climate project:

GS12062 / CTL Landfill Gas Project / 10 tCO₂

Date: 15/07/2026



John Buckley
Managing Director, Carbon Footprint Ltd
www.carbonfootprint.com