

Which initiatives are we aligning with to support our ambition?



RE100 is a global initiative bringing together the world’s most influential businesses committed to 100% renewable electricity. Led by the Climate Group and in partnership with CDP, their mission is to accelerate change towards zero carbon grids at scale. RE100 member companies are already driving enough renewable electricity demand to power a medium sized country.



EV100 is a global initiative led by the Climate Group and in partnership with CDP to bring together forward-looking companies committed to accelerating the transition to electric vehicles, who commit to transition their fleets to EV and install EV charging for staff and customers by 2030.



The Science-Based Target initiative (SBTi) is a partnership between CDP, the United Nations Global Compact (UNGC), World Resources Institute (WRI) and the Worldwide Fund for Nature (WWF). SBTi facilitates a third-party validation process which assesses whether corporate climate targets are in line with the emissions reductions required by climate science.