

Map of One Kendall Square Complex

Quanterix Corporation
One Kendall Square
Building 1400W, 2nd Floor
Cambridge, MA 02139

