

memo

CSR活動の一つとして、
農業体験に参加しました！
皆様、ご協力いただき
ありがとうございました！
おおやま

農業体験 オーガニックファーム みなみあそ

みずの
うまれる
さと

トマトの脇芽摘み
トマトの収穫

6
Jun

7
Jul

トマトの収穫

10
Oct

ニンニクの植え付け
桃の剪定

5
May

トマトの脇芽摘み
桃の摘果
ニンニクの根切り

