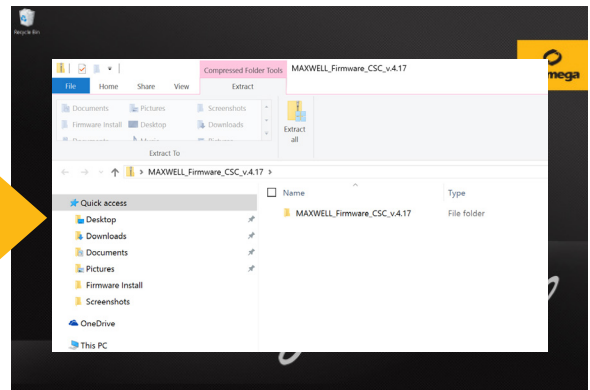
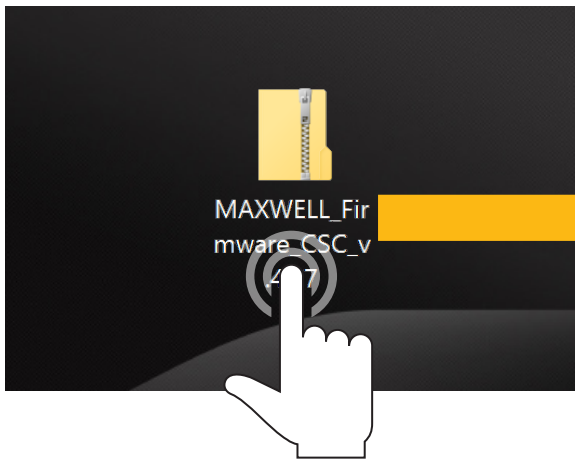


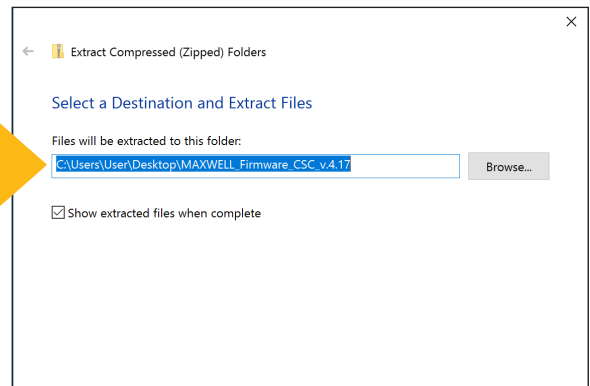
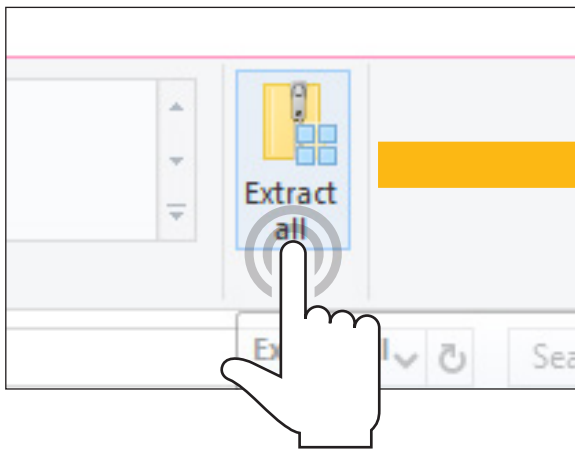
1

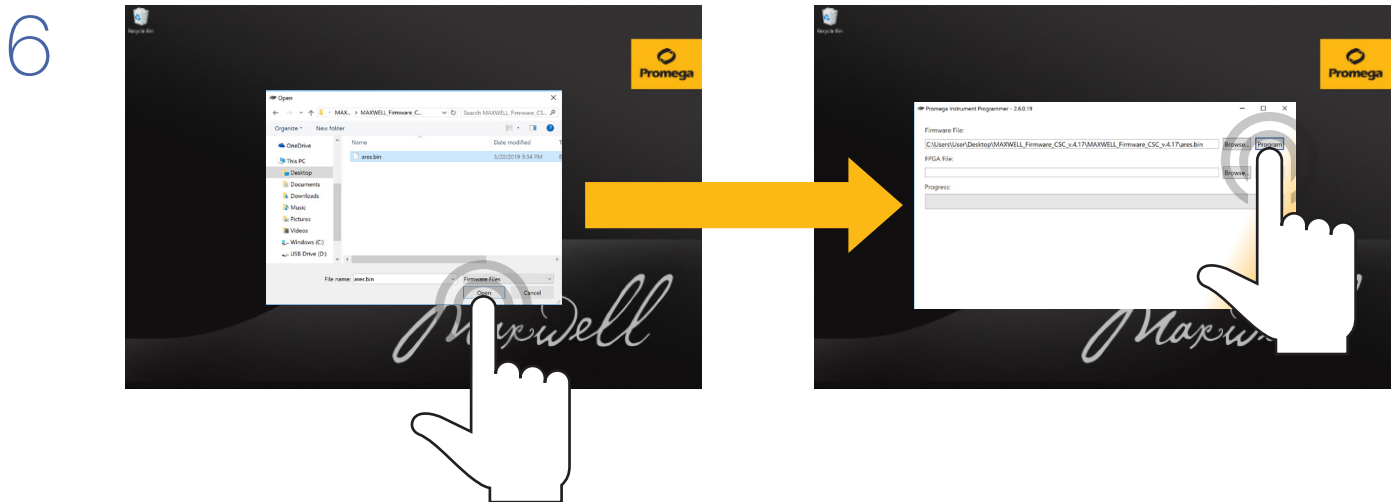
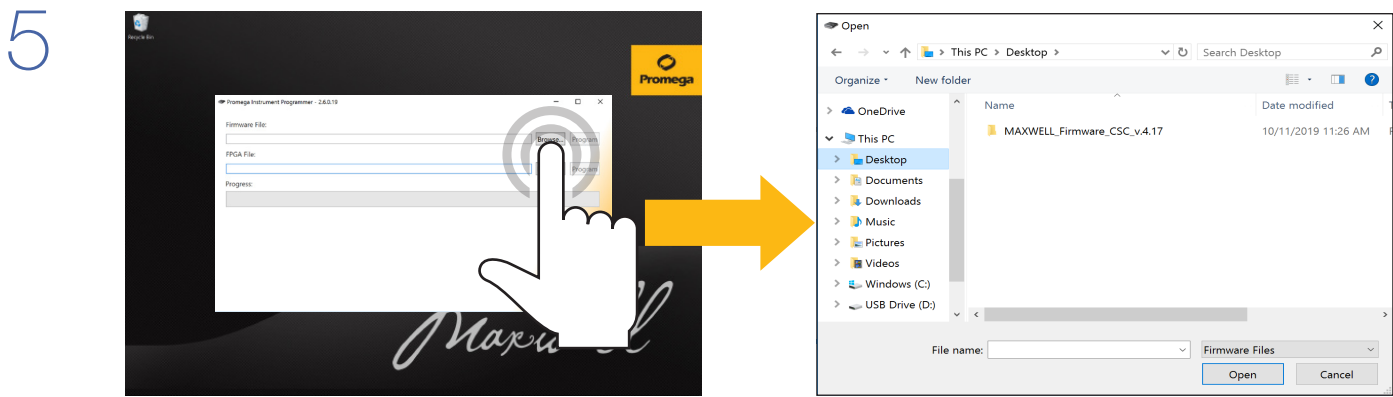


2



3





7

