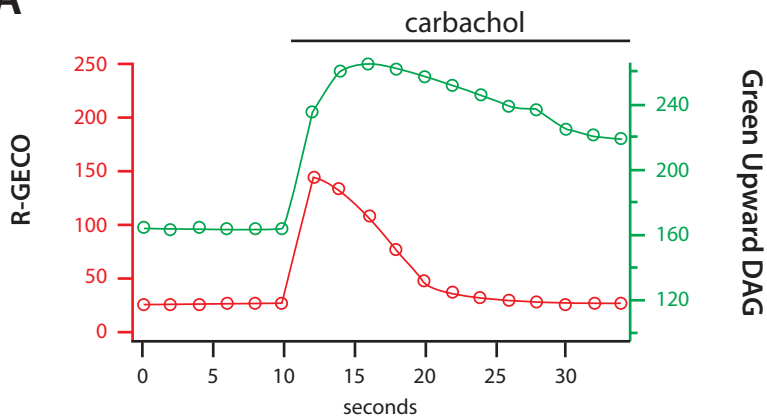


Multiplex Detection

A



B

