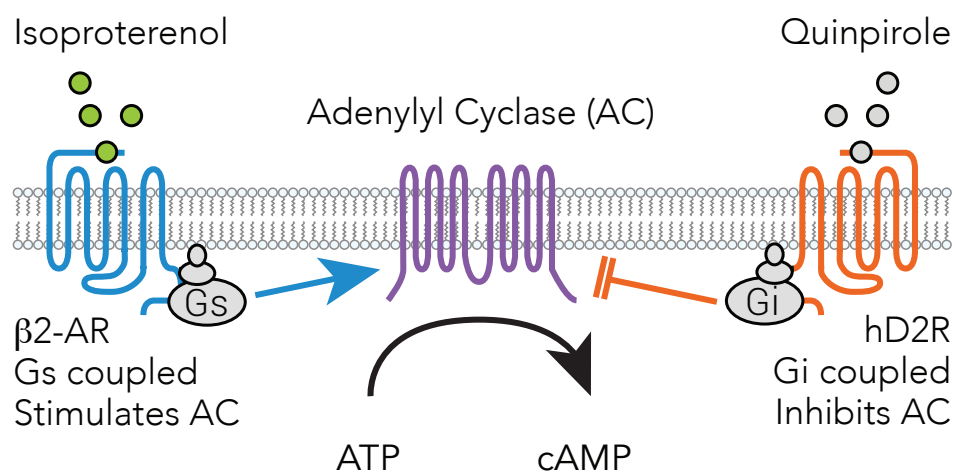
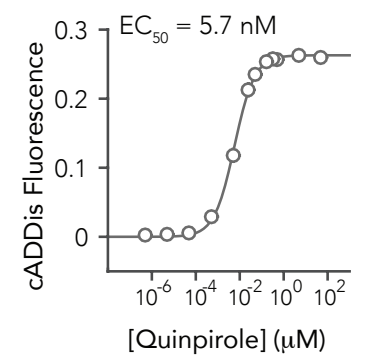
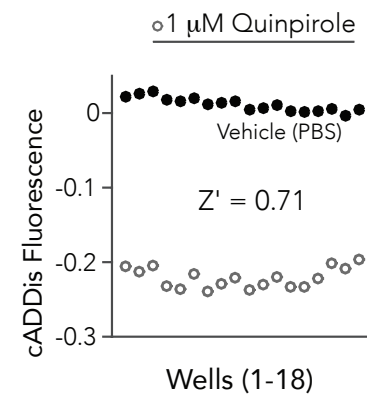
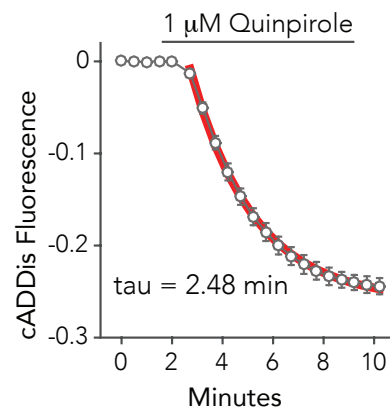
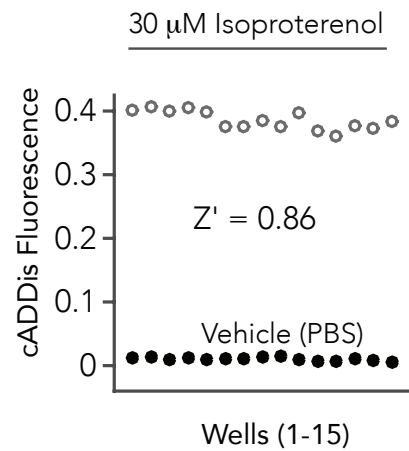
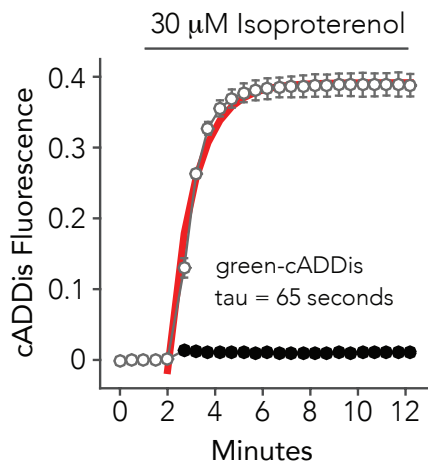


# G<sub>s</sub>



# G<sub>i</sub>



# G<sub>q</sub>

