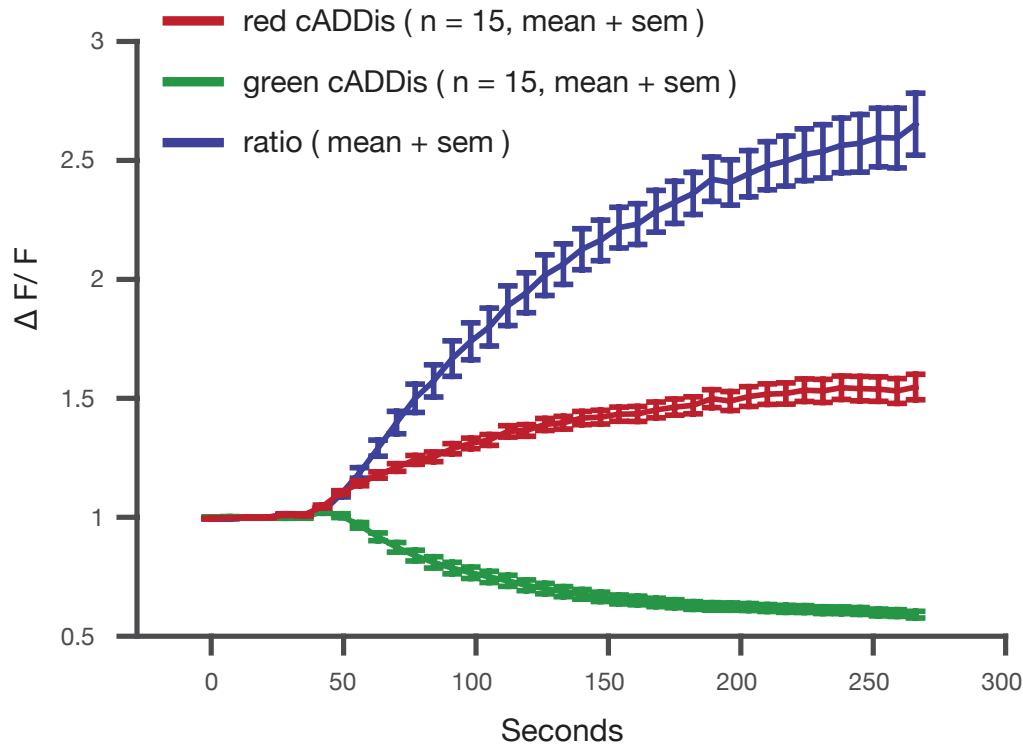
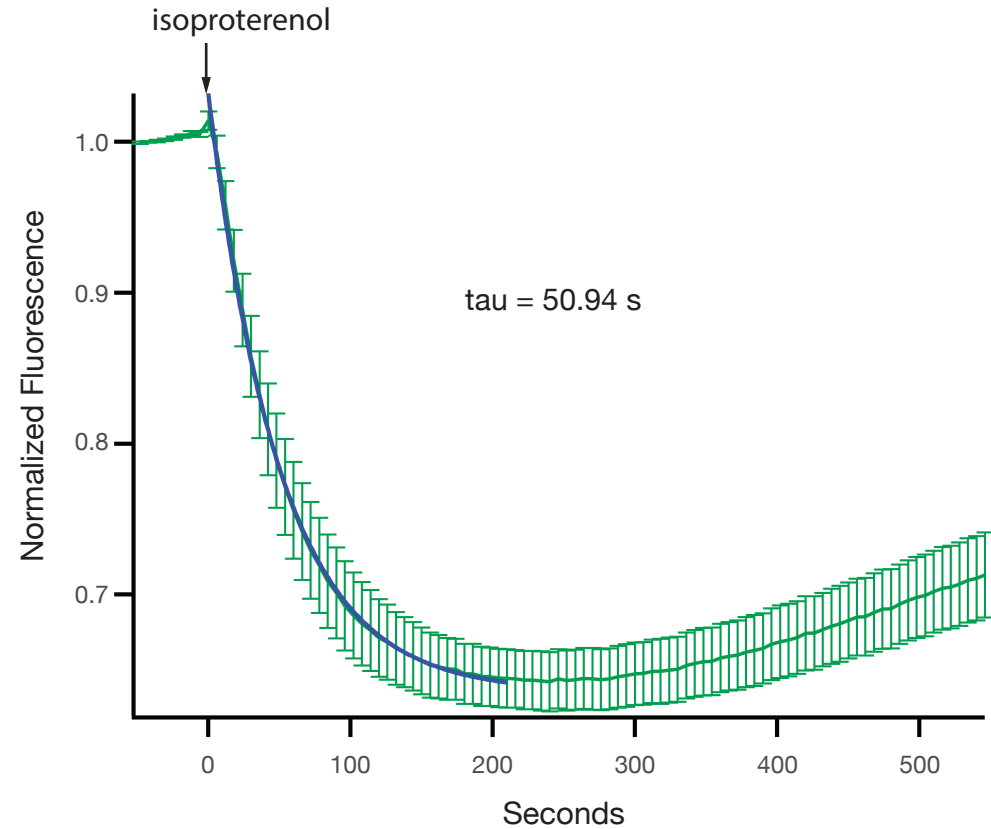


Ratiometric cAMP measurement



Red fluorescent cADDis sensor for cAMP combined with the green fluorescent cADDis sensor is a ratiometric measurement with a > 2.5 fold increase

G_s signaling detection in living cells



The downward green fluorescent cAMP sensor, cADDis, starts bright and decreases when cAMP increases.