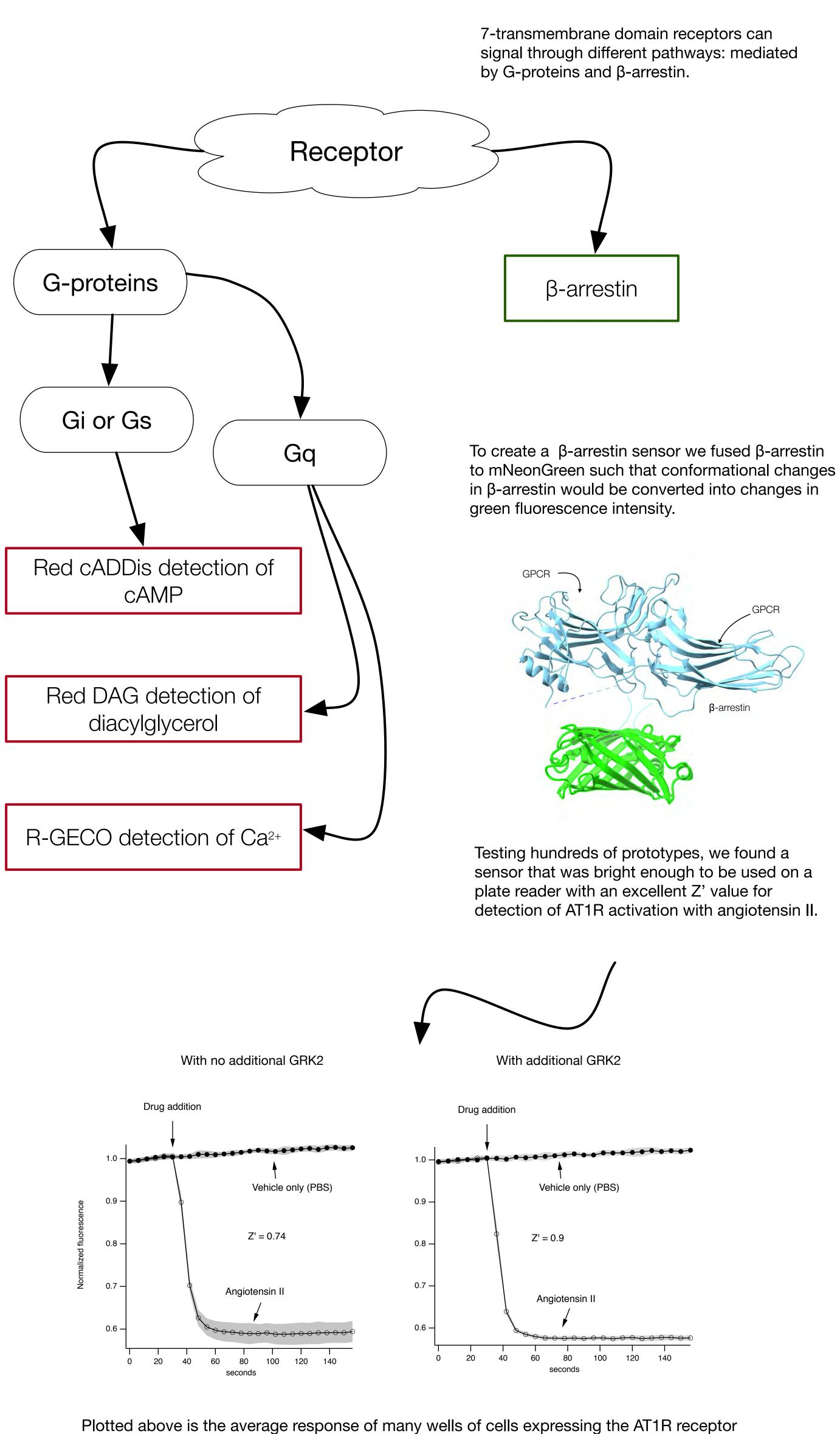


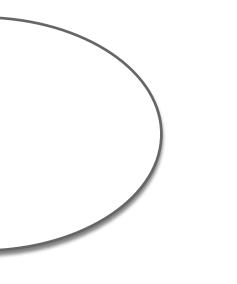
## We created a green fluorescent $\beta$ -arrestin sensor to pair with our red sensors for Gprotein signaling in living cells.

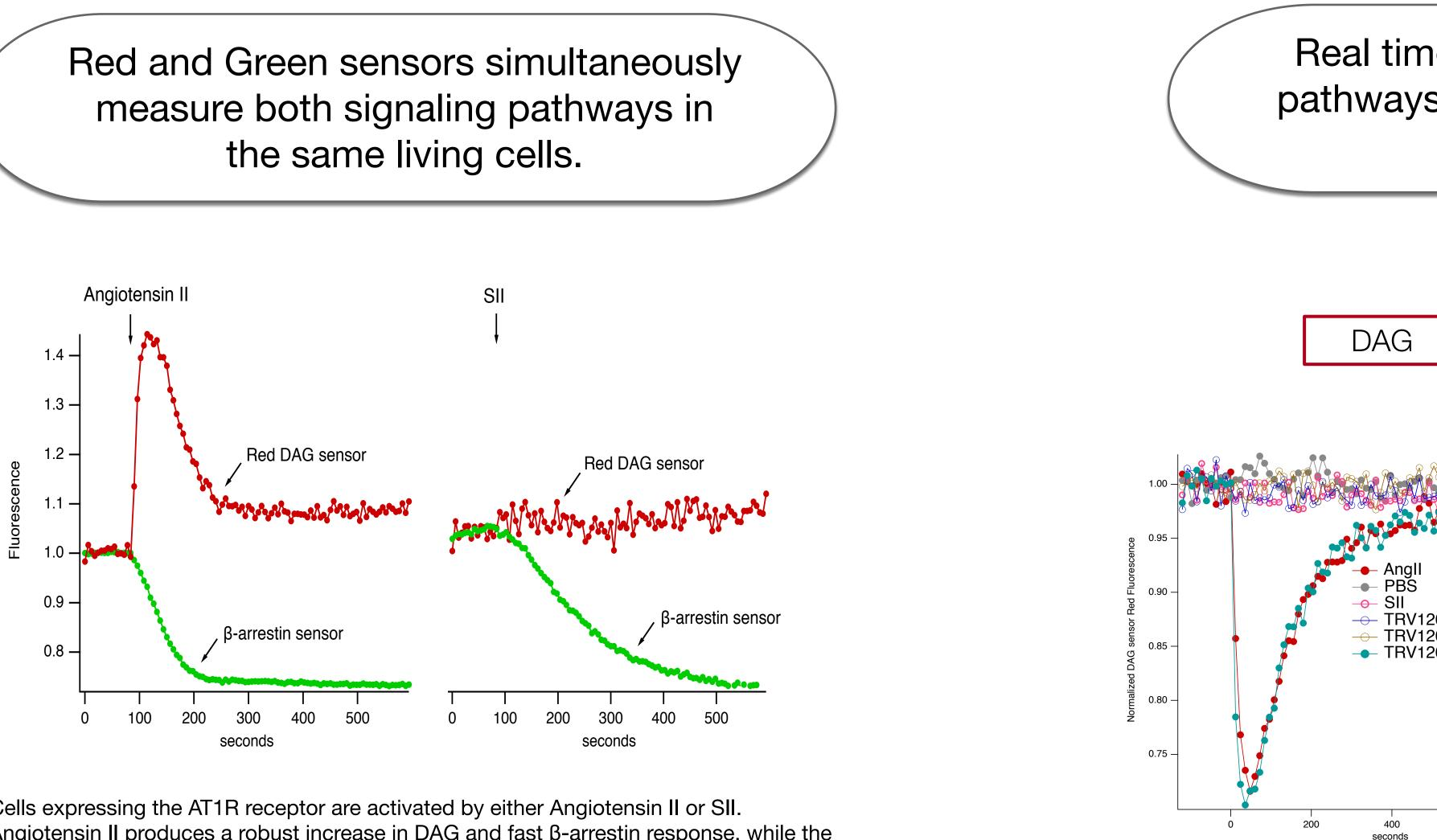


as well as the  $\beta$ -arrestin sensor. The standard deviation is illustrated in gray. The Z' value for detecting Angiotensin II activation (30 uM) was 0.74. Additional GRK2 reduced the variability and improved the Z'.

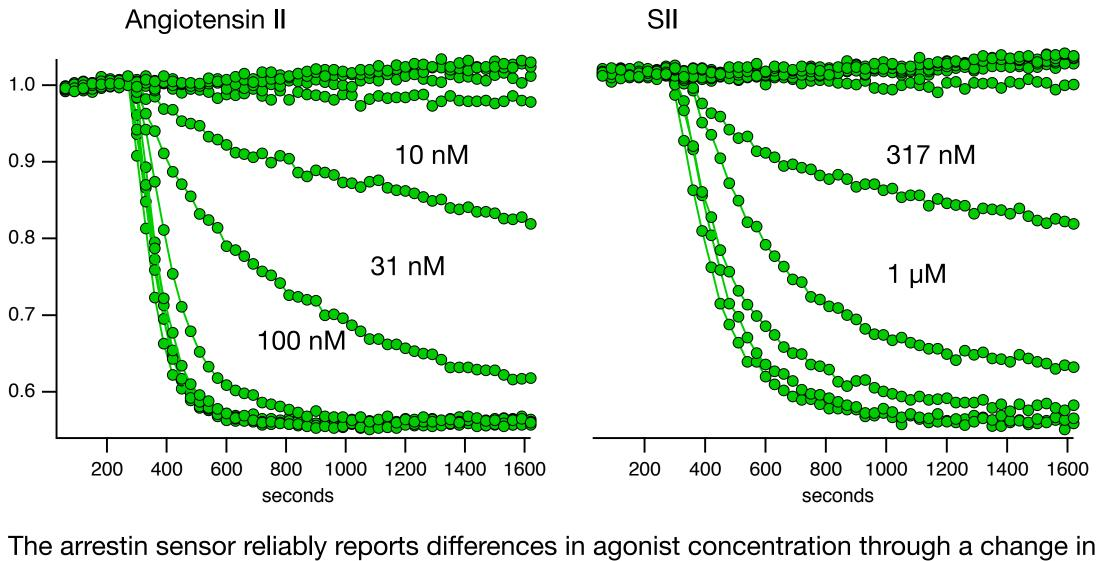
## Coloring $\beta$ -arrestin signaling green, and G-protein signaling red, a new live cell assay for quantitatively measuring agonist bias at seven transmembrane domain receptors.

Paul Tewson<sup>1</sup>, Sam Hoare<sup>2</sup>, Kevin Harlen<sup>1</sup>, Scott Martinka<sup>1</sup>, Anne Marie Quinn<sup>1</sup>, Thomas E. Hughes<sup>1</sup> (Montana Molecular <sup>1</sup>, Pharmechanics<sup>2</sup>)

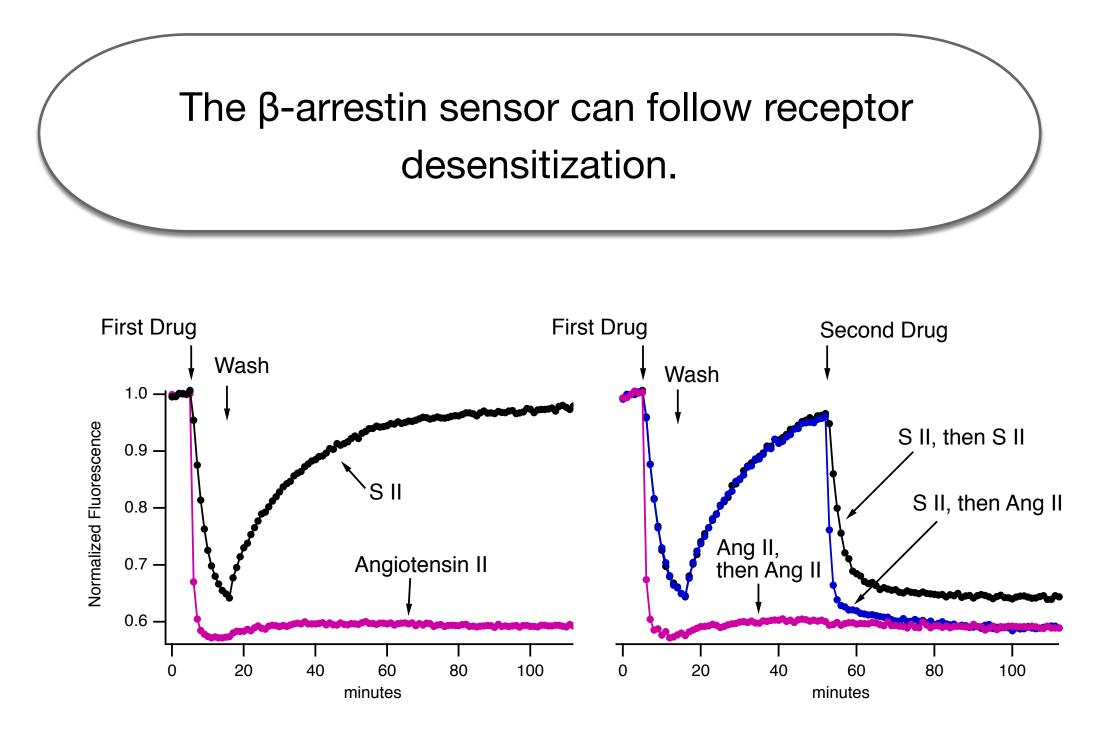




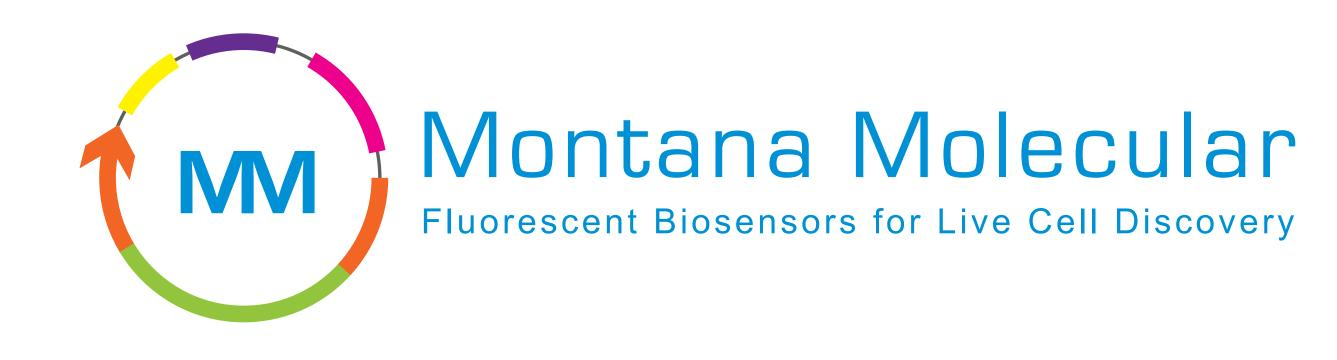
Cells expressing the AT1R receptor are activated by either Angiotensin II or SII. Angiotensin II produces a robust increase in DAG and fast  $\beta$ -arrestin response, while the  $\beta$ -arrestin biased ligand SII produces no detectable DAG response and a slow  $\beta$ -arrestin response.

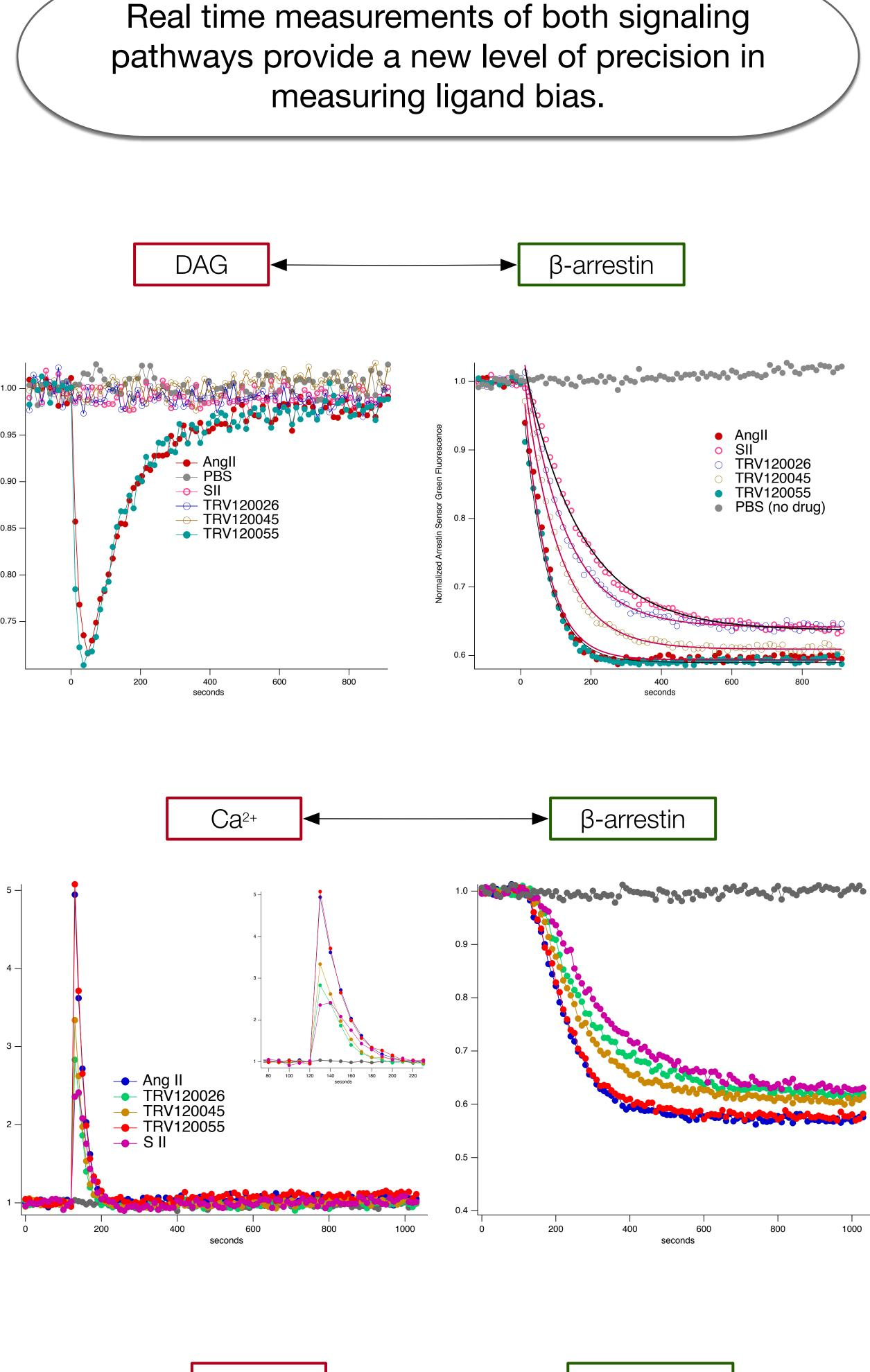


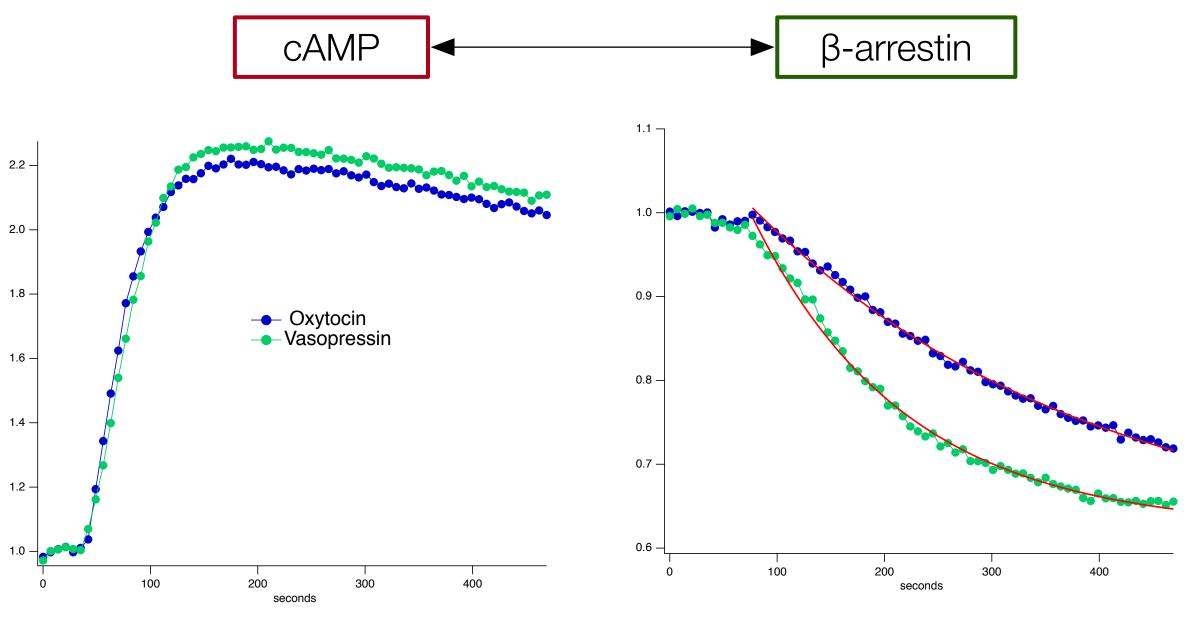
kinetics as well as the amplitude.



To explore receptor desensitization we first washed out the drug. Cells expressing the angiotensin receptor (AT1R) and activated with Angiotensin II did not return to baseline. However if they were treated with SII, they did recover. A second application then reactivated the  $\beta$ -arrestin sensor.







The human vasopressin receptor (human AVPR2) signals through cAMP and  $\beta$ -arrestin. Oxytocin and Vasopressin both activate the receptor and elevate cAMP. The  $\beta$ -arrestin responses are consistently different.