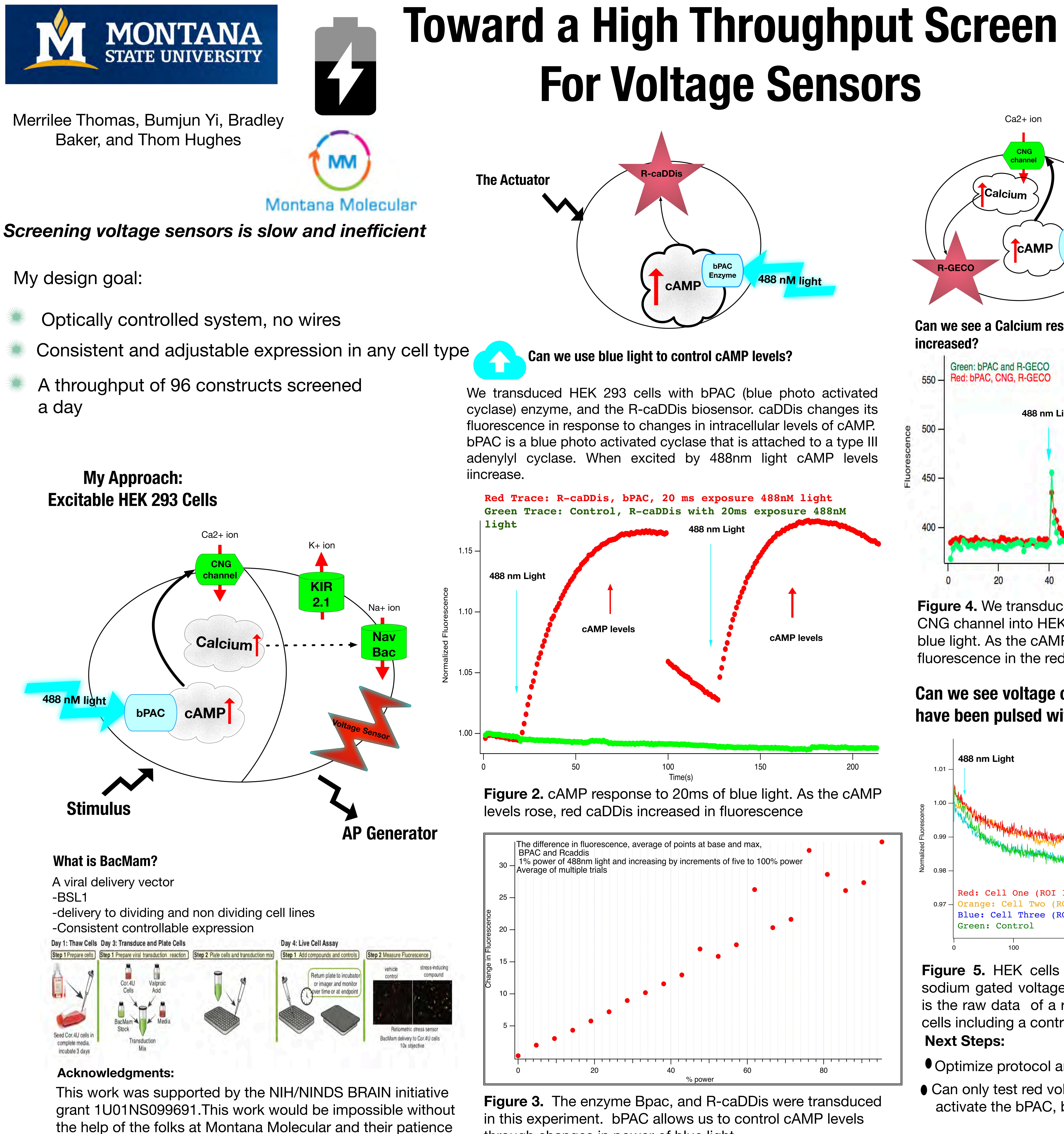


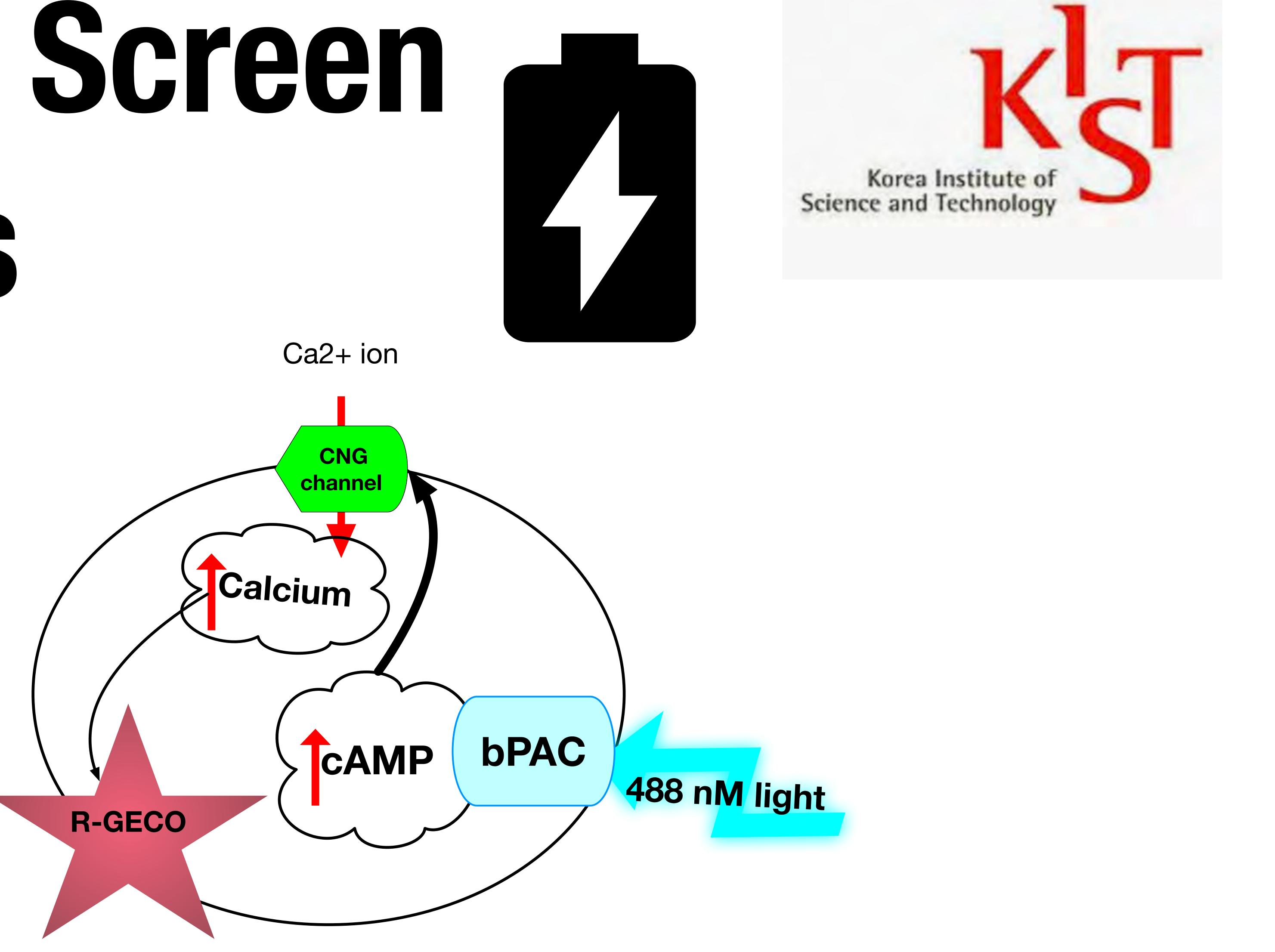
Baker, and Thom Hughes

- a day

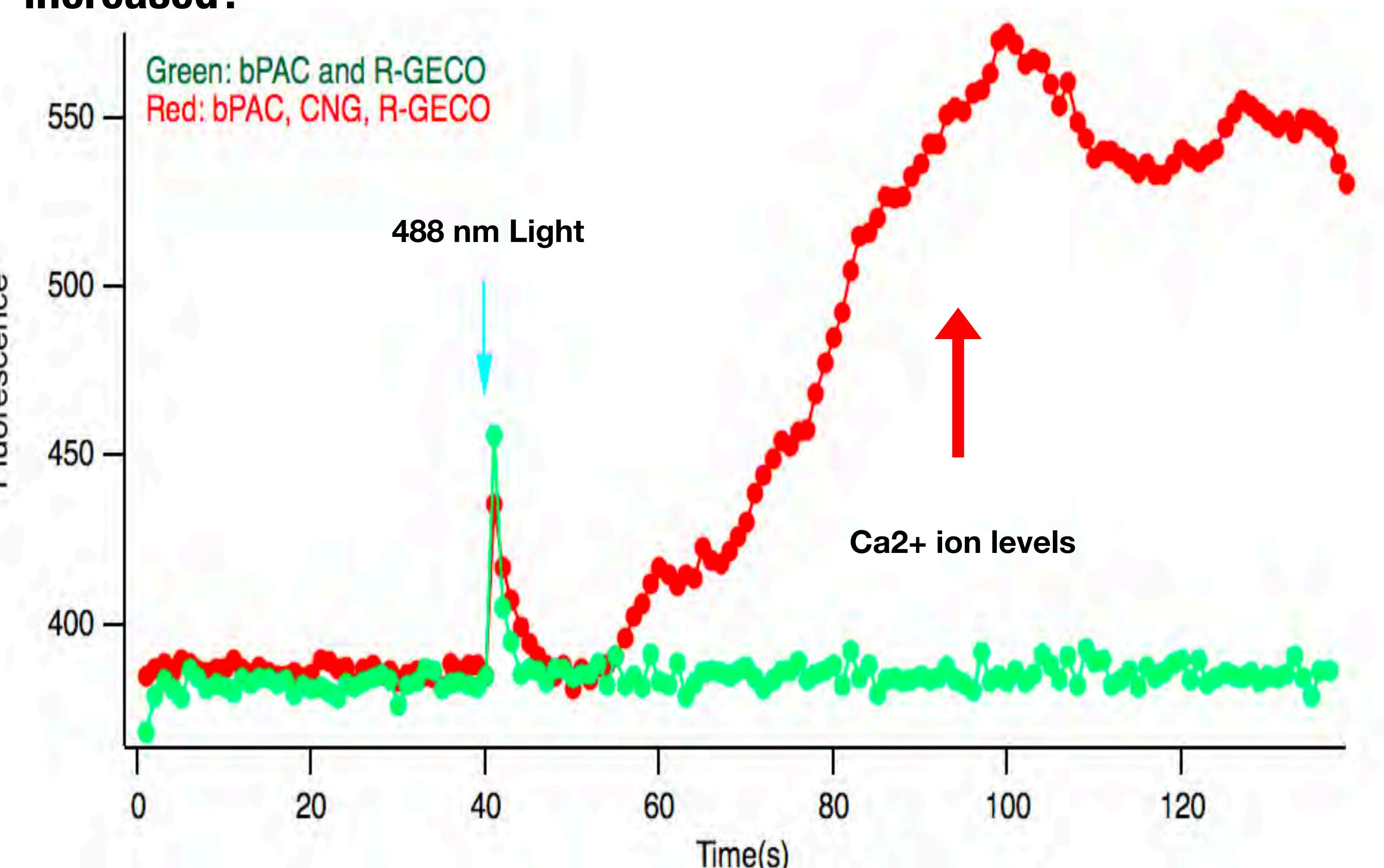
and never ending supply of HEK 293 cells.



through changes in power of blue light.



increased?



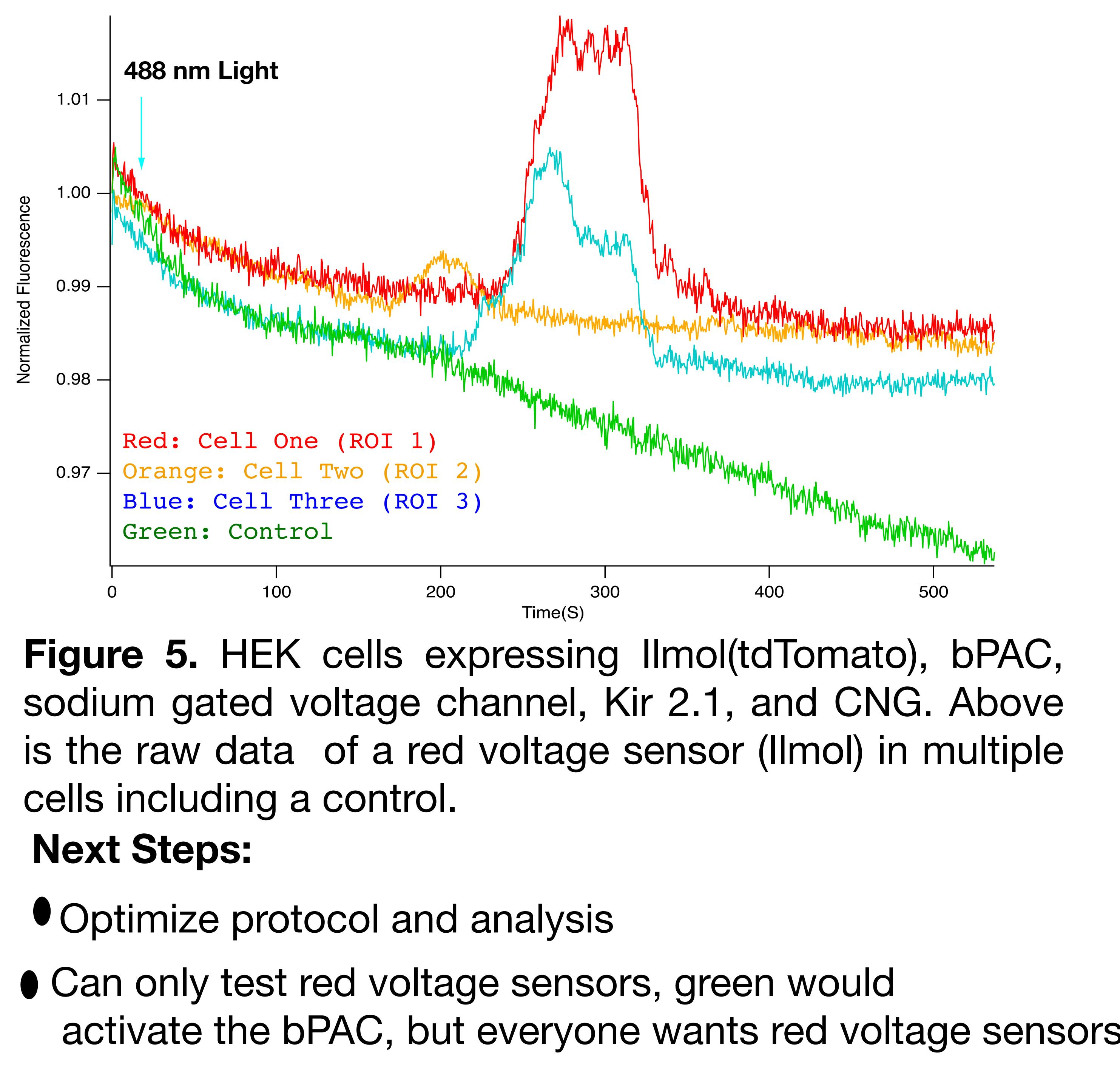




Figure 4. We transduced bPac Enzyme, R-Geco, and the CNG channel into HEK cells. A dose response to 20ms of blue light. As the cAMP levels rose, there was a increase in fluorescence in the red caDDis sensor.

## Can we see voltage changes in HEK cells that have been pulsed with blue light?