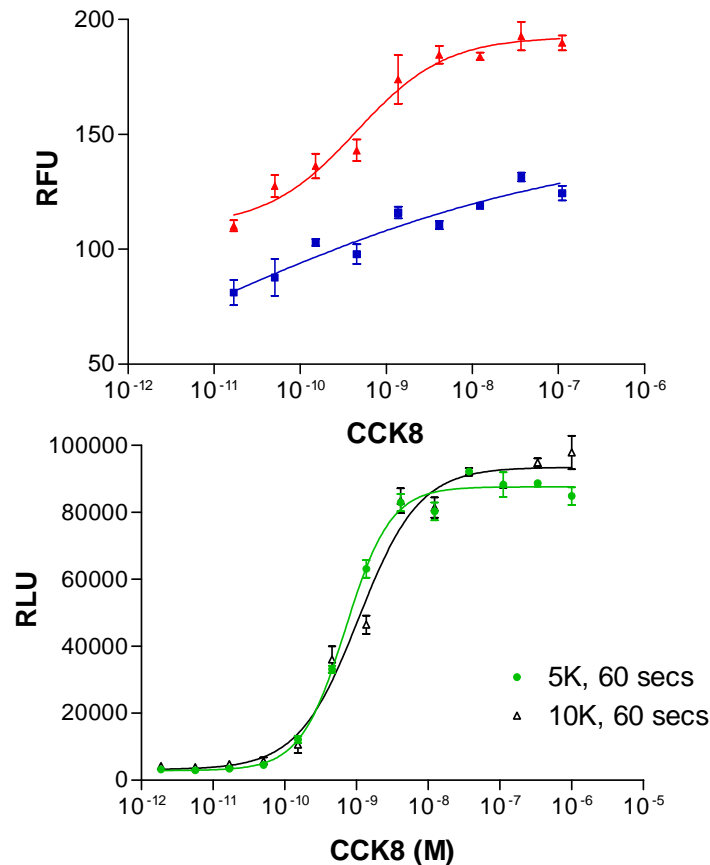
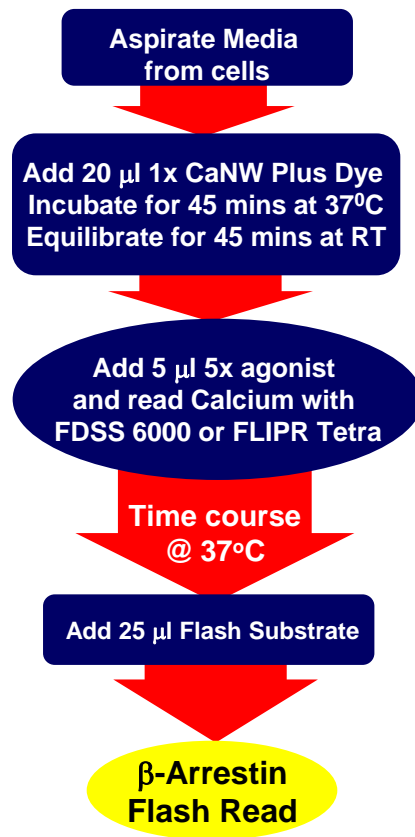


Sequential Measurement of Calcium and PathHunter Arrestin in the Same Well



Sequential Analysis of Calcium and Arrestin in the Same Well. Using an instrument with onboard liquid handling, we were able to demonstrate that the same plate treated with compounds can be sequentially analyzed first for calcium mobilization and then for arrestin signaling using our dual-purpose PathHunter cell lines. The upper panel shows the calcium measurement in CCK8 stimulated CCKAR cells (5k red, 10k cells blue), and the bottom panel shows the arrestin response in the same cells.