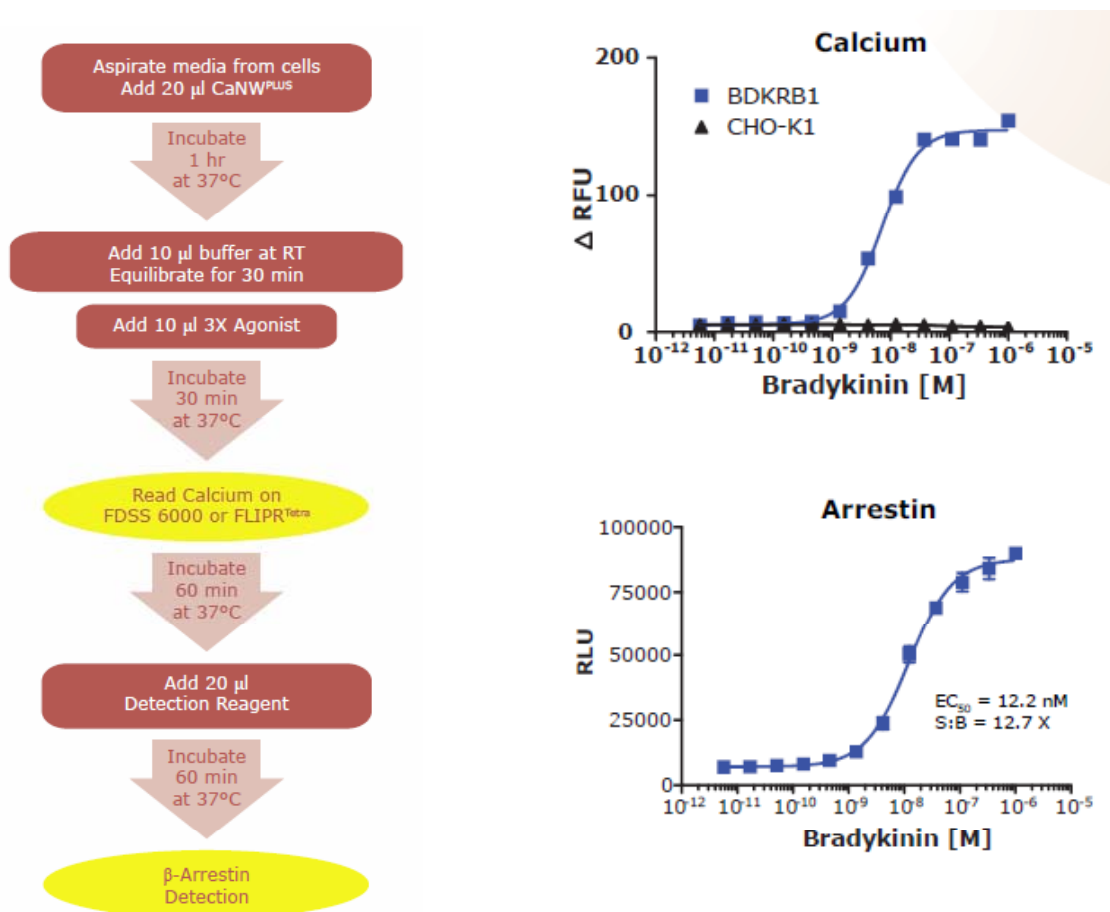


## Sequential Measurement of Calcium and Arrestin Recruitment in the Same Well



**Sequential Analysis of Calcium and Arrestin in the Same Well.** Using an instrument with onboard liquid handling, we were able to demonstrate the same plate treated with compounds can be sequentially analyzed first for calcium mobilization using the Calcium NoWash<sup>PLUS</sup> kit (90-0091L) and then for arrestin recruitment using the PathHunter Detection Reagents (93-0001) and our dual purpose PathHunter Bradykinin B1 Receptor (BDKRB1) cell line.