



Figure 2. EasyScreen™ β-Arrestin Assays enable selective confirmation of hits from a screen

EasyScreen CHO-K1 cells expressing human OPRD1 receptor were assayed for Arrestin signaling and second messenger activation in the presence of DADLE, a control agonist for OPRD1, and 3 known agonists for alternate GPCRs. **(A)** Only the target-specific DADLE compound elicited a measurable Arrestin response. **(B)** Treatment of cells with DADLE resulted in the expected cAMP response since OPRD1 couples to the cAMP pathway. A non-specific PAR agonist, TFLLR-NH2 also gave a robust response due to the activation of off-target, endogenous GPCRs. **(C)** DADLE treatment did not increase cytosolic calcium levels while the 3 other compounds resulted in robust increases in calcium via signaling through endogenous GPCRs on CHO-K1 cells.