



Discover_{Rx}

Figure 2. (A) In the PathHunter OPRD1 cell line exposure to DADLE, a known agonist to OPRD1 results in a robust arrestin response. As expected, the PathHunter OPRK1 cell line fails to respond. (B) In the heterodimer cell line, activation of the untagged OPRD1 results in transactivation of OPRK1 as shown by recruitment of arrestin and increased enzyme activity.