



### Figure 1. EasyScreen $\beta$ -Arrestin GPCR Assays result in excellent reproducibility

CHO-K1 cells expressing Murine CMKLR1 (mCMKLR1) were stimulated with 50 nM Chemerin, a known agonist, for 90 min prior to the addition of Detection Reagents. mCMKLR1 cells were also treated with vehicle alone (PBS/BSA). Eighty replicates were used for each treatment (%CV for vehicle = 7; %CV for agonist = 6). The Signal to Background was 13.0;  $Z' > 0.8$ .