



### Figure 3. Selectivity and Ligand Rank Ordering

(A) A series of known chemokine agonists were tested against select cAMP Hunter eXpress targets. CXCR2, CXCR3, and CXCR4 maintain agonist selectivity demonstrating that these assays are uniquely suited for profiling and/or target prioritization. (B) Cells expressing  $\beta$ 2-Adrenergic receptor (ADRB2) were treated with increasing concentrations of four known agonists for 30 minutes prior to the addition of Detection Reagents. Isoproterenol and Salmeterol are partial agonists.