



Discover_{Rx}

Figure 2. Performance comparison between live cells and frozen eXpress cells

(A) Correlation between EC_{50} values of control agonists obtained using live cells and PathHunter™ eXpress cells. Data is presented for 100 receptors. EC_{50} values above the line indicate higher sensitivity in the eXpress format compared to live cells. **(B)** CHO-K1 PathHunter™ eXpress cells and live cells expressing the Somatostatin 2 (SSTR2) receptor were stimulated with increasing concentrations of SST-28, a known agonist, for 90 minutes prior to the addition of detection reagents. PathHunter™ eXpress cells show equivalent pharmacology to live cells.