

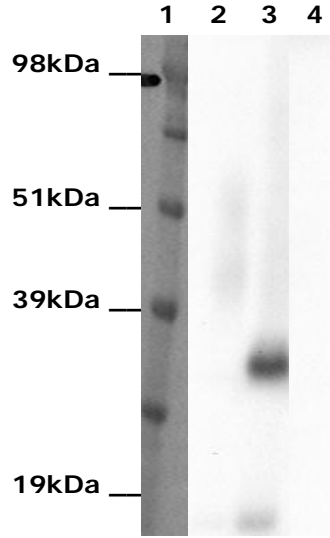
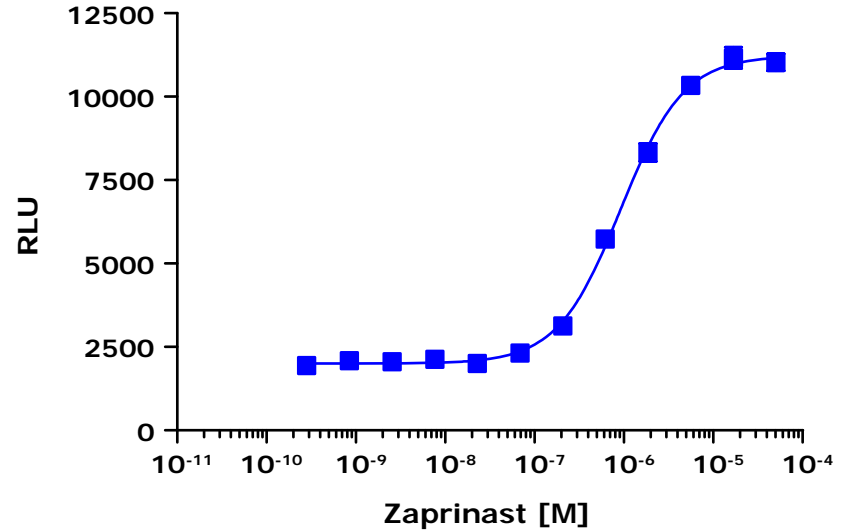
A.**B.****Discover_x**

Figure 1. Easily deorphanize any orphan GPCR using validated cell lines

(A) Expression of the GPR35 oGPCR receptor was confirmed by Eastern blot, as shown by the presence of a 34 KD band (lane 3). **(B)** Dose response curve for Zaprinst, a “surrogate” ligand for GPR35 showing a robust response to agonist.