

Detection of "Super" Agonists

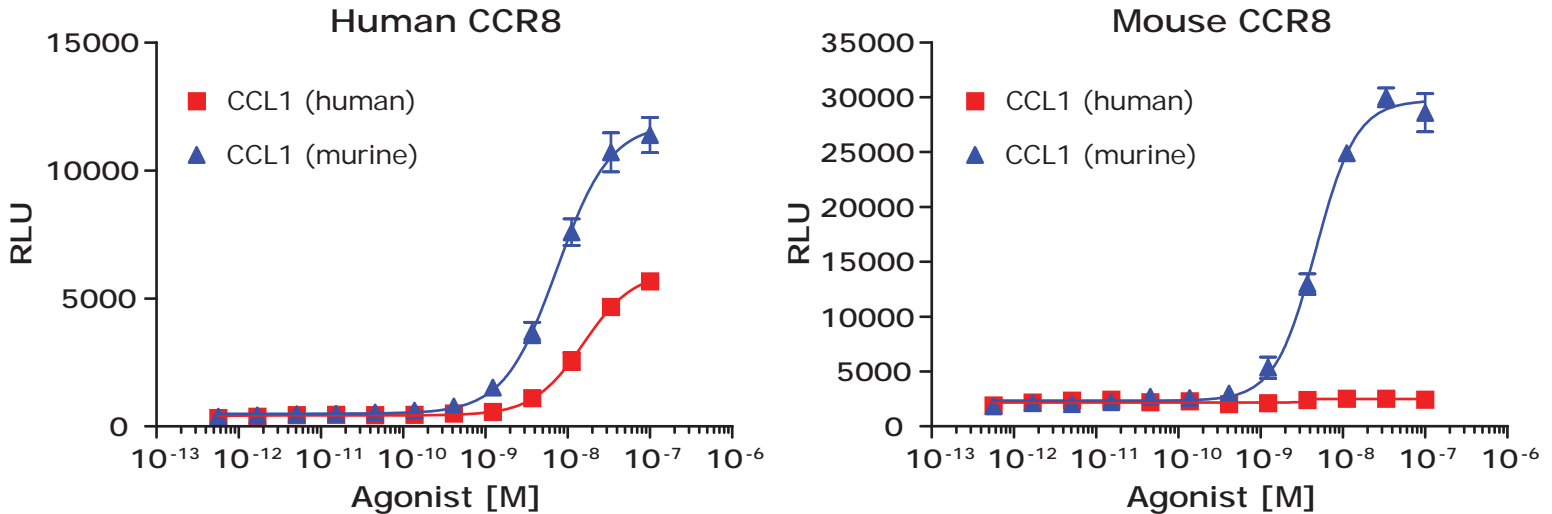


Figure 2. PathHunter™ β -Arrestin cells expressing human Chemokine (C-C motif) Receptor 8 (CCR8) were treated with CCL1 ligand. The efficacy of *murine* CCL1 for the *human* CCR8 receptor is much higher than the efficacy of the *human* CCL1. Thus, the *murine* CCL1 behaves as a "super" agonist at the *human* CCR8 receptor. At the same time, the *human* CCL1 shows little, if any, efficacy for the *murine* CCR8. A similar effect is seen with CCR9.