

Detection of Species Selective Agonists

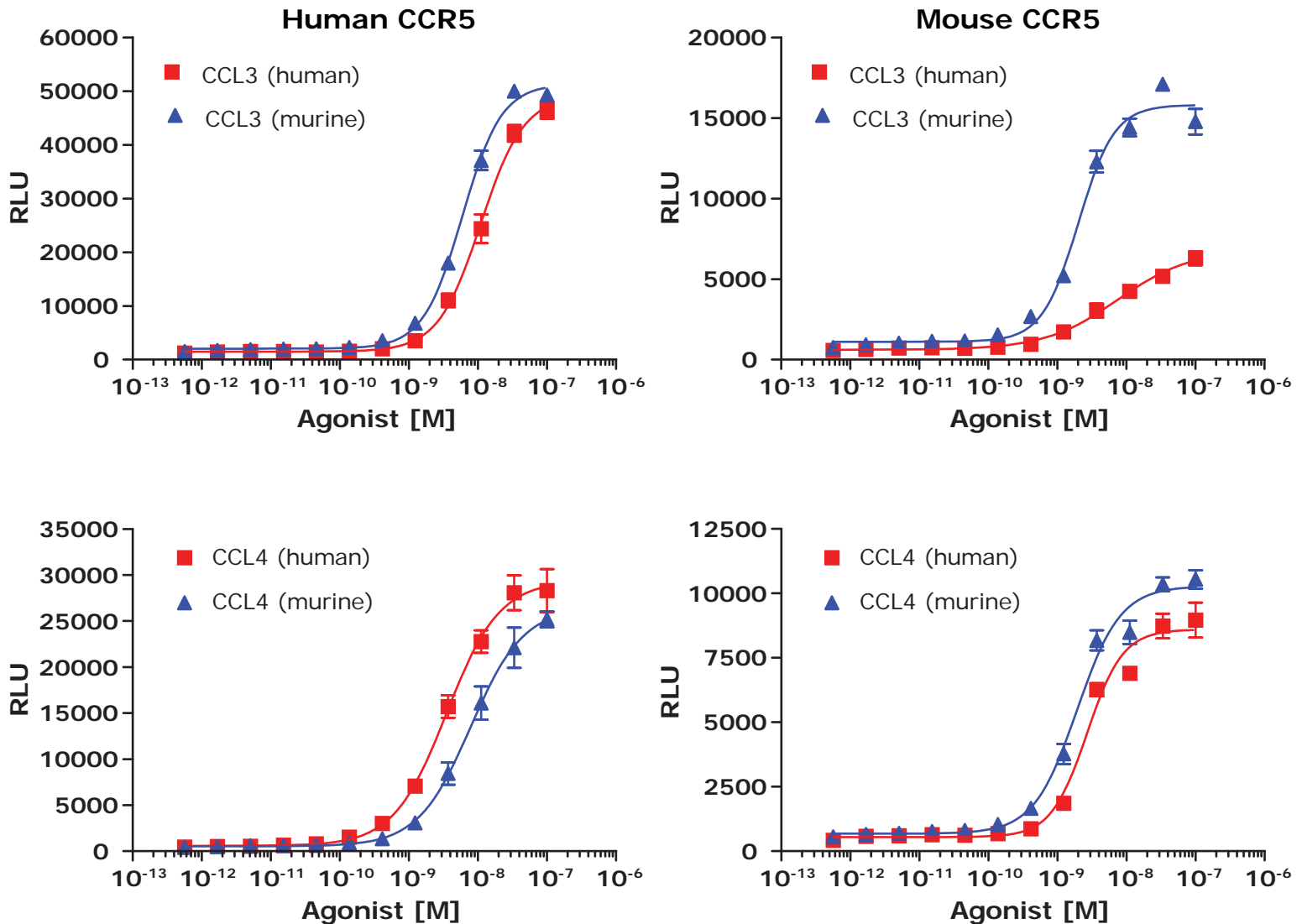


Figure 1. PathHunter™ β -Arrestin cells expressing human and murine Chemokine (C-C motif) Receptor 5 (CCR5) were treated with CCL3 and CCL4 agonists. *Human CCL3* is a partial agonist for *murine CCR5*, but both species of ligand are full agonists for the *human CCR5*. *CCL4* ligand showed full agonist activity on both the *human* and *murine CCR5* cell lines.