



## Figure 2. Agonist and antagonist dose response curves

cAMP Hunter eXpress CHO-K1 cells expressing the Melatonin 1B (MTNR1B) receptor were stimulated with increasing concentrations of Iodomelatonin, a known agonist, for 30 minutes prior to addition of Detection Reagents (Left Panel). MTNR1B cells were also treated with increasing concentrations of Luzindole, a known antagonist, in the presence of Iodomelatonin (Right Panel).