



Figure 2. PathHunter® Total GPCR Internalization Assays

The PathHunter® Total GPCR Internalization cell line overexpressing the b2 Adrenergic receptor (ADRB2) was treated with a panel of known agonists. Agonists and partial agonists were readily identified and expected pharmacology and rank ordering was observed.