



Figure 1. PathHunter® Activated GPCR Internalization Assays

The PathHunter® Activated GPCR Internalization cell line overexpressing the Opioid δ Receptor (OPRD1) was treated with a panel of compounds. SNC-80 (purple curve), a known strongly internalizing compound and functional antagonist of the OPRD1 receptor is clearly identified as a super agonist in the internalization assay.