



**Discover<sub>Rx</sub>**

**Figure 1. PathHunter™  $\beta$ -Arrestin assay for monitor functional GPCR heterodimers**

The PathHunter  $\beta$ -Arrestin assay only generates signal when the ProLink tagged GPCR binds to the Arrestin-EA fusion protein. By using PathHunter Arrestin clones as the starting material, a second untagged GPCR can be introduced into the cell line and the effects of the untagged GPCR on the ProLink tagged GPCR can be measured.