

How does it work?

PathHunter® Signal Detection

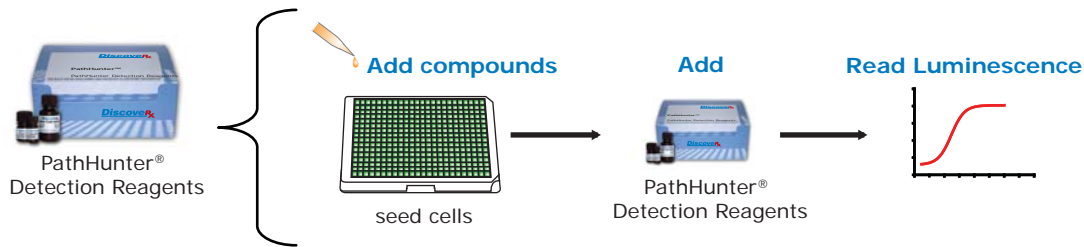


Figure 2. The PathHunter® System include engineered clonal cell lines validated for stable GPCR receptor tagged with ProLink™ and a β -Arrestin fused with enzyme acceptor (EA); as well as unique PathHunter® Certified Reagents especially optimized to yield high quality, reproducible results. These system utilize a simple, one-step addition protocol and standard chemiluminescent detection and are ideally suited for 96-, 384-, 1536-well compound screening.