



Figure 1. Confirm pharmacology of your compound regardless of whether it is an agonist, antagonist, inverse agonist or allosteric modulator using a simple, homogeneous protocol combined with chemiluminescent detection.