

# NanoBRET™ Target Engagement Assay for CSNK1G2

## Protocol

HEK293 cells were transfected with NanoLuc®-CSNK1G2 Fusion Vector and seeded into the wells of 96-well plates. The cells were incubated with test compound for 2 hours following the addition of the NanoBRET™ tracer reagent. The BRET signal was measured on a luminometer after dispensing the NanoBRET™ Nano-Glo® Substrate and Extracellular NanoLuc® Inhibitor into the wells. IC50 values were calculated by fitting the data to the following equation.

$$\text{BRET ratio} = \text{Min} + (\text{Max} - \text{Min}) \times \frac{1}{(1 + (\text{IC50}/[\text{Compound}])^{\text{Slope}})}$$

