



Catalog No. PXC-PPI-2

## Split Luc HEK293 I $\kappa$ B $\alpha$ / $\beta$ -TRCP Complex Cell Line

### Product Summary

The N-terminal and C-terminal fragments of unique split click beetle luciferase are fused to I $\kappa$ B $\alpha$  and  $\beta$ -TRCP respectively. The formation of I $\kappa$ B $\alpha$ / $\beta$ -TRCP complex brings the N-terminal luciferase into proximity with the C-terminal luciferase, and bioluminescence activity is recovered. To detect this reaction, Split Glo Cell Assay Reagent (Catalog No. PXR-SG001) is required.

### Description

Accession Number	I $\kappa$ B $\alpha$ : NM_020529 $\beta$ -TRCP: NM_001256856
Description	I $\kappa$ B $\alpha$ : NF $\kappa$ B inhibitor alpha (NFKBIA) $\beta$ -TRCP: beta-transducin repeat containing E3 ubiquitin protein ligase (BTRC), transcript variant 3
Cell Line	HEK293
Species	Human
Storage	Liquid Nitrogen

### Function

Control Agonist	TNF $\alpha$
Control Antagonist	IKK inhibitor
Assay Plate	96 well plate
Cell number / well	6 x 10 <sup>4</sup>
Assay wells	3
Incubation time (hour)	1
Incubation temperature (°C)	37
Antagonist concentration ( $\mu$ M)	Various
S/B Ratio (w/wo inhibitor)	18.8

