

## Split Luc HEK293 AGTRL1 $\beta$ -Arrestin-2 Cell Line

### Product Summary

The N-terminal and C-terminal fragments of unique split click beetle luciferase are fused to  $\beta$ -arrestin and GPCR, respectively. Binding of a ligand to GPCR triggers phosphorylation of GPCR, thereby inducing its interaction with  $\beta$ -arrestin. This interaction 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

Receptor Family	Apelin
Target GPCR	AGTRL1
Coupling	Gi
Accession Number	NM_005161
Description	Apelin receptor
$\beta$ -Arrestin Isoform	$\beta$ -Arrestin-2
Cell Line	HEK293
Species	Human
Storage	Liquid Nitrogen

### Function

Control Agonist	[Pyl1]-apelin-13
Assay Plate	96 well plate
Cell number / well	$2 \times 10^4$
Assay wells	4
Incubation time (min)	20
Incubation temperature ( $^{\circ}$ C)	37
Agonist concentration ( $\mu$ M)	1
S/B Ratio	90.5

