

Split Luc HEK293 NPY1R β -Arrestin-2 Cell Line

Product Summary

The N-terminal and C-terminal fragments of unique split click beetle luciferase are fused to β -arrestin and GPCR, respectively. Binding of a ligand to GPCR triggers phosphorylation of GPCR, thereby inducing its interaction with β -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	Neuropeptide Y
Target GPCR	NPY1R
Coupling	Gi
Accession Number	NM_000909
Description	Neuropeptide Y receptor Y1
β -Arrestin Isoform	β -Arrestin-2
Cell Line	HEK293
Species	Human
Storage	Liquid Nitrogen

Function

Control Agonist	Neuropeptide Y
Assay Plate	96 well plate
Cell number / well	2×10^4
Assay wells	3
Incubation time (min)	60
Incubation temperature ($^{\circ}$ C)	37
Agonist concentration (μ M)	1
S/B Ratio	53.1

