

Split Luc HEK293 CCKBR 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, 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	Cholecystokinin
Target GPCR	CCKBR
Coupling	Gq
Accession Number	NM_17685
Description	Cholecystokinin B receptor
β -Arrestin Isoform	β -Arrestin-2
Cell Line	HEK293
Species	Human
Storage	Liquid Nitrogen

Function

Control Agonist	Gastrin-1
Assay Plate	96 well plate
Cell number / well	2×10^4
Assay wells	4
Incubation time (min)	90
Incubation temperature ($^{\circ}$ C)	37
Agonist concentration (μ M)	1
S/B Ratio	46.0

