

Catalog No. PXC-014-1

Split Luc HEK293 SSTR2 β -Arrestin-1 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 | Somatostatin |
| Target GPCR | SSTR2 |
| Coupling | Gi |
| Accession Number | NM_001050 |
| Description | Somatostatin receptor 2 |
| β -Arrestin Isoform | β -Arrestin-1 |
| Cell Line | HEK293 |
| Species | Human |
| Storage | Liquid Nitrogen |

Function

| | |
|--|-----------------|
| Control Agonist | Somatostatin |
| Assay Plate | 96 well plate |
| Cell number / well | 2×10^4 |
| Assay wells | 4 |
| Incubation time (min) | 20 |
| Incubation temperature ($^{\circ}$ C) | 37 |
| Agonist concentration (μ M) | 1 |
| S/B Ratio | 166.8 |

