

Catalog No. PXC-PPI-1

Split Luc HEK293 Hif-1 Complex Cell Line

Product Summary

The N-terminal and C-terminal fragments of unique split click beetle luciferase are fused to Hif-1 α and Hif-1 β , respectively. The formation of Hif-1 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	Hif-1 α : NM_001530 Hif-1 β : NM_001668
Description	Hif-1 α : hypoxia inducible factor 1, alpha Hif-1 β : hydrocarbon receptor nuclear translocator (ARNT)
Cell Line	HEK293
Species	Human
Storage	Liquid Nitrogen

Function

Control Antagonist	Acriflavine
Assay Plate	96 well plate
Cell number / well	2×10^4
Assay wells	3
Incubation time (hour)	24
Incubation temperature (°C)	37
Antagonist concentration (μ M)	10
S/B Ratio (w/wo inhibitor)	6.1

