

Catalog No. PXC-PPI-3

Split Luc HEK293 ARIA/PTEN Complex Cell Line

Product Summary

The N-terminal and C-terminal fragments of unique split click beetle luciferase are fused to ARIA and PTEN, respectively. The formation of ARIA/PTEN complex brings the N-terminal luciferase into proximity with the C-terminal luciferase, and bioluminescence activity is recovered. To detect this reaction, D-Luciferin (Catalog No. CAY-14681-1G) is required.

Description

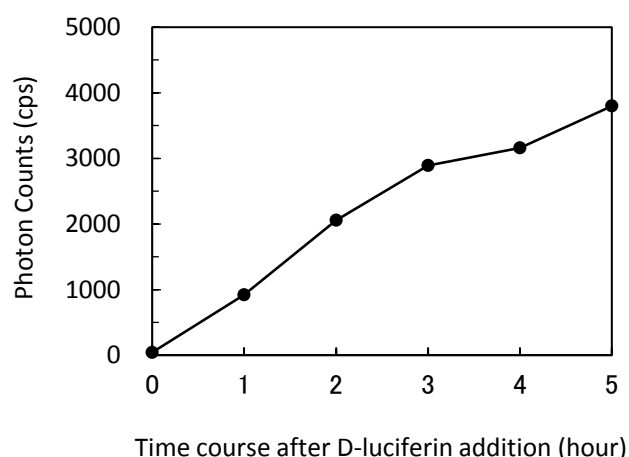
Accession Number	ARIA: DQ462572 PTEN: U96180
Description	ARIA: endothelial cell-specific molecule-2 (ECSM2) PTEN: protein tyrosine phosphatase (TEP1)
Cell Line	HEK293
Species	Human
Storage	Liquid Nitrogen
Remarks	ARIA and PTEN is constitutively associated. For inhibition assay, the effect of the antagonist can be evaluated both before and after luciferase addition. The signal is expected to be linear up to 10 hour after adding D-Luciferin*.

***Split Glow Cell Assay Reagent is not suitable to detect the signal.**

Function

ARIA/PTEN assay

Assay Plate	96 well plate
Cell number / well	6×10^4
Assay wells	1
Incubation temperature (°C)	Room Temp. (25)



1. Harvest the cells by treating with Trypsin-EDTA. Pellet the cells (180 x g for 2 minutes at room temperature) and wash with fresh complete medium.
2. Resuspend cells in complete medium and adjust the density to 6×10^5 cells/mL.
3. Add 100 μ L/well cell suspension to a 96 well plate (6×10^4 cells per well) and incubate for 48 hours in a humidified CO₂ atmosphere maintained at 37°C.
4. Remove the medium and add gently 2mM D-Luciferin (in HBSS) (100 μ L/well).
5. Incubate up to 10 hours at room temperature to get appropriate signal and measure luminescence*.

Please cover the bottom of the plate before adding Reagent. Transparent bottom is not suitable for luminescence measurements.