



Secrete-Pair™ Dual Luminescence Assay Kit

Overview

Secrete-Pair™ Dual Luminescence Assay Kit is designed to analyze the activities of Gaussia Luciferase (GLuc) and Secreted Alkaline Phosphatase (SEAP) of a dual-reporter system side-by-side using the same sample from the cell culture medium. Both GLuc and SEAP are secreted reporter proteins. Samples can be easily obtained from cell culture medium without lysis of the cells.

Two buffer conditions are provided in the kit for GLuc assays depending on the applications. Buffer GL-S contains a stabilizer and can be used for stabilized activity by overcoming the quick decay of the GLuc signal. When higher sensitivity is required for detecting low expression of GLuc, Buffer GL-H can be used for higher enzyme activity.

Secrete-Pair measures dual reporter signals and allows transfection normalization. The normalized GLuc activities can be compared across samples free of the impact of transfection variation.

Advantages

Live cell assays

- Secreted GLuc and SEAP
- Lysis of the cells is not necessary

Robust and flexible conditions

- Buffer for stable activity extends the half-life of light emission to approximately 30 minutes
- Buffer for higher sensitivity can be used to detect low GLuc expression

Dual-reporter detection

- Detects GLuc and SEAP
- Enables transfection-normalization for true cross-sample comparison

High-throughput compatible

- Quick and easy assay format
- High sample number compatible

To Order

Secrete-Pair™ kits

Catalog#	Product	Description	Price
LF031	Secrete-Pair™ Dual Luminescence Assay Kit (100 rxns)	Detects Gaussia luciferase (GLuc) and secreted alkaline phosphatase (SEAP)	\$139
LF032	Secrete-Pair™ Dual Luminescence Assay Kit (300 rxns)		\$379
LF033	Secrete-Pair™ Dual Luminescence Assay Kit (1000 rxns)		\$959
LF061	Secrete-Pair™ Gaussia Luciferase Assay Kit (100 rxns)	Detects Gaussia luciferase (GLuc)	\$77
LF062	Secrete-Pair™ Gaussia Luciferase Assay Kit (1000 rxns)		\$399

Cloning vectors

Catalog#	Vector name	Description	Promoter	Selection marker
ZX101	pEZX-GN01	Gaussia luciferase (GLuc) reporter cloning vector	N/A	Puromycin
ZX102	pEZX-GN03		miniCMV	Puromycin
ZX103	pEZX-GA01	Gaussia luciferase (GLuc) and secreted alkaline phosphatase (SEAP) reporter cloning vector	N/A	Puromycin
ZX104	pEZX-GA02		SV40	Neomycin
ZX105	pEZX-GA03		miniCMV	Puromycin
ZX106	pEZX-LvGN01	Gaussia luciferase (GLuc) reporter cloning vector (lentiviral)	N/A	Puromycin
ZX107	pEZX-LvGA01	Gaussia luciferase (GLuc) and secreted alkaline phosphatase (SEAP) reporter cloning vector (lentiviral)	N/A	Puromycin

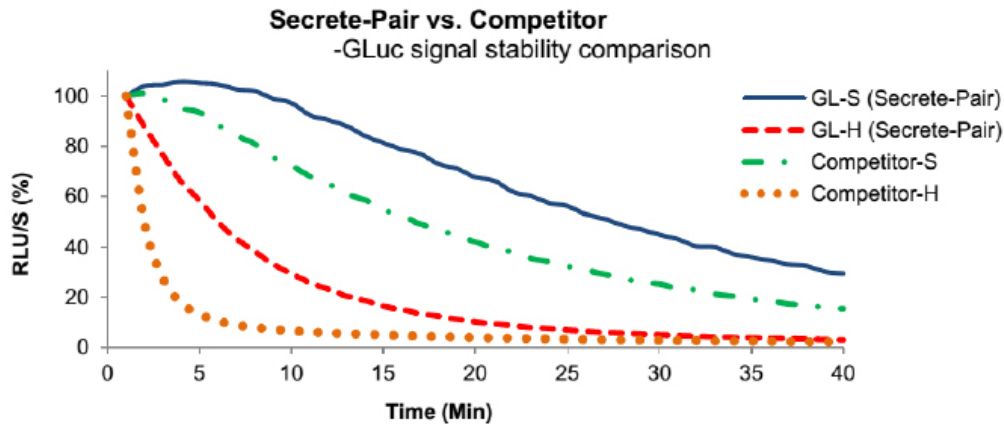


Figure 1. Comparison of GLuc signal stability in different buffer systems from Secretre-Pair and a competitor Gaussia luciferase assay kit. Cell culture medium was collected from cells transfected with the humanized wild type GLuc reporter clones. 10 μ l of the medium was used in each assay. Two buffer systems of each kit were tested and the assays were performed according to the manufacturer protocols. The percentage of signal retained (Y axis) is used as an indicator for signal stability. For both kits, the GLuc activities in buffers with a stabilizer (-S) are much more stable than those in buffers without a stabilizer (-H). However, when compared side-by-side, Secretre-Pair buffer systems provide more stable GLuc signal than the competitor kit.

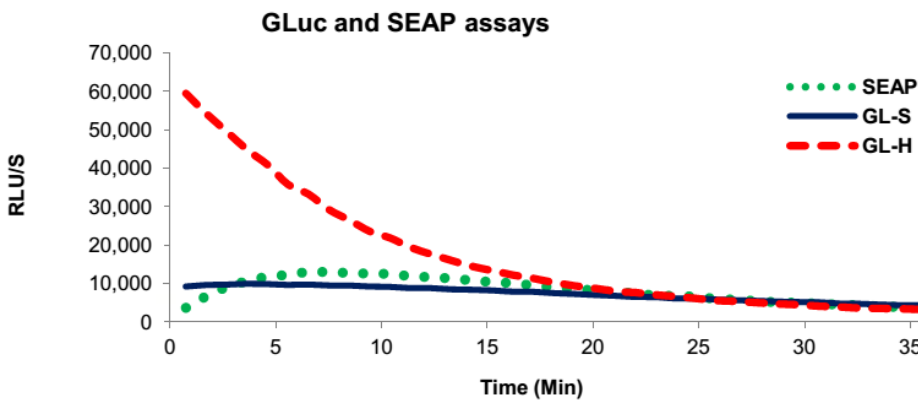
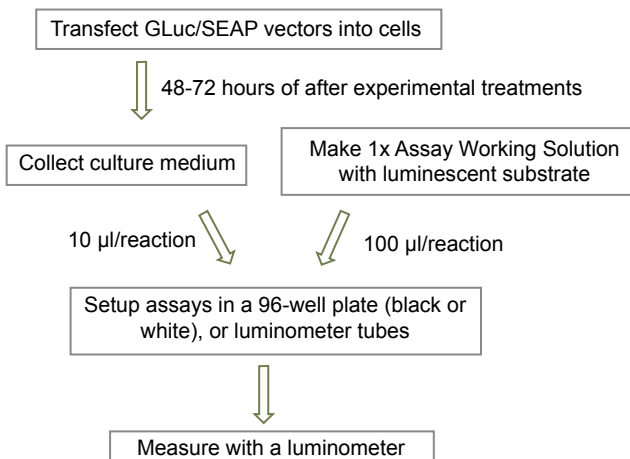


Figure 2. GLuc and SEAP assays. Cell culture medium was collected from cells transfected with GLuc-SEAP dual-reporter clone. 10 μ l of the medium was used in each assay. At the beginning, the GLuc activity in Buffer GL-H is about 4-6 times higher than that in Buffer GL-S. Then it quickly decays. The GLuc activity in Buffer GL-S, however, is much more stable.

Protocol Overview



Related products

- GLuc-ON™ Promoter Reporter Clones
- GLuc-ON™ SEAP Expression Clone
- miTarget™ miRNA 3'UTR Target Clones
- GLuc-ON™ Transcriptional Response Element (TRE) clones

GeneCopoeia, Inc.

9620 Medical Center Drive, Suite 101
Rockville, MD 20850, USA

Email inquiry@genecopoeia.com
Tel +1 (301) 762-0888
Toll free +1 (866) 360-9531
Fax +1 (301) 762-3888
Website www.genecopoeia.com

SDL-031318