

Duo-Luciferase Vectors

Cat. No.	Product	Description	Promoter	Reporter Luc	Control Luc	Marker
ZX001	pEZX-FR01	Firefly (FLuc) + Renilla luciferase (RLuc) cloning vector	None	Firefly (FLuc)	Renilla (RLuc)	Puro
ZX002	pEZX-FR02		SV40			None
ZX003	pEZX-FR03		miniCMV			Puro

Luc-Pair™ Firefly or Renilla Luciferase Assay Kits (Single Luciferase Assay Kits)

	Cat. No.	Luciferase	Product	Price
Ultra Sensitivity	LF007	Firefly (FLuc)	Luc-Pair™ Firefly Luciferase HS Assay Kit (100 rxns)	\$77
	LF008		Luc-Pair™ Firefly Luciferase HS Assay Kit (300 rxns)	\$125
	LF009		Luc-Pair™ Firefly Luciferase HS Assay Kit (1000 rxns)	\$375
	LF010	Renilla (RLuc)	Luc-Pair™ Renilla Luciferase HS Assay Kit (100 rxns)	\$77
	LF011		Luc-Pair™ Renilla Luciferase HS Assay Kit (300 rxns)	\$125
	LF012		Luc-Pair™ Renilla Luciferase HS Assay Kit (1000 rxns)	\$375
High-throughput Enhanced Stability	LF016	Firefly (FLuc)	Luc-Pair™ Firefly Luciferase HT Assay Kit (10 ml)	\$80
	LF017		Luc-Pair™ Firefly Luciferase HT Assay Kit (30 ml)	\$130
	LF018		Luc-Pair™ Firefly Luciferase HT Assay Kit (100 ml)	\$390
	LF019	Renilla (RLuc)	Luc-Pair™ Renilla Luciferase HT Assay Kit (10 ml)	\$80
	LF020		Luc-Pair™ Renilla Luciferase HT Assay Kit (30 ml)	\$130
	LF021		Luc-Pair™ Renilla Luciferase HT Assay Kit (100 ml)	\$390

Contact us:  
GeneCopoeia, Inc.  
9620 Medical Center Dr. 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



Related Products

- Secrete-Pair™ Dual Luminescence Assay Kits
- Secrete Pair™ Gaussia Luciferase Assay Kit
- GLuc-ON™ Promoter Reporter Clones
- miTarget™ miRNA 3' UTR Target Clones
- GLuc-ON™ Transcription Response Element (TRE) Clones
- Luciferase+GFP Dual-Labeled Caner Cell Lines

LPL-031318

• Luc-Pair™ Duo-Luciferase Assay Kits •

Ultra SensitivityEnhanced Stability

Features

- » Duo reporter system: firefly luciferase + Renilla luciferase (as internal control)
- » Ultrasensitive detection of bioluminescence with low background and wide dynamic range compared to fluorescent reporters
- » Gene regulation studies based on the expression of luciferase gene fused to the regulatory elements of interest (promoters, enhancers or 5' or 3' UTRs)
- » All Luc-Pair Assay kits are compatible with GeneCopoeia and Promega FLuc and RLuc vectors

Ultra Sensitivity	Luc-Pair™ Duo-Luciferase HS Assay Kit
Enhanced Stability	Luc-Pair™ Duo-Luciferase Assay Kit 2.0
High-throughput Enhanced Stability	Luc-Pair™ Duo-Luciferase HT Assay Kit

Luc-Pair™ Duo-Luciferase HS Assay Kit

The Luc-Pair™ Duo-Luciferase HS Assay Kit works great for many different eukaryotic cell types (e.g. mammalian, invertebrates, etc.) using micro-plate or single-tube luminometers. It is ideal for many applications, such as measuring the inhibitory effect of a miRNA on a target sequence.



Cat. No.	Product	Price
LF004	Luc-Pair™ Duo-Luciferase HS Assay Kit (100 rxns)	\$149
LF005	Luc-Pair™ Duo-Luciferase HS Assay Kit (300 rxns)	\$357
LF006	Luc-Pair™ Duo-Luciferase HS Assay Kit (1000 rxns)	\$979

Luc-Pair™ Duo-Luciferase HS Assay Kit compared with Promega kit

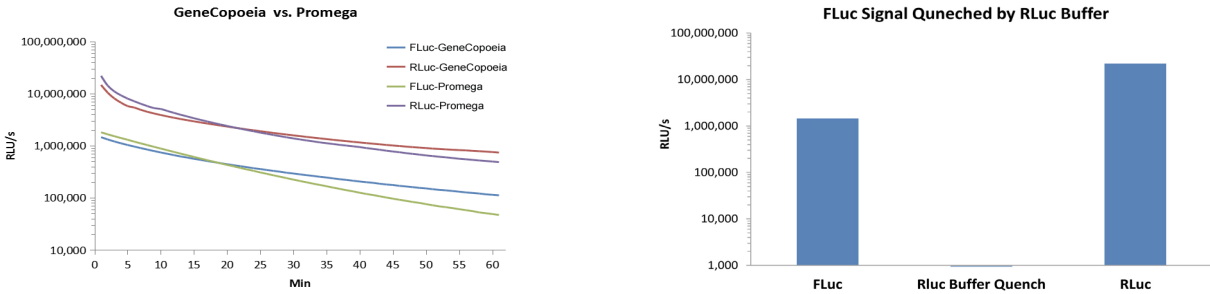


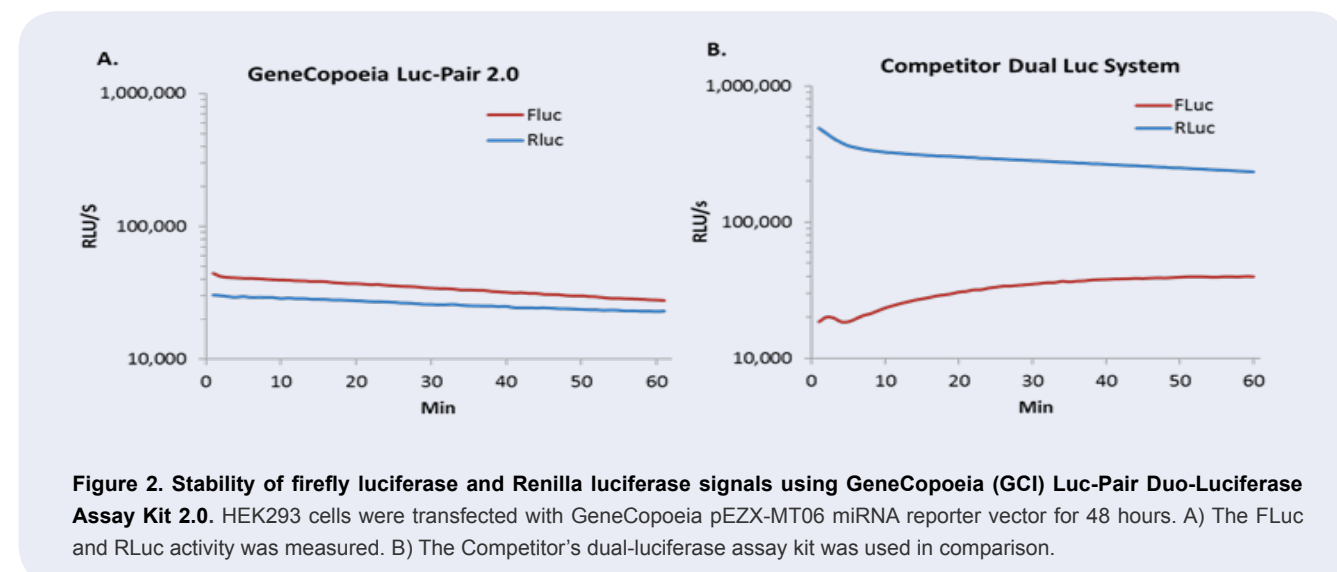
Figure 1. Activity of Firefly luciferase and Renilla luciferase signals using GeneCopoeia Luc-Pair™ Duo-Luciferase HS Assay vs Promega dual-luciferase assay kit. Firefly luciferase activity is quenched with the addition of Luc-H Buffer II.

## Luc-Pair™ Duo-Luciferase Assay Kit 2.0 (Enhanced Stability)

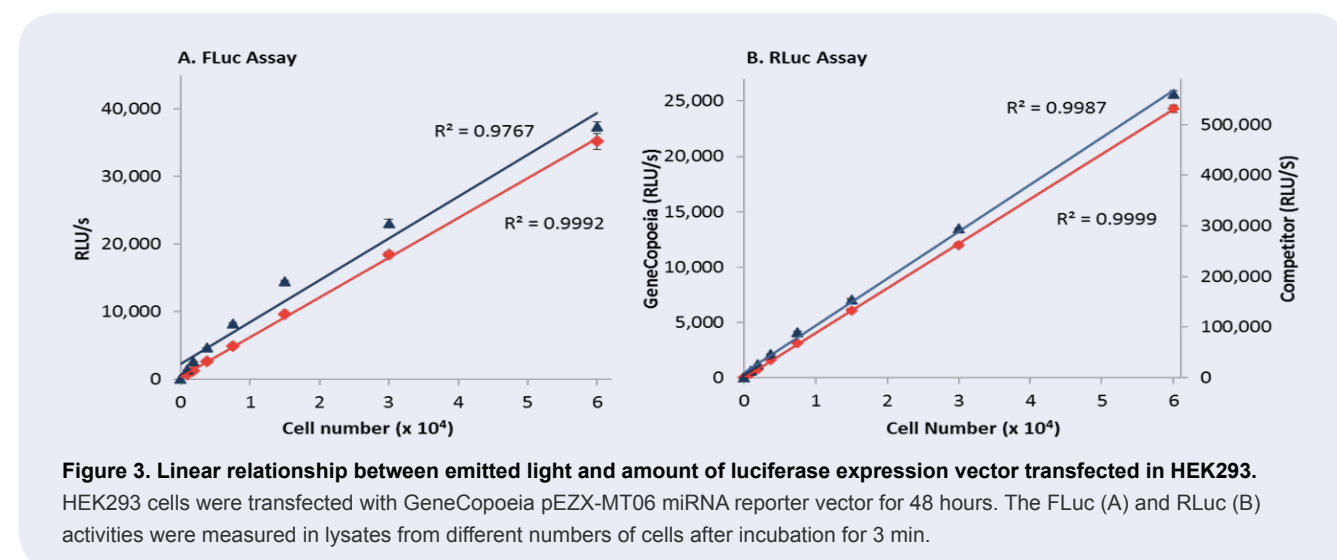
The Luc-Pair™ Duo-Luciferase Assay Kit 2.0 provides a more flexible system in the same convenient 96-well plate format as before for measuring Firefly and *Renilla* luciferases sequentially. Due to its enhanced stability, the Luc-Pair™ Duo-Luciferase Assay Kit 2.0 is the ideal solution for measuring the inhibitory effect of a miRNA on a target sequence.

Cat. No.	Product	Price
LF001	Luc-Pair™ Duo-Luciferase Assay Kit 2.0 (100 rxns)	\$149
LF002	Luc-Pair™ Duo-Luciferase Assay Kit 2.0 (300 rxns)	\$357
LF003	Luc-Pair™ Duo-Luciferase Assay Kit 2.0 (1000 rxns)	\$979

## Luc-Pair™ Duo-Luciferase Assay Kit 2.0 more stable than Promega kit



## Side-by-Side Comparison of Luc-Pair™ 2.0 vs. Promega kit

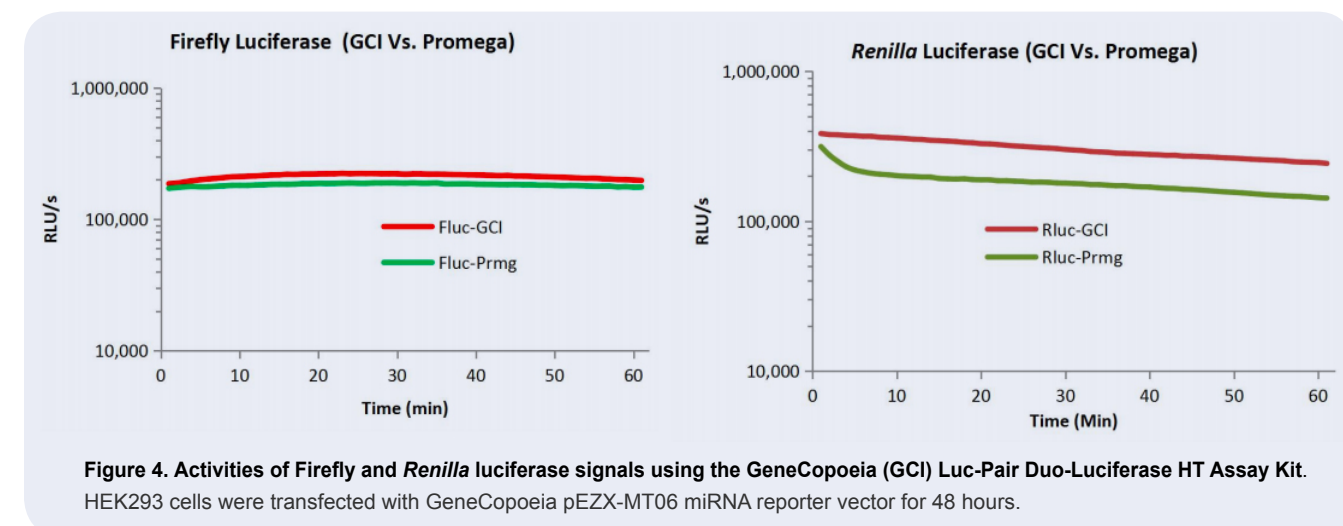


## Luc-Pair™ Duo-Luciferase HT Assay Kit

GeneCopoeia's Luc-Pair™ Duo-Luciferase HT Assay Kit is a convenient system for high-throughput detection of Firefly luciferase (FLuc) and *Renilla* luciferase (RLuc) activities in 96- or 384-well plates and works great for many different eukaryotic cell types (e.g. mammalian, invertebrates, etc.). The Luc-Pair™ Duo-Luciferase HT Assay reagents **can be added directly to cells in growth medium** without washing or preconditioning, and are optimized for use with many different media containing 0–10% serum (e.g. DMEM, RPMI1640, etc.). Firefly luciferase luminescence is produced by one reagent, while a second reagent simultaneously quenches the Firefly luciferase and produces *Renilla* luciferase luminescence.

Cat. No.	Product	Price
LF013	Luc-Pair™ Duo-Luciferase HT Assay Kit (10 ml)	\$156
LF014	Luc-Pair™ Duo-Luciferase HT Assay Kit (30 ml)	\$375
LF015	Luc-Pair™ Duo-Luciferase HT Assay Kit (100 ml)	\$1022

## Luc-Pair™ Duo-Luciferase Assay HT Kit more stable than Promega kit



## Luc-Pair™ Duo-Luciferase Assay HT Kit stability in various culture media

