

Premade Labeled Cancer Cell Lines

Luciferase+GFP Dual Labeled Cancer Cell Lines

GeneCopoeia offers premade stable cancer cell lines with dual labels (luciferase + GFP). They are ready to use and express high levels of luciferase and GFP.

Applications

- Study tumor cell properties using biochemical and genetic methods *in vitro*
- Monitor tumor growth, progression, and metastasis *in vivo* in response to a variety of environmental stimuli, drugs, and other therapeutic treatments.

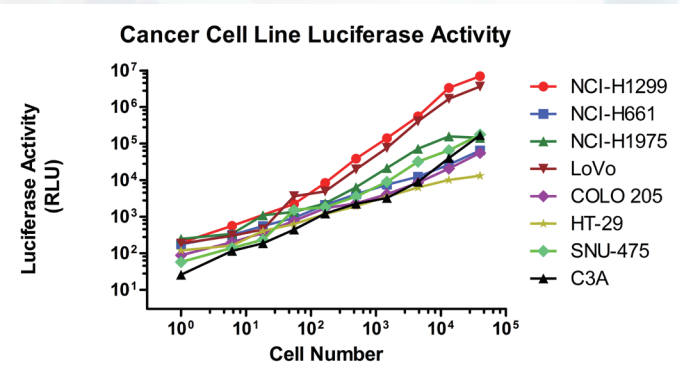
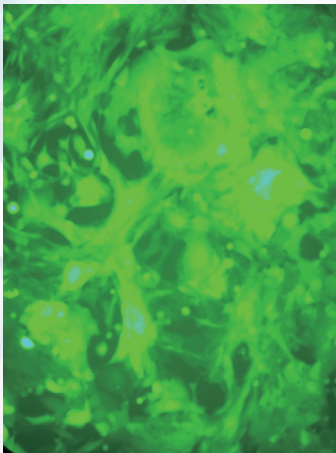


Figure 1. Luciferase activity of dual labeled cancer cell lines

Figure 2. SCL-C03-HLG dual-labeled cancer cell line expressing GFP

GFP fluorescence after exposure for 100 mS at 95% cell confluence using a Nikon Fluorescence Microscope



Detection Marker: Firefly Luciferase and GFP

Cat No.	Cancer Cell Line	Tissue
SCL-C01-HLG	NCI-H1299	Lung
SCL-C02-HLG	NCI-H661	Lung
SCL-C03-HLG	NCI-H1975	Lung
SCL-C04-HLG	LoVo	Colon
SCL-C05-HLG	COLO 205	Colon
SCL-C06-HLG	HT-29	Colon
SCL-C07-HLG	SNU-475	Liver
SCL-C08-HLG	C3A	Liver
SCL-C09-HLG	PLC/PRF/5	Liver
SCL-C010-HLG	Panc 10.05	Pancreas
SCL-C011-HLG	CFPAC-1	Pancreas
SCL-C012-HLG	BxPC-3	Pancreas
SCL-C013-HLG	KATOIII	Stomach
SCL-C014-HLG	NCI-N87	Stomach
SCL-C015-HLG	AGS	Stomach
SCL-C016-HLG	HCC70	Breast
SCL-C017-HLG	MCF7	Breast

Premade Labeled Cancer Cell Lines

GFP Labeled Cancer Cell Lines

GeneCopoeia offers premade stable cancer cell lines labeled with GFP. They are ready to use and express high levels of GFP.

Applications

- Study tumor cell properties using biochemical and genetic methods *in vitro*
- Monitor tumor growth, progression, and metastasis *in vivo* in response to a variety of environmental stimuli, drugs, and other therapeutic treatments.

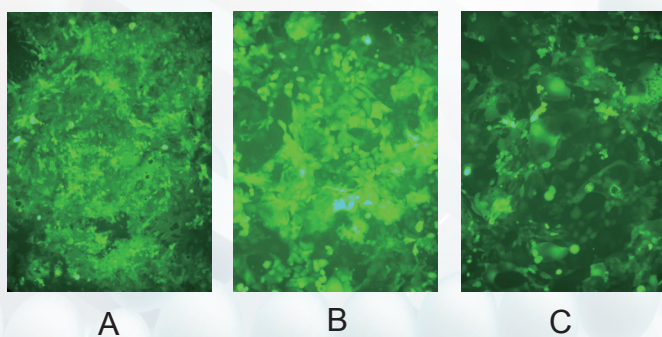


Figure 3. GFP labeled cancer cell lines

- A:** SCL-C08-GFP (C3A Liver Cancer Cell Line)
B: SCL-C09-GFP (PLC/PRF/5 Liver Cancer Cell Line)
C: SCL-C10-GFP (Panc 10.05 Pancreat Cancer Cell Line)

GFP fluorescence after exposure for 100 mS

For more informations about GeneCopoeia™ **Mammalian Stable Cell Line Development Services** or premade labeled cancer cell lines, contact us at inquiry@genecopoeia.com or 1-866-360-9531 (toll-free)

Detection Marker: eGFP

Cat No.	Cancer Cell Line	Tissue
SCL-C01-GFP	NCI-H1299	Lung
SCL-C02-GFP	NCI-H661	Lung
SCL-C03-GFP	NCI-H1975	Lung
SCL-C04-GFP	LoVo	Colon
SCL-C05-GFP	COLO 205	Colon
SCL-C06-GFP	HT-29	Colon
SCL-C07-GFP	SNU-475	Liver
SCL-C08-GFP	C3A	Liver
SCL-C09-GFP	PLC/PRF/5	Liver
SCL-C010-GFP	Panc 10.05	Pancreas
SCL-C011-GFP	CFPAC-1	Pancreas
SCL-C012-GFP	BxPC-3	Pancreas
SCL-C013-GFP	KATOIII	Stomach
SCL-C014-GFP	NCI-N87	Stomach
SCL-C015-GFP	AGS	Stomach
SCL-C016-GFP	HCC70	Brest
SCL-C017-GFP	MCF7	Brest
SCL-C018-GFP	MDA-MB-231	Brest

STA2-091514