



OmicsLink™ Expression-Ready ORF Clones

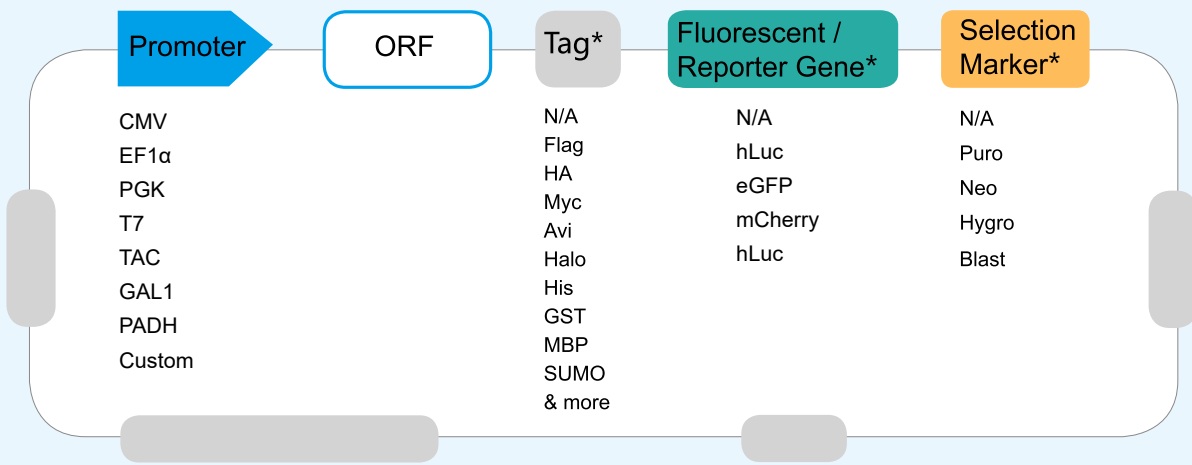
GeneCopoeia offers the largest collection of ORF clones, with genome-wide coverage for human, mouse, and rat. All ORFs can be readily expressed in combination with a variety of fusion tags (fluorescence, antibody, solubility, purification or no tag), promoters and selection markers, for easy expression in different host systems.

Advantages:

- 200+ vector types for mammalian, lentiviral, and more
- Next-day shipping of 140,000 Nextday-ORF™ clones
- Starting from \$69 for selected clones
- Custom services also available



OmicsLink™ ORF Clone Collections



Custom Services:

- ORF mutant clones
- ORF cDNA lentivirus and AAV
- Stable cell lines with ORF clones
- Custom gene synthesis services

Clone Sets (Human Gene or Custom Design)

Set type	# of Genes	# of Clones
CDs	341	430
Cytokines and cytokine receptors	347	420
Drug target genes	5022	6144
GPCRs	526	583
Innate immune proteins	565	702
Ion channels	297	366
Kinases	707	940
Nuclear hormone receptors	56	82
Phosphatases	213	261
Protease inhibitors	111	134
Secretory proteins	1497	1788

OmicsLink™ Expression-Ready ORF Clones

Selected Mammalian Expression Vectors

Vector	promoter	Tag/Reporter	Selection Marker
M02 ^{Next day}	CMV	-	Neo
M35 ^{Next day}	CMV	C-Flag	Neo
M98 ^{Next day}	CMV	C-eGFP (monomeric)	Neo
M61	CMV	IRES2-eGFP	Neo
M56	CMV	C-mCherry	Neo
M29	CMV	N-eGFP	Neo
M03	CMV	C-eGFP	Neo
M06	CMV	N-3×HA	Neo
M11	CMV	N-Flag	Neo
M09	CMV	C-Myc	Neo
M12	CMV	N-3×Flag	Neo
M55	CMV	N-mCherry	Neo
M14	CMV	C-3×Flag	Neo
M07	CMV	C-3×HA	Neo
M77	CMV	C-His	Neo
M01	CMV	N-His	Neo
M43	CMV	N-Myc	Neo
M46	CMV	C-Flag-IRES-eGFP	Neo
M51	CMV	C-His-IRES-eGFP	Neo
M16	CMV	C-eYFP	Neo
M10	CMV	C-Myc-His	Neo
M45	CMV	C-3×HA-IRES-eGFP	Neo
M72	CMV	C-Myc-IRES2-eGFP	Neo
M15	CMV	N-eYFP	Neo
M17	CMV	C-Avi	Neo
M48	CMV	N-Avi-IRES-Biotin ligase	Neo
M50	CMV	C-Halo/Tev protease	Neo
M08	CMV	C-3×HA-His	Neo
M49	CMV	N-Halo/Tev protease	Neo
M05	CMV	N-Avi	Neo
M04	CMV	N-GST/EK protease	Neo
M33	CMV	C-eCFP	Neo
M32	CMV	N-eCFP	Neo

Selected Lentiviral Expression Vectors

Vector	promoter	Tag/Reporter	Selection Marker
Lv105 ^{Next day}	CMV	-	Puro
Lv152	CMV	-	Hygro
Lv242 ^{Next day}	CMV	C-Avi-Flag	Puro
Lv205	CMV	IRES2-eGFP	IRES-Puro
Lv120	CMV	C-3×HA	Puro
Lv102	CMV	N-Flag	Puro
Lv203	CMV	C-Flag-SV40-eGFP	IRES-Puro
Lv241	CMV	C-Flag	Puro
Lv201	CMV	SV40-eGFP	IRES-Puro
Lv207	CMV	SV40-eGFP	IRES-Hygro
Lv130	CMV	C-mCherry	Puro
Lv213	CMV	IRES2-mCherry	IRES-Puro
Lv103	CMV	N-eGFP	Puro
Lv216	CMV	C-Flag-SV40-mCherry	Puro
Lv125	CMV	C-Myc	Puro
Lv111	CMV	N-mCherry	Puro
Lv217	CMV	SV40-hLUC	IRES-Puro
Lv129	CMV	C-Halo	Puro
Lv211	CMV	C-Halo	IRES-Hygro
Lv206	CMV	SV40-mCherry	IRES-Puro
Lv109	CMV	N-eCFP	Puro
Lv128	CMV	C-His	Puro
Lv110	CMV	N-Halo	Puro
Lv123	CMV	C-eYFP	Puro
Lv126	CMV	C-Myc-His	Puro
Lv202	CMV	C-Myc-SV40-eGFP	IRES-Puro
Lv204	CMV	C-Myc-SV40-mCherry	IRES-Puro
Lv104	CMV	N-eYFP	Puro
Lv108	CMV	N-Avi	Puro
Lv127	CMV	C-eCFP	Puro

Not all available tags, reporter genes and selection markers are listed here. Please visit our website www.genecopoeia.com for a complete list.

Contact us:

GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,
Rockville, MD 20850 USA



Email
Tel
Fax
Website

inquiry@genecopoeia.com
+1 (301) 762-0888
+1 (301) 762-3888
www.genecopoeia.com