

# Complete Safe Harbor System

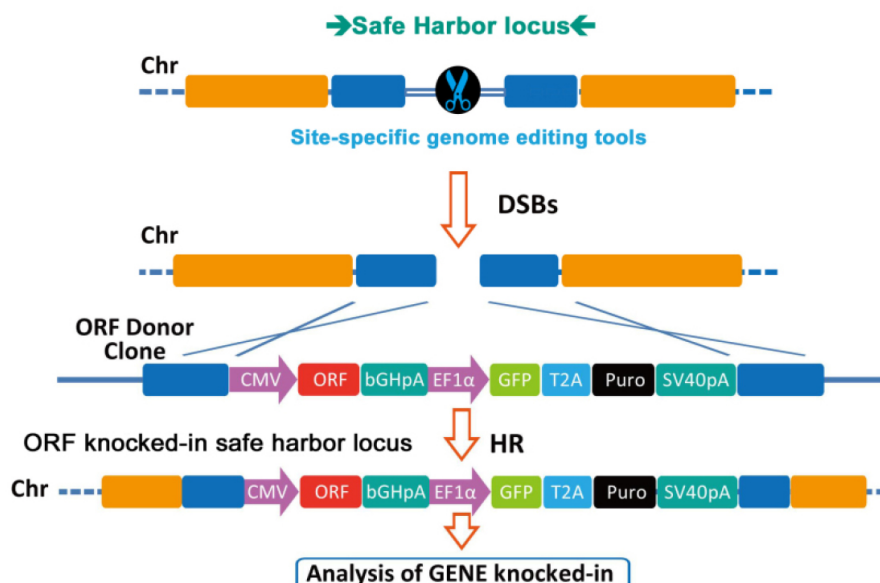
● Efficient CRISPR- or TALEN- mediated gene knock-in for transgenesis ●

GeneCopoeia's Safe Harbor system is the ideal solution for expressing transgenes in cells and animals. Transgene insertion at the AAVS1 and ROSA26 sites in humans and mice, respectively, permits consistent and stable expression with no known adverse effects on fitness.

GeneCopoeia now offers a large variety of individual Safe Harbor kits with knock-in donor cloning vectors, ORF knock-in donors, CRISPR Cas9 nuclease knock-in donors, and Cas9 Safe Harbor stable cell lines, providing a convenient means to carry out genome editing applications with high efficiency.

## Advantages

- ◎ **Safe integration.** Ensures consistent, stable transgene expression with no known adverse effects on cell viability or fitness.
- ◎ **Comprehensive system.** CRISPR or TALEN-based kits come with a choice of donor cloning vectors, premade knock-in clones, or made-to-order knock-in clones, as well as verification PCR primers.
- ◎ **Compatible knock-in ORFs.** More than 20,000 human and more than 15,000 mouse sequence-verified ORFs in custom-built knock-in donor clones.



## Do-it-yourself knock-in kits: With donor cloning vector

Product	Catalog#	Price
Human AAVS1 safe harbor gene knock-in kit (with donor)	SH004, SH029	\$1395
Mouse ROSA26 safe harbor gene knock-in kit (with donor)	SH054, SH079	\$1395

## ORF Knock-in: Pre-made Safe Harbor donor clones carrying ORFs

Product	Catalog#	Price
Human AAVS1 safe harbor gene knock-in kit (without donor)	SH000, SH025	\$1195
Mouse ROSA26 safe harbor gene knock-in kit (without donor)	SH050, SH075	\$1195
AAVS1/ROSA26 knock-in ORF donor clone	Varies	From \$495

## Cas9 knock-in: Express the CRISPR Cas9 nuclease from Safe Harbor sites

Product	Catalog#	Price
<b>Kits</b>		
Human AAVS1 Cas9 safe harbor gene knock-in kit	SH012, SH013, SH037, SH038, SH039	\$1595
Mouse ROSA26 Cas9 safe harbor gene knock-in kit	SH062, SH063, SH064, SH087, SH088, SH089	\$1595
<b>Stable cell lines</b>		
Human cell line NCI-H1299 stably expressing CRISPR Cas9, single clone (CMV / Puro / AAVS1)	SL501	\$1395
Human cell line HEK293T stably expressing CRISPR Cas9, single clone (CMV / Puro / AAVS1)	SL502	\$945
Human cell line HeLa stably expressing CRISPR Cas9, single clone (CBh/ Hygro / AAVS1)	SL503	\$1395
Mouse cell line Neuro2a stably expressing CRISPR Cas9, single clone (CBh/ Hygro / ROSA26)	SL509	\$1395
Visit our website, <a href="http://www.genecopoeia.com/product/cas9-cell-line">www.genecopoeia.com/product/cas9-cell-line</a> , to see a complete list.		
Custom order	GeneCopoeia also offers custom-built Cas9-expressing stable cell line services.	

### Contact us:

#### GeneCopoeia, Inc.

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

Email	<a href="mailto:inquiry@genecopoeia.com">inquiry@genecopoeia.com</a>
Tel	+1 (301) 762-0888
Toll free	+1 (866) 360-9531
Fax	+1 (301) 762-3888
Website	<a href="http://www.genecopoeia.com">www.genecopoeia.com</a>