

# Mammalian Stable Cell Line

## Development Services



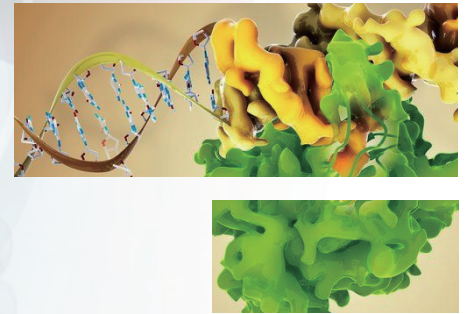
We now offer state-of-the-art services for establishing stable cell lines that **meet your specific research needs.**



### Stable cell lines with ORF clones

- Cell-based functional assays
- Protein production for biochemical assays
- Protein production for crystal or NMR structural determination
- Fusion tagging for pulldown/immunoprecipitation
- Fusion tagging for live cell imaging
- Monoclonal antibody production
- Drug target analysis

#### Applications



### Stable cell lines expressing shRNAs

- Expression profiling to identify candidate downstream genes in a network or pathway
- Determination of gene function in cells
- Drug target validation

#### Applications

### Stable cell lines constructed with CRISPR or TALEN

- Gene knockout
- Introduction of point mutations or defined insertions/deletions
- Correction of disease mutations back to wild type
- Gene tagging
- Replacing promoter or gene sequence
- Transgene knockin at human and mouse safe harbor sites or at a customer-specified locus

#### Applications

# Mammalian Stable Cell Line Services

## Advantages

### Stable cell lines

ORF	shRNA	Genome editing
-----	-------	----------------

State-of-the-art facility for stable cell line development and QC following industrial standards.

√	√	√
---	---	---

Largest in-house collection of sequence-verified and expression-ready ORF, shRNA clones and lentiviral particles reduces turnaround time and ensures the highest quality.

√	√	N/A
---	---	-----

Large selection of promoters including inducible promoters for toxic gene expression.

√	√	N/A
---	---	-----

Large selection of different cell based assays for kinases, GPCR, ion channel, growth factor, and other protein-protein or protein-compound reactions.

√	√	N/A
---	---	-----

Large selection of vectors with different fusion tags such as Flag<sup>®</sup>, HA, V5, His, and GFP to meet your needs.

√	N/A	N/A
---	-----	-----

DHFR or GS gene amplification stable cell lines for protein production and process development of Fed-batch cell culture.

√	N/A	N/A
---	-----	-----

MAR option for increased open-chromatin site integration.

√	N/A	N/A
---	-----	-----

≥ 70% knockdown efficiency in shRNA applications.

N/A	√	N/A
-----	---	-----

Complete solutions for CRISPR- or TALEN-based targeted genomic modifications.

N/A	N/A	√
-----	-----	---

We also provide **premade labeled cancer cell lines** either GFP labeled or dual labeled (luciferase + GFP). Contact us for more details!

### GeneCopoeia, Inc.

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

Website: [www.genecopoeia.com](http://www.genecopoeia.com)

For **more information**, or to **request a quote** for services, contact us at

[inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)

or **1-866-360-9531 (toll free)**

STA1-091514