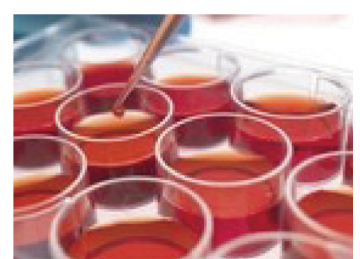


# Mammalian Stable Cell Line

## 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

Applications

- 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



### Stable cell lines expressing shRNAs

Applications

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

### Stable cell lines constructed with CRISPR or TALEN

Applications

- 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

# Mammalian Stable Cell Line Services

## Advantages

- › 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.
- › Large selection of promoters including inducible promoters for toxic gene expression.
- › Large selection of different cell based assays for kinases, GPCR, ion channel, growth factor, and other protein-protein or protein- compound reactions.
- › Large selection of vectors with different fusion tags such as Flag®, HA, V5, His, and GFP.
- › DHFR or GS gene amplification stable cell lines for protein production and process development of Fed-batch cell culture.
- › MAR option for increased open-chromatin site integration.
- › ≥ 70% knockdown efficiency in shRNA applications.
- › Complete solutions for CRISPR- or TALEN-based targeted genomic modifications.

We also provide premade stable cancer cell lines either GFP labeled, dual labeled (luciferase + GFP) or Cas9 stable cell lines. 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 )

STA1-031518