

GeneCopoeia ORF cDNA Clones

Including OmicsLink™, ORFEXPRESS™ Gateway® PLUS, and ORFeome clones

Product format	Storage and Handling
Glycerol stock	<ol style="list-style-type: none"> 1. Upon receiving, mix the glycerol stock well before streaking it onto selective plates to ensure that selective makers are not lost. Then freeze the glycerol stocks immediately at -70°C to -80°C for long term storage. Note: Plates of streaked bacteria can be sealed with Parafilm® and stored upside-down at 4°C for several weeks. 2. Glycerol stocks can potentially contain mutants arising from cultures used to prepare them, or can deteriorate during storage. Inoculate liquid cultures from healthy single colonies picked from a freshly streaked selective plate to ensure that cells growing in the culture are all descended from a single founder cell and have the same genetic makeup. 3. Avoid repeated thawing and re-freezing of glycerol stocks as this can reduce the viability of bacteria. Note: Scraping from the top of the frozen glycerol stock often provides enough cells to streak on a selective plate and can help avoid fully thawing the stock. Alternatively, removing some culture from the middle of stock as it unthaws can provide more cells.
Purified plasmid	Store at 4°C for short term storage or -80°C for long term storage.

Plasmid Preparation

For lentiviral clones, please see GeneCopoeia [Plasmid Prep Protocol for Lentiviral Clones](#).

For non-viral clones, please follow instruction manuals for plasmid prep kits from manufacturers, such as Qiagen.

Related Products from GeneCopoeia

[EndoFectin™ Transfection Reagents](#)

[GeneCopoeia Chemically Competent Cells](#)

[OmicsLink™ Anti-tag Antibodies](#)

[Lentiviral Products and Services](#)

[All-in-One™ qPCR mix and primers](#)

[OmicsLink™ shRNA Clones](#)

[GeneCopoeia microRNA Products](#)

[GLuc-ON™ Promoter Clones](#)

[Custom Gene Synthesis Services](#)

[CoolCutter™ SUMO Protease](#)

Warranty, cancellation, Quality assurance

All ORF cDNA clones (the Product) are subject to the following policy. Customers are referred to as Buyers.

Changes or Cancellation

Any changes or cancellation must be done by 5 p.m. EST on the day the order is placed. Changes or cancellation to the order after 5 p.m. EST (prior to shipment of the Product ordered) is allowed only upon mutual agreement between GeneCopoeia and the Buyer and is subjected to a cancellation fee 20% of the purchase order. Once shipment of the Product has been made by GeneCopoeia, cancellation will not be permitted.

Quality Assurance

GeneCopoeia quality assurance policy ensures that all **OmicsLink™ ORF cDNA Expression Clones are free of artificial frame-shifting variations** including deletions(s) and additions(s) of base(s) and translation termination mutation(s) (non-stop codon to stop codon as the result of point mutation). **This policy does not cover** point variations due to possible naturally occurring polymorphisms and/or mutations that may be introduced by PCR in very rare cases. If Buyer requests a specific version of a gene from different cDNA libraries, GeneCopoeia will make its best effort to obtain that particular version but without guarantee of success. For this recloning effort, Buyer is to pay a fee that will be fairly negotiated between Buyer and GeneCopoeia. **Splicing variants are replaceable** subject to availabilities and a modest fee to cover the cloning cost. In the event that requested ORF clone(s) of specific splicing variant(s) are not available, GeneCopoeia will make its best effort to clone it and Buyer will pay a fee that will be fairly negotiated between Buyer and GeneCopoeia.

Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund or credit. This limited warranty shall not extend to anyone other than the original Buyer of the Product. Notice of nonconforming products must be made to GeneCopoeia within 1 year of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product or a refund or credit limited to the actual purchase price. Shipping charges will not be credited. GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

Limited Use License

A limited use license is granted to the Buyer of the Product. The Product shall be used by the Buyer for internal research purposes only and should not be transferred (in any way) to any third party. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale (includes but is not limited to the transferring of full-length ORF(s) from GeneCopoeia ORF clones to other expression vectors), or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us.

GeneCopoeia, Inc.

9620 Medical Center Drive, #101
Rockville, MD 20850
Tel: 301-762-0888; Fax: 301-762-3888
Email: support@genecopoeia.com
Web: www.genecopoeia.com