



Lentifect™ Lentiviral Particles: Cancellation, Quality Assurance, Limited Use License and Warranty

All lentiviral particles (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 the initial titer in TU/mL as determined by qRT-PCR for the lentiviral particles purchased will be as indicated in the accompanying datasheet. However, GeneCopoeia does not guarantee the resulting transduction efficiency for specific target cells as the transducibility may vary from one cell type to another. The following statements are pursuant to Lentifect™ shRNA lentiviral particles only. The quality assurance guarantee of knockdown efficiency for the OmicsLink™ shRNA Expression Clones does not extend to the Lentifect™ shRNA lentiviral particles. If knockdown efficiency is of considerable concern, we recommend purchasing your shRNA of interest as an OmicsLink™ shRNA Expression Clone first to validate the knockdown efficiency before proceeding with packaging into Lentifect™ particles.

Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. Notice of nonconforming or damaged products must be made to GeneCopoeia within 30 days from the day of receipt of the Product. 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 Buyer with a refund or credit. This limited warranty shall not extend to anyone other than the original Buyer 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, 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
T: 301-762-0888/F: 301-762-3888
Email: support@genecopoeia.com