

Tear Away License for Gaussia Luciferase Users

By opening this package or by using the materials enclosed within; Recipient is legally bound and accepts the following terms and conditions:

Non-Commercial entities: This product is covered by US Patent Nos. 6,232,107, 6,436,682, 6,780,974, 7,045,599, 7,109,315 7,238,497 World Patent WO1999/049019 entitled "LUCIFERASES, FLUORESCENT PROTEINS, NUCLEIC ACIDS ENCODING THE LUCIFERASES AND FLUORESCENT PROTEINS AND THE USE THEREOF IN DIAGNOSTICS" and other patents that are the legal property assigned to Prolume Ltd / NanoLight Technologies, POB 2746 Pinetop, Arizona, USA 85935. This product is licensed only to the purchasing laboratory-research group. Recipient agrees <u>not to transfer</u> this plasmid or derivatives of this vector, or product to any other laboratory, person or research group, even if within the same institution.

Recipient agrees not to alter or make any changes to the nucleotide coding sequence or secretory coding sequence of the luciferase(s) or fluorescent protein(s) contained within without prior written permission from Prolume Ltd / NanoLight Technologies (www.nanolight.com). Recipient agrees not to file for any patent rights or to make any inventions claiming any portion of the luciferase(s) or fluorescent protein(s) within this material without prior written permission from Prolume Ltd/NanoLight Technologies.

Commercial Recipients defined herein as ANY recipient using Prolume/NanoLight materials for any commercial purpose whatsoever, not limited to and including the search for new drugs, or new drug discovery methods, research, develop, derive products, engage In the sale or license of any products, or make inventions by use of the luciferase and fluorescent protein materials enclosed, fully agree to the terms mentioned above for Non-Commercial entities; AND ADDITIONALLY agree to and agree to be bound to use the materials FOR EVALUATION PURPOSES ONLY. Commercial Recipient agrees to obtain a valid license from Prolume/NanoLight Technologies for any use of these or any related materials beyond 180 days of receipt, or destroy all constructs and progeny containing or derived from the enclosed materials beyond 180 days from receipt. Multiple product purchases does not grant recipient continuous license extensions.

GeneCopoeia, Inc.
9620 Medical Center Drive
Suite 101
Rockville, MD 20878
USA
+1 (301) 762-0888
inquiry@genecopoeia.com