



www.nanolight.com

www.luciferin.com

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 Patents 6,232,107, 6,436,682, 6,780,974, 7,045,599, 7,109,315, 7,238,497 and world counterparts 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 Technology. This product is licensed only to the purchasing laboratory-research group. Recipient agrees not to transfer this plasmid or derivatives of this vector, or product to any other laboratory, person or research group, even 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.

Universities, Non-profit Foundations, & Commercial For-profit Entities If product and/or materials derived therefrom are being used in any commercial way you are defined as a "Commercial Recipient" meaning: using in search of new drugs, or new drug discovery methods, research to invent, patent, or for any gain, develop, derive products, engage in the sale or license of any products, or make inventions by use of the luciferases and/or fluorescent protein materials enclosed, fully agree to the terms mentioned above for Non-Commercial entities; AND ADDITIONALLY agree to and are hereby 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 materials beyond 180 days of receipt, or destroy all derivatives from/and the original enclosed materials 180 days from receipt.**