

RapidReporter® Limited-use Label License

This License Agreement applies to the purchased RapidReporter® materials: This license applies to the materials contained in the attached package (the “**materials**”). The use of such materials is subject to the terms and conditions of this Label License. Any use of the materials other than in accordance with these terms and conditions will amount to a breach of the license for which you as the recipient may be liable, including for patent infringement. The materials comprising the product in the attached package are sold under license from GeneStream Pty Ltd and are covered by patent families WO02072844 and WO1999/049019; including U.S. Patent Nos. 6,232,107; 6,436,682; 6,780,974 and various other patents worldwide.

Recipient agrees to be bound by these terms and conditions: By purchase of this product and upon opening of this package or by use of the materials enclosed within, the recipient organization and each of its employees, agents or contractors having access to such materials (the “**Recipient**”) of this product agrees to be bound by and accepts the following terms and conditions for the use of these materials.

Before using the materials For-profit Entities (as defined below) must sign these terms to acknowledge acceptance and fax or e-mail to GeneStream Pty Ltd at the contact details specified below.

Limited use permitted: GeneStream Pty Ltd’s RapidReporter (whether supplied as-is or incorporated into another product) must only be used with the reporter gene contained therein and with reagents provided by GeneStream Pty Ltd or its licensed distributor, Active Motif. Use with any other reporter gene or reagent will amount to a breach of this label license, in which case no warranty whatsoever is given by GeneStream Pty Ltd or any licensee or distributor and the user of RapidReporter hereby indemnifies GeneStream and its distributors and licensees in respect of such use or application. In the event of such breach, GeneStream Pty Ltd will be entitled to immediately terminate this license and require delivery-up or destruction of all RapidReporter material in the licensee’s possession or control and all information or materials related to use of the RapidReporter material and require a certificate of confirmation that all such materials have been delivered to GeneStream or destroyed.

Non-profit Entities: This product is licensed only to the purchasing laboratory/research group. Recipient hereby agrees not to transfer the materials or derivatives of the materials including any vector, nucleotide sequence or protein to any other laboratory, person or research group, even if within the same institution or organization.

Recipient hereby agrees not to alter nor make any changes to the nucleotide coding sequence of the materials, or derived from the materials, without prior written permission from GeneStream Pty Ltd.

Recipient agrees not to file for any patent rights or file claims to any inventions disclosing information about any portion of the reporter gene (from transcription start site to transcription stop site) contained within the materials without prior written permission from GeneStream Pty Ltd.

A “**Non-profit Entity**” means (1) a university or other institution of higher education located in any country; (2) an entity of the type described in section 501(c)(3) of the Internal Revenue Code of 1954 [26 U.S.C.

501(c)(3)] and exempt from taxation under section 501(a) of the Internal Revenue Code [26 U.S.C. 501(a)]; or (3) any Non-profit Entity located in a foreign country that would qualify as a Non-profit Entity under condition (2) of this section if it were located in the United States.

Any entity that is not a Non-profit Entity is considered a For-profit Entity (“**For-profit Entity**”). A Non-profit Entity performing research for a For-profit Entity or organization will, for the purposes of this label license, be considered to be a For-profit Entity.

For-profit Entities: For-profit Entities receiving this product agree to be bound by all the terms and conditions referred to above AND ADDITIONALLY agree to and are hereby bound to use the materials **FOR EVALUATION PURPOSES ONLY**. For-profit Entities purchasing this product shall have six (6) months to evaluate the product in their research. Thereafter, such For-profit Entities shall either contact GeneStream Pty Ltd and take a license for their continued use of the product or destroy the product and all derivatives thereof and notify GeneStream Pty Ltd of such destruction.

Not for use in high-throughput drug screening: Irrespective of the status of an organization as “Non-profit” or “For-profit” under this label license, the product must not in any circumstances be used in high-throughput drug screening without the prior written consent of GeneStream Pty Ltd. For the purposes of this label license, “high-throughput drug screening” means testing more than 100,000 compounds in a year or more than 100,000 samples in a single experiment.

Any Recipient that does not accept these label license terms shall return the unopened package and materials to Active Motif for a full refund.

Recipient organization name:	
Purchase order or sales order number:	
Signed by: (for and on behalf of the Recipient organization)	
Printed name:	
Address:	
Phone:	
E-mail:	

**Complete the above table and then fax to GeneStream Pty Ltd. at +61 8 9245 8586,
or scan and e-mail to: info@genestream.com.au.**