

**Silencer® GAPDH siRNA,  
In Vivo Ready**

Store at or below –20°C.  
*Do not store in a frost-free freezer.*

**Catalog # (P/N):** 4404025

**Amount:** 250 nmol

**Appearance:** Powder

**Target Information:**

Gene Symbol: GAPD  
Full Gene Name: Glyceraldehyde-3-phosphate dehydrogenase  
Organism(s): Human, Mouse, and Rat  
RefSeq Number(s): NM\_002046 (human), NM\_008084.2 (mouse), and NM\_017008 (rat)  
Entrez Gene ID(s): 2597 (human), 14433 (mouse), and 24383 (rat)

**Format:** Annealed

**Purity:** In Vivo Ready (HPLC purified)

**Storage Conditions:** Store at or below –20 °C. *Do not store in a frost-free freezer.* (Dried oligonucleotides are shipped at ambient temperature.)

**Safety Information:** Read the Safety Data Sheet, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

---

**USER INFORMATION**

---

**Product Description:**

*Silencer®* GAPDH siRNA, In Vivo Ready, provides a positive control for experiments involving *Silencer* siRNA delivery to animals.

In Vivo Ready siRNAs are high quality siRNAs that are purified especially for introduction into animals. Each siRNA strand is individually purified by HPLC, desalted, and annealed with its complementary strand. In Vivo Ready siRNAs are then further purified utilizing a process that removes excess salt, via a semi-permeable membrane. The result is a highly pure siRNA with minimal salt content, suitable for in vivo applications. In Vivo Ready siRNAs are then filtered through a 0.2 µm pre-sterilized filter and tested for the presence of endotoxin.

At a concentration of 50 µM in de-ionized water, In Vivo Ready siRNA contains <5.0 mM Na<sup>+</sup>, <0.06 mM K<sup>+</sup>, and <0.02 mM Mg<sup>2+</sup>.

*Silencer* GAPDH siRNA, In Vivo Ready, has the same sequence that has been shown to reduce the expression of GAPDH by 70-95% in several common human, mouse and rat cell lines, and is purified to rigorous In Vivo Ready specifications.

**Handling Instructions:**

RNA oligonucleotides are susceptible to degradation by exogenous ribonucleases introduced during handling. Wear gloves when handling this product. Use RNase-free reagents, tubes, and barrier pipette tips.

Upon receipt, store dried oligonucleotides at 4°C, or in a non-frost-free freezer at or below –20°C (dried oligonucleotides are shipped at ambient temperature). For long-term storage, store at or below –20°C in a non-frost-free freezer.

Use standard biological sterile techniques when handling In Vivo Ready siRNA that will be administered to animals.

**Using In Vivo Ready siRNA with  
InvivoFectamine® 2.0 Reagent:**

InvivoFectamine® 2.0 Reagent (P/N 1388501) is a proprietary, animal-origin-free, lipid-based transfection reagent for systemic, in vivo siRNA delivery to mouse liver tissue. InvivoFectamine 2.0 Reagent is ideally suited for use with Ambion *In Vivo* siRNA, with high in vivo transfection efficiency in liver following tail-vein injection. Low volume delivery of siRNA using low pressure, combined with the low toxicity of InvivoFectamine 2.0 Reagent, avoids a stress response in the animal.

**Resuspension of siRNA for use with InvivoFectamine® 2.0 Reagent**

1. Briefly centrifuge the tube to ensure that the dried siRNA is at the bottom of the tube.
2. Resuspend the siRNA in nuclease-free sterile water and vortex to thoroughly resuspend.

For best results, prepare the siRNA stock solution at the highest concentration that is workable for your experiments. Dilute the siRNA stock as needed for immediate use.

The working siRNA concentration for use with InvivoFectamine 2.0 is 200 µM (~3 mg/mL).

A calculator for suspension of dry oligonucleotides is available at:  
[www4.appliedbiosystems.com/techlib/append/oligo\\_dilution.html](http://www4.appliedbiosystems.com/techlib/append/oligo_dilution.html)

**Storage of resuspended siRNA:** Store at or below  $-20^{\circ}\text{C}$ . Stock solutions of siRNA at concentrations  $\geq 2\text{ }\mu\text{M}$  can undergo up to 50 freeze-thaw cycles without significant degradation. Storage in a frost-free freezer is not recommended, however. Long-term storage at  $-70^{\circ}\text{C}$  has traditionally been recommended, but siRNA stock solutions at concentrations  $\geq 2\text{ }\mu\text{M}$  can be stored at  $-20^{\circ}\text{C}$  for extended periods (up to 1 year).

#### **Preparation of InvivoFectamine® 2.0-siRNA complexes and in vivo delivery**

Follow the instructions provided with InvivoFectamine 2.0 Reagent, available at the web catalog page at [www.invitrogen.com](http://www.invitrogen.com) (search for InvivoFectamine 2.0). An siRNA dose of 7 mg/kg is recommended as a starting point for experiments. This dose corresponds to 200  $\mu\text{L}$  of a 0.7 mg/mL solution injected into a 20-g mouse.

#### **Using In Vivo Ready siRNA with Other In Vivo Delivery Strategies:**

##### **Resuspension of In Vivo Ready siRNA**

Follow the instructions for resuspension with use with InvivoFectamine 2.0 Reagent. At step 2, resuspend the siRNA in sterile water or a sterile buffer appropriate for your application. Common examples are provided below.

##### **Suggested Buffers for Systemic Delivery**

- Sterile, phosphate buffered saline (PBS)
- Sterile saline (0.9% NaCl), or variants containing sugars such as mannitol or glucose (5–15%)
- Ringer's solution: 147 mM NaCl, 4 mM KCl, 1.13 mM  $\text{CaCl}_2$

##### **Suggested Non-irritating Buffers for Central Nervous System Delivery**

- Sterile saline (0.9% NaCl)
- Isotonic buffer (100 mM potassium acetate, 30 mM HEPES-KOH, 2 mM magnesium acetate, 26 mM NaCl, pH 7.4)

Store resuspended siRNA as described under "Resuspension of siRNA for use with InvivoFectamine 2.0 Reagent."

##### **Suggested dosing for a 20-g mouse**

A typical volume for systemic delivery is  $\sim 200\text{ }\mu\text{L}$  (for low pressure injection) at concentrations of 50–500  $\mu\text{M}$ . This combination corresponds to an siRNA dosing range of 5–50 mg/kg for a 20-g mouse.

Dose (mg/kg)	nmol siRNA/Dose	Concentration for 200 $\mu\text{L}$ Dose ( $\mu\text{M}$ *)
1.0	1.5	7.4
5.0	7.4	37.0
10.0	14.8	74.0
20.0	29.6	148.0
50.0	74.0	370.0

\* In Vivo Ready siRNAs are soluble in aqueous solution at concentrations up to 1.5 mM.

##### **For more information**

Visit these web resources for more information on buffers and injection routes:  
[www4.appliedbiosystems.com/RNAi/invivo](http://www4.appliedbiosystems.com/RNAi/invivo)

[www.invitrogen.com](http://www.invitrogen.com) > **Products & Services** > **Applications** > **RNAi, Epigenetics & Gene Regulation** > **RNAi** > **in vivo RNAi** > **in vivo RNAi Protocols**

---

#### **RELATED PRODUCTS**

---

##### **Anti-GAPDH, Mouse Monoclonal 6C5**

P/N AM4300

Ideal for detecting knockdown of GAPDH at the protein level by Western blot or immunofluorescence.

##### **Ambion® In Vivo Pre-designed and Validated siRNAs**

P/N Various (see [www4.appliedbiosystems.com/geneassist](http://www4.appliedbiosystems.com/geneassist))

An all-new class of modified siRNAs with superior serum stability for in vivo applications, designed with the *Silencer*® Select algorithm. Ambion® In Vivo siRNAs are non-toxic in vivo (mouse) and non-immunogenic (cell-based assays), while exhibiting potency and specificity equivalent to or better than *Silencer* Select siRNAs.

##### **InvivoFectamine® 2.0 Reagent**

P/N 1377-505, 1377-501

A proprietary lipid-based reagent designed to efficiently deliver siRNA to mouse liver tissue in an easy-to-complex preparation. See [www.invitrogen.com](http://www.invitrogen.com).

##### **TaqMan® Gene Expression Assays**

[www.allgenes.com](http://www.allgenes.com) or [www4.appliedbiosystems.com/geneassist](http://www4.appliedbiosystems.com/geneassist)

A comprehensive collection of over 700,000 probe and primer sets for quantitative gene expression analysis using real-time PCR. Search the GeneAssist™ Atlas at [www4.appliedbiosystems.com/geneassist](http://www4.appliedbiosystems.com/geneassist) to find suggested TaqMan Gene Expression Assays corresponding to your siRNA targets of interest.

---

## QUALITY CONTROL

---

<b>Identity:</b>	The mass of a sample of each single-stranded RNA oligonucleotide is analyzed using MALDI-TOF mass spectrometry and compared to the calculated mass.
<b>Purity:</b>	Analytical HPLC of a sample of the final purified single-stranded RNA oligonucleotides is used to confirm $\geq 95\%$ purity.
<b>Annealing:</b>	A sample of the annealed siRNA is analyzed by nondenaturing gel electrophoresis.
<b>Endotoxin:</b>	A <i>Limulus</i> Amebocyte Lysate (LAL) assay of a sample of the siRNA is used to confirm $<10$ Endotoxin Units (EU) per nmol.

---

## OTHER INFORMATION

---

**Safety Data Sheets:** Safety Data Sheets (SDSs; previously known as MSDSs) for any chemical product supplied by Applied Biosystems or Ambion are available 24 hours a day. At [www.appliedbiosystems.com](http://www.appliedbiosystems.com), select Support, then SDS/MSDS. Search by chemical name, product name, product part number, or SDS/MSDS part number. Right-click to print or download the SDS of interest. At [www.ambion.com](http://www.ambion.com), go to the web catalog page for the product of interest. Select SDS/MSDS, then right-click to print or download. Or, e-mail (MSDS\_Inquiry\_CCRM@appliedbiosystems.com), telephone (650-554-2756; USA), or fax (650-554-2252; USA) your request, specifying the catalog or part number(s) and the name of the product(s). We will e-mail the associated SDSs unless you request fax or postal delivery. Requests for postal delivery require 1-2 weeks for processing.

**Warranty and Liability:** *For research use only. Not intended for any animal or human therapeutic or diagnostic use.*

Ambion, a part of Life Technologies Corporation, is committed to delivering superior product quality and performance, supported by industry-leading global service and technical support teams. Warranty information for the accompanying consumable product is available at [www.ambion.com/info/warranty](http://www.ambion.com/info/warranty) in "Limited Warranty for Consumables," which is subject to the exclusions, conditions, exceptions, and limitations set forth under the caption "EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS" in the full warranty statement. Please contact Ambion if you have any questions about our warranties or would like information about post-warranty support.

Information in this document is subject to change without notice. Ambion, a part of Life Technologies Corporation, assumes no responsibility for any errors that may appear in this document.

Ambion disclaims all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. In no event shall Ambion be liable, whether in contract, tort, warranty, or under any statute or on any other basis for special, incidental, indirect, punitive, multiple or consequential damages in connection with or arising from this document, including but not limited to the use thereof.

**Trademarks, Patents, and Licensing:**

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. TaqMan is a registered trademark of Roche Molecular Systems.

### Limited Use Label License

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact [outlicensing@lifetech.com](mailto:outlicensing@lifetech.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

This product is licensed under European Patents 1144623, 1214945 and foreign equivalents from Alnylam Pharmaceuticals, Inc., Cambridge, USA and is provided only for use in academic and commercial research including in vitro and in vivo identification and validation of drug targets (but excluding the evaluation or characterization of this product as the potential basis for a siRNA-based drug) and not for any other commercial purposes. Information about licenses for commercial use (including discovery and development of siRNA-based drugs) is available from Alnylam Pharmaceuticals, Inc., 300 Third Street, Cambridge, MA 02142, USA.

This product is covered by U.S. Patent Nos. 6,506,559; 7,056,704 and 7,078,196.

**Purchaser notification:** Unless expressly stated, the purchase of this product conveys no license or grant of any intellectual property rights, whether express, implied, by estoppel or otherwise.

© 2010 Life Technologies Corporation. All rights reserved. 4412542C