

SAIVI™ Alexa Fluor® 680 and Alexa Fluor® 750 Injectable Contrast Agents *human serum transferrin*

Table 1. Contents and storage information.

Material	Amount	Concentration	Storage	Stability
Human serum transferrin conjugates	1 mL, 12.5 nmol per vial	1.25 nmol transferrin/100 µL (recommended volume per injected dose) in azide-free phosphate buffered saline (PBS), pH 7.2; passed through a 0.2 µm sterile filter	<ul style="list-style-type: none"> • 2–6°C • Do not freeze • Protect from light 	When stored as directed, product is stable for at least 3 months.
Approximate Fluorescence Excitation and Emission, in nm : 679/702, for SAIVI™ Alexa Fluor® 680 agent; 749/775, for SAIVI™ Alexa Fluor® 750 agent.				

Introduction

Transferrin is a monomeric serum glycoprotein (~80,000 daltons) that binds to receptors that are overexpressed in many types of rapidly proliferating tumor cells.¹ Transferrin can accumulate intracellularly in tumors through receptor-mediated endocytosis, and may also accumulate in tumor tissue via the enhanced permeability and retention (EPR) effect attributed to pathological alterations of tumor vasculature. The high blood vessel density, increased permeability, and ineffective lymphatic drainage common to some tumors can contribute to this effect;^{1,2} similar accumulation may occur in association with vascular irregularities found in inflammatory processes such as rheumatoid arthritis.³

Molecular Probes provides researchers with human serum transferrin conjugated to near-infrared–fluorescent Alexa Fluor® 680 and Alexa Fluor® 750 dyes (Figure 1) for use as injectable contrast agents in small animal *in vivo* imaging. SAIVI™ injectable contrast agents have been optimized for emission intensity and tested by *in vivo* imaging after injection in disease models established in mice.

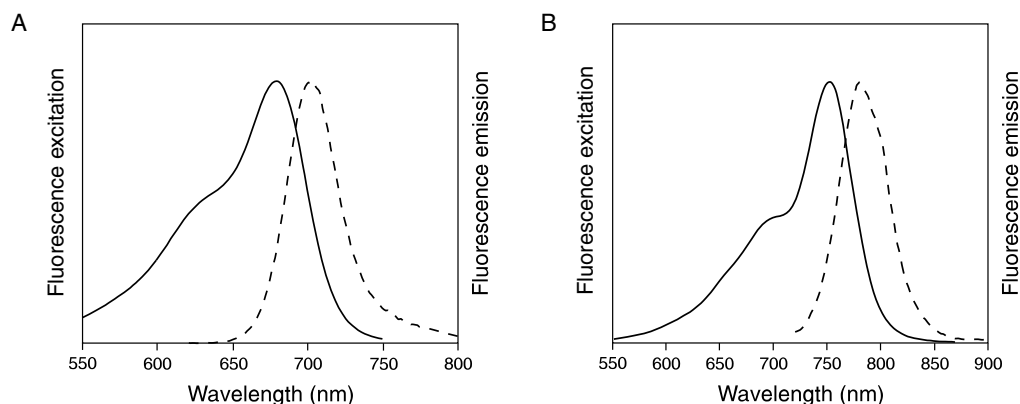


Figure 1. Fluorescence excitation and emission spectra of Alexa Fluor® 680 dye (A) and Alexa Fluor® 750 dye (B).

Guidelines for Use

Using Imaging Agents

Allow the injectable contrast agent to equilibrate to room temperature before use.

The recommended procedure for *in vivo* imaging with SAIVI™ Alexa Fluor® 680 and Alexa Fluor® 750 imaging agents is administration via tail-vein injection and imaging 10 min–24 hr after injection, depending upon the experiment. In our experience, substantial clearance of circulating agent (mainly through the kidneys) will occur over ~24 hr; maximal signal may occur within the first few hours of injection. We recommend imaging frequently in initial experiments to determine the appropriate time course for each type of experiment.

These imaging agents have been used to visualize xenograft subcutaneous tumors and to characterize vascular changes in rodent models of inflammatory disease, using appropriate near-infrared imaging equipment. We have observed evolution of signal to background in xenografted solid tumors over time periods ranging from 1 to 48 hr.

Properties

The human holotransferrin (transferrin saturated with iron) used to prepare these transferrin conjugates is determined to be >98% pure by SDS-polyacrylamide gel electrophoresis.

Warning

These transferrin conjugates contain human transferrin. The venous blood from which the human transferrin is isolated has been tested for the presence of Hepatitis B Surface Antigen (HBsAg) and for HIV (HTLV III) antibody and was found to be negative for both. However, in accordance with good laboratory procedure, these products must be handled as if they are capable of transmitting hepatitis or other infectious agents.

References

1. Cancer Res 46, 6387 (1986);
2. Photochem Photobiol 72, 234 (2000);
3. Arthritis Rheum 50, 961 (2004).

Product List

Current prices may be obtained from our website or from our Customer Service Department.

Cat #	Product Name	Unit Size
S34790	SAIVI™ Alexa Fluor® 680 injectable contrast agent *human serum transferrin*	1 mL
S34791	SAIVI™ Alexa Fluor® 750 injectable contrast agent *human serum transferrin*	1 mL

Contact Information

Molecular Probes, Inc.

29851 Willow Creek Road
Eugene, OR 97402
Phone: (541) 465-8300
Fax: (541) 335-0504

Customer Service:

6:00 am to 4:30 pm (Pacific Time)
Phone: (541) 335-0338
Fax: (541) 335-0305
probesorder@invitrogen.com

Toll-Free Ordering for USA:

Order Phone: (800) 438-2209
Order Fax: (800) 438-0228

Technical Service:

8:00 am to 4:00 pm (Pacific Time)
Phone: (541) 335-0353
Toll-Free (800) 438-2209
Fax: (541) 335-0238
probestech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd.
3 Fountain Drive
Inchinnan Business Park
Paisley PA4 9RF, UK
Phone: +44 (0) 141 814 6100
Fax: +44 (0) 141 814 6260
Email: euroinfo@invitrogen.com
Technical Services: eurotech@invitrogen.com

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Assistance Department in Eugene, Oregon.

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Limited Use Label License No. 223: Labeling and Detection Technology

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, Tel: (541) 465-8300. Fax: (541) 335-0354.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2006, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.