

SAIVI™ Alexa Fluor® 680 and Alexa Fluor® 750 Injectable Contrast Agents *bovine serum albumin*

Table 1. Contents and storage information.

Material	Amount	Concentration	Storage	Stability
Bovine serum albumin conjugates	1 mL, 15 nmol per vial	1.5 nmol albumin/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

Serum albumin, a monomeric protein that constitutes approximately two-thirds of the protein mass of serum, transports an assortment of molecules including fatty acids, bile acids, eicosanoids, vitamins, hormones, ions, toxins, and drugs. Serum albumin is widely used as a macromolecular carrier for targeting tumors with anticancer drugs, photosensitizers, and cytostatic and contrast-enhancing agents.¹ Serum albumin conjugates have also been used for intraoperative sentinel lymph node mapping.² Pooling of this macromolecule in areas of perturbed vasculature may occur by means of the enhanced permeability and retention effect (EPR effect) attributed to pathological alterations of tumor vasculature. High blood vessel density, increased permeability, and ineffective lymphatic drainage can contribute to this effect.³ Such accumulation may occur in association with vascular irregularities found in tumors and in inflammatory processes such as rheumatoid arthritis.^{4,5}

Molecular Probes provides researchers with bovine serum albumin (BSA) 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 as reagents for the *in vivo* imaging of regions of blood pooling and vascularization in an inflammatory disease model in mice.

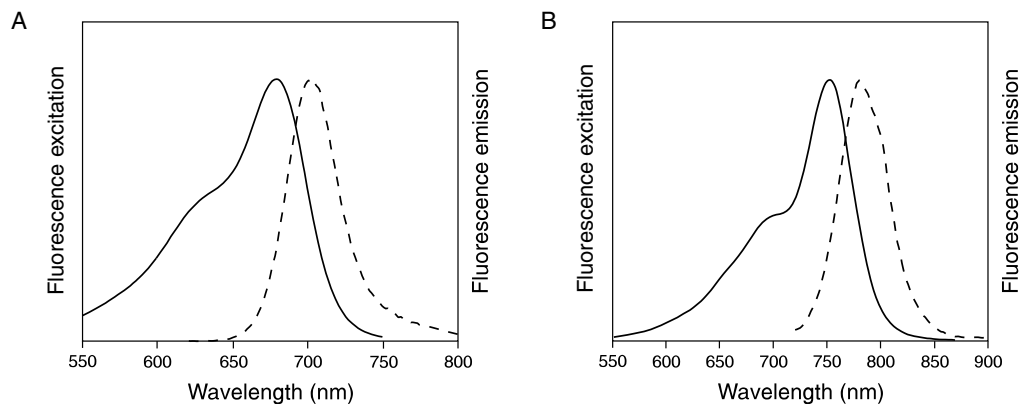


Figure 1. Fluorescence excitation and emission spectra of Alexa Fluor® 680 dye (A) and Alexa Fluor® 750 dye (B).

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 BSA imaging agents is administration via tail-vein injection and imaging 10 min–24 hr after injection, depending upon the experiment. We recommend imaging frequently in initial experiments to determine the appropriate time course for each type of experiment.

These imaging agents have been successfully used to characterize regions of inflammation in a mouse model of collagen-induced arthritis, using appropriate near-infrared imaging equipment. We have observed rapid evolution of signal to background at the site of vascular alterations within the first hour following injection; pooling of circulating agent (mainly in the intestines and bladder) was observed within a few hours and persisted throughout a seven-day imaging period. No signs of acute or long-term toxicity or discomfort were observed in test animals.

Properties

The molecular weight of bovine serum albumin is approximately 66,000 daltons.

References

1. *All About Albumin: Biochemistry, Genetics, and Medical Applications*, Academic Press, New York (1996);
2. *Molecular Imaging* 4, 172 (2005);
3. *Cancer Res* 46, 6387 (1986);
4. *Photochem Photobiol* 72, 234 (2000);
5. *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
S34788	SAIVI™ Alexa Fluor® 680 injectable contrast agent *bovine serum albumin*	1 mL
S34789	SAIVI™ Alexa Fluor® 750 injectable contrast agent *bovine serum albumin*	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.