

FluoSpheres™ Fluorescent Microspheres

Properties and modifications

Pub. No. MAN0001849 Rev. B.0



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

Invitrogen™ FluoSpheres™ Fluorescent Microspheres are manufactured using high-quality, ultraclean polystyrene microspheres. These microspheres are loaded with Invitrogen™ proprietary dyes, resulting in bright fluorescent microspheres. With our special staining methods, all of the fluorescent dye molecules are contained inside each polystyrene microsphere, instead of merely on the bead's surface. The protective environment within the bead shields the dye from many of the environmental effects that cause quenching or photobleaching of exposed fluorophores. The method we employ also ensures the narrow distribution of fluorescence intensity and size. The stability, uniformity, and reproducibility of the FluoSpheres™ microspheres, as well as the extensive selection of colors available, make them the preferred tools for research assays that use fluorescence. In addition, fluorescent microspheres are potentially more sensitive than colorimetric methods in most of the major microsphere-based test systems presently in use, including microsphere-agglutination tests, filter-separation tests, particle-capture ELISA methods, and two-particle sandwich techniques. Precautions are taken throughout the manufacturing process to ensure that the microparticles are kept free of contaminating agents. The final product is sold as a suspension in ultrapure water, in most cases containing 2-mM azide or 0.02% thimerosal as a preservative.

Surface properties

FluoSpheres™ are available with 3 different surface functional groups, making them compatible with many conjugation strategies. Our fluorescent dyes have a negligible effect on surface properties of the polystyrene beads or on their protein adsorption. Anionic FluoSpheres™ with sulfate- or carboxylate-modified surface groups have been used most frequently in biological applications due to their broader availability and tendency to bind less to negatively charged cell surfaces. Cationic microspheres with amine-modified surface groups have been used less frequently, but may have some distinct advantages for certain applications, since they are stable in alkaline pH conditions and high concentrations of multivalent anions such as calcium and magnesium. We caution, however, that the surface properties have an important role in the functional utility of the microspheres; we cannot guarantee the suitability of a particular bead type for all applications.

Carboxylate-modified FluoSpheres™

The carboxylate-modified microsphere products are made by grafting polymers containing carboxylic acid groups to sulfate microspheres. The resulting microsphere has a highly charged, relatively hydrophilic, slightly porous layer that is only a few angstroms thick, which does not change the seed particle size significantly. Carboxylate-modified microspheres will adsorb proteins and other biomolecules, but not as strongly as hydrophobic microspheres. Carboxylate-modified microspheres are often superior for applications in biological systems because they are more highly charged, which reduces their attraction to cells. It is also easier to further reduce nonspecific binding by the introduction of additives such as bovine serum albumin (BSA) or dextrans. A further potential advantage of carboxylate-modified microspheres is that they can be covalently coupled to proteins, nucleic acids, and other biomolecules. Covalent coupling requires more effort than passive adsorption, but can result in conjugates with greater specific activity and products that remain active longer. Covalent coupling to carboxylate-modified microspheres is the method of choice for conjugating low molecular weight peptides and oligonucleotides. Carbodiimide-mediated coupling of proteins to carboxylate-modified microspheres is discussed in more detail in “Covalent coupling of proteins to carboxylate-modified microspheres” on page 4. Their pendent carboxyl groups also make these microspheres suitable for covalent coupling of amine-containing biomolecules using water-soluble carbodiimide reagents such as EDAC.

Amine-modified FluoSpheres™

Amine-modified microsphere products are prepared by further chemical modification of carboxylate-modified microspheres to give hydrophilic particles with positively charged amine groups. The charge density is high, permitting their use in high ionic strength buffers. Amine-modified microspheres contain aliphatic amine surface groups that can be coupled to a wide variety of amine-reactive molecules, including succinimidyl esters and isothiocyanates of haptens and drugs or carboxylic acids of proteins, using a water-soluble carbodiimide. The amine surface groups can also be reacted with SPDP (Cat. No. [S1531](#)) to yield (after reduction) microspheres with sulfhydryl groups.

Sulfate FluoSpheres™

The FluoSpheres™ with sulfate surface groups are relatively hydrophobic particles that will passively adsorb almost any protein, including BSA, IgG, and avidin or streptavidin. The microsphere suspensions are stable at up to about 0.2 M univalent electrolyte concentrations, but will readily agglomerate in the presence of low concentrations of divalent cations unless stabilized by a hydrophilic coating. Sulfate microspheres ($pK_a < 2$) are stable at acidic pH above their approximate pK_a . Even though they have charged surfaces, the hydrophobic microspheres will bind strongly to any molecule that has hydrophobic character, including proteins, nucleic acids, and many small biomolecules such as drugs and hormones. The hydrophobic microsphere products are usually suitable for applications in systems that are free of biologicals and need no further modifications. In biological systems, including immunoassay applications, the microspheres can be easily coated with various proteins or polysaccharides that will greatly reduce their capacity to adsorb biomolecules nonspecifically. Specific, stable adsorption of proteins such as avidin, streptavidin, and antibodies is accomplished simply by mixing the microspheres and the protein together and then separating the microsphere-bound protein from the unbound protein. Refer to “Passive adsorption of proteins to hydrophobic microspheres” on page 3 for details.

Optimization of buffers

The type and density of the microsphere surface charges dictate the best buffer system choice for use in experiments.

- Avoid cationic buffers, such as Tris, when using sulfate- and carboxylate-modified anionic microspheres.
- Avoid borate, citrate, or phosphate buffers when using amine-modified cationic microspheres.

The ionic strength of the buffer should be kept as low as possible, especially when the microspheres are very small or have a low charge density. Since anionic microspheres are very sensitive to low concentrations of multivalent cations, calcium and magnesium salts should be avoided if at all possible. Cationic microspheres are not sensitive to these ions and may be best for applications in which high concentrations of these ions are anticipated. Because of their hydrophobic character, microsphere particles are great scavengers, therefore, the water used for preparation of buffers should be as pure as possible. Either doubly distilled water or high-purity ion-exchanged water is strongly recommended. In general, the smaller the particle size, the more critical are these requirements, since very small microspheres have fewer charge groups for stabilization. The buffer pH can be important when using carboxylate-modified or amine-modified microspheres. The carboxylate-modified microspheres should be used at a pH greater than ~5.0, while the amine-modified microspheres require a pH of less than ~9.0. If these conditions are not followed, the charge groups on these particles may be neutralized, leading to agglomeration. If agglomeration does occur as a result of incorrect pH, the particles can usually be redispersed by adjusting the pH to the correct range followed by gentle sonication.

Controlling nonspecific binding of microspheres

Nonspecific binding is a common problem that is encountered in working with microspheres and is often the major reason for abandoning an otherwise well-conceived experiment with microsphere particles. Microspheres are generally hydrophobic, and although various modifications tend to make them less hydrophobic, these particles are polystyrene-based and therefore always retain some hydrophobic characteristics. In biological systems, most of the nonspecific binding problems are a result of hydrophobic interactions; however, some of the problems may also be caused by charge-based interactions (for example, a positively charged molecule attracted to a negatively charged microsphere surface). The best way to minimize these nonspecific binding events is to coat the microsphere with a large macromolecule such as a protein or a polysaccharide, which reduces nonspecific binding by blocking the hydrophobic or charged binding sites on the microsphere surface. Although many types of coating agents can be used, the most frequently employed are bovine serum albumin (BSA), egg albumin, and whole serum. Egg albumin should be avoided in systems that employ biotin–avidin binding. When using the hydrophobic microspheres, suspending the particles in a 1% solution of the protein-based coating agent is typically sufficient to allow the particles to be completely and stably coated.

Dextrans can be used as coating agents in place of or in addition to proteins. Unlike proteins, the hydrophilic dextrans bind reversibly to microspheres. They form a layer at the surface of the particles and make them more hydrophilic, thus reducing nonspecific interactions. If dextrans are used as coating agents, 40,000 MW dextran at a 2% weight/volume ratio is recommended. If the particle is a hydrophilic carboxylate-modified microsphere, the coating agent may not bind strongly enough to the particles and may fail to prevent nonspecific binding. In this case, covalent coupling of a coating agent such as BSA is recommended. In this method, specific binding proteins, such as immunoglobulin or avidin, can be mixed with BSA and simultaneously coupled covalently, resulting in a specifically active microsphere with a covalently bound BSA coating. As a last resort, or in situations where the use of detergents is acceptable, nonionic surfactants such as polyoxyethylenes (Triton™ X-100 or Tween™ 80) can be coated onto the microsphere at concentrations ranging from 0.01–0.1% (the exact amount to be determined by experimentation).

Our BlockAid™ Blocking Solution (Cat. No. [B10710](#)), a protein-based blocking solution, is designed for use with our streptavidin-, NeutrAvidin™ -, biotin-, and protein A-labeled FluoSpheres™ microspheres. In flow cytometry applications, we find BlockAid™ Blocking Solution to be superior to other commercially available blocking solutions and to a number of “home-made” blocking solutions described in the scientific literature. We expect BlockAid™ Blocking Solution to be useful for preventing the nonspecific binding of protein-coated or other macromolecule-coated microspheres in a variety of flow cytometry and microscopy applications. BlockAid™ Blocking Solution is available in a 50-mL unit size.

Keeping FluoSpheres™ in a monodisperse state

Microsphere particles, which are hydrophobic by nature, will always tend to agglomerate. In aqueous suspensions, the surface charge on the particles prevents agglomeration. Surfactant-free microspheres do not have detergents to aid in dispersion, so these preparations are slightly more sensitive to conditions that can lead to agglomeration. These conditions include: 1) high concentration of particles; 2) high electrolyte concentration; and 3) neutralization of surface charge groups. To minimize these adverse conditions, it is recommended to keep the microsphere suspensions dilute. Recommended particle concentrations to be used when coating microsphere particles with proteins (both passive adsorption and covalent coupling) is 0.5–1.0% solids. Reaction buffers and storage buffers of relatively low ionic strength (100 mM or less) are best. The use of multivalent cations should be especially avoided with anionic microsphere particles. Finally the pH should be maintained so that all of the surface charges on the microsphere particles are fully ionized. If agglomeration does occur, the particles can frequently be rescued by either diluting the microsphere suspension, adjusting the pH, or reducing the ionic strength and then redispersing the suspension by means of a bath sonicator. The use of a bath sonicator greatly aids in working with microsphere suspensions, and it is strongly recommended that this device be utilized if possible. Routine sonication of microsphere preparations is advised before each use, especially in critical applications where a high degree of monodispersity is required. In the case of very small particles (less than 0.1 μm), the sonicated suspension can be briefly centrifuged at high speed (12,000 rpm) to further remove agglomerates from the suspension (the monodisperse particles will remain in suspension under these conditions).

Passive adsorption of proteins to hydrophobic microspheres

Passive adsorption of proteins and other molecules with hydrophobic domains to microspheres is the simplest method of coating, since no chemical reactions are necessary. The coated microsphere product can often be purified from unbound ligand by a simple centrifugation and washing procedure. Passive adsorption should be used only with the hydrophobic microspheres (sulfate surface groups). The hydrophobic interactions that bind macromolecules to the microsphere particles are essentially independent of pH; however, pH and charge can influence the conformation of protein molecules and thus facilitate their binding. Virtually every protein studied to date has been shown to bind to hydrophobic microspheres, and, in general, proteins have been shown to bind most efficiently at a pH that is near their isoelectric point.

The microsphere particle concentration most suitable for adsorption is in the range of 0.5–1% solids. At this relatively low concentration of particles, the aggregation caused by protein bridging is minimized. The ligand to be attached should be added to the dilute suspension of microsphere particles at a concentration of 20–50 μg/mL of final suspension in a buffer with an ionic strength of less than 100 mM. The suspension is stirred, shaken, or rocked gently for a period of a few minutes to a day or more at room temperature. While the physical adsorption is very fast and is complete in just a few seconds, protein-dependent conformational changes can take an hour or two for completion. It should be noted that partial coverage of the microsphere particles with a ligand usually produces a system with greater binding specificity than one where full surface coverage is achieved.

To avoid nonspecific adsorption of further proteins when coated microspheres are used in applications such as a diagnostic test, the microspheres can be back-coated with albumin, gelatin, or other macromolecules (0.5–1 mg/mL). These substances will fill in any remaining hydrophobic areas on the particles. If the coated microsphere particles are used in applications that tolerate detergents, a nonionic surfactant such as Tween™ 20 or Triton™ X-100 can be added to increase hydrophilicity of the particles.

The coated microsphere particles can be separated from unbound ligand by centrifugation and washing if the particle diameter is greater than about 0.2 µm. The smaller the particles, the greater the centrifugation force and time that will be required to cause them to sediment. Some care should be taken to avoid excessive centrifugation force, however; otherwise, the particles may be packed together too tightly and will overcome the repulsive forces between the particles. After the supernatant is carefully removed, the particles are resuspended in washing buffer by vortexing or sonication. In the case of microsphere particles with diameters of less than about 0.2 µm, some type of filtration process will be necessary for separation of unbound ligand. If the ligand is small, ordinary dialysis tubing (12,000–14,000 MW cut-off) can be used; otherwise, cellulose ester dialysis tubing with a MW cut-off of 300,000 daltons can be used for most proteins, including IgG. When dialyzing the particles, the buffer should be the same as that used for the adsorption process, and at least five changes of buffer should be made. Other separation processes such as gel filtration can also be used, although in our experience, the microsphere particles tend to stick nonspecifically to some types of gels.

Covalent coupling of proteins to carboxylate-modified microspheres

There are many procedures published in the literature that describe covalent coupling of proteins and other macromolecules to carboxylate-modified microspheres. Almost all of these use a water-soluble carbodiimide (EDAC) to activate the surface carboxyl groups on the microsphere particles. The following procedure is a simple one-step method we have used with excellent results in our laboratory for coupling avidin, streptavidin, BSA, and goat anti-mouse to our carboxylate-modified FluoSpheres™ microspheres. The reaction can be easily scaled up or down to fit individual needs.

1. Prepare 100 mL of 50 mM phosphate buffer, pH 7.4, containing 0.9% NaCl (50 mM PBS).
2. Prepare 100 mL of 50 mM MES buffer, pH 6.0.
3. Dissolve 10–25 mg of protein (e.g., avidin, streptavidin, IgG, BSA, etc.) at 2–5 mg/mL in MES buffer in a glass centrifuge tube.
4. Add 5 mL of a 2% aqueous suspension of carboxylate-modified microsphere. Incubate at room temperature for 15 minutes.

Note: Agglomeration of the microsphere particles may be observed at this point in the procedure. This agglomeration can be caused by bridging of the particles by protein, neutralization of the charged carboxyl groups, or both. Adjusting the pH to 6.5 and sonication of the mixture in a bath sonicator usually will redisperse the particles. If the particles do not redisperse with these treatments, try a lower concentration of particles and reagents (begin with a 50% reduction in concentration).

5. Add 40 mg of EDAC (1-ethyl-3-(3-dimethylaminopropyl)-carbodiimide; Cat. No. [E2247](#)). Mix by vortexing (see Note in step 4).
6. Adjust the pH to 6.5 ± 0.2 with dilute NaOH. Incubate the reaction mixture on a rocker or orbital shaker for 2 hours at room temperature or overnight, if desired (see Note in step 4).
7. Add glycine to give a final concentration of 100 mM to quench the reaction. Incubate 30 minutes at room temperature.
8. Centrifuge to separate the protein-labeled microsphere particles from unreacted protein. The time and speed of the centrifugation will vary with the diameter of the microsphere particles. As a guideline, 0.5-µm particles and smaller should be centrifuged at $25,000 \times g$ for 30–60 minutes. Particles 1.0 µm and greater can usually be sedimented at $3,000\text{--}5,000 \times g$ for 20 minutes. It may not be possible to centrifuge 20 nm and 40 nm particles without extended spin times. We therefore recommend that dialysis be used with beads of this size.

Note: Particles with a diameter <0.2 µm can be separated from unbound ligand by dialysis using 100,000–300,000 MW cut-off cellulose ester dialysis tubing (Spectra/Por™). We have confirmed that this method effectively removes BSA, avidin, and IgG from microsphere particles with average diameters ranging from 0.03 to 1.0 µm.

9. Resuspend the pellet in 50 mM PBS by gentle vortexing or by use of a bath sonicator. Centrifuge as described in step 8.
10. Repeat step 9 twice for a total of 3 washes.
11. Resuspend the protein-conjugated microspheres in 5 mL of 50 mM PBS. Other buffers compatible with the microspheres (see “Optimization of buffers” on page 2) are also suitable. If desired, the microspheres can be resuspended in a final buffer containing 1% BSA. The BSA will adsorb to the remaining hydrophobic sites on the microspheres and help to provide a more stable suspension that may be less prone to nonspecific interactions with other proteins.
12. Add 2-mM sodium azide and store the microspheres at 4°C. DO NOT FREEZE.

Ordering information

FluoSpheres™ Fluorescent Microspheres are supplied as aqueous suspensions containing 2% solids, except for the 0.04- μm microspheres, which are supplied as aqueous suspensions containing 5% solids. All sizes fall within a narrow range. Sizes indicated are nominal and may vary from batch to batch. Actual sizes, as determined by electron microscopy, are specified on the product labels.

Table 1 Summary of FluoSpheres™ Fluorescent Microspheres

Excitation/emission (nm) ^[1]	0.02 μm	0.04 μm	0.1 μm	0.2 μm	0.5 μm	1.0 μm	2.0 μm	4.0 μm
FluoSpheres™ Carboxylate-Modified Fluorescent Microspheres								
Blue (350/440 nm)	—	—	F8797	—	—	F8815	—	—
Blue (365/415 nm)	F8781	—	—	F8805	—	F8814	F8824	—
Orange (365/610 nm)	—	—	—	F20881	—	—	—	—
Yellow-green (505/515 nm)	F8787 F8888 ^[2]	F10720 ^[3] F8795	F8803 F8888 ^[2]	F8811 F8888 ^[2]	F8813 F8888 ^[2]	F8823 F8888 ^[2]	F8827 F8888 ^[2]	—
Nile red (535/575 nm)	F8784	—	—	—	—	F8819	F8825	—
Orange (540/560 nm)	—	F10720 ^[3] F8792	F8800	F8809	—	F8820	—	—
Red-orange (565/580 nm)	—	F8794	—	—	—	—	—	—
Red (580/605 nm)	F8786	F10720 ^[3] F8793	F8801	F8810	F8812	F8821	F8826	—
Crimson (625/645 nm)	F8782	—	—	F8806	—	F8816	—	—
Dark red (660/680 nm)	F8783	F10720 ^[3] F8789	—	F8807	—	—	—	—
Infrared (715/755 nm)	—	—	F8799	—	—	—	—	—
FluoSpheres™ Sulfate Fluorescent Microspheres								
Yellow-green (505/515 nm)	F8845	—	—	F8848	—	F8852	F8853	F8859
Red (580/605 nm)	—	—	—	—	—	F8851	—	F8858
FluoSpheres™ Amine-Modified Fluorescent Microspheres								
Yellow-green (505/515 nm)	—	—	—	F8764	—	F8765	—	—
Red (580/605 nm)	—	—	—	F8763	—	—	—	—

^[1] Approximate fluorescence excitation and emission in nm are indicated in parentheses.

^[2] The FluoSpheres™ Size Kit includes 1 mL each of the 0.02-, 0.1-, 0.2-, 0.5-, 1.0-, and 2.0- μm sizes.

^[3] The FluoSpheres™ Sample Kit includes 1 mL each of the 505/515-, 540/560-, 580/605-, and 660/680-nm excitation/emission microspheres.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Revision history: Pub. No. MAN0001849 B.0

Revision	Date	Description
B.0	19 December 2022	The microspheres table was updated.
A.0	11 July 2022	The format and content were updated. The version numbering was reset to A.0 in conformance with internal document control.
1.0	2 June 2004	New document for the FluoSpheres™ Fluorescent Microspheres.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.