

Biotech: Diagnostics | Immunodiagnostics

# Protein and cell analysis reagents and instrumentation

For rapid development of diagnostics

A scientist in a white lab coat, safety glasses, and a white face mask is working in a laboratory. The scientist is wearing blue gloves and is focused on a task. The background is a blurred laboratory setting with various pieces of equipment and a large window. The overall tone is professional and scientific.

## Supporting your success in the development of new diagnostics

Your work to develop innovative diagnostics that deliver answers as quickly and precisely as possible can lead to greatly improved health outcomes. Effective tests empower health care providers with the ability to assess risk and identify disease early, monitor the efficacy of treatments, and provide reliable prognoses. We have listened to your feedback, and we know you face a myriad of challenges on the journey from discovery to commercialization.

Looking at your needs, our capabilities to help you navigate these challenges are far superior. When you partner with Thermo Fisher Scientific to help enable development of a new diagnostic, your success is our wholehearted purpose. Take advantage of the industry's broadest range of high-quality and cost-effective solutions; connect with innovative tools and support across your entire workflow; leverage our unique ability to scale solutions to your needs as you grow; and discover the advantage of having our company's unrivaled resources and expertise alongside you at every step, helping you compete and succeed in bringing your diagnostic solutions to market.





IL-6

IFN

TNF- $\alpha$



# Contents

<b>Introduction</b>	<b>7</b>		
<hr/>			
<b>Manufacturing capabilities</b>	<b>11</b>	<b>Linking mechanisms</b>	<b>47</b>
<hr/>			
<b>Capture surfaces</b>	<b>15</b>	Crosslinkers	47
Latex beads	15	PEGylation reagents	54
Coated multiwell plates	18	Biotinylation reagents	56
Custom plate coating service	22	Biotin quantitation kits	60
Biotin-binding proteins	23	Pierce premium grade reagents	61
<hr/>		Fluorescent antibody and protein labeling kits	63
<b>Antibodies and detection probes</b>	<b>25</b>	Fluorescent dyes	65
Primary antibodies	26	Enzyme labeling kits	68
Recombinant primary antibodies	27	<hr/>	
Secondary antibodies	28	<b>Nonspecific binding</b>	<b>71</b>
Recombinant secondary antibodies	29	Blocking buffers	71
Fluorescent and enzyme-conjugated secondary antibodies	29	Wash buffers	75
SARS-CoV-2 antibodies	30	Detergents	76
Antibody affinity purification buffers	31	Protein stabilizers	77
Custom antibody development and production services	32	<hr/>	
ELISA products	35	<b>Detection substrates</b>	<b>79</b>
ELISA kits	35	Chemiluminescent substrates	80
Antibody pair kits	36	Colorimetric substrates	85
Biotin-binding protein conjugates	37	Fluorescent substrates	86
<hr/>		<hr/>	
<b>Protein biology</b>	<b>44</b>	<b>Instruments</b>	<b>89</b>
Gel electrophoresis	44	Attune Flow Cytometers	89
Protein standards and ladders	44	EVOS M7000 Imaging System	94
Mass spectrometry	45	Plate readers and washers	96
Tandem mass tags (TMT)	45	Multiskan SkyHigh Microplate Spectrophotometer	98
<hr/>		Qubit Flex Fluorometer	99
		<hr/>	



# Introduction

## Who understands your challenges? We do.

Immunodiagnostic assay development and commercialization are challenging and time-consuming. Selecting the best combination of tools and technologies to create a high-performing, competitively positioned assay is demanding. We understand this challenge from our work with leading immunodiagnostics companies who face many of the same difficulties you do. Working with a broad array of companies gives us a unique perspective on this immunodiagnostics market, allowing us to see what differentiates the merely good companies from the industry leaders.

We see companies juggling costs, assay development time, and assay quality, which forces them to focus on certain aspects while sacrificing others. This balancing act can result in failure to fully maximize proprietary technology or delays to market, but assay development doesn't have to be this way.

Then there are the other obstacles that you might face. What if you hit a reliability issue with the raw materials used in an assay already commercially available? That could result in the loss of market share and millions of dollars in revenue while the assay is redesigned. Or what happens if someone mistakenly performs an assay based on a raw material (for example, an antibody) that can't be used for diagnostics? You have to start over—and unfortunately, we see this all the time.

And all you want is to create the best assays quickly and cost-effectively. Does it have to be this hard?

### The best solution: partnership

One of the best ways to meet your goals is to partner with an experienced supplier. Think about it. You leverage your proprietary technology and in-house knowledge with the experience of a supplier who has worked with many of the leading immunodiagnostics companies and understands raw materials and new technologies. Using our experience for your benefit can be very powerful. It is the difference we see that separates the most successful companies because it enables them to break free of the cost, quality, and time paradigm. These are the companies who are developing high-quality assays quickly and cost-effectively.

### Why choose Thermo Fisher Scientific?

By choosing us, you gain an ally with a dedicated diagnostics partnering business that is always available for your projects. We have one of the largest ranges of technologies for immunodiagnostics development, with integrated solutions that you can leverage. Our customized solutions enable you to get the best possible assay, and our experienced teams can help troubleshoot when you hit snags in development or post-launch. By partnering early in the development process, we can help you save time and produce an optimal assay.



# How to partner with us

Our Global OEM and Licensing team supports your assay development needs, providing access to our experienced team and providing the rights to commercially use our comprehensive tool kit that includes many different integrated product solutions. These tools will not only help improve the performance of your immunoassay, but can minimize development costs and lot-to-lot assay variation. We offer leading reagent products and technologies with superior levels of performance, together with the high quality and lot-to-lot consistency expected by our customers.

We will provide samples from our broad product portfolio that best meet your requirements. Our immunodiagnostic development portfolio delivers a unique, integrated solution with all of the products necessary for developing an immunoassay, including:

- **Capture surfaces**—latex particles and coated microplates that are the base support for the immunoassay
- **Antibodies and detection probes**—monoclonal, polyclonal, primary, secondary, and recombinant antibodies that bind the analyte of interest, link the analyte to the capture surface for detection; we offer a comprehensive portfolio of over 2,000 ELISA products

- **Linking mechanisms**—crosslinkers, PEGylation, biotinylation, and labeling reagents that connect capture surfaces to antibodies, and antibodies to enzymes and dyes used with detection technologies.
- **Blocking buffers and detergents**—a wide selection to minimize nonspecific binding in the immunoassay, improving assay sensitivity and dynamic range while minimizing false positives.
- **Detection substrates**—enzyme-triggered colorimetric, fluorescent, and chemiluminescent substrates.
- **Instrumentation and equipment**—an unmatched offering of innovative technologies including the Invitrogen™ Attune™ NxT Flow Cytometer, Invitrogen™ EVOS™ M7000 Imaging System, Invitrogen™ Qubit™ Flex Fluorometer, and Gibco™ plate readers and washers

If our off-the-shelf products do not meet your desired specifications, they can be further customized to suit your exact requirements. Our business development manager will determine if a license is required, and can offer special pricing for large-volume orders.

**We are the single supplier to help you accelerate assay development and kit manufacturing.**



### Evaluate workflow needs

Experienced representatives with technology and market experience

### Select products from our portfolio

Extensive collection of technologies and tools for use in immunodiagnostics development

### Customize product design as needed

- Product knowledge
- Process control

### Licensing and supply agreements

>5,000 patents and licenses

### Manufacture products

- cGMP/QSR
- ISO 9001 and/or ISO 13485
- Private labeling
- Custom packaging
- Global distribution







# Manufacturing capabilities

## World-class manufacturing

Many of our products are manufactured in ISO 9001– and/or ISO 13485–certified facilities. Our resources include bioprocess, a HEPA-filtered cleanroom, chemical synthesis laboratories, and QC testing suites to help ensure consistent product performance. Our manufacturing suite is designed to process bulk production batches and customized formulations. For large volume and custom projects, we provide:

### Reagent formulation

- Formulated in a class 100,000–equivalent cleanroom
- Large-scale liquid formulation (up to 5,000 L)
- Hazardous formulation labs
- Custom formulations with in-process testing
- Chromatography media preparation, and column filling
- Buffer dry-blends

### Expert organic chemical synthesis

- Synthetic organic chemistry (milligrams to >300 kg)
- Specialties include synthesis of:
  - Crosslinking and biotinylation reagents
  - GC derivatization reagents
  - Activated supports for chromatography media
  - Enzyme substrates
  - Fluorescent dyes
- Class 1 division 1 production facility with 10–300 gallon Pfaudler reactors
- Chemical purification
- Custom synthesis

## Bioprocessing

- Large-scale fermentation of fine enzymes
- Cytokine standards and quantitation kits
- Protein chemical modification
- Protein immobilization onto solid supports

## Custom services

- Monoclonal and polyclonal antibody production and purification
- Custom hybridoma development
- Custom antibody conjugation to enzymes, biotin, and dyes including Alexa Fluor and DyLight fluorescent dyes
- Custom peptide synthesis
- Automated high-speed 96- and 384-well microplate coating

## Custom packaging and labeling

- Automated high-speed liquid filling/capping/labeling lines
- Aseptic filtration
- Controlled-atmosphere fills for dry materials
- Lyophilization
- Chromatography column filling
- Various packaging options from single-use packages to large-volume carboys
- Hazardous chemical and reagent handling
- Private labeling and kit production

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Quality assurance

Our controlled processes are designed for product quality and consistency, and we provide regulatory documentation and support.

### Analytical testing methods

We manufacture each product lot to high quality specifications using validated methods. Customizable quality specifications can be established for any large volume or custom product, with accompanying certificates of analysis. Our testing methods include:

### Purity

- Quantitative <sup>1</sup>H NMR for accurate purity measurement
- LC-MS
- Infrared spectroscopy
- HPLC
- Melting point analysis
- Solvent panel by GC headspace
- Water content by Karl Fischer titration
- ICP-MS for metal content
- Benzoyl and peroxide levels for detergents

### Consistency

- Appearance
- Particle size analysis
- Solubility
- Concentration and pH

### Activity

- Immunoassay
- Enzyme activity
- Cell-based assays
- Binding capacity

### Regulatory documentation and support

In addition to standard material safety data sheets, we can provide documentation for large-volume and custom products to help meet your regulatory needs. Our logistics team is experienced in global distribution and our quality assurance team can work directly with your regulatory staff to help manage change notifications and site audits.

- Custom certificate of analysis
- Certificate of country origin
- Certificate of animal origin
- OSHA and department of transportation (DOT) hazardous material handling and shipping
- Import and export compliance expertise for chemicals and instruments
- Change control notification
- Site audits welcome





# Capture surfaces

Capture surfaces are the solid support for an immunoassay. We offer a range of capture surfaces including latex beads and surface-coated multiwell plates. We can also work with you to provide custom supports for your application.

## Latex beads

### UltraClean surfactant-free microspheres

We offer a wide selection of Invitrogen™ UltraClean™ surfactant-free microspheres (latex beads) for research and commercial applications (Table 1). Latex beads are colloidal particles typically made of polystyrene (Figure 1). They are available as standard products in many different sizes and surface functionalities or can be tailored to your specifications. Latex beads can be easily modified with proteins such as antibodies or streptavidin via passive adsorption or covalent linkage.

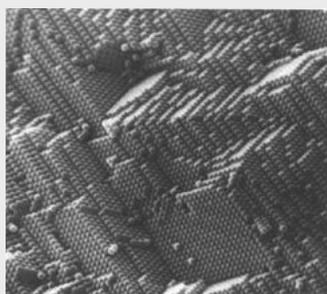


Figure 1. 250 nm polystyrene latex particles.

Table 1. Latex beads applications.

<b>Immunoassays</b>	<ul style="list-style-type: none"><li>• Agglutination tests (lateral flow) (Figure 2)</li><li>• Sandwich assays (ELISA)</li><li>• Particle capture</li><li>• Contrast agents</li></ul>
<b>Flow cytometry</b>	<ul style="list-style-type: none"><li>• Instrument calibration</li><li>• Assay performance</li><li>• Particle capture assays (multiplexing)</li></ul>
<b>Microscopy</b>	<ul style="list-style-type: none"><li>• Instrument calibration</li><li>• Assay performance</li></ul>
<b>Fluid flow</b>	<ul style="list-style-type: none"><li>• Blood flow determination</li><li>• Microfluids</li><li>• Water flow</li><li>• Air flow (flow of airborne particles)</li></ul>
<b>Cell biology</b>	<ul style="list-style-type: none"><li>• Tracing</li><li>• Cell differentiation</li><li>• Cell migration</li></ul>
<b>HTS and HCS</b>	<ul style="list-style-type: none"><li>• Instrument calibration (excitation, emission, focus, etc.)</li></ul>

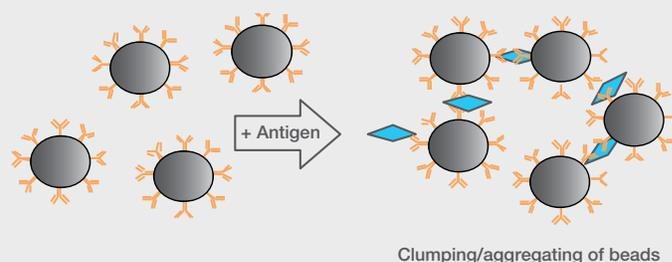


Figure 2. Agglutination assay. Antibody-coated particles assemble via antigens. Detection is accomplished by visualization on a plate (size 0.7–1  $\mu\text{m}$ ), lateral flow assay using colored particles captured in a membrane (0.1–0.3  $\mu\text{m}$ ), or by optical density (turbidometric) measurement (<0.15  $\mu\text{m}$ ).

For more information on our latex beads, go to [thermofisher.com/latexbeads](https://www.thermofisher.com/latexbeads)

## UltraClean high-activity latex beads

Invitrogen™ UltraClean™ high-activity latex beads are hydrophobic and surfactant-free. They are stabilized against aggregation by covalently linked charge groups (sulfate, carboxyl, amidine) and have ~95% of their surface available for passive adsorption of proteins (Table 2).

**Table 2. Latex beads are available in a wide range of surface functional types.**

<b>Physical</b>	Hydrophobic and hydrophilic	
<b>Chemical</b>	<b>Functional group</b>	<b>Type</b>
	-SO <sub>4</sub> <sup>2-</sup>	Strong acid
	-COOH	Weak acid
	-(NH)(NH <sub>2</sub> )	Strong base
	-NH <sub>2</sub>	Weak base
	-CHO	Aldehyde
	-CH <sub>2</sub> Cl	Chloromethyl
<b>Sizes</b>	~20 nm to 15 μm size range, surface-dependent	

### Amidine

Amidinated positively charged hydrophobic latex is particularly suitable for the preparation of latex intermediates. Amidine latex beads should be used in low- to neutral-pH environments and are available in a range of sizes and surface charge densities (70–1,000 Å<sup>2</sup> per charged group).

### Carboxyl

Carboxyl charge-stabilized hydrophobic latex beads are available in a range of sizes and surface charge densities (70 Å<sup>2</sup> per charged group down to 3,000 Å<sup>2</sup> per charged group). These beads may be used either for physical adsorption of antigens or antibodies, or for covalent coupling of components to the particle surface.

### Carboxyl/sulfate

These hydrophobic polystyrene latex beads possess carboxyl and sulfate groups in comparable numbers. As a consequence, the total effective charge is pH-dependent. These beads have been designed for applications in which the reactivity of the carboxyl group combined with the charge-stabilizing characteristics of the sulfate groups is beneficial. They are available in a range of proportions of surface charge groups to one another and particle size.

### Sulfate

These latex beads are stabilized by sulfate charges. Depending upon manufacturing conditions and particle size, the surface charge density of sulfate groups ranges from about one charged group for every 200 Å<sup>2</sup> of particle surface down to one group for every 2,000 Å<sup>2</sup> of surface.

The pK<sub>a</sub> of the sulfate group is <2; consequently, these particles are stable in acidic media and may be used in media of physiological ionic strength. Sulfate latex beads are suitable for calibration of particle size analysis equipment and appropriate for immunoassays that rely upon physical adsorption of antigens or antibodies.

## UltraClean superactive latex beads

Invitrogen™ UltraClean™ superactive latex beads are hydrophilic and contain a very high density of functional groups for covalent coupling of proteins. The superactive layer is a three-dimensional layer, which increases the colloidal stability of the particles and provides a “soft landing” for proteins during interactions with the beads. There is less distortion of the protein structure than if it were physically adsorbed to a rigid surface.

### Carboxylate-modified latex (CML)

Carboxylate-modified latex beads are produced by copolymerizing carboxylic acid-containing polymers. The result is a hydrophilic and somewhat “fluffy” surface layer. The charge density of the CML particles ranges from about 10 Å<sup>2</sup> to 100 Å<sup>2</sup> per charged group.

### Chloromethyl

The chloromethyl latex bead has a high density of chloromethyl groups attached to the styrene monomeric unit. These surface functional groups react directly with amino groups in antibodies, antigens, or other ligands under mild aqueous conditions to yield a stable covalent product by a one-step process. The hydrophobic particles are stabilized by negatively charged sulfate groups. This type of particle can be used at both high- and low-pH conditions.

### Aldehyde/sulfate

These hydrophilic super-active latex beads contain an abundance of aldehyde groups grafted to the surface of the polymer particle. Typical aldehyde density is ~50 Å<sup>2</sup>/group. The high density of aldehyde groups enables facile coupling of proteins and other materials to the latex particles in a one-step process. These particles are ideal candidates for a variety of applications in diagnostic assay development.

### Aldehyde/amidine

These particles are similar to aldehyde/sulfate latex beads but contain a positively charged amidine functional group to provide colloidal stability.

### Aliphatic amine

The aliphatic amine latex bead contains a high density of amine groups and can be used to covalently couple proteins. Location of the amine at the end of the spacer arm minimizes steric hindrance, thereby improving the kinetics of latex agglutination reactions. The particles are stabilized by the positively charged amine groups under low- to neutral-pH conditions. Care should be taken not to use them at high pH.

## Specialty latexes

A variety of specialty latexes and crosslinked particles, including NIST-traceable microspheres, are also available.

For more information on our latex beads, go to [thermofisher.com/latexbeads](https://www.thermofisher.com/latexbeads)

## Coated multiwell plates

We offer a wide selection of high-performance, surface-coated plates (precoated and preblocked polystyrene 96-well and 384-well microplates) in clear, white, and black for use with standard or fluorescence plate readers (Table 3). The choice of plate color depends on the detection chemistry. Clear polystyrene flat-bottom plates are used for colorimetric assays, while black or white opaque plates are used for fluorescence and chemiluminescence applications. Each lot is functionally tested to help ensure minimal variance between wells and between plates. Custom-coated multiwell plates are also available.

**Table 3. Thermo Scientific™ Pierce™ coated polystyrene microplates.**

Microplate coating	Application
Protein A, G, or A/G	For binding antibodies via their Fc regions
Protein L	For binding Fab antibody fragments and single-chain variable fragments (scFvs) through the kappa light chain
Secondary antibodies	For binding antibodies, as an alternative to Protein A, G, or L
NeutrAvidin protein or streptavidin	For binding biotinylated proteins, peptides, or nucleic acids; also available in black or white opaque microplates
Biotin	For binding avidin, streptavidin, or Thermo Scientific™ Pierce™ NeutrAvidin™ biotin-binding protein
Ni <sup>2+</sup> or glutathione	For binding recombinantly expressed proteins containing polyhistidine or glutathione S-transferase (GST)
Maleic anhydride	For binding large or small amine-containing molecules
Maleimide-activated	For binding sulfhydryl-containing molecules
Anti-GST	For capturing proteins expressing glutathione S-transferase



For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## NeutrAvidin and Streptavidin Coated Plates

Thermo Scientific™ Pierce™ Streptavidin and NeutrAvidin™ Coated Plates are preblocked, ready-to-use plates for binding biotinylated antibodies or nucleic acid probes. The plates are available in standard binding capacity, high binding capacity, and high-sensitivity formats (Table 4).

**Table 4. Comparison of Pierce NeutrAvidin and Streptavidin Coated Plates.** Detection ranges were determined using black plates of each product type and Thermo Scientific™ QuantaBlu™ Fluorogenic Peroxidase Substrate Kit.

	High sensitivity (HS)	High binding capacity (HBC)	Standard binding capacity (SBC)
<b>Application</b>	Detect low concentrations of biotinylated molecules	Detect high concentrations of biotinylated molecules	General ELISA screening applications
<b>Biotinylated protein minimum size</b>	>26 kDa	>8 kDa	>8 kDa
<b>Detection range, NeutrAvidin plates</b>	5–125 ng/mL	15–2,500 ng/mL	15–300 ng/mL
<b>Detection range, streptavidin plates</b>	5–300 ng/mL	62–10,000 ng/mL	31–1,250 ng/mL

## Pierce Biotin Coated Plates

Thermo Scientific™ Pierce™ Biotin Coated Plates can be used in immunoassays with NeutrAvidin, streptavidin, avidin, or other biotin-binding proteins. The plates are preblocked to help reduce nonspecific binding.

## Pierce Protein A, G, A/G, and L Coated Plates

Thermo Scientific™ Pierce™ Protein A, G, A/G, and L Coated Plates provide alternatives to direct, passive adsorption methods for immobilizing antibodies for ELISA and other plate-based assay techniques. These plates are uniformly and stably coated with one of four popular immunoglobulin-binding proteins (protein A, protein G, protein A/G, or protein L). They bind to the Fc region of antibodies, promoting optimal orientation for maximum antigen capture (Figure 3). Their consistent coating helps ensure minimal variation.

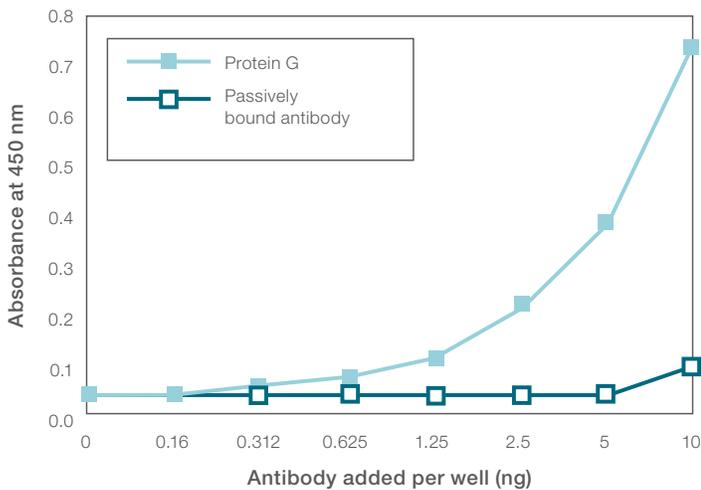


Figure 3. Properly oriented antibodies retain higher activity.

## Pierce Anti-GST Coated Plates

Thermo Scientific™ Pierce™ Anti-GST Coated Plates are polystyrene microplates coated with mouse monoclonal anti-GST antibody and preblocked for immediate use. Unlike glutathione-coated plates, these anti-GST plates effectively bind both native and denatured forms of GST. In most cases, prepurification of cell lysates is not necessary before screening and analysis of recombinant GST-tagged protein expression by ELISA using the plates.

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Pierce Glutathione, Nickel, and Copper Coated Plates

Thermo Scientific™ Pierce™ Glutathione, Nickel, and Copper Coated Plates are used to capture and detect fusion proteins; glutathione-coated plates capture and detect glutathione S-transferase (GST) fusion proteins, whereas nickel and copper bind His-tagged fusion proteins (Figures 4 and 5).

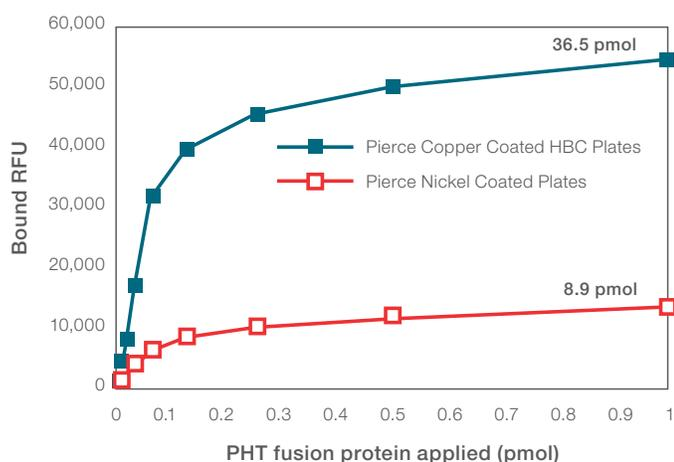


Figure 4. Binding comparison of a histidine-tagged fluorescent fusion protein to standard Pierce Nickel Coated and Copper Coated High Binding Capacity (HBC) Plates. Pierce Copper Coated HBC Plates exhibit a 4-fold greater capacity for binding purified polyhistidine-tagged protein when assayed using a 100 µL volume. Incubation time was 2 hours for binding.

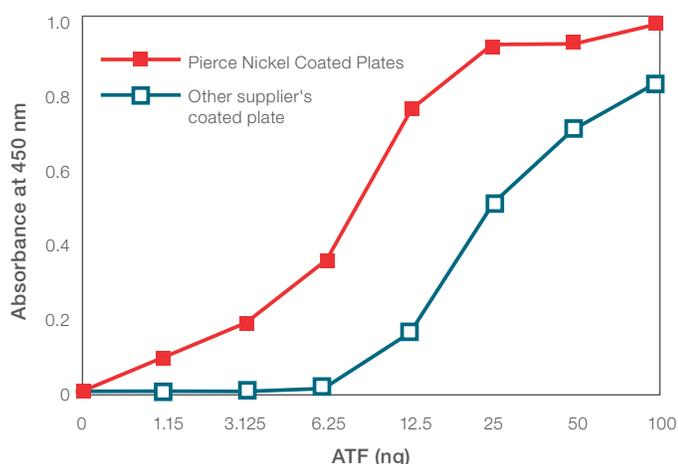


Figure 5. Binding comparison of histidine-tagged ATF fusion protein to Pierce Nickel Coated Plates vs. another supplier's plate.

## Pierce Amine and Sulfhydryl-Binding Plates

Thermo Scientific™ Pierce™ Maleic Anhydride Activated Plates allow covalent attachment of proteins and other primary amine-containing compounds to microplate wells. These plates are useful for immobilizing compounds that do not readily bind to plain polystyrene plates.

Thermo Scientific™ Pierce™ Maleimide Activated Plates are ideal for binding sulfhydryl-containing molecules that are difficult to coat onto polystyrene plates, such as peptides that contain a terminal cysteine. Our coated plates are an especially useful tool for assessing specific anti-hapten antibody titers during antibody production.

## Pierce Antibody Coated Plates

Thermo Scientific™ Pierce™ Antibody Coated Plates are useful for binding assays when available antibodies are in low quantities, denatured, or become inactive upon direct adsorption to polystyrene plates. Because these plates are IgG-specific, purified antibodies are not required. These plates may be used for direct, indirect, competitive, or sandwich assays.

- Anti-mouse plates: binding capacity:  
~7 pmol IgG/well
- Anti-rabbit plates: binding capacity:  
~12 pmol IgG/well

For more information on our coated plates, go to [thermofisher.com/coatedplates](https://thermofisher.com/coatedplates)

## Custom plate coating service

We have developed novel chemistries and coatings for a number of applications, and have the capacity to process up to 5,000 microplates per day. Put our experience to work for you to develop the right combination of ligand, plate, and blocker for your application. We offer customized quality assurance tests to help ensure you receive reliable products that consistently meet your needs. Additionally, we can provide five test plates before we ship any lot of Thermo Scientific™ Pierce™ custom coated plates to you.

**Table 5. Plate coating options.** Choose any plate, ligand, and buffer combination.

Plate type (96- or 384-well)	Coating ligand	Blocking buffer
Clear plates	Antibodies	Thermo Scientific™ SuperBlock™ Blocker
White plates	Peptides	Thermo Scientific™ StartingBlock™ Blocker
Black plates	Fusion proteins	Protein-free blocker
Clear bottom, black plates	Metal chelates	Purified casein
Clear bottom, white plates	Biological polymers	BSA
Filter plates	Your own ligand	Serum
Your own plate		Your own blocker

## Custom microplate packaging options

We can provide the appropriate packaging for your custom plates based on your usage. For example, your plates can be packaged for large screening applications (e.g., 25 plates/pack, ready for stacking) or for inclusion in a kit for resale (e.g., single-pouch packages).

### Partner with us to get the coated plate of your choice

#### Do you need:

- Coated 96- or 384-well plates, slides, or other coated surfaces?
- A coated plate using a certain type of plate or a specific supplier's plate?
- A specific surface chemistry that is not shown here?

#### We can help

Go to [thermofisher.com](https://www.thermofisher.com) to learn about our plate coating service. Outside the US, contact your local distributor or branch office.

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Biotin-binding proteins

Our family of biotin-binding proteins includes streptavidin, avidin, and NeutrAvidin protein (Table 6). Each protein binds four biotins per molecule with high affinity and selectivity. These proteins are ideal for solid phase modification for immunoassays.

- Streptavidin is most commonly used—it is unglycosylated and exhibits low levels of nonspecific binding
- Avidin is a highly cationic glycoprotein with an isoelectric point of about 10—it can cause nonspecific background in some applications due to its positively charged residues and oligosaccharide components
- NeutrAvidin protein has been processed to remove the carbohydrate and lower its isoelectric point, resulting in reduced nonspecific background

**Table 6. Comparison of biotin-binding proteins.**

	Avidin	Streptavidin	NeutrAvidin protein
<b>Molecular mass</b>	67 kDa	53 kDa	60 kDa
<b>Biotin-binding sites</b>	4	4	4
<b>Isoelectric point (pI)</b>	10	6.8–7.5	6.3
<b>Specificity</b>	Low	High	Highest
<b>Affinity for biotin (<math>K_d</math>)</b>	$10^{-15}$ M	$10^{-15}$ M	$10^{-15}$ M
<b>Nonspecific binding</b>	High	Low	Lowest

### Pierce streptavidin protein

Streptavidin is a biotin-binding protein that was originally isolated from *Streptomyces avidinii*. In contrast to avidin, streptavidin has no carbohydrate and has a mildly acidic pI of 5. Pierce streptavidin products use a recombinant form of streptavidin having a mass of 53 kDa and a near-neutral pI. Streptavidin is a tetrameric protein, with each subunit binding one molecule of biotin with affinity similar to that of avidin. Guanidinium chloride will dissociate avidin and streptavidin into subunits, but streptavidin is more resistant to dissociation.

### NeutrAvidin biotin-binding protein

Thermo Scientific™ NeutrAvidin™ biotin-binding protein is a form of avidin that has been processed to remove carbohydrate and to lower its isoelectric point, which can substantially decrease background due to nonspecific binding. The method used to deglycosylate the avidin retains its specific binding.

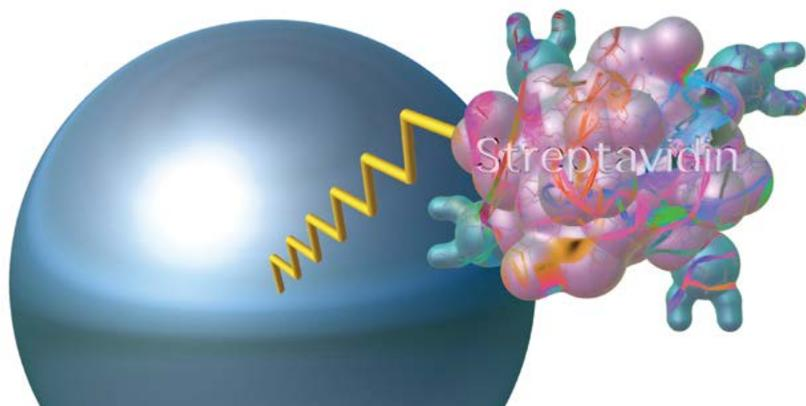
### Pierce avidin protein

Thermo Scientific™ Pierce™ Avidin is a purified avidin protein from chicken egg white that effectively binds biotin, a small molecule that is frequently used to tag antibodies and other probes for immunodetection methods.

#### Highlights:

- **Native avidin**—glycoprotein from chicken egg white, 67 kDa, pI 10
- **Soluble**—glycosylation and high isoelectric point renders avidin more soluble than streptavidin
- **Affordable**—significantly less expensive than recombinant streptavidin
- **ABC staining**—avidin is usually preferred over streptavidin for avidin–biotin complex (ABC) staining methods in immunohistochemistry (IHC)

Avidin is a glycoprotein found in the egg white and tissues of birds, reptiles, and amphibians. This protein contains four identical subunits having a combined molecular mass of 67–68 kDa. Each subunit binds one molecule of biotin, and studies have shown that tryptophan and lysine are involved in the binding site for biotin. The sequencing of the subunit indicates it consists of 128 amino acids. Avidin has an isoelectric point of 10–10.5 and is very soluble in water and salt solutions. Avidin is stable over a wide range of pH values and temperatures. Extensive chemical modification has little effect on the activity of avidin, making it useful for detection and protein purification.





# Antibodies and detection probes

Our extensive line of antibodies includes polyclonal and monoclonal antibodies and Invitrogen™ recombinant rabbit antibodies that can be used to probe, stain, locate, purify, and detect virtually any protein or posttranslational modification. Our offering of antibodies also includes secondary antibodies conjugated to the wide array of fluorescent and enzymatic labels, including Invitrogen™ Alexa Fluor™ and Thermo Scientific™ DyLight™ fluorescent dyes for high-sensitivity detection. We offer a comprehensive portfolio of thousands of ELISA products, ranging from antibody pairs to sensitive and accurate ELISA kits for over 2,000 targets. We also provide enzyme- and fluorescent dye-labeled streptavidin and avidin conjugates.

## Over 50,000 products:

- Monoclonal antibodies—mouse, rat, rabbit
- Polyclonal antibodies—goat, rabbit
- Conjugates for flow cytometry and cellular imaging
- Cleavage site-specific antibodies
- Antibodies to posttranslational modifications, including phosphorylation, acetylation, methylation, and more

## Customization options:

- Alexa Fluor fluorescent dyes
- DyLight fluorescent dyes
- Invitrogen™ fluorescent dyes
  - Pacific Blue™, Pacific Green™, and Pacific Orange™ dyes
  - pHrodo™ Red and pHrodo™ Green dyes
- Common enzymes such as horseradish peroxidase (HRP) and alkaline phosphatase (AP)
- Biotin
- Invitrogen™ Qdot™ nanocrystals
- Traditional fluorophores and labels



## Primary antibodies

We have a diverse collection of thousands of highly specific primary antibodies to targets for use in a variety of assays. These antibodies are validated using a combination of tests, including western blotting, immunocytochemistry, immunohistochemistry, ELISAs, turbidimetric, lateral flow, and flow cytometry assays.

Our portfolio includes antibodies directed against cluster of differentiation (CD) markers, cancer markers, immunology targets, stem cell markers, cell signaling targets, cytokines and chemokines, cell organelles and cell structure targets, cell junctions, and adhesion molecules, as well as other key targets involved in a variety of cellular processes.

### Highlights:

- **Validated**—extensive validation allows for highly specific and consistent antibodies
- **Robust**—they perform in many applications, including immunofluorescence, immunohistochemistry, and flow cytometry
- **Selection**—large menu of unique specificities
- **Quality**—most antibodies are produced under ISO 9001
- **Convenient antibody search tool**—use our antibody search tool at [thermofisher.com/antibodies](https://www.thermofisher.com/antibodies) to find the primary antibody that suits your application

### Applications:

- ELISA
- Immunocytochemistry
- Immunofluorescence
- Immunohistochemistry
- Immunoprecipitation
- Agglutination
- Immunodiffusion
- Competition assay
- Flow cytometry
- ChIP assay
- Cytotoxicity assay
- Electron microscopy
- FACS
- FRET
- Fluorescent quenching
- Functional assay
- Infection
- Gel shift assay
- Hemagglutination assay
- Immunoradiometric assay
- Inhibition assay
- Neutralization
- Radioimmunoassay
- Western blot

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Recombinant primary antibodies

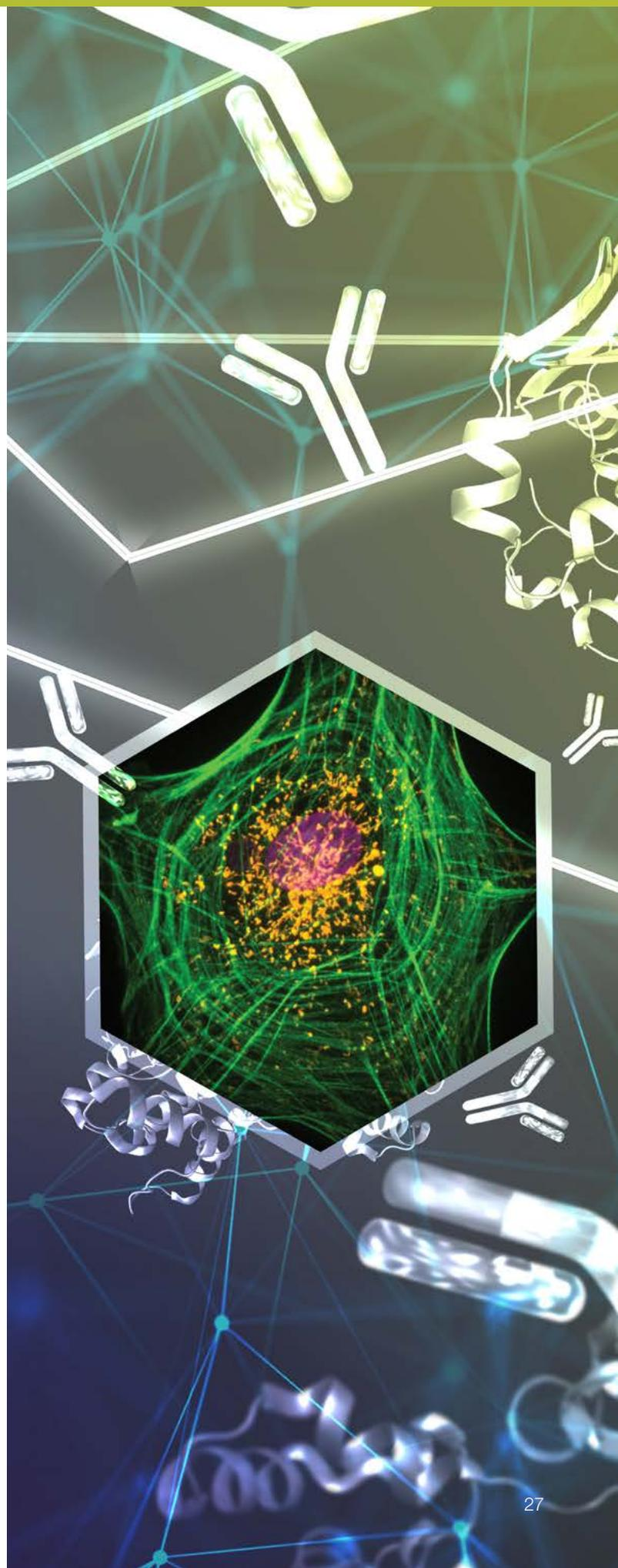
Invitrogen™ recombinant rabbit monoclonal antibodies are developed by immunizing animals, screening for functionality, and then cloning the immunogen-specific antibody genes into high-level mammalian expression vectors. The monoclonal antibodies resemble rabbit monoclonals isolated from serum or produced by hybridomas, but they demonstrate greater specificity and sensitivity.

### Monoclonals: highly consistent performance

Monoclonal antibodies offer peak specificity for antibody tools; however, due to the nature of the production process required, lot-to-lot variation and cell-line drift are both potential issues.

As a consequence, the performance of a traditional monoclonal antibody can change from lot to lot, requiring that you revalidate each lot before committing to valuable samples and time.

Because recombinant monoclonal antibodies are derived from cloned DNA sequences of the heavy and light antibody chains, they are not susceptible to cell-line drift or lot-to-lot variation, thus allowing for peak specificity and performance.



## Secondary antibodies

Secondary antibodies are used for the indirect detection of target antigens. Secondary antibodies can offer significant advantages such as increased sensitivity through the signal amplification and greater flexibility in labeling and detection options.

We have developed an extensive selection of high-quality conjugated and unconjugated secondary antibodies that can be used for fluorescent, colorimetric, and chemiluminescent detection of primary antibodies in a wide range of applications, such as cell imaging and flow cytometry (Table 7). Our Alexa Fluor dye- and DyLight dye-conjugated secondary antibodies provide exceptionally bright and photostable conjugates. All of our secondary antibodies are designed to perform against intended species, across all applications.

A wide selection of highly cited research antibodies is available, many conjugated to a broad range of dyes and enzymes (Table 7):

- Alexa Fluor and DyLight fluorescent dyes
- Classic fluorescent dyes, such as FITC, R-phycoerythrin (RPE), and allophycocyanin (APC)
- Biotin
- Enzyme conjugates, such as HRP and AP

**Table 7. Overview of available secondary antibodies and conjugate options.**

Reactivity	Alexa Fluor dyes	DyLight dyes	Other fluorescent conjugates	Enzyme- and biotin-labeled conjugates	Target Ig class
Anti-chicken	Alexa Fluor 350	DyLight 350	FITC	Enzyme labeled with HRP	IgG
Anti-goat	Alexa Fluor 405	DyLight 405	TRITC	Enzyme labeled with AP	IgM
Anti-mouse	Alexa Fluor 488	DyLight 488	Rhodamine	Biotin	IgA
Anti-human	Alexa Fluor 532	DyLight 550	Texas Red dye	... and more	IgG2a
Anti-rabbit	Alexa Fluor 546	DyLight 594	Texas Red-X dye		IgG1
Anti-rat	Alexa Fluor 568	DyLight 633	RPE		... and more
... and more	Alexa Fluor 680	DyLight 650	APC		
	Alexa Fluor 647	DyLight 680	Qdot probes		
	Alexa Fluor 750	DyLight 755	Pacific dyes		
		DyLight 800	... and more		
		DyLight 800-4xPEG			

For more information on our wide selection of secondary antibodies and conjugates, go to [thermofisher.com/secondary-antibody-selection](https://www.thermofisher.com/secondary-antibody-selection)

## Recombinant secondary antibodies

Invitrogen™ recombinant secondary antibodies represent a breakthrough in recombinant antibody technology designed to provide precise and accurate detection of mouse, rabbit, and goat primary antibodies in a variety of applications. The proprietary screening and production process yields specific mixtures of recombinant goat or rabbit secondary antibodies that bind with the epitope precision of monoclonal antibodies, while also achieving the multi-epitope coverage (e.g., H+L) and sensitivity of polyclonal antibodies. Each of our secondary antibodies is formulated and optimized to help achieve excellent results in ELISA, western blotting, and cell imaging.

### Highlights:

- Designed to eliminate cross-reactivity in detection of primary antibodies (Figure 6)
- Developed as recombinant monoclonal antibodies to enable precise and accurate detection
- Formulated to recognize both heavy- and light-chain epitopes (H+L) of target IgG molecules
- Selected and optimized for use with cell imaging, ELISA, and western blotting applications
- Offered in four types: goat anti-mouse (GAM), goat anti-rabbit (GAR), rabbit anti-mouse (RAM), and rabbit anti-goat (RAG) antibodies
- Available unconjugated and with biotin, HRP, or selected Alexa Fluor dye conjugates

## Fluorescent and enzyme-conjugated secondary antibodies

Fluorescent dye–conjugated secondary antibodies provide a much-needed tool for identifying proteins in many applications including fluorescence cell imaging, western blotting, and immunohistochemistry. The advantages of using a fluorescently labeled secondary antibody include brighter signal, multiplexing capabilities, and ease of use. We offer a wide selection of fluorescent dye–conjugated secondary antibodies for your research, with over 60 different fluorescent dyes and dye combinations, and nanocrystals in a wide range of specific target and host species. In addition, we offer HRP- and AP-labeled antibodies.

### Highlights:

- Available with Alexa Fluor dyes, DyLight dyes, Qdot nanocrystals, biotin, HRP, AP, and more
- Affinity-purified; some antibodies are available in cross-adsorbed formats to minimize cross-reactivity

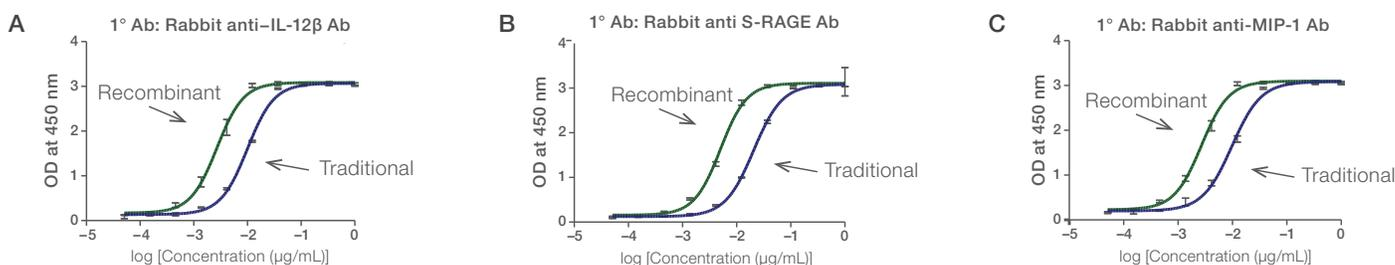
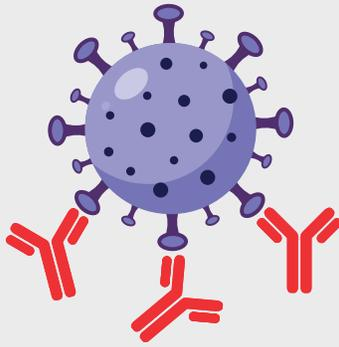


Figure 6. Data demonstrating performance of Invitrogen recombinant goat anti–rabbit IgG (H+L) secondary antibody and traditional polyclonal goat anti–rabbit IgG (H+L) secondary antibody in an ELISA assay. Indirect ELISA to detect recombinant human IL-12β (A), S-RAGE (B), and MIP-1β (C), using the respective primary antibodies and serial dilutions (3-fold starting at 1 µg/mL) of secondary antibody to test performance. An HRP-conjugated rabbit anti-goat (H+L) secondary antibody was used as a detector (1:2,500). A nonlinear regression analysis was performed. EC<sub>50</sub> values for the Invitrogen recombinant antibody were 0.007 (IL-12β), 0.003 (S-RAGE), and 0.005 (MIP-1β) µg/mL, and for the traditional polyclonal antibody were 0.01 (IL-12β), 0.01 (S-RAGE), and 0.02 (MIP-1β) µg/mL.

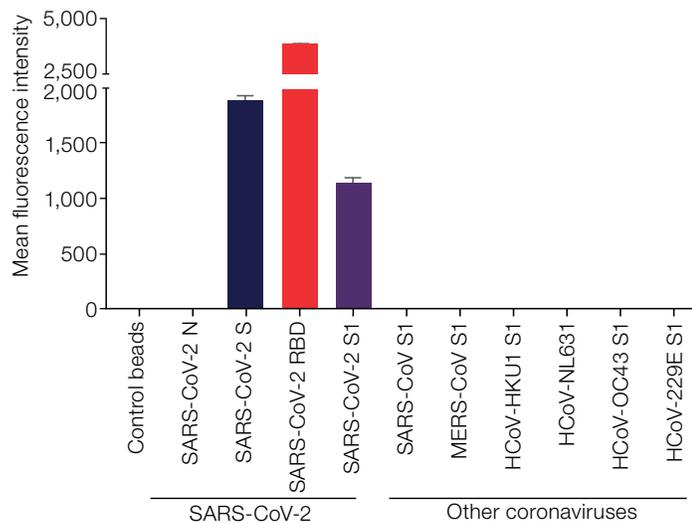
For more information, go to [thermofisher.com/superclonal](https://www.thermofisher.com/superclonal)



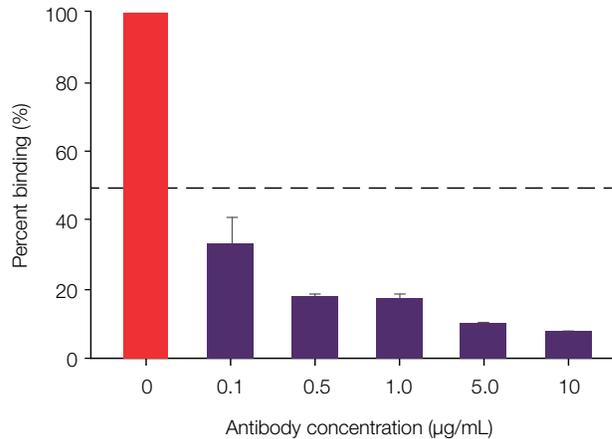
### Introducing antibodies to support SARS-CoV-2 research

Learn more about our new recombinant antibodies to SARS-CoV-2, which are specific and neutralizing

These antibodies were developed using complementarity-determining regions (CDRs) computationally derived from published antibody–virus interactions, cloned into our proprietary rabbit and human antibody backbones, and expressed in our animal-free production system.



**Figure 7. Specific for the SARS-CoV-2 spike protein receptor-binding domain (RBD)** No interaction with SARS-CoV, responsible for the 2002–2004 health crisis, or other coronaviruses. Specifically binds to the RBD of the S1 subunit of the spike protein of SARS-CoV-2



**Figure 8. Blocks binding to the ACE2 receptor.** Blocks the binding of the SARS-CoV-2 spike protein RBD to human ACE2 in ELISA-based assays.

Learn more about our entire SARS-CoV-2 antibody portfolio at [thermofisher.com/sarscov2abs](https://thermofisher.com/sarscov2abs)

## Antibody affinity purification buffers

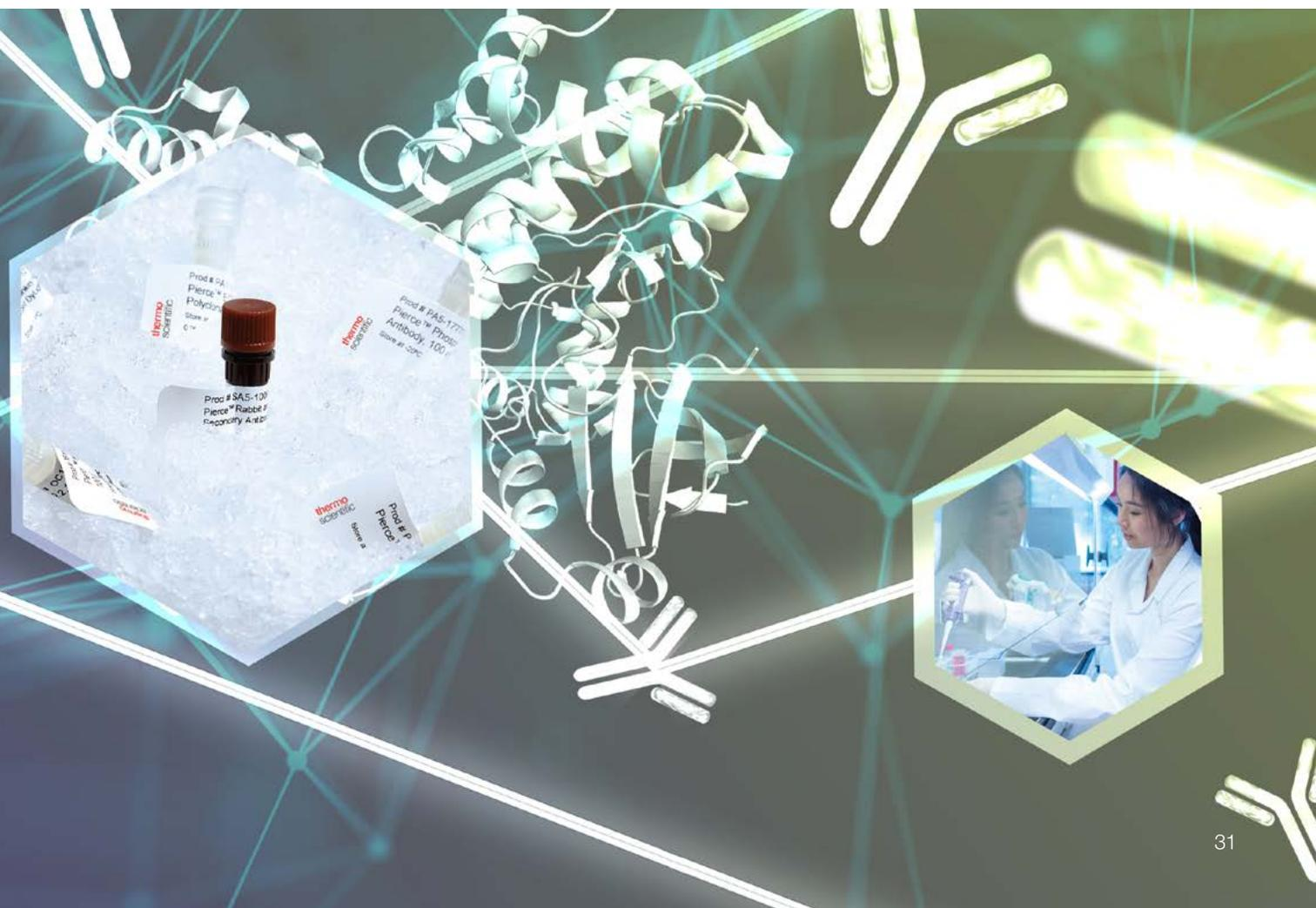
Thermo Scientific™ Pierce™ Binding and Elution Buffers ensure high-yield, nondenaturing antibody purification with IgG affinity purification supports such as Protein A and Protein G agarose beads.

### Highlights:

- **Ready to use**—carefully formulated and prefiltered 1X solutions
- **Three binding buffers**—optimized for Protein A, Protein G, or Protein A/G
- **Two elution buffer options**—acidic (pH 2.0) or high-salt (near-neutral)
- **Versatile**—elution buffers are appropriate for antibody–antigen affinity purification

Products available for each antibody purification method:

- Protein A IgG Binding Buffer: pH 8.0; phosphate-based; contains glycerol and EDTA
- Protein A/G IgG Binding Buffer: pH 8.0; phosphate-based; contains EDTA
- Protein G IgG Binding Buffer: pH 5.0; contains 0.02% sodium azide
- IgG Elution Buffer: pH 2.0; amine-based
- Gentle Ag/Ab Binding Buffer: pH 8.0; phosphate-free; contains 0.05% sodium azide
- Gentle Ag/Ab Elution Buffer: pH 6.6; high salt (ionic strength)



## Custom antibody development and production services

### Full-service, customized production of monoclonal and polyclonal antibodies

Our custom antibody development service leverages our experience in making more than 18,500 antibodies to peptides and recombinant proteins (Table 8). Our proprietary antigen design tools, including Thermo Scientific™ Antigen Profiler software, produce robust antibodies that perform well in your targeted assays.

We specialize in antigen design and generation of custom peptide antibodies and monospecific peptide antibodies to highly discrete epitopes. Our detailed knowledge of antigen-determining factors allows us to produce custom antibodies with superior specificity, affinity, and assay utility.

#### Highlights:

- **Peptide design, synthesis, and conjugation**—we will help you design the best antigenic peptide using our exclusive Antigen Profiler system; then we'll synthesize the peptide for you and conjugate it to the carrier protein of your choice
- **Fusion protein expression and purification**—we offer bacterial, mammalian, and insect protein expression services as part of our antibody development platform; we'll make your protein and purify it to prepare it for immunization and ELISA testing
- **Polyclonal or monoclonal antibody protocols**—choose one of six species options (eight immunization schedules) for polyclonal antibody production, or one of three package sizes (number and strains of mice) for monoclonal antibody production
- **Recombinant rabbit monoclonal antibody production services**—develop recombinant rabbit monoclonal antibodies to your target; screening can either be completed by our development scientists or in your labs
- **Biomarker antibody development**—choose from several standardized packages for producing polyclonal or monoclonal antibodies to specific biomarkers of interest
- **Screening and titering analysis**—we screen and characterize the antisera or hybridoma supernatants produced
- **Purification specificities and deliverables**—we purify all antibodies; in addition, we offer many specialized purification options for obtaining monospecific antibodies, such as against phosphorylated, acetylated, or other posttranslational modification states

For more information on our custom antibodies, go to [thermofisher.com/customabs](https://thermofisher.com/customabs)

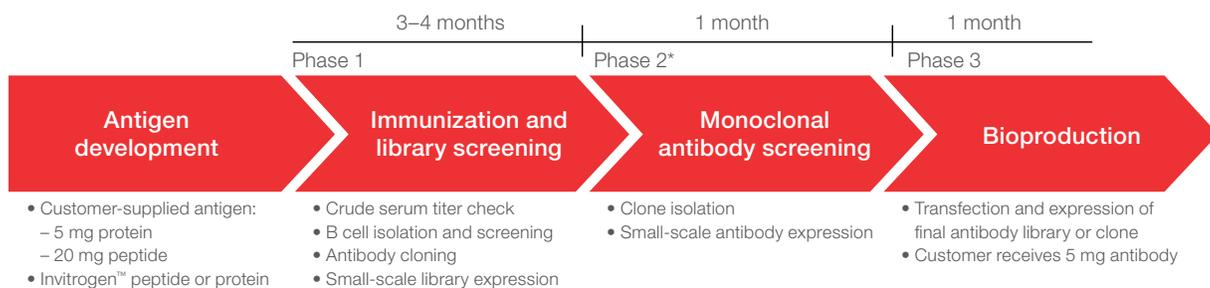
**Table 8. Summary of custom polyclonal and monoclonal antibody development and production service capabilities and options.**

Polyclonal antibody production				Monoclonal antibody production			
	Recombinant protein antibody	Standard anti-peptide antibody	Modification-specific antibody*	Modification-specific antibody*	Standard anti-peptide antibody	Recombinant protein antibody	
Antigen design with Antigen Profiler	✓	✓	✓	✓	✓	✓	Antigen design with Antigen Profiler
Peptide antigen synthesis	NA	✓	✓	✓	✓	NA	Peptide antigen synthesis
Depleting peptide synthesis	NA	NA	✓	✓	NA	NA	Control peptide synthesis
Protein expression (vs. you supply)	Optional	NA	NA	NA	NA	Optional	Protein expression (vs. you supply)
Animal immunization (8 protocol options)	✓	✓	✓	✓	✓	✓	Animal immunization (3 protocol options)
Bleeds, serum prep, ELISA screening, and titration	✓	✓	✓	✓	✓	✓	Hybridoma fusion, cloning, screening, and ELISA titration
Affinity purification and depletion (several options)	Optional	Optional	✓	Optional	Optional	Optional	Culture, production, and purification (several options)
Specific assay validation or antibody labeling	Optional	Optional	Optional	Optional	Optional	Optional	Specific assay validation or antibody labeling

\* Modification-specific antibodies are phosphospecific antibodies or monospecific antibodies against peptides with other posttranslational modifications or specific states (phosphorylation, sumoylation, myristoylation, acetylation, polymorphisms, drug binding, ubiquitination, glycosylation, isoforms, splice variants, ligand binding, prenylation, protein cleavage neoepitopes, mutations, species cross-reactivity).

## Recombinant rabbit monoclonal antibody development and production services

Recombinant antibodies are rabbit monoclonal antibodies that are developed by immunizing animals, screening for functionality, and then cloning the immunogen-specific antibody genes into high-level mammalian expression vectors.



\* Phase 2 is only performed in the monoclonal protocol. During development, the protocol includes phases 1 and 3 only.

The recombinant antibodies are then produced by large-scale cell culture and purified with protein A.

Choose the Invitrogen™ custom antibody service to obtain antibodies with the following features:

- Greater specificity and sensitivity compared to standard antibodies
- Lot-to-lot consistency due to recombinant technology
- Animal origin-free antibodies, expressed in a mammalian cell system

### Custom conjugation of antibodies or proteins

We offer different fluorescent labels for primary antibodies, secondary antibodies, anti-dye and anti-hapten antibodies, and streptavidin, including proprietary labels such as Alexa Fluor and DyLight dyes, Qdot nanocrystals, Invitrogen™ Texas Red™-X, and Invitrogen™ Pacific Blue™ dyes, which span the visible light spectrum from deep blue to near-infrared emission, generic fluorophores such as fluorescein and tetramethylrhodamine, RPE, APC, and peridinin chlorophyll protein complex (PerCP). We also offer conjugates of biotin, 2,4-dinitrophenyl (DNP), digoxigenin, dansyl, and other haptens.

We have more than 30 years of experience in preparing custom antibody conjugates that are optimized to meet your application-specific assay needs and match your instrument and optical specifications. The type of conjugation chemistry and linkers used can affect the functional outcome of your antibody or protein. Our expertise and our broad range of technologies enable you to get superior results. We can couple antibodies in a site-specific way and to labels like gold particles, polystyrene microspheres and nanospheres, and magnetic particles.

We are also highly experienced with conjugating antibodies and streptavidin to enzymes including AP, HRP, beta-galactosidase, and others.



Our custom conjugation service is efficient and confidential, and we stand behind the quality of our work. We can conjugate your antibody or protein to many labels, including:

- Alexa Fluor dyes
- DyLight dyes
- Qdot nanocrystals
- Pacific Blue, Pacific Green, and Pacific Orange dyes
- Traditional fluorophores such as FITC, TRITC, and Texas Red dyes
- Biotin
- RPE, APC, and Alexa Fluor tandem dyes
- HRP, AP, and other enzymes

Additional testing is available upon request and includes:

- Activity validation/bioassay testing (Biacore or other applications such as IF, FACS, etc.)
- Purity testing and/or mass spectrometry (MS) analysis
- Endotoxin testing and removal
- Special purification

\* See Terms and Conditions at [thermofisher.com/performanceguaranteed](https://www.thermofisher.com/performanceguaranteed)

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## ELISA products

### Over 2,000 ELISA products, ranging from antibody pairs to ready-to-use kits

The enzyme-linked immunosorbent assay (ELISA) is a benchmark for the quantitation of proteins, using a solid-phase enzyme immunoassay (EIA) to detect and measure protein targets in various sample types. ELISAs are designed to provide rapid and consistent results that are relatively easy to analyze. We offer a comprehensive portfolio for research areas including immunology, inflammation, neurobiology, and cancer.

### ELISA kits

Our ELISA kits help provide accurate, sensitive, and consistent quantitative results. Each target protein is tested in biologically relevant models and are calibrated to NIBSC (National Institute for Biological Standards and Control) standards, if available. In addition, kits are validated using common sample types, including serum, plasma, and cell culture supernatant. Cell lysates are used to validate kits that detect signaling proteins or phosphorylation.

These ELISA kits must meet rigorous quality control specifications and are manufactured in an ISO facility to help ensure excellent quality and reproducibility.

### Advantages of our ELISA kits:

- Broad menu of over 800 targets
- Optimized for sensitive, accurate, and consistent performance
- Thorough instructions to complete protocol in 2.5 to 4 hours (varies by kit)
- Validated for typical sample types (e.g., serum, plasma, supernatant, lysates)

### Protein targets:

- Cytokines
- Chemokines
- Interleukins
- Inflammation targets
- Signaling proteins
- Receptors
- Neurobiology markers
- Phosphorylated proteins
- Growth factors
- Adhesion molecules

### Ready-to-use ELISA kits typically include:

- Antibody-coated 96-well plate
- Standards
- Primary detection antibody (typically biotinylated)
- Secondary detection reagent (usually streptavidin-HRP)
- Diluent buffers
- Wash buffers
- Substrate and stop solutions
- Plate covers

For more information on our ELISA kits, go to [thermofisher.com/elisa](https://www.thermofisher.com/elisa)

## Antibody pair kits

Antibody pair kits contain matched, pretitered, and fully optimized capture (coating) and detection antibodies. These kits enable you to build your own ELISA or any other assay platform that utilizes a matched antibody pair.

These matched pair kits are designed to accurately quantify cytokines, chemokines, growth factors, signaling pathway targets, and proteins associated with immunology, inflammation, cancer, cardiovascular, and neurodegenerative disease research.

### Advantages of the antibody pair kits for ELISA:

- **Quality**—reliable antibodies, prematched antibody pairs and proven detection reagents
- **Ease of use**—simplified protocol and optimized reagents
- **Flexibility**—multiple detection technologies (fluorescence, absorbance, or chemiluminescence) possible
- **Cost savings**—more economical than complete, ready-to-use ELISA kits that contain precoated plates

### Build and customize your own immunoassays

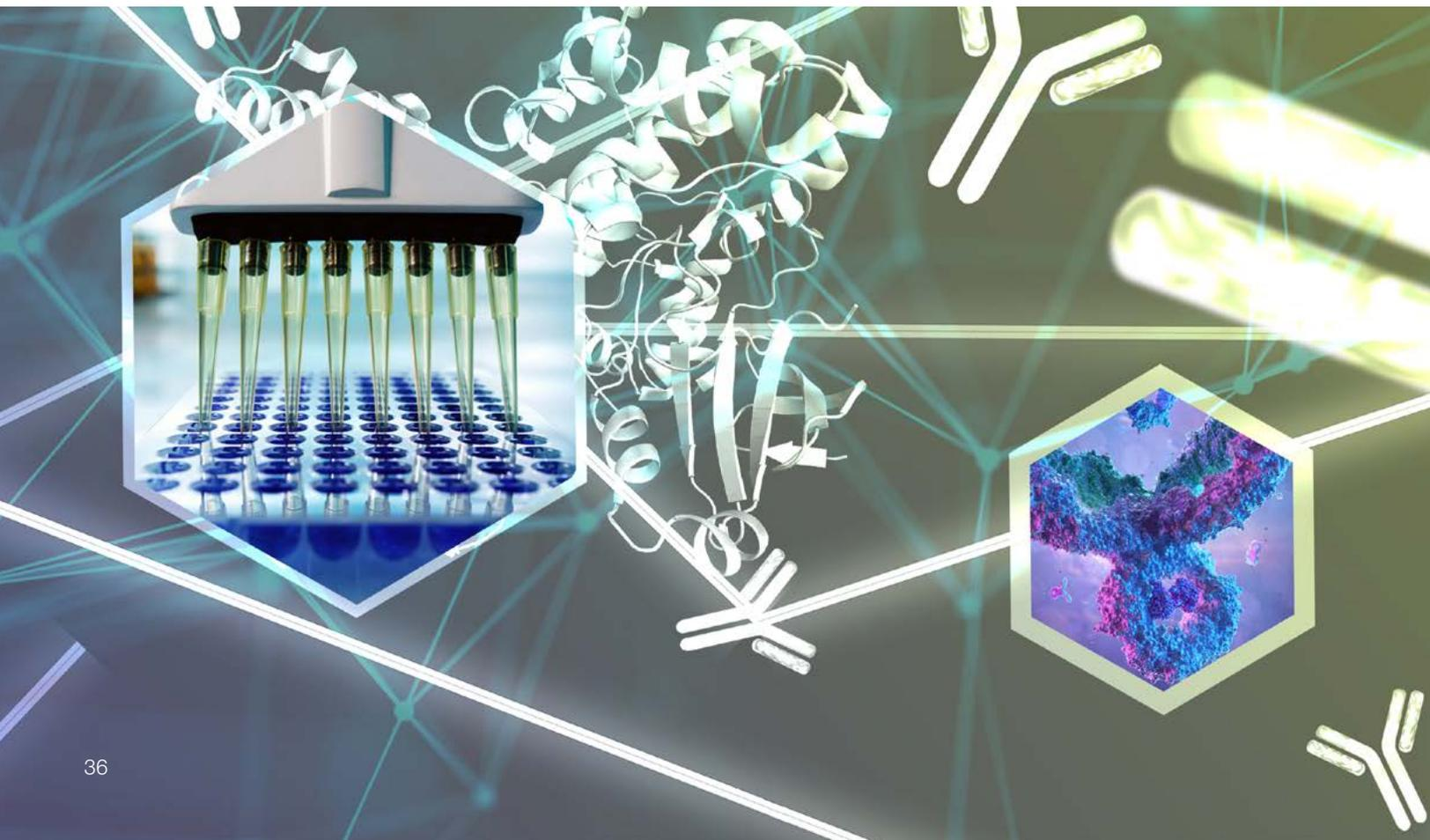
The kits are designed for use with a variety of sample types such as serum, plasma, cell culture supernatant, cell lysate, tissue homogenate, urine, and cerebrospinal fluid (CSF).

For convenience, we also offer a buffer set for antibody pairs that contains premade, easy-to-use buffers and solutions that are optimized for use with the antibody pair kits.

Each antibody pair kit supplies sufficient reagents for 10 ELISA plates (5 plates for intracellular targets) and includes:

- Capture antibody
- Detector antibody
- Recombinant standard
- HRP conjugates

For more information on our antibody pair kits, go to [thermofisher.com/antibodypairs](https://www.thermofisher.com/antibodypairs)



## Biotin-binding protein conjugates

### Enzyme-labeled streptavidin and avidin conjugates

#### Pierce High Sensitivity Streptavidin-HRP

The Thermo Scientific™ Pierce™ High Sensitivity Streptavidin-HRP conjugate is an exclusive peroxidase-conjugated biotin-binding protein that provides signal amplification and exceptional storage stability.

#### Highlights:

- **High sensitivity**—detect low levels of target without background; obtain high signal-to-noise ratios (Figure 9)
- **Cost-effective**—use less conjugate in western blotting and ELISA applications and still obtain excellent results
- **Flexible**—compatible with typical chemiluminescent, fluorescent, and colorimetric peroxidase substrates
- **Convenient**—ready-to-use stabilized liquid format means there is no waiting for thawing and no need to aliquot

This specially manufactured variety of HRP-conjugated streptavidin is designed to meet the needs of today's scientists for more sensitive detection in immunoassay applications. The conjugate is suitable for use with chemiluminescent, chemifluorescent, or colorimetric substrates. Each High Sensitivity HRP conjugate is packaged in an easy-to-use stabilized solution for convenient storage at 4°C for at least 1 year.

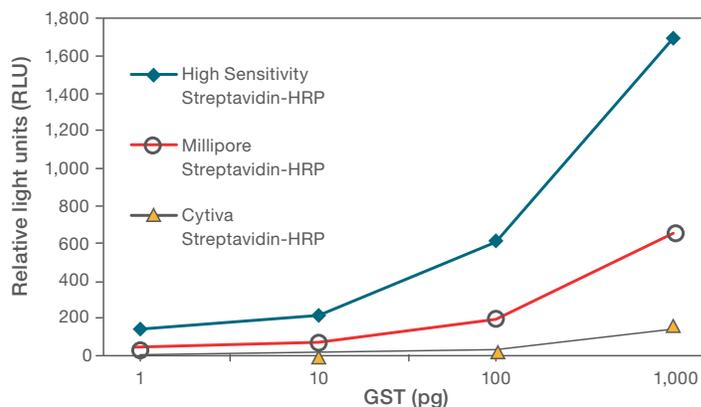


Figure 9. Pierce High Sensitivity Streptavidin-HRP conjugate enables low-level target detection with high signal-to-noise ratios. Recombinant Thermo Scientific™ Pierce™ GST (glutathione S-transferase) was serially diluted (0–10,000 pg/mL) with Thermo Scientific™ StartingBlock™ (PBS) Blocking Buffer. Each dilution (100  $\mu$ L) was added to a 96-well Thermo Scientific™ Pierce™ glutathione-coated plate in four replicates, including a negative control. The plate was incubated for 60 minutes at room temperature and washed three times with Pierce™ PBS Tween™ 20 detergent. Biotinylated anti-GST (100  $\mu$ L at 250 ng/mL; Santa Cruz) was added to all wells. The plate was incubated for 30–60 minutes at room temperature and washed three times. Streptavidin-HRP conjugates were diluted in blocking buffer as per manufacturer's directions. Pierce High Sensitivity Streptavidin-HRP conjugate was diluted 1:10,000, and conjugates from other suppliers were diluted 1:1,000. The conjugate solutions (100  $\mu$ L) were added to the plate and incubated for 60 minutes (light protected) at room temperature. The plate was washed five times and 150  $\mu$ L/well of Thermo Scientific™ SuperSignal™ ELISA Pico Substrate was added. Signal intensity was measured using a luminometer.

### Pierce Streptavidin Poly-HRP conjugate

The Thermo Scientific™ Pierce™ Streptavidin Poly-HRP conjugate is a biotin-binding protein conjugated with polymers of HRP, enabling signal amplification and detection of biotinylated antibodies for IHC and other methods. Pierce Streptavidin Poly-HRP conjugate is designed to deliver the highest sensitivity and low background in immunoassays where sample volume is limited or when the target molecule is present at low levels. Streptavidin Poly-HRP conjugate is purified to remove unconjugated streptavidin molecules that reduce signal intensity by competing for binding sites with HRP-conjugated molecules. In addition, the conjugate is devoid of HRP monomers that can cause background signal.

#### Highlights:

- **High sensitivity**—detect low-abundance targets (low picogram to femtogram range) with high signal-to-noise ratios
- **Robust**—consistent manufacturing and purification to minimize low and unconjugated molecules for minimal background and high sensitivity
- **Flexible**—compatible with chromogenic, fluorogenic, and chemiluminescent substrates
- **Versatile**—compatible with ELISA, western blotting, and IHC
- **Easy to use**—can be directly substituted into immunoassays and other detection assays
- **Convenient**—ready-to-use, stabilized liquid format stored at 4°C
- **Cost-effective**—requires less conjugate per assay than standard HRP conjugates

### Pierce Streptavidin, HRP conjugate

The Thermo Scientific™ Pierce™ HRP-conjugated streptavidin (Cat. No. 21124) includes streptavidin in a purified form, conjugated to peroxidase for substrate-based detection.

The Pierce streptavidin HRP conjugate enables detection of biotinylated antibodies and other probes in a variety of standard assay methods, including western blotting, ELISA, IHC, and fluorescence imaging. The conjugate is supplied as a lyophilized powder in phosphate-based buffers for immediate reconstitution with water.

### Streptavidin, alkaline phosphatase conjugate

The Invitrogen™ streptavidin alkaline phosphatase conjugate can be used to detect biotin in a signal amplification scheme in conjunction with chromogenic or fluorogenic substrates.

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

### Pierce High Sensitivity NeutrAvidin HRP conjugate

The Thermo Scientific™ Pierce™ High Sensitivity NeutrAvidin™ HRP conjugate is a specially prepared peroxidase-conjugated form of the biotin-binding protein avidin that provides signal amplification similar to poly-HRP conjugate and exceptional storage stability. This specially manufactured variety of HRP-conjugated avidin is designed to meet the needs of today's scientists for more sensitive detection in ELISA, western blotting, and IHC applications (Figure 10). The conjugate is suitable for use with chemiluminescent, chemifluorescent, or colorimetric substrates. Each High Sensitivity HRP conjugate is packaged in an easy-to-use, stabilized solution for convenient storage at 4°C for at least 1 year.

NeutrAvidin protein is deglycosylated native avidin from egg white. Removal of the excess carbohydrate by a proprietary process yields a protein with a more neutral isoelectric point and less nonspecific binding. NeutrAvidin protein exhibits high assay specificity and sensitivity with high signal-to-noise ratios. Purified and conjugated forms of NeutrAvidin protein provide exceptional performance in western blot, ELISA, and IHC applications that require biotin-binding probes.

#### Highlights:

- **NeutrAvidin protein**—a specially deglycosylated form of avidin that provides highly specific, low-background binding of biotinylated antibodies in many applications
- **High sensitivity**—detect low levels of target without background; obtain high signal-to-noise ratios (Figure 10)
- **Cost-effective**—use less conjugate in western blotting and ELISA applications and still obtain excellent results
- **Flexible**—compatible with typical chemiluminescent, fluorescent, and colorimetric peroxidase substrates
- **Convenient**—ready-to-use, stabilized liquid format, which means there is no waiting for thawing and no need to aliquot

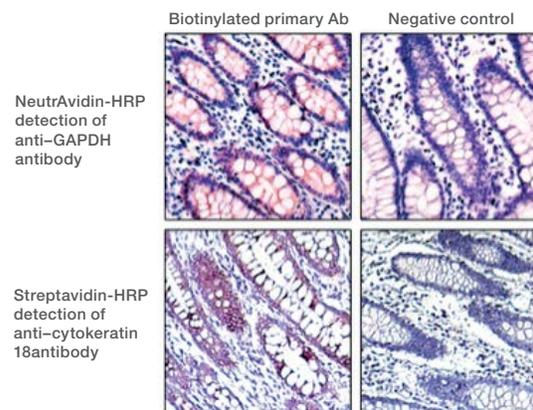


Figure 10. Excellent IHC staining of GAPDH and cytokeatin 18 in human colon carcinoma tissues with Pierce High Sensitivity HRP conjugates. Formalin-fixed, paraffin-embedded (FFPE) human colon carcinoma tissues were stained using a metal-enhanced DAB substrate (Thermo Scientific™ Pierce™ Immunohistochemistry Peroxidase Detection Kit). The tissues were incubated with either biotinylated anti-GAPDH or rabbit anti-cytokeatin 18 antibody followed by biotinylated anti-rabbit IgG (left panels) or blocking buffer only (right panels). The sections were subsequently incubated with Pierce High Sensitivity NeutrAvidin HRP (top panels) or Streptavidin HRP conjugate (bottom panels). Tissues were counterstained using the Harris-modified hematoxylin solution (blue staining in all panels). GAPDH and cytokeatin 18, stained using Pierce NeutrAvidin HRP and Streptavidin HRP conjugates, appear brown (left panels) while the negative control panels shows no brown staining in the tissue.

### Pierce NeutrAvidin HRP conjugate

Thermo Scientific™ Pierce™ NeutrAvidin™ HRP conjugate is a specially prepared form of the biotin-binding protein avidin, conjugated to peroxidase, that decreases background in western blotting and ELISA applications.

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

### NeutrAvidin Alkaline Phosphatase Conjugate

Thermo Scientific™ alkaline phosphatase–conjugated NeutrAvidin protein is a specially prepared form of avidin that decreases background in biotin binding. Conjugated NeutrAvidin protein offers exceptional performance in western blot, ELISA, and IHC applications that require biotin-binding probes. NeutrAvidin protein exhibits high assay specificity and sensitivity with high signal-to-noise ratios.

#### Highlights:

- **Near-neutral isoelectric point**—pI 6.3, more neutral than native avidin
- **Nearly devoid of glycosylation**—decreased possibility of lectin binding compared to native avidin
- **No RYD recognition sequence**—no known off-target binding domains like those contained in streptavidin
- **Affordable**—significantly less expensive than streptavidin

### Pierce Horseradish Peroxidase Conjugated Avidin

Thermo Scientific™ Pierce™ Horseradish Peroxidase Conjugated Avidin is used in biotin-based detection methods, including western blotting or ELISA detection with appropriate enzyme substrates. Avidin from chicken egg white is a glycoprotein that effectively binds biotin, which is frequently used to tag antibodies and other probes for immunodetection methods.

#### Highlights:

- **Native avidin**—glycoprotein from chicken egg white, 67 kDa, pI 10
- **Soluble**—glycosylation and isoelectric point render avidin more soluble than streptavidin
- **Affordable**—significantly less expensive than recombinant streptavidin
- **ABC staining**—avidin is usually preferred over streptavidin for avidin–biotin complex (ABC) staining methods in IHC
- **HRP conjugate**—for membrane- and plate-based assays

### Fluorescent streptavidin and avidin conjugates

#### Alexa Fluor fluorescent streptavidin, avidin, and NeutrAvidin conjugates

We offer a broad range of Invitrogen™ Alexa Fluor™ fluorescent streptavidin, avidin, and NeutrAvidin conjugates (Table 9). Benefits of the Alexa Fluor Dyes and their conjugates include exceptionally bright conjugates, excellent photostability, and a broad range of emission colors (Figure 11).

Table 9. Alexa Fluor fluorescent streptavidin conjugates.

Streptavidin conjugates	Ex (nm)*	Em (nm)*
Alexa Fluor 350 streptavidin	346	442
Alexa Fluor 405 streptavidin	402	421
Alexa Fluor 430 streptavidin	434	539
Alexa Fluor 488 streptavidin	495	519
Alexa Fluor 500 streptavidin	503	525
Alexa Fluor 514 streptavidin	518	540
Alexa Fluor 532 streptavidin	530	554
Alexa Fluor 555 streptavidin	555	565
Alexa Fluor 546 streptavidin	556	573
Alexa Fluor 568 streptavidin	578	603
Alexa Fluor 594 streptavidin	590	617
Alexa Fluor 610 streptavidin	612	628
Alexa Fluor 633 streptavidin	632	647
Alexa Fluor 635 streptavidin	633	647
Alexa Fluor 647 streptavidin	650	668
Alexa Fluor 660 streptavidin	663	690
Alexa Fluor 680 streptavidin	679	702
Alexa Fluor 700 streptavidin	702	723
Alexa Fluor 750 streptavidin	749	775
<b>NeutrAvidin and avidin conjugates</b>		
Alexa Fluor 350 NeutrAvidin	346	442
Alexa Fluor 488 avidin	495	519

\* Approximate fluorescence excitation (Ex) and emission (Em) maxima for conjugates, in nm.

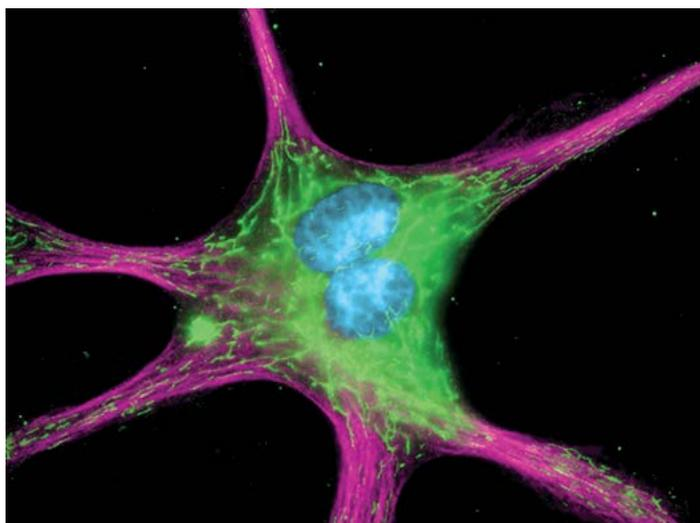


Figure 11. The cytoskeleton of a fixed and permeabilized bovine pulmonary artery endothelial (BPAE) cell. Tubulin was detected using mouse monoclonal anti- $\alpha$ -tubulin antibody and visualized with Alexa Fluor 647 goat anti-mouse IgG antibody (pseudocolored magenta). Endogenous biotin in mitochondria was labeled with green-fluorescent Alexa Fluor 488 streptavidin; DNA was stained with blue-fluorescent DAPI.

### Alexa Fluor streptavidin phycobiliprotein conjugates

Invitrogen™ Alexa Fluor™ streptavidin phycobiliprotein conjugates are fluorescent streptavidin conjugates with tandem dyes RPE (R-phycoerythrin) and APC (allophycocyanin) labeled with long-wavelength Alexa Fluor dyes (Table 10). These tandem conjugates are comprised of a donor phycobiliprotein-RPE conjugate or APC conjugated to longer wavelength-emitting fluorescence acceptors (Table 10). By the process of fluorescence resonance energy transfer (FRET), an energy transfer cascade is established wherein most of the light absorbed by the donor RPE (or APC) results in fluorescence of the acceptor dye. This process can be quite efficient, resulting in almost total transfer of energy to the acceptor dye.

Table 10. Tandem Alexa Fluor streptavidin phycobiliprotein conjugates.

Tandem RPE conjugates		
Streptavidin conjugates	Ex (nm)*	Em (nm)*
Alexa Fluor 610–RPE streptavidin	496, 546, 565	630
Alexa Fluor 647–RPE streptavidin	496, 546, 565	668
Alexa Fluor 680–RPE streptavidin	496, 546, 565	702
Alexa Fluor 750–RPE streptavidin	496, 546, 565	771
APC and tandem APC conjugates		
Streptavidin conjugates	Ex (nm)*	Em (nm)*
Allophycocyanin (APC) streptavidin	650	660
Alexa Fluor 680–APC streptavidin	650	702
Alexa Fluor 700–APC streptavidin	650	723
Alexa Fluor 750–APC streptavidin	650	775

\* Approximate fluorescence excitation (Ex) and emission (Em) maxima for conjugates, in nm.

### Qdot streptavidin conjugates

Invitrogen™ Qdot™ streptavidin conjugates combine the highly specific binding properties of streptavidin with the exceptional photostability of Qdot nanocrystals. The large surface area afforded by the Qdot nanocrystal allows simultaneous conjugation of multiple streptavidin molecules to a single fluorophore. Advantages conferred by this approach include increased avidity for targets, the potential for cooperative binding in some cases, and the use of efficient signal amplification methodologies. For example, combining biotin-functionalized products with the streptavidin labels allows for successive enhancements in signal via “sandwiching” (streptavidin/biotin/streptavidin, etc.) following an initial labeling step. Unlike the Qdot streptavidin conjugates, the Invitrogen™ Qdot™ ITK™ streptavidin conjugates have the streptavidin covalently attached to the inner amphiphilic coating without a polyethylene glycol (PEG) linker (Table 11).

Table 11. Qdot streptavidin conjugates.

Streptavidin conjugates	Dye
Qdot 525 Streptavidin Conjugate	Qdot 525
Qdot 525 ITK Streptavidin Conjugate Kit	Qdot 525
Qdot 545 ITK Streptavidin Conjugate Kit	Qdot 545
Qdot 565 Streptavidin Conjugate	Qdot 565
Qdot 565 ITK Streptavidin Conjugate Kit	Qdot 565
Qdot 585 Streptavidin Conjugate	Qdot 585
Qdot 585 ITK Streptavidin Conjugate Kit	Qdot 585
Qdot 605 Streptavidin Conjugate	Qdot 605
Qdot 605 ITK Streptavidin Conjugate Kit	Qdot 605
Qdot 625 Streptavidin Conjugate	Qdot 625
Qdot 655 Streptavidin Conjugate	Qdot 655
Qdot 655 ITK Streptavidin Conjugate Kit	Qdot 655
Qdot 705 Streptavidin Conjugate	Qdot 705
Qdot 705 ITK Streptavidin Conjugate Kit	Qdot 705
Qdot 800 Streptavidin Conjugate	Qdot 800
Qdot 800 ITK Streptavidin Conjugate Kit	Qdot 800



# Protein biology

## Gel electrophoresis

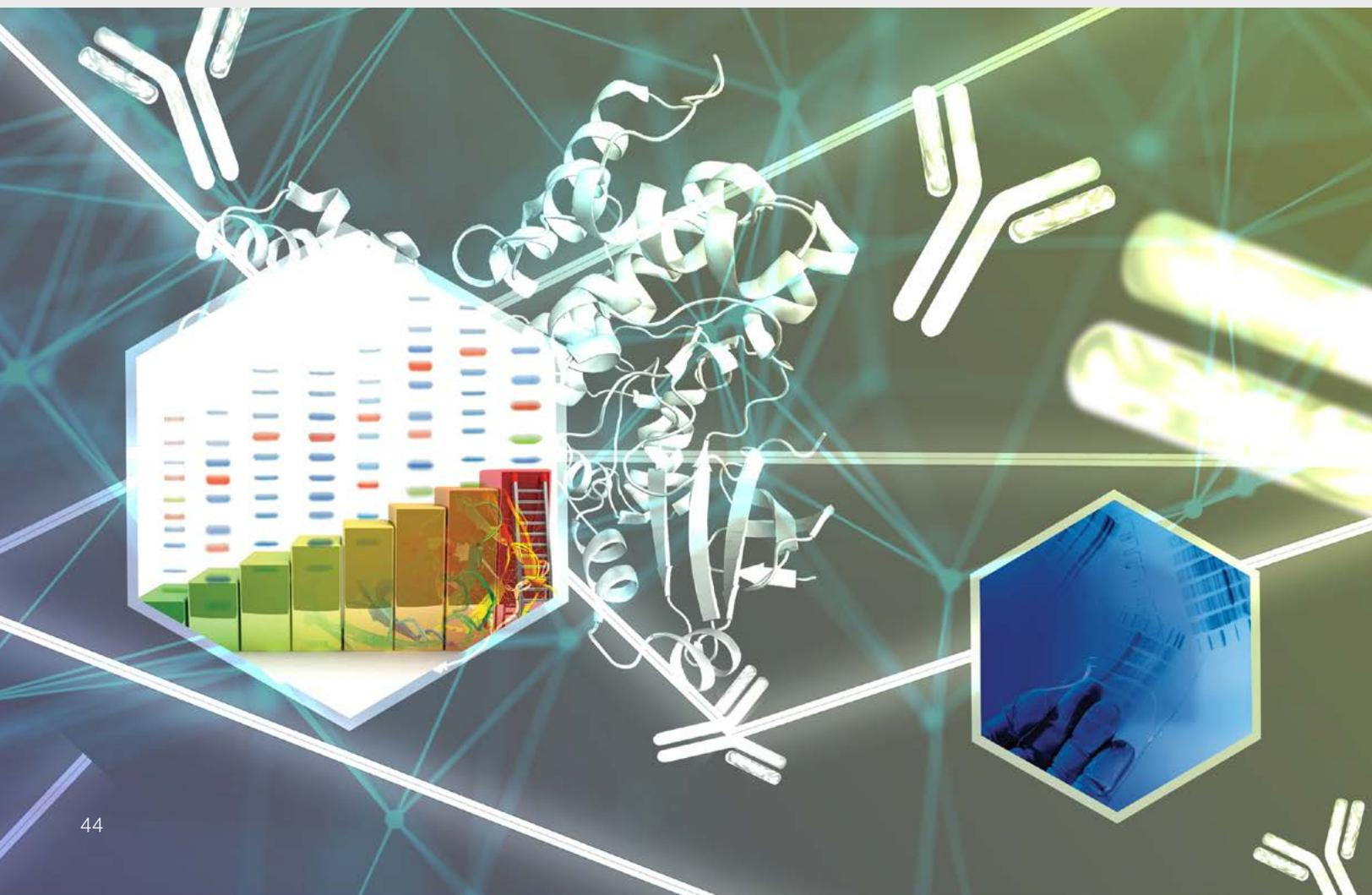
### Protein standards and ladders

Choose from a variety of protein standards and ladders (molecular weight markers) for **protein electrophoresis** and **western blotting applications**, including SDS-PAGE and isoelectric focusing (IEF). All standards and ladders are supplied in ready-to-use formats, eliminating the need to reduce, pre-mix, or add loading dyes. These protein standards are consistent from lot to lot and are strictly quality controlled on appropriate gels to help ensure consistent band migration and intensity. All prestained protein ladders are compatible for use in western blotting.

### PageRuler Prestained Protein Ladder

Thermo Scientific™ PageRuler™ Prestained Protein Ladder is a mixture of 10 blue-, orange-, and green-stained proteins, ranging from 10 to 180 kDa, for use as size standards in protein electrophoresis (SDS-PAGE) and western blotting. The ladder is supplied in a ready-to-use format for direct loading onto gels—no need to heat, reduce, or add sample buffer prior to use.

Find out more about our diverse array of protein ladders on [thermofisher.com/proteinladders](https://www.thermofisher.com/proteinladders)



# Mass spectrometry

## Tandem mass tags

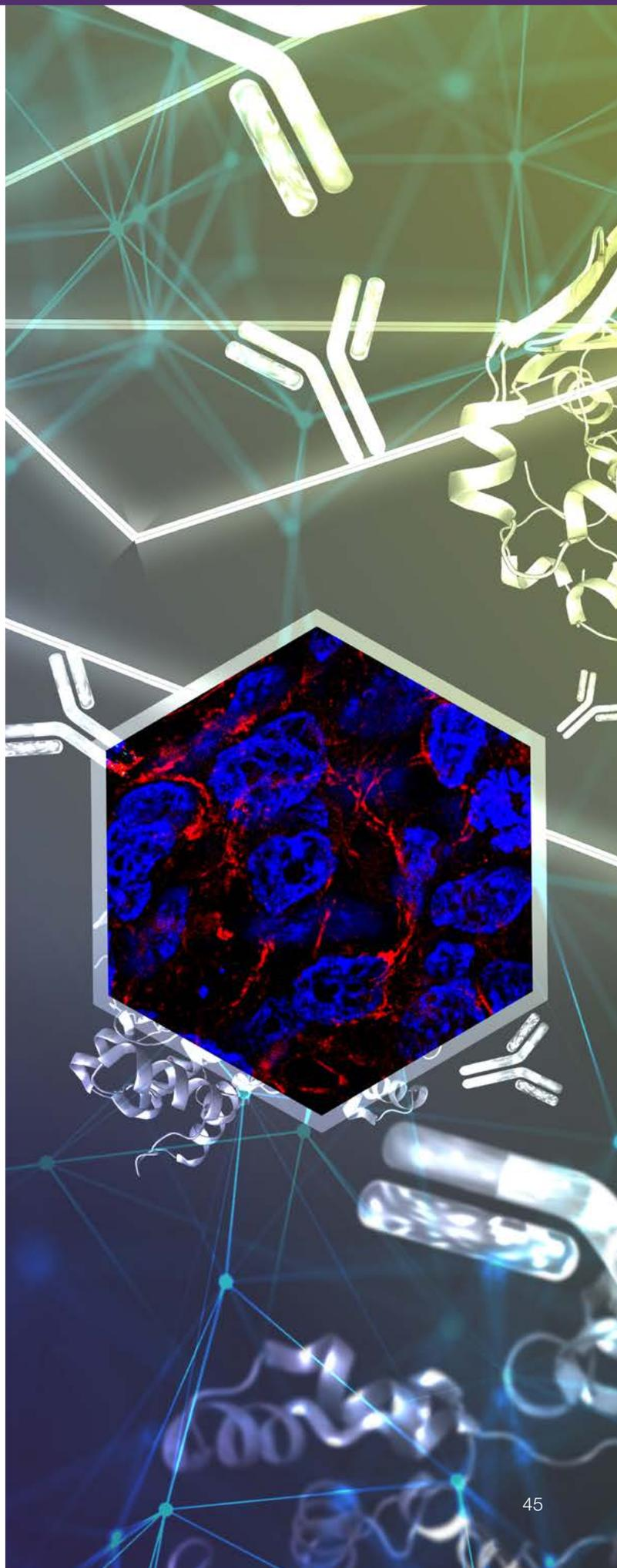
Thermo Scientific™ Tandem Mass Tag (TMT) reagents are designed to enable identification and quantitation of proteins in different samples using tandem mass spectrometry (MS). All mass tagging reagents within a set have the same nominal mass (i.e., are isobaric) and chemical structure composed of an amine-reactive NHS ester group, a spacer arm (mass normalizer), and a mass reporter. Commercial use of TMT Reagents as a service requires a license.

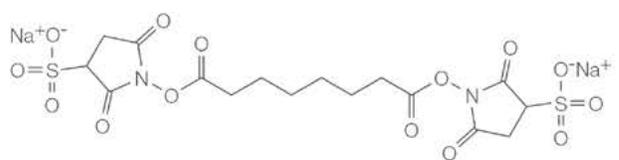
## New! TMTpro 16plex Label Reagent Set

Thermo Scientific™ TMTpro™ 16plex label reagents are the next generation of tandem mass tags, designed to increase the level of sample multiplexing without compromising on protein identification and quantitation. The new TMTpro label reagents are similar in design to TMT label reagents in that they are both isobaric and amine-reactive, but TMTpro reagents differ in structure, having a longer spacer region and isobutyl proline mass reporter region. The TMTpro 16plex reagents have the same labeling efficiency and peptide/protein ID rates and provide the same level of quantitative precision as the original TMT reagents, with the benefit of improved quantitative accuracy with larger sample sets.

### Highlights:

- **Multiplex**—concurrent MS analysis of up to 16 samples derived from cells, tissues, or biological fluids
- **Robust**—increased multiplex capability results in fewer missing quantitative values among samples and higher confidence among replicates
- **Efficient**—amine-reactive, NHS ester-activated reagents ensure efficient labeling of all peptides regardless of protein sequence or proteolytic enzyme specificity
- **Compatible**—optimized for use with high resolution Thermo Scientific MS/MS platforms, such as the Thermo Scientific™ Q Exactive™ and Orbitrap Fusion™ Tribrid™ instrument series, including the Orbitrap Eclipse™ Tribrid™ and Orbitrap Exploris™ 480 mass spectrometers with data analysis fully supported by Thermo Scientific™ Proteome Discoverer™ Software.





**BS<sup>3</sup>**

Bis(sulfosuccinimidyl)suberate

MW 572.43

Spacer arm 11.4 Å

## Linking mechanisms

### Crosslinkers

Crosslinking reagents are used in many techniques to assist in immobilization of a ligand to a solid phase, and also for the conjugation of a detection antibody to an enzyme or fluorescent dye. We offer crosslinkers, PEGylation reagents, biotinylation kits and reagents, and antibody and enzyme labeling kits. In addition, we provide custom conjugation of antibodies and proteins.

Crosslinking is the process of chemically joining two or more molecules by a covalent bond. Crosslinking reagents, or crosslinkers, are molecules that contain two or more reactive ends capable of chemically attaching to specific functional groups (primary amines, sulfhydryls, etc.) on proteins or other molecules.

Crosslinkers are defined by their reactivities:

**Homobifunctional crosslinkers** have identical reactive groups at either end of a spacer arm. Generally, they must be used in one-step reaction procedures to randomly “fix” or polymerize molecules containing like functional groups. This is useful for capturing a snapshot of all protein interactions, but it cannot provide the precision needed for other types of crosslinking applications.

**Heterobifunctional crosslinkers** possess different reactive groups at either end. These reagents not only allow for single-step conjugation of molecules that have the respective target functional groups, but they also allow for sequential (two-step) conjugations that minimize undesirable polymerization or self-conjugation. In sequential procedures, heterobifunctional reagents are reacted with one protein using the most labile group of the crosslinker first. After removing excess unreacted crosslinker, the modified first protein is added to a solution containing the second protein where reaction through the second reactive group of the crosslinker occurs.

Heterobifunctional crosslinkers are the preferred choice for antibody-enzyme or other protein-to-protein crosslinking applications. Unwanted self-conjugation inherent when using homobifunctional NHS-ester reagents or glutaraldehyde can be avoided by reagents such as SMCC or sulfo-SMCC. In separate reactions, one protein can be reacted to the amine-specific end of this reagent while the other protein is treated with a reducing agent or sulfhydryl-addition reagent to expose or create sulfhydryl groups. Finally, after removing excess unreacted reagents, the two proteins can be mixed to allow the sulfhydryl-reactive groups (e.g., maleimide) of the first protein to conjugate with the sulfhydryl groups of the second protein (Figure 12).

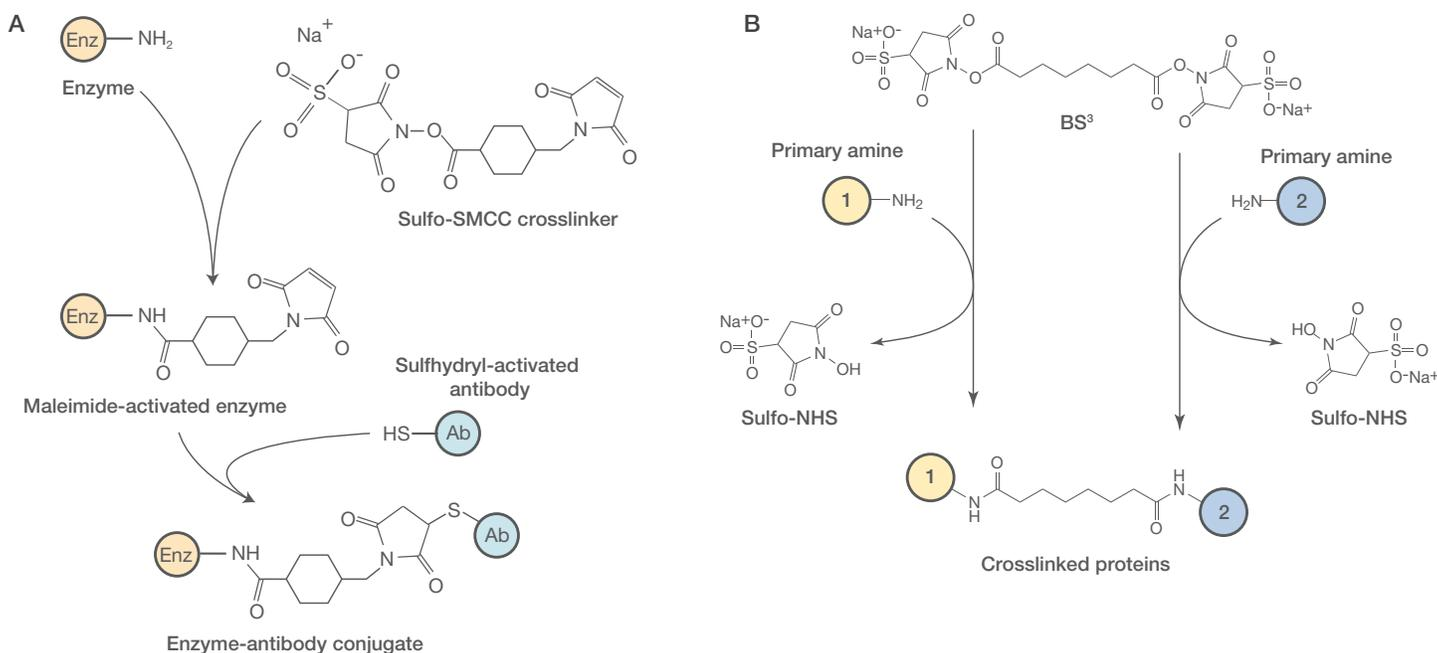


Figure 12. Example reactions for producing detection conjugates. **(A)** Enzyme–antibody conjugation with heterobifunctional crosslinker Sulfo-SMCC. **(B)** Use of homobifunctional crosslinker BS<sup>3</sup>.

## Pierce crosslinkers

Many types of Thermo Scientific™ Pierce™ crosslinkers are available for protein, peptide, and other macromolecular immobilization and conjugation needs. Both homobifunctional and heterobifunctional crosslinkers are offered with a variety of spacer-arm lengths, solubility, and cleaving characteristics.

Our wide selection of crosslinking reagents now includes those that contain discrete-length polyethylene glycol (PEG) spacers. These PEG groups increase reagent and conjugate solubility, minimize toxic and immunological effects compared to non-PEG spacers, and provide several options for accommodating specific crosslinking distances (Figure 13).

### PEG crosslinkers offer:

- Increased water solubility
- Decreased aggregation
- Better accessibility

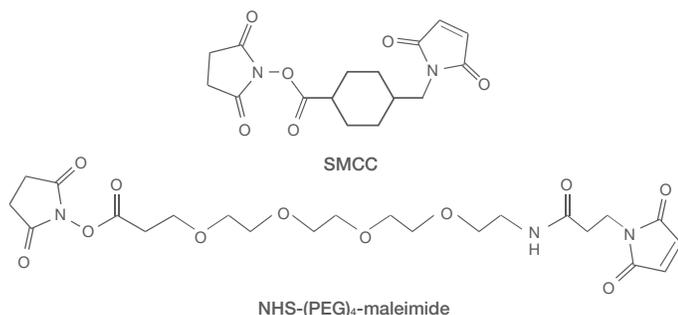
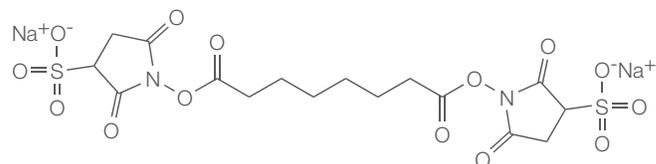


Figure 13. Chemical structures of SMCC and PEG alternative NHS-(PEG)<sub>4</sub>-maleimide.

Pierce crosslinkers are manufactured in our organic synthesis lab at scales ranging from milligram to hundreds of kilograms. Available quality tests include visual inspection, structural confirmation by infrared spectrometry (IR), purity by HPLC or quantitative NMR (qNMR), solvent analysis by GC headspace, water content by Karl Fischer titration, and elemental analysis. Custom specifications can also be set.

## Pierce BS<sup>3</sup> crosslinker



**BS<sup>3</sup>**

Bis(sulfosuccinimidyl)suberate  
MW 572.43  
Spacer arm 11.4 Å

Thermo Scientific™ Pierce™ Premium Grade BS<sup>3</sup> (Sulfo-DSS) crosslinker is bis(sulfosuccinimidyl)suberate, an amine-to-amine crosslinker that is homobifunctional, water-soluble, noncleavable, and membrane-impermeant.

BS<sup>3</sup> contains an amine-reactive *N*-hydroxysulfosuccinimide (NHS) ester at each end of an 8-carbon spacer arm. NHS esters react with primary amines at pH 7–9 to form stable amide bonds, along with release of the *N*-hydroxysulfosuccinimide leaving group. Proteins, including antibodies, generally have several primary amines in the side chain of lysine (K) residues and the N-terminus of each polypeptide that are available as targets for NHS-ester crosslinking reagents.

### Features of Pierce BS<sup>3</sup> crosslinker:

- Reactive groups: sulfo-NHS ester (both ends)
- Reactive with amino groups (primary amines)
- Amine-reactive sulfo-NHS ester reacts rapidly with any primary amine-containing molecule
- Water-soluble, compared to Thermo Scientific™ Pierce™ DSS
- Membrane-impermeant, allowing for cell surface labeling
- High-purity, crystalline reagent can be used to create high-purity crosslinked conjugates

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

Because it contains the hydrophilic sulfonyl moiety, Pierce BS<sup>3</sup> crosslinker is soluble up to ~100 mM in water and many commonly used buffers, thus avoiding the use of organic solvents, which may perturb protein structure. Pierce DSS, the non-water-soluble analog of BS<sup>3</sup>, is also available for applications that require a less hydrophilic crosslinker (e.g., to effect intracellular crosslinking). DSS and BS<sup>3</sup> have essentially identical crosslinking activity toward primary amines.

**Properties of Pierce BS<sup>3</sup>:**

- Alternative names: Sulfo-DSS
- Molecular formula: C<sub>16</sub>H<sub>18</sub>N<sub>2</sub>O<sub>14</sub>S<sub>2</sub>Na<sub>2</sub>
- Molecular weight: 572.43
- Spacer-arm length: 11.4 Å (8 atoms)
- CAS number: 82436-77-9
- Reactive groups: sulfo-NHS esters react with primary amines at pH 7–9

**Applications:**

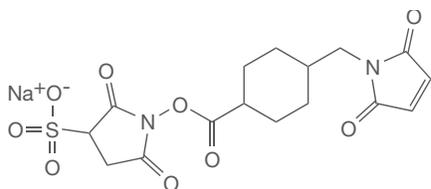
- Protein crosslinking to create bioconjugates via single-step reactions
- Crosslink cell-surface proteins prior to cell lysis and immunoprecipitation
- Immobilize proteins onto amine-coated surfaces
- Identify receptor-ligand interactions by chemical crosslinking
- “Fix” protein interactions to allow identification of weak or transient protein interactions
- Identify near-neighbor protein interactions

**Pierce BS<sup>3</sup> specifications:**

We manufacture Pierce BS<sup>3</sup> crosslinker to high specifications to produce specific bioconjugates, to help ensure the integrity of your data, and to offer you a high degree of consistency. Each lot of Pierce BS<sup>3</sup> crosslinker is tested to meet the following minimum specifications:

- Purity: >93% by quantitative NMR (the highest standard for crosslinker purity)
- Solubility: >5.8 mg/mL in DI water, clear solution with no insoluble material

## Pierce Sulfo-SMCC crosslinker



### Sulfo-SMCC

Sulfosuccinimidyl 4-(N-maleimidomethyl)cyclohexane-1-carboxylate  
MW 436.37  
Spacer arm 8.3 Å

Thermo Scientific™ Pierce™ Sulfo-SMCC is a water-soluble, amine-to-sulfhydryl crosslinker that contains NHS-ester and maleimide reactive groups at opposite ends of a medium-length cyclohexane spacer arm (8.3 Å).

#### Features of Pierce Sulfo-SMCC crosslinker:

- Reactive groups: sulfo-NHS ester and maleimide
- Reactive with amino and sulfhydryl groups
- Water-soluble (compared to SMCC), so crosslinking can be done in physiologic solutions
- High-purity, crystalline reagent can be used to create high-purity, maleimide-activated derivatives
- Cyclohexane bridge confers added stability to the maleimide group making Sulfo-SMCC the ideal crosslinking agent for maleimide activation of proteins. Maleimide groups are stable for 64 hours in 0.1 M sodium phosphate buffer, pH 7 at 4°C

Sulfosuccinimidyl-4-(N-maleimidomethyl)cyclohexane-1-carboxylate (Sulfo-SMCC) is a noncleavable and membrane-impermeant crosslinker. It contains an amine-reactive NHS ester and a sulfhydryl-reactive maleimide group. NHS esters react with primary amines at pH 7–9 to form stable amide bonds. Maleimides react with sulfhydryl groups at pH 6.5–7.5 to form stable thioether bonds. The maleimide groups of Sulfo-SMCC and SMCC are unusually stable up to pH 7.5 because of the cyclohexane bridge in the spacer arm. Because it contains the hydrophilic sulfonyl moiety, Pierce Sulfo-SMCC crosslinker is soluble up to ~10 mM in water and many commonly used buffers, thus avoiding the use of organic solvents, which may perturb protein structure.

#### Applications:

- Enzyme labeling of antibodies—both enzyme and antibody specificity can be preserved
- Create specific bioconjugates via one- or two-step crosslinking reactions
- Create sulfhydryl-reactive, maleimide-activated carrier proteins for coupling haptens

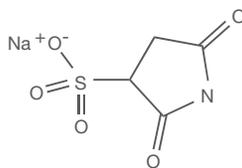
#### Specifications:

Pierce Sulfo-SMCC crosslinker is manufactured to high specifications to produce specific bioconjugates and maleimide-activated proteins, to help ensure the integrity of your data and to offer you a high degree of crosslinking consistency. Each lot of Pierce Sulfo-SMCC crosslinker is tested to meet the following minimum specifications.

- Identity: IR scan shows only peaks characteristic of the structure and functional groups of Sulfo-SMCC crosslinker
- Purity: ≥90% by quantitative NMR (the highest standard for crosslinker purity)
- Solubility: ≥10 mg/mL in DI water

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Pierce Sulfo-NHS reagent



Sulfo-NHS Acetate  
M.W. 259.17

Thermo Scientific™ Pierce™ Sulfo-NHS (*N*-hydroxysulfo-succinimide) is a chemical modification reagent for converting carboxyl groups to amine-reactive NHS esters for bioconjugation, crosslinking, labeling, and immobilization methods.

### Features of Pierce Sulfo-NHS reagent:

- Efficiency of EDC-mediated coupling is increased in the presence of NHS or sulfo-NHS
- Amine-reactive NHS esters or sulfo-NHS esters can be made with any carboxyl-containing molecule
- Sulfo-NHS derivatives are usually readily water-soluble (can be added directly to physiologic buffers) and membrane-impermeant (can be used for cell surface labeling)
- High-purity, crystalline sulfo-NHS can be used to create high-quality activated derivatives

Sulfo-NHS enables control and modification of carbodiimide crosslinking reactions involving activation of carboxylates ( $-\text{COOH}$ ) for conjugation with primary amines ( $-\text{NH}_2$ ).

Derivatives are easily synthesized by mixing the sulfo-NHS with a carboxyl-containing molecule and a dehydrating agent such as the carbodiimide EDC (EDAC). The method is the basis for generating many types of protein labeling reagents, including amine-reactive fluorescent dyes, biotin affinity tags, and PEGylation compounds.

### Applications:

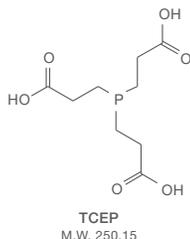
- Improve efficiency of EDC coupling reactions
- Convert carboxyls to amine-reactive sulfo-NHS esters
- Crosslink proteins to carboxyl-coated beads or surfaces more efficiently
- Activate nanoparticles with amine-reactive sulfo-NHS esters

### Specifications for Pierce Sulfo-NHS reagent:

We manufacture Sulfo-NHS to high specifications to produce specific bioconjugates, to help ensure the integrity of your data, and to offer you a high degree of consistency. Each lot of Pierce sulfo-NHS reagent is tested to meet the following minimum specifications:

- **Purity**—greater than 95% by quantitative NMR (the highest standard for crosslinker purity); average lot purity is greater than 99%
- **Solubility**—sample dissolves at 2 mg/mL in deionized water to yield a clear, colorless solution
- **Identity**—IR scan shows only peaks characteristic of sulfo-NHS

## Bond-Breaker TCEP Solution, Neutral pH



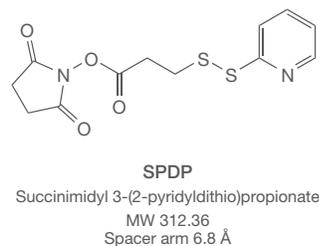
Thermo Scientific™ Bond-Breaker™ TCEP Solution, Neutral pH, is a stable, 0.5 M solution of the thiol-free, phosphine-based TCEP compound, useful as a 10X stock for addition to SDS-PAGE sample loading buffers to reduce protein disulfide bonds.

### Features of Bond-Breaker TCEP Solution, Neutral pH:

- Unlike dithiothreitol (DTT) and beta-mercaptoethanol (beta-ME), TCEP is odor-free
- Efficient—5 to 50 mM TCEP is designed to thoroughly reduce most peptide or protein disulfide bonds within a few minutes (i.e., just as effective as DTT)
- Enables specific and complete reduction of even the most stable water-soluble alkyl disulfides
- Reduces protein disulfides at room temperature and pH 5, typically in less than 5 minutes
- Resistant to air oxidation; nonvolatile and nonreactive toward other protein functional groups
- Reduces peptides and proteins over a broad range of pH, salt, detergent, and temperature conditions
- Removal of the reducing agent is not necessary before most applications (e.g., histidine-tagged protein purification, maleimide conjugations), because TCEP does not contain sulfhydryl groups

Bond-Breaker TCEP Solution is a potent, odorless, thiol-free reducing agent with broad application to protein modification and other research involving reduction of disulfide bonds. It is an effective and convenient replacement for beta-ME or DTT in SDS-PAGE sample buffers. The neutral pH of this reagent provides sharp bands and avoids exposing proteins to the strong acid of TCEP-HCl, which can result in hydrolysis and carbohydrate modification.

## Pierce SPDP crosslinker



Thermo Scientific™ Pierce™ SPDP (succinimidyl 3-(2-pyridyldithio)propionate) is a short-chain crosslinker for amine-to-sulfhydryl conjugation via NHS-ester and pyridyldithiol reactive groups that form cleavable (reducible) disulfide bonds with cysteine sulfhydryl groups.

### Features of Pierce SPDP crosslinker:

- Reactive groups: NHS ester and pyridyldithiol
- Reactive with amino and sulfhydryl groups
- Releases a detectable by-product when reacted with a free sulfhydryl group; by measuring the release of pyridine-2-thione at 343 nm, the reaction can be easily monitored
- Disulfide bond in the spacer arm is readily cleaved by 10–50 mM DTT or TCEP at pH 8.5
- Spacer arm is also easily cleaved using reducing SDS-PAGE sample loading buffer
- Cleavable crosslinker allows separation of crosslinked products
- Water-insoluble (dissolve first in DMF or DMSO)
- SPDP crosslinker is membrane-permeant, so crosslinking can be done inside the cells

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

**Applications:**

- Reversible crosslinking of proteins
- Adding sulfhydryl groups to proteins by attachment to lysine residues
- Protein crosslinking to create antibody-enzyme and other bioconjugates via a two-step reaction
- Attaching haptens to carrier proteins for antibody production
- Preparing immunotoxin conjugates

**Specifications:**

We manufacture Pierce SPDP crosslinker to high specifications to produce specific bioconjugates to help ensure the integrity of your data and to offer you a high degree of consistency. Each lot of SPDP is tested to ensure it is >95% pure by quantitative NMR (the highest standard for crosslinker purity).

We offer three analogs of SPDP: the standard version (SPDP), a derivative with a longer spacer arm (LC-SPDP), and a sulfonated water-soluble variety (sulfo-LC-SPDP). SPDP (*N*-succinimidyl 3-(2-pyridyldithio)-propionate) and LC-SPDP (succinimidyl 6-(3-[2-pyridyldithio]-propionamido)hexanoate) are heterobifunctional, thiol-cleavable, and membrane-permeant crosslinkers. They each contain an amine-reactive NHS ester that will react with lysine residues to form a stable amide bond. The other end of the spacer arm is terminated in the pyridyl disulfide group that will react with sulfhydryls to form a reversible disulfide bond.

**Pierce EDC crosslinker**

Thermo Scientific™ Pierce™ EDC (1-ethyl-3-(3-dimethylamino-propyl)carbodiimide hydrochloride) is a water-soluble carbodiimide crosslinker that activates carboxyl groups for spontaneous reaction with primary amines, enabling peptide immobilization and hapten–carrier protein conjugation.

EDC (or EDAC) is a zero-length crosslinking agent used to couple carboxyl groups to primary amines. This crosslinker has been used in diverse applications such as forming amide bonds in peptide synthesis, attaching haptens to carrier proteins to form immunogens, labeling nucleic acids through 5' phosphate groups, and creating amine-reactive NHS-esters of biomolecules. EDC reacts with a carboxyl to form an amine-reactive *O*-acylisourea intermediate. If this intermediate does not encounter an amine, it will hydrolyze and regenerate the carboxyl group. In the presence of sulfo-NHS, EDC can be used to convert carboxyl groups to amine-reactive sulfo-NHS esters. This is accomplished by mixing the EDC with a carboxyl-containing molecule and adding sulfo-NHS.

**Features of Pierce EDC crosslinker:**

- Reactive group: carbodiimide
- Reaction target: activates carboxyl groups to conjugate to amino groups (primary amines)
- React EDC alone with target groups or include NHS or sulfo-NHS to increase reaction efficiency or to stabilize active intermediate for later reaction to amines
- Forms neutral amide bonds between carboxyls and amines
- Water-soluble reagent can be added directly to reactions in aqueous, physiological buffers
- Soluble reaction byproducts are easily removed by washing with water or dilute acid
- High-purity, crystalline reagent can be used to create high-quality, activated derivatives

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

### Properties of Pierce EDC crosslinker:

- Molecular formula:  $C_8H_{17}N_3 \cdot HCl$
- Molecular weight: 191.7
- Spacer arm length: 0.0 Å
- CAS Number: 25952-53-8
- Reactive group: carbodiimide
- Reactivity: forms active intermediate with carboxyl groups at pH 4.7–6.0 (optimal), then intermediate reacts with primary amines

### Applications:

- Conjugate carboxyl and amino groups among peptides and proteins
- Couple haptens to immunogenic carrier proteins (e.g., attach a peptide to keyhole limpet hemocyanin (KLH))
- Immobilize peptide antigens to affinity-purify antibodies
- Create NHS-activated, amine-reactive labeling compounds
- Crosslink proteins to carboxyl-coated beads or surfaces
- Activate nanoparticles with amine-reactive sulfo-NHS esters
- DNA labeling through 5' phosphate groups

Also see Thermo Scientific™ Pierce™ Premium Grade EDC, page 62.

For information on all of our crosslinkers, go to [thermofisher.com/crosslinkers](https://www.thermofisher.com/crosslinkers)

## PEGylation reagents

We offer activated linear and branched derivatives of PEG for the PEG-modification of peptides and proteins via primary amines and sulfhydryl groups. These reagents are typically used to increase solubility, prolong stability, and reduce immunogenicity. Certain experimental systems and assay platforms depend on the ability to alter the mass, solubility, or other properties of proteins, immunogens, reaction vessels, and other materials.

### Amine-reactive PEGylation reagents

Covalent modification with PEG groups requires PEG compounds that contain a reactive or targetable functional group at one end. The simplest method for PEGylating proteins, which are rich in surface primary amines, is to use a PEG compound that contains an NHS ester group at one end.

- MS(PEG)<sub>n</sub> reagents are available in various PEG lengths (n = 4, 8, 12, and 24)
- TMS(PEG)<sub>n</sub> is a branched form of this reagent, containing three methyl-(PEG)<sub>12</sub> arms attached to a (PEG)<sub>4</sub>-NHS ester stem

### Sulfhydryl-reactive PEGylation reagents

MM(PEG)<sub>n</sub> compounds are linear reagents for PEGylating sulfhydryl groups. At the end, opposite the inert methyl group(s), is a maleimide moiety, which reacts to form stable thioether bonds with sulfhydryl groups.

- MM(PEG)<sub>n</sub> is available in two PEG lengths (n = 12 and 24)

### PEGylated amino acids and amine compounds

MA(PEG)<sub>n</sub> and CA(PEG)<sub>n</sub> are polyethylene glycol compounds of discrete length (n = 4, 8, 12, and 24) containing methyl-and-amine or carboxyl-and-amine ends. Although these functional groups are not spontaneously reactive, they are easily targeted by various crosslinking and immobilization reagents for the construction of surface chemistries.

## Pierce MS(PEG)<sub>n</sub> reagents

Thermo Scientific™ Pierce™ MS(PEG)<sub>n</sub> reagents are methyl-terminated, polyethylene glycol compounds (n = 4, 8, 12, and 24) activated as NHS esters for covalent PEGylation of primary amines on proteins (e.g., lysines) or assay surfaces.

### Features of MS(PEG)<sub>4</sub>:

- NHS-activated for efficient PEGylation of primary amines at pH 7–9; reaction of NHS-ester group results in formation of stable, irreversible amide bonds
- Fully characterized PEGylation reagents with defined PEG chain lengths; molecules of discrete molecular weight for consistency of performance in protein-modification applications
- Provided as a series of 4, 8, 12, and 24 ethylene glycol units, enabling modification procedures to be optimized for a specific application while retaining all the benefits associated with protein PEGylation
- PEG spacer provides unique advantages, including increased stability, reduced tendency toward aggregation, and reduced immunogenicity
- Easy-to-follow instructions help increase the likelihood of a successful outcome

MS(PEG)<sub>n</sub> is the abbreviation for a set of compounds having PEG spacers with methyl (–CH<sub>3</sub>) and amine-reactive NHS-ester groups at opposite ends. The unbranched, hydrophilic, discrete-length molecules have the form methyl-(PEG)<sub>n</sub>-NHS ester, where the subscript “n” denotes 4, 8, 12, or 24 ethylene glycol units. The NHS ester is spontaneously reactive with primary amines (–NH<sub>2</sub>), providing for efficient PEGylation of proteins, peptides, and other amine-containing molecules or surfaces.

### PEGylation applications:

- PEGylate amine surfaces
- Add inert mass to proteins, immunogens, drug compounds, and probes
- Improve solubility (decrease aggregation) of proteins or peptides without affecting function
- Protect proteins from proteolysis

### Advantages of discrete-length methyl-PEG-NHS ester compounds

These reagents are specially synthesized as homogeneous compounds of discrete chain length and defined molecular weight. As such, they enable precise control and optimization of surface protein modification experiments. By contrast, typical preparations of PEG compounds are heterogeneous mixtures composed of multiple chain lengths and a range of molecular weights.

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Why PEGylate a protein or peptide?

PEG-containing reagents have been used to modify proteins to provide specific advantages. Protein PEGylation can improve the stability of the modified protein, protect it from proteolytic digestion, increase its half-life in biological applications, mask it from causing an immunogenic response, decrease its antigenicity or potential toxicity, improve its solubility, diminish the potential for aggregation, and minimize interference for both *in vitro* and *in vivo* applications. Polyethylene glycol, also called polyethylene oxide (PEO), has these effects because it is nontoxic, non-immunogenic, hydrophilic, water-soluble, and highly flexible.

For more information on our PEGylation reagents, go to [thermofisher.com/pegylation](http://thermofisher.com/pegylation)

## Biotinylation reagents

The biotin–avidin interaction is commonly exploited to detect and/or purify proteins because of the high specificity that these two molecules have for each other. Biotinylation is the process of attaching biotin to proteins and other macromolecules. Biotinylation reagents are available for targeting specific functional groups or residues, including primary amines, sulfhydryls, carboxyls, and carbohydrates (Figure 14). Photoreactive biotin compounds that react nonspecifically upon exposure to ultraviolet (UV) light are also available and expand the scope of the molecules that may be biotinylated.

The variety of biotinylation reagents with different functional group specificities is extremely useful, allowing one to choose a reagent that does not inactivate the target macromolecule.

- Primary amine groups ( $-\text{NH}_2$ ) are the most commonly targeted functional groups for biotinylation because of the abundance of lysine side chains and N-terminal amines
- Sulfhydryl groups ( $-\text{SH}$ ), which are found in exposed cysteine residues, are the second most common targets for biotinylation
- Carbohydrate residues containing *cis*-diols can be oxidized to create active aldehydes ( $-\text{CHO}$ )

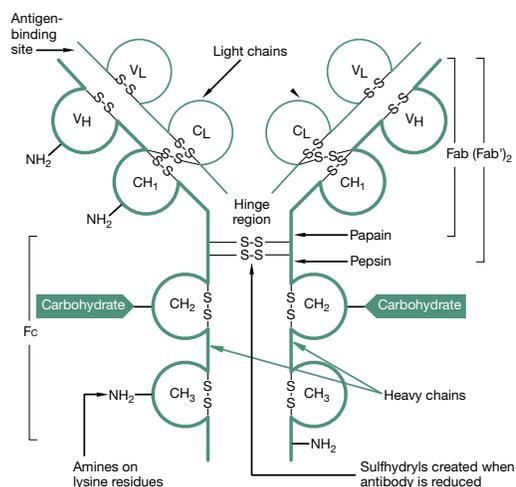


Figure 14. Functional groups available on an antibody for labeling.

## EZ-Link biotinylation reagents

Thermo Scientific™ EZ-Link™ biotinylation reagents and kits are used to chemically tag and label antibodies, proteins, and peptides with biotin at particular amino acid functional groups to create labeled antibody or other probes for streptavidin affinity binding, purification, and detection.

Our wide selection of reagents and kits includes traditional biotin-labeling reagents and PEG forms of biotinylation reagents. Compared to traditional reagents, PEG biotinylation reagents offer:

- Long-chain biotin alternatives
- Water-soluble, hydrophilic spacers
- PEG-biotinylated antibodies are more soluble after biotinylation
- Increased signal in ELISA applications (Figure 15)
- No aggregation issues (Figure 16)

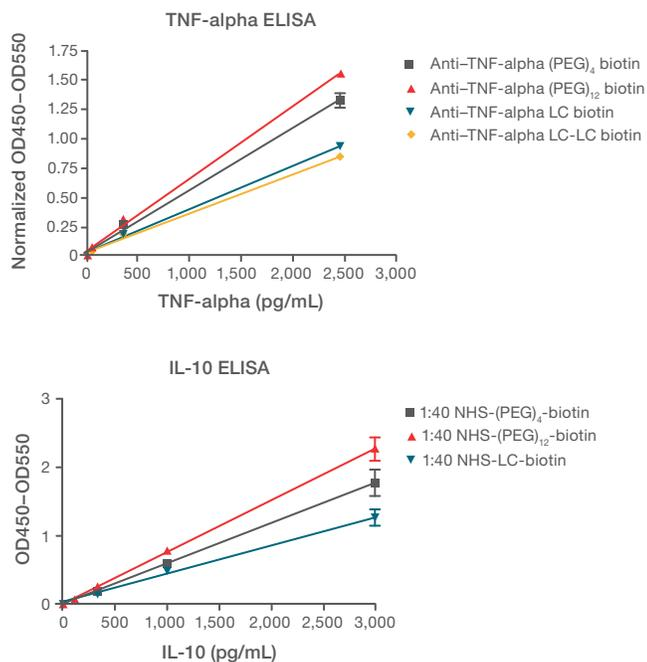


Figure 15. Increased signal with PEG biotinylation reagents.

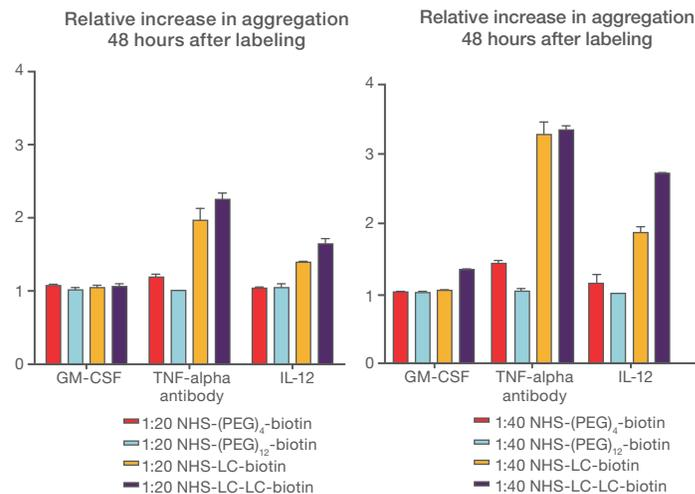


Figure 16. Reduction of aggregation with PEG biotinylation reagents.

## EZ-Link NHS-PEG<sub>4</sub>-Biotin

Thermo Scientific™ EZ-Link™ NHS-PEG<sub>4</sub>-Biotin is a long (29.0 Å), PEGylated, water-soluble reagent for simple and efficient biotin labeling of antibodies, proteins, and other primary amine-containing macromolecules. The NHS ester group reacts specifically and efficiently with lysine and N-terminal amino groups to form stable amide bonds. The hydrophilic polyethylene glycol (PEG) spacer arm imparts water solubility that is transferred to the biotinylated molecule, thus reducing aggregation of labeled proteins stored in solution. The PEG spacer arm also gives the reagent a long and flexible connection to minimize steric hindrance for binding to avidin molecules.

### Highlights:

- **Protein labeling**—biotinylate antibodies or other proteins for detection or purification using streptavidin probes or resins
- **Amine-reactive**—reacts with primary amines (–NH<sub>2</sub>), such as the side-chain of lysines (K) or the amino-termini of polypeptides
- **Hydrophilic**—spacer arm contains a hydrophilic, 4-unit, PEG group
- **Enhances solubility**—PEG spacer imparts water solubility to the biotinylated molecule, helping to prevent aggregation of biotinylated antibodies stored in solution
- **Irreversible**—forms permanent amide bonds; spacer arm cannot be cleaved
- **Long reach**—spacer arm (total length added to target) is 29 Å; this reduces steric hindrance when binding to avidin molecules

## EZ-Link Sulfo-NHS-LC-Biotin

Thermo Scientific™ EZ-Link™ Sulfo-NHS-LC-Biotin is an intermediate-length, water-soluble biotinylation reagent for labeling antibodies, proteins, and other molecules that have primary amines.

### Highlights:

- **Protein labeling**—biotinylate antibodies to facilitate immobilization, purification or detection using streptavidin resins or probes
- **Cell surface labeling**—biotinylates only surface proteins of whole cells because the negatively charged reagent does not permeate cell membranes
- **Amine-reactive**—reacts with primary amines (–NH<sub>2</sub>), such as lysine side-chains, or the amino-termini of polypeptides
- **Soluble**—charged sulfo-NHS group increases reagent water solubility compared to ordinary NHS-ester compounds
- **Irreversible**—forms permanent amide bonds; spacer arm cannot be cleaved
- **Medium length**—spacer arm (total length added to target) is 22.4 Å; provides excellent balance between adequate length (minimal steric hindrance for biotin binding) and modest mass

Sulfo-NHS-LC-biotin is one of three very similar EZ-Link reagents that are water-soluble and noncleavable and enable simple and efficient biotinylation of antibodies, proteins, and any other primary amine-containing macromolecules in solution. Specific labeling of cell surface proteins is another common application for these uniquely water-soluble and membrane-impermeant reagents. Differing only in their spacer arm lengths, the three sulfo-NHS-ester reagents offer the possibility of optimizing labeling and detection experiments where steric hindrance of biotin binding is an important factor (Table 12).

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

Table 12. EZ-Link biotinylation reagents.

Amine biotinylation reagents						
Product categories	Product	Product features	Spacer arm (Å)	Water-soluble?	Cleavable?	
NHS-PEG <sub>4</sub> biotinylation kits	• EZ-Link NHS-PEG <sub>4</sub> Biotinylation Kit	Maximize the solubility and stability of antibodies and other proteins	29	Yes	No	
	• EZ-Link Micro NHS-PEG <sub>4</sub> Biotinylation Kit					
Sulfo-NHS-LC biotinylation kits	• EZ-Link Sulfo-NHS-LC Biotinylation Kit	Complete kit with an optimized procedure for labeling a protein and determining how much biotin has been attached	22.4	Yes	No	
	• EZ-Link Micro Sulfo-NHS-LC Biotinylation Kit					
NHS-PEG solid-phase biotinylation kits	• EZ-Link NHS-PEG Solid-Phase Biotinylation Kit—1 mL Column • EZ-Link NHS-PEG Solid-Phase Biotinylation Kit	A column-based approach to antibody biotinylation and clean-up	NA	NA	NA	

Carbonyl biotinylation reagents						
Product categories	Product	Product features	Spacer arm (Å)	Water-soluble?	Cleavable?	Membrane-permeant?
Hydrazide-biotin reagents	• EZ-Link Hydrazide-Biotin	Short and medium-length reagents for labeling glycoproteins and other carbohydrate-containing compounds having oxidizable sugars or aldehydes	15.7	No	No	Yes
	• EZ-Link Biotin-LC-Hydrazide		24.7	No	No	Yes
Hydrazide-PEG <sub>4</sub> -biotin	• EZ-Link Hydrazide-PEG <sub>4</sub> -Biotin	A carbohydrate biotinylation reagent with a long hydrophilic spacer arm	31.3	Yes	No	No

Sulfhydryl biotinylation reagents						
Product categories	Product	Product features	Spacer arm (Å)	Water-soluble?	Cleavable?	Membrane-permeant?
Maleimide-PEG <sub>11</sub> -biotin	• EZ-Link Maleimide PEG <sub>11</sub> -Biotin	A column-based approach to antibody biotinylation and clean-up	59.1	Yes	No	No
Maleimide-PEG <sub>2</sub> -biotin	• EZ-Link Maleimide PEG <sub>2</sub> -Biotin	A simple, water-soluble biotinylation reagent specific for sulfhydryls	29.1	Yes	No	No
Maleimide solid-phase biotinylation kits	• EZ-Link Maleimide-PEG Solid-Phase Biotinylation Kit—Prepacked Column	A column-based approach to antibody biotinylation and clean-up	NA	NA	NA	NA
	• EZ-Link Maleimide-PEG Solid-Phase Biotinylation Kit—Mini-Spin Columns					

Carboxyl biotinylation reagents						
Product categories	Product	Product features	Spacer arm (Å)	Water-soluble?	Cleavable?	Membrane-permeant?
Amine-PEG-biotin reagents	• EZ-Link Amine-PEG <sub>2</sub> -Biotin	Short, hydrophilic reagents for biotinylation of carboxyls via EDC	20.4	Yes	No	No
	• EZ-Link Amine-PEG <sub>3</sub> -Biotin		22.9	Yes	No	No
	• EZ-Link Amine-PEG <sub>11</sub> -Biotin	Long, hydrophilic reagent for biotinylation of carboxyls via EDC	53.2	Yes	No	No

For information on all of our biotinylation reagents, go to [thermofisher.com/biotinylation](https://www.thermofisher.com/biotinylation)

## Biotin quantitation kits

These simple and accurate 4-hydroxyazobenzene-2-carboxylic acid (HABA) dye assay and fluorescence detection kits are used to measure the amount of biotin attached to proteins or other biomolecules following biotinylation, allowing you to determine the molar ratio of biotin to labeled antibody molecule in an easy, efficient manner (Table 13).

Table 13. Biotin quantitation kits and reagents.

Product	Product features	Detection method
Pierce Fluorescence Biotin Quantitation Kit	Fluorescent assay reagent and HABA dye for rapid and sensitive determination of biotinylation levels in labeled antibodies and other proteins with minimal sample quantities	Fluorescence
Pierce Biotin Quantitation Kit	A simple, colorimetric method for determining biotinylation efficiency	Colorimetric
Pierce HABA	HABA dye reagent and biotinylated protein standard for convenient colorimetric determination of biotinylation levels in antibodies and other proteins	Colorimetric

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)



## Pierce premium grade reagents

Our Thermo Scientific™ Pierce™ premium grade reagents are high-quality formulations of selected chemical modification reagents, specially characterized for applications where product integrity and risk minimization are paramount. The current menu of reagents offered in this premium format includes several popular crosslinkers and biotinylation reagents, and a reducing agent. The reagents are ideal for projects that require long-term, consistent performance and an extra level of troubleshooting support afforded by robust chemical characterization and documentation. Available Pierce premium grade reagents include:

- Pierce Premium Grade Sulfo-NHS
- Pierce Premium Grade EDC
- Pierce Premium Grade Sulfo-NHS-LC-Biotin
- Pierce Premium Grade Sulfo-NHS-SS-Biotin
- Pierce Premium Grade DSP
- Pierce Premium Grade BS<sup>3</sup>
- Pierce Premium Grade Sulfo-SMCC
- Pierce Premium Grade SPDP
- Pierce Premium Grade TCEP-HCl

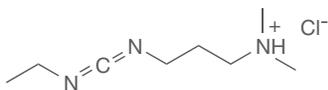
### Highlights:

- **High quality**—identity and purity confirmed by several tests, including quantitative NMR
- **Product integrity**—enhanced level of testing and characterization compared to standard grade (Table 14)
- **Lot retention**—ample supply of past lots retained to help ensure future process testing
- **Change management**—Change Control Notification (CCN) service
- **Consistent manufacturing**—batch-specific manufacturing documentation review

Table 14. Comparison of premium grade and standard grade crosslinking reagents.

Test parameter	Test method	Premium grade	Standard grade
Purity	Quantitative NMR using an internal standard	Yes	Yes
Visual	Color assessment	Yes	Yes
Solubility	Sample dissolves at specified concentration in deionized water to yield a clear, colorless solution	Yes	Yes
Identity	Infrared spectroscopy	Yes	Yes
Mass identity	Mass spectrometry	Yes	NA
Water content	Karl Fischer titration	Yes	NA
Trace metals	Inductively coupled plasma mass spectrometry (ICP-MS)	Yes	NA
Elemental analysis	Combustion analysis (values reported for C, H, N, O, and S)	Yes	NA
Residual solvent analysis	Headspace gas chromatography	Upon request	NA

## Pierce Premium Grade EDC crosslinker



### EDC

1-Ethyl-3-(3-dimethylaminopropyl)carbodiimide • HCl  
MW 191.70  
Spacer arm 0.0 Å

Thermo Scientific™ Pierce™ Premium Grade EDC (1-ethyl-3-(3-dimethylaminopropyl)carbodiimide•HCl) crosslinker is our highest-quality formulation of this popular carbodiimide crosslinker, specially characterized for applications where product integrity and risk minimization are paramount.

### Features of Pierce Premium Grade EDC crosslinker:

- Carbodiimide crosslinker for conjugation between carboxylates and amines
- Identity and purity confirmed by several tests for exceptional quality
- High level of testing and characterization compared to standard grade for high product integrity
- Ample supply of past lots retained to help ensure future process testing
- Change control notification (CCN) service
- Batch-specific manufacturing documentation review

Compared to the standard grade product, Pierce Premium Grade EDC crosslinker provides more clearly defined quality and product support by including (a) increased analytical testing and product characterization, (b) greater batch-specific information and quality assurance review, (c) extensive lot sample retention, and (d) change control notification. EDC is a water-soluble carbodiimide crosslinker that activates carboxyl groups for spontaneous reaction with primary amines, enabling peptide immobilization and hapten-carrier protein conjugation. In the presence of sulfo-NHS, EDC can be used to convert carboxyl groups to amine-reactive sulfo-NHS esters. This is accomplished by mixing the EDC with a carboxyl containing molecule and adding sulfo-NHS.

## Specifications of Pierce Premium Grade

### EDC crosslinker:

We manufacture crosslinking reagents to offer high overall product integrity, consistency, and performance for the intended research applications.

- Visual: white crystalline powder, free of foreign matter
- Identity: IR scan shows only peaks characteristic of EDC
- Water content: Karl Fischer titration, ≤2%
- Trace metals: all metals, including lead, ≤50 ppm
- Elemental analysis: reported values for C, H, and N

### Applications:

- Conjugate carboxyl and amino groups among peptides and proteins
- Couple haptens to immunogenic carrier proteins (e.g., attach a peptide to KLH)
- Immobilize peptide antigens to affinity-purify antibodies
- Create NHS-activated, amine-reactive labeling compounds
- Crosslink proteins to carboxyl-coated beads or surfaces
- Activate nanoparticles with amine-reactive, sulfo-NHS esters

For information on all of the Pierce Premium Grade Reagents, go to

[thermofisher.com/premiumreagents](https://thermofisher.com/premiumreagents)

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Fluorescent antibody and protein labeling kits

We offer a number of antibody and protein labeling kits for the direct attachment of a broad range of fluorescent dyes and labels including Alexa Fluor dyes, DyLight dyes, Qdot labels, R-phycoerythrin, APC tandem dyes, and biotin, at scales from less than 10 µg up to 1 mg of IgG antibody. Direct labeling of antibodies allows you to use more than one same-species antibody in a single staining experiment. You can use traditional labeling chemistries optimized for your application or site-specific labeling using Invitrogen™ SiteClick™ or Zenon™ labeling technology.

### Kits for fluorescent site-specific antibody labeling

**SiteClick labeling technology** specifically attaches the label to the heavy chains of an IgG antibody helping to ensure that the antigen-binding domains remain available for binding to your antigen target. This site selectivity is achieved by targeting the carbohydrate domains present on essentially all IgG antibodies regardless of isotype and host species.

**Zenon labeling technology** provides a versatile, easy-to-use system for labeling human, mouse IgG1, IgG2a, and IgG2b antibodies and rabbit IgG antibodies. Zenon fragments are specifically designed to target and bind to the Fc portion of the primary antibody only, giving a rapid, noncovalent method of quickly labeling small quantities of primary antibody. Zenon labeling technology is simple, efficient, and scalable; the entire labeling procedure takes only 10 minutes.

### Kits for conventional antibody and protein labeling

**Invitrogen™ APEX™ antibody labeling kits** provide an efficient, convenient, and reliable method for covalently attaching fluorescent labels to very small amounts of IgG antibody in the presence of any contaminants, including stabilizing proteins and amine-containing buffers.

**Thermo Scientific™ DyLight™ antibody labeling kits** are designed for fast, efficient labeling of antibodies with DyLight dyes.

**Invitrogen™ Alexa Fluor™ antibody labeling kits** enable direct labeling of monoclonal and polyclonal antibodies.

**Invitrogen™ microscale protein labeling kits** conveniently label small amounts of purified protein or antibodies (MW 12,000–150,000) with a bright, photostable Alexa Fluor dye or biotin.

Our **protein labeling kits** provide a nearly effortless way to stably label larger amounts of IgG antibody with a fluorescent dye or hapten in two hours.

**Invitrogen™ Alexa Fluor™ SAIVI™ Rapid Antibody Labeling Kits** are uniquely designed to provide the rigorous requirements to control the degree of labeling (DOL) using our near-infrared (NIR) Alexa Fluor dyes.

For more information on all of our antibody and protein labeling kits, go to [thermofisher.com/antibodylabeling](https://www.thermofisher.com/antibodylabeling)

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## SiteClick antibody labeling system

The Invitrogen™ SiteClick™ antibody labeling system allows simple and gentle site-selective attachment of detection molecules to heavy-chain *N*-linked glycans far from the antigen-binding domain, enabling excellent reproducibility from labeling to labeling and from antibody to antibody (Figures 17 and 18). A number of different detection molecules can be site-selectively attached to the heavy-chain glycans, including phycobiliproteins (e.g., RPE), Qdot probes, fluorescent dyes, metal-chelating compounds, and other small molecules such as biotin, allowing multiplex analysis with antibodies from the same species.

### Highlights:

- Highly efficient, site-specific, reproducible labeling chemistry that preserves antigen-binding domains
- Qdot labels for fluorescence microscopy and flow cytometry (Figure 19)
- RPE labels for flow cytometry
- Alexa Fluor labels using Invitrogen™ Click-iT™ Alexa Fluor™ DIBO Alkynes

### How the SiteClick antibody labeling system works

In general, IgG antibodies contain two *N*-linked glycans attached to specific conserved asparagine residues located in the antibody heavy-chain Fc domain. These sugar chains are structurally quite homogeneous, and the terminal sequences of the glycan branches are highly consistent. Most of the antibody glycan branches terminate with galactose-*N*-acetylglucosamine (Gal-GlcNAc) or with *N*-acetylglucosamine (GlcNAc). Removal of the terminal Gal residue with beta-galactosidase unmasks the majority of terminal GlcNAc labeling sites for the subsequent enzymatic beta-galactosyl transferase (GalT) reaction. After cleavage of terminal Gal residues by beta-galactosidase, each *N*-linked glycan will contain, on average, two terminal GlcNAc residues per heavy chain (four terminal GlcNAc per antibody), which can be labeled with a DIBO alkyne (Figure 17).

The SiteClick method is compatible with antibodies from a number of different species including, but not limited to, human, rabbit, mouse, rat, goat, hamster, and chicken. Additionally, SiteClick labeling is effective with several antibody classes such as IgG, IgM, and IgY; note that chicken IgY antibodies have six heavy-chain glycans instead of two, and therefore can be labeled to a higher extent.

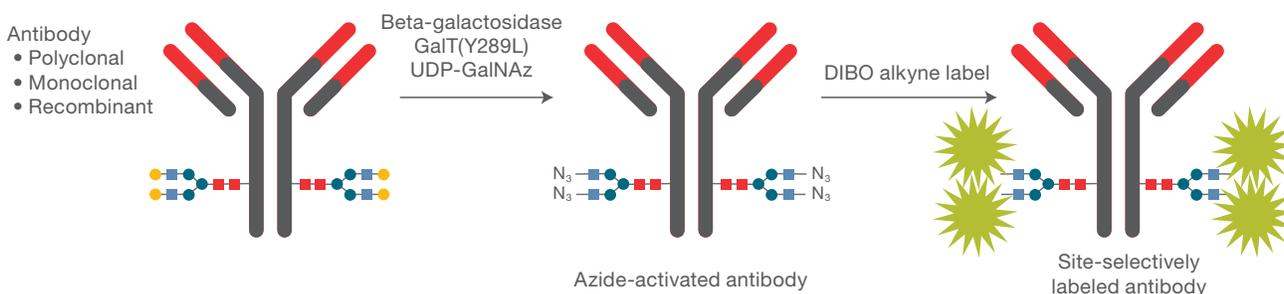


Figure 17. The SiteClick antibody labeling system. The first step in the SiteClick antibody labeling process involves removal of terminal galactose residues from the heavy-chain *N*-linked glycans using beta-galactosidase, exposing essentially all possible modifiable GlcNAc residues. Second, the free terminal GlcNAc residues are activated with azide tags by enzymatic attachment of GalNAz to the terminal GlcNAc residues using the GalT(Y289L) enzyme. In the third step, the azide residues are reacted with the dibenzocyclooctyne (DIBO)-functionalized probe of choice (e.g., Alexa Fluor 488 DIBO Alkyne). The average degree of labeling is 3–3.5 labels per antibody.

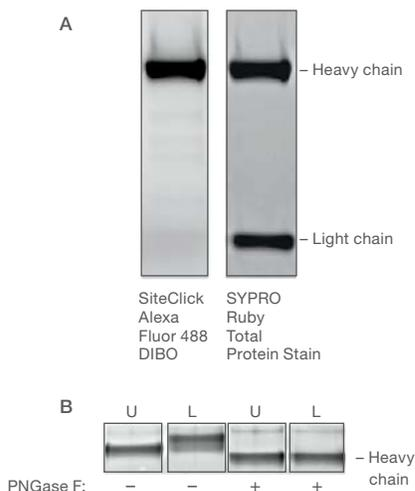


Figure 18. Site selectivity of SiteClick antibody labeling. **(A)** A beta-tubulin monoclonal antibody was azide-activated using the GalT(Y289L) enzyme, labeled with DIBO-functionalized Alexa Fluor 488 dye, and then analyzed by gel electrophoresis (left panel). Once imaged, the same gel was post-stained with Invitrogen™ SYPRO™ Ruby Total Protein Gel Stain (right panel), showing that only the heavy chain of the antibody is click-labeled with the Alexa Fluor 488 DIBO alkyne. **(B)** Anti-beta-tubulin mouse monoclonal antibodies were labeled with a small-molecule DIBO-PET chelating agent (L) or left unlabeled (U) (left two lanes) and then post-treated with PNGase F, which selectively cleaves *N*-linked glycans from asparagine residues (right two lanes). After treatment with PNGase F, both chelator-labeled (L) and unlabeled (U) species are shifted to the same lower molecular weight, confirming the site-selective labeling of heavy-chain *N*-linked glycans.

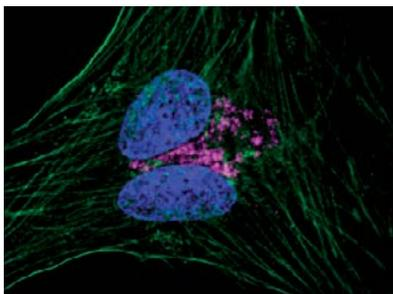


Figure 19. Immunocytochemistry with antibodies labeled by the SiteClick method. HeLa cells were fixed, permeabilized, and incubated with 10 nM Qdot 655 nanocrystal-conjugated anti-golgin 97 antibody (magenta). This conjugate was generated using the SiteClick Qdot 655 Antibody Labeling Kit and mouse monoclonal anti-golgin 97 antibody (CDF3). After antibody incubation, the cells were counterstained with Invitrogen™ NucBlue™ Live (blue) and Invitrogen™ ActinGreen™ 488 (green) ReadyProbes™ reagents prior to imaging.

## Fluorescent dyes

As one of the world's leading developers of fluorescence-based technology, we offer many fluorescent labels for primary antibodies, secondary antibodies, anti-dye and anti-hapten antibodies, and streptavidin, including proprietary labels such as Alexa Fluor, DyLight, Texas Red-X, and Pacific Blue dyes.

### Alexa Fluor dyes

The Alexa Fluor dyes (Table 15) produce exceptionally bright and photostable conjugates. These dyes are available in several reactive forms for labeling biopolymers and derivatizing low-molecular weight molecules. All Alexa Fluor dyes are available as an amine-reactive succinimidyl ester, and many are available in thiol, aldehyde, and carboxylic acid reactive forms.

#### Highlights:

- Compatible with all common excitation sources and instruments
- Bright and unusually photostable fluorescence of their bioconjugates
- Good water solubility, which makes the reactive dyes easy to conjugate and the conjugates resistant to precipitation and aggregation
- Insensitivity of their absorption and emission spectra to pH over a broad range
- Well-differentiated spectra, providing many options for multicolor detection and fluorescence resonance energy transfer (FRET); spans the near-UV, visible, and near-IR spectra (Figure 20)
- High quantum yields and long fluorescence lifetimes
- Extremely high FRET efficiency, with calculated  $R_0$  values of up to 84 Å between pairs of Alexa Fluor dyes and up to 77 Å between Alexa Fluor dyes and some nonfluorescent quenchers

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

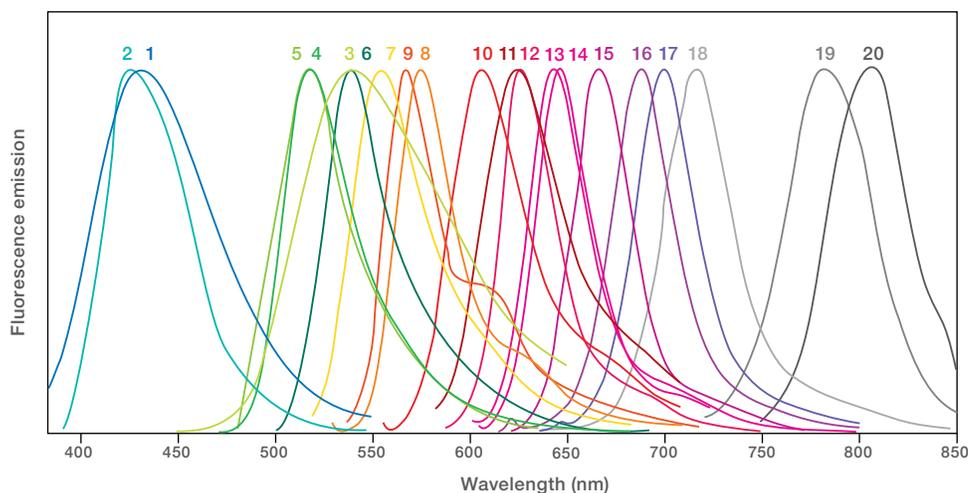
Table 15. Properties of Alexa Fluor dyes.

Alexa Fluor dye	Absorption max. (nm)	Emission max. (nm)	Emission color*	Extinction coefficient**
Alexa Fluor 350	346	442	Blue	19,000
Alexa Fluor 405	401	421	Blue	34,000
Alexa Fluor 430	433	541	Green/yellow	16,000
Alexa Fluor 488	496	519	Green	71,000
Alexa Fluor 532	532	553	Yellow	81,000
Alexa Fluor 546	556	573	Orange	104,000
Alexa Fluor 555	555	565	Orange	150,000
Alexa Fluor 568	578	603	Orange/red	91,000
Alexa Fluor 594	590	617	Red	73,000
Alexa Fluor 610	612	628	Red	138,000
Alexa Fluor 633	632	647	Far-red	239,000
Alexa Fluor 635	633	647	Far-red	140,000
Alexa Fluor 647	650	665	Near-IR†	239,000
Alexa Fluor 660	663	690	Near-IR†	132,000
Alexa Fluor 680	679	702	Near-IR†	184,000
Alexa Fluor 700	702	723	Near-IR†	192,000
Alexa Fluor 750	749	775	Near-IR†	240,000
Alexa Fluor 790	784	814	Near-IR†	270,000

\* Typical emission color seen through the eyepiece of a conventional fluorescence microscope with appropriate filters.

\*\* Extinction coefficient at  $I_{max}$  in  $M^{-1} cm^{-1}$ .

† Human vision is insensitive to light beyond ~650 nm; it is not possible to directly view near-IR fluorescent dyes.



1. Alexa Fluor 350 dye
2. Alexa Fluor 405 dye
3. Alexa Fluor 430 dye
4. Alexa Fluor 488 dye
5. Alexa Fluor 500 dye
6. Alexa Fluor 514 dye
7. Alexa Fluor 532 dye
8. Alexa Fluor 546 dye
9. Alexa Fluor 555 dye
10. Alexa Fluor 568 dye
11. Alexa Fluor 594 dye
12. Alexa Fluor 610 dye
13. Alexa Fluor 633 dye
14. Alexa Fluor 635 dye
15. Alexa Fluor 647 dye
16. Alexa Fluor 660 dye
17. Alexa Fluor 680 dye
18. Alexa Fluor 700 dye
19. Alexa Fluor 750 dye
20. Alexa Fluor 790 dye

Figure 20. Emission spectra for Alexa Fluor dyes.

## DyLight PEGylated dyes

The Thermo Scientific™ DyLight™ PEGylated dyes are derivatives of our high-performance DyLight dyes that can be used to fluorescently label antibodies and other proteins (Tables 16 and 17). Conjugates made with DyLight PEGylated dyes can be used as molecular probes for cellular imaging and other fluorescence detection methods. DyLight PEGylated dyes contain 2 to 4 nontoxic polyethylene glycol (PEG) chains. The PEG chains enhance fluorescence, reduce nonspecific binding of conjugates, improve solubility of the dyes and labeled molecules in aqueous solution, aid in cell permeability, and improve retention, especially in tumors [1]. The near-infrared (NIR) to far-red fluorescence properties of some of the PEGylated dyes such as DyLight 680, 755, and 800 dyes make them useful in biological, chemical, and pharmaceutical applications, including *in vivo* imaging (Figure 21).

The DyLight PEGylated amine- and sulfhydryl-reactive dyes provide superior fluorescence characteristics (e.g., brightness, photostability, pH insensitivity, and water solubility). In addition, the high fluorescence intensity of DyLight PEGylated dyes provides outstanding sensitivity and requires less conjugate in most applications.

### Highlights:

- **Soluble**—PEGylated dyes and labeled molecules improve solubility in aqueous solution, aid in cell permeability, and improve retention
- **High fluorescence intensity**—significantly brighter fluorescence than many other dyes
- **Efficient labeling methods**—well-characterized chemistry and optimized protocols enable reliable, high-quality labeling
- **Reduced nonspecific binding**

**Table 16. Properties of the PEGylated DyLight NHS-ester fluorescent dyes.**

DyLight™ Fluor	Ex/Em*	$\epsilon^{**}$	MW	Spectrally similar dyes
DyLight 550-2xPEG	557/571	150,000	1,102	Alexa Fluor 555, Cy3, DyLight 550, CF555
DyLight 650-4xPEG	658/681	250,000	1,425	Alexa Fluor 647, Cy5, DyLight 650, DyLight 650, CF647
DyLight 680-4xPEG	684/706	180,000	1,729	Alexa Fluor 680, Cy5.5, DyLight 680, CF680, IRDye 680
DyLight 755-4xPEG	754/777	220,000	1,451	Alexa Fluor 750, DyLight 755, CF750
DyLight 800-4xPEG	783/797	270,000	1,685	Alexa Fluor 790, Cy7, DyLight 800, CF790, IRDye 800

\* Excitation and emission maxima in nanometers.

\*\* Molar extinction coefficient ( $M^{-1} cm^{-1}$ ).

**Table 17. Properties of the PEGylated DyLight maleimide fluorescent dyes.**

DyLight Fluor	Ex/Em*	$\epsilon^{**}$	MW	Spectrally similar dyes
DyLight 550-2xPEG	557/571	150,000	1,127	Alexa Fluor 555, Cy3, DyLight 550, CF555
DyLight 650-4xPEG	656/675	250,000	1,450	Alexa Fluor 647, Cy5, DyLight 650, CF647
DyLight 680-4xPEG	684/706	180,000	1,754	Alexa Fluor 680, Cy5.5, DyLight 680, CF680, IRDye 680
DyLight 755-4xPEG	757/778	220,000	1,476	Alexa Fluor 750, DyLight 755, CF750
DyLight 800-4xPEG	784/798	270,000	1,710	Alexa Fluor 790, Cy7, DyLight 800, CF790, IRDye 800

\* Excitation and emission maxima in nanometers.

\*\* Molar extinction coefficient ( $M^{-1} cm^{-1}$ ).

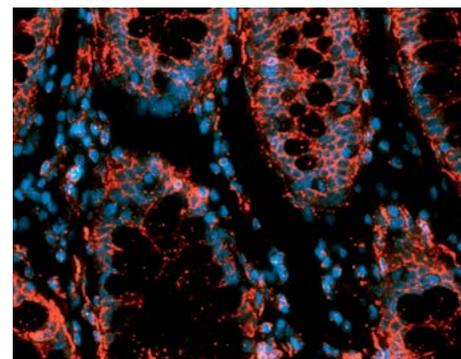


Figure 21. DyLight 550-2xPEG dye performance in IHC. Paraffin-embedded normal colon tissue was fluorescently probed for the EGF receptor with rabbit anti-EGFR antibody and DyLight 550-2xPEG dye-conjugated goat anti-rabbit secondary antibody. Nuclei were also counterstained with Hoechst 33342 stain.

### Reference

1. Knop K et al. (2010) Poly(ethylene glycol) in drug delivery: Pros and cons as well as potential alternatives. *Angew Chem Int Ed* 49:6288–6308.

## Enzyme labeling kits

### EZ-Link HRP enzyme labeling kits and reagents

Thermo Scientific™ EZ-Link™ preactivated HRP makes it easy to convert almost any protein to a detection reagent. HRP is the enzyme most commonly used for immunoassay detection systems. This enzyme catalyzes reactions with substrates to form soluble color responses or colored precipitates, or to generate the chemical emission of light (chemiluminescence). Enzyme conjugates make stable assay reagents and can be stored for long periods at  $-20^{\circ}\text{C}$ .

### Pierce Maleimide Activated Horseradish Peroxidase (HRP)

Thermo Scientific™ EZ-Link™ Maleimide Activated Horseradish Peroxidase is for preparation of HRP conjugates with proteins, peptides, or other ligands that contain sulfhydryl groups, such as reduced cysteines.

#### Highlights:

- **Activated HRP**—HRP modified with maleimide groups for conjugation to sulfhydryl molecules
- **Sulfhydryl-reactive**—maleimide groups conjugate to reduced thiols ( $-\text{SH}$ ), as in the side chain of cysteine residues
- **High-activity HRP**—enzyme activity is greater than 240 units/mg; lyophilized, activated enzyme is stable for at least 12 months at  $4^{\circ}\text{C}$

### EZ-Link Plus Activated Peroxidase

Thermo Scientific™ EZ-Link™ Plus Activated Peroxidase is an amine-reactive form of HRP that provides coupling efficiencies of greater than 95% with antibodies and other proteins.

#### Highlights:

- **Activated HRP**—periodate-treated, aldehyde-activated horseradish peroxidase, ready for conjugation to antibodies and other proteins at sites of primary amines (e.g., lysines)
- **Permanent conjugation**—reacts efficiently (95%) with primary amines to form covalent amide bonds upon treatment with sodium cyanoborohydride (included in kit)
- **High-activity HRP**—enzyme activity is 120 to 200 units/mg; lyophilized, activated enzyme is stable for at least 12 months at  $-20^{\circ}\text{C}$
- **Convenient quantities**—each 1 mg quantity of activated enzyme is sufficient for reaction with 1 mg of IgG to produce about 0.5 mL of conjugate
- **Customizable**—vary the molar ratios, reaction buffer and pH, and other parameters to achieve conjugates with different levels of HRP incorporation and activity

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Pierce Horseradish Peroxidase

Thermo Scientific™ Pierce™ Horseradish Peroxidase is purified HRP enzyme for use in activity assays and conjugation to antibodies for ELISA, western blot, and immunohistochemistry applications.

### Highlights:

- **Purified form**—lyophilized salt-free powder; ready to dissolve and use
- **High specific activity**—typically greater than 300 units/mg (lot-specific value reported)
- **Compared to AP**—HRP is smaller (40 kDa) than alkaline phosphatase (AP; 140 kDa) and has higher specific enzyme activity than both AP and beta-galactosidase (beta-Gal)
- **Many options**—numerous substrate solutions and assay techniques are available for HRP

This purified HRP is supplied lyophilized as a salt-free powder for reconstitution and use in protein research methods. The main application for HRP in molecular biology and protein research is as a reporter system for immunoassays and other probe-based assay techniques such as ELISA, western blotting, EMSA, and Southern blotting. The enzyme is usually conjugated to specific secondary antibodies or streptavidin, and its activity is detected with a color-forming (or light-generating) substrate.

One unit catalyzes the production of 1 mg of purpurogallin from pyrogallol in 20 seconds at 20°C and pH 6.0.

## Pierce Alkaline Phosphatase

This purified calf intestinal alkaline phosphatase (CIP) is supplied in Tris buffer and 50% glycerol. It can be conjugated to specific primary or secondary antibodies and its activity detected with a color-forming (or light-generating) substrate.

### Highlights:

- **Purified form**—ready to dilute and conjugate
- **Concentrated**—approximately 20 mg/mL (lot-specific value reported)
- **High specific activity**—typically greater than 1,600 units/mg (lot-specific value reported)
- **Tris-buffered solution**—supplied in 5 mM Tris, 5 mM magnesium chloride, and 0.1 mM zinc chloride, pH ~7.0 in 50% glycerol

One unit equals the amount of protein required to hydrolyze one micromole of *p*-nitrophenyl phosphate (PNPP) per minute at 25°C in a glycine buffer, pH 9.6.



# Nonspecific binding

We offer an extensive line of high-quality buffers and detergents designed for use in a variety of immunoassay procedures. Our detergents undergo extensive processing to minimize levels of oxidants and carbonyls.

## Blocking buffers

Designed to reduce nonspecific binding, Thermo Scientific™ Pierce™ blocking buffers can be used in protein detection systems such as ELISA, western blotting, lateral flow devices, and protein arrays. We have a variety of blocking buffer formulations with different blocking and buffering agents (Table 18).

Table 18. Choose the right blocking buffer for your experiment.

Product	ELISA	WB	IHC	Nucleic acid detection
SuperBlock Blocking Buffers	✓	✓	✓*	
StartingBlock Blocking Buffers	✓	✓	✓	
Blocker Casein Blocking Buffers	✓	✓	✓	✓
Blocker BSA Blocking Buffers	✓	✓	✓	✓
Blocker BLOTTO Blocking Buffer	✓	✓	✓	
SEA BLOCK Blocking Buffer	✓	✓		
Pierce Protein-Free Blocking Buffers	✓	✓		
I-Block™ Protein-Based Blocking Reagent	✓	✓		

\* Ideal for IHC with antibody or avidin/biotin probes (blocker is serum and biotin-free).

## SuperBlock Blocking Buffer

Highlights of Thermo Scientific™ SuperBlock™ Blocking Buffer include:

- **Fast**—block membranes typically in 5 to 10 minutes and ELISA plates in 2 minutes
- **Flexible**—biotin-free for use with streptavidin systems
- **Convenient**—optimized PBS or TBS solutions available in multiple package sizes (100 mL, 1 L, 5 L, and dry-blend pouches)
- **Low background**—the nonserum protein solution yields a high signal-to-noise ratio
- **Stable**—store buffer at 4°C for 1 year; store blocked plates dry for up to 12 months

## StartingBlock Blocking Buffer

Highlights of Thermo Scientific™ StartingBlock™ Blocking Buffer include:

- **Compatible with many detection systems**—western blot, ELISA, and IHC with antibody or avidin/biotin probes (blocker is serum- and biotin-free)
- **Short blocking times**—less than 15 minutes for nitrocellulose or PVDF membranes, even faster for polystyrene microplate wells
- **Strip and reprobe without reblocking**—blots stay blocked even after stripping with Thermo Scientific™ Restore™ Stripping Buffer
- **High signal-to-noise ratios**—biotin- and serum protein-free formulation contributes to achieving signal-to-noise ratios from 10:1 to 20:1
- **Convenient formats**—ready-to-use 1X formulations in PBS and TBS, with and without Tween 20 detergent

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Blocker Casein

Highlights of Thermo Scientific™ Blocker™ Casein include:

- **Purified casein**—single-protein blocking buffer provides fewer chances of cross-reaction with assay components than serum or milk solutions
- **Easy to use**—1% solutions of Hammarsten-grade casein are ready to use; can be diluted further as needed
- **Flexible**—available in PBS and TBS formulations to suit a variety of applications
- **Safe**—stable, thimerosal-free formulations

## Blocker BSA

Highlights of Thermo Scientific™ Blocker™ BSA include:

- **Purified protein**—10% solution of high-quality bovine serum albumin; a single purified protein provides fewer chances of cross-reaction with assay components than serum or milk solutions
- **Convenient**—concentrated formulation helps save storage space and can be diluted easily, enabling optimal blocking results for specific applications
- **Easy to use**—no waiting for powder to dissolve with this ready-to-dilute liquid concentrate

## Blocker BLOTTO Blocking Buffer

Highlights of Thermo Scientific™ Blocker™ BLOTTO Blocking Buffer include:

- **Popular**—nonfat milk has been used for many years in a variety of protein methods, although it is not recommended for avidin-based techniques because it contains some endogenous biotin
- **Convenient**—supplied as a ready-to-use 1X TBS solution; can be diluted as needed
- **Easy to use**—formulated with anti-foaming agent and thimerosal-free preservative
- **Flexible**—may be used for multiple applications, including as a diluent for antibodies

## SEA BLOCK Blocking Buffer

Highlights of Thermo Scientific™ SEA BLOCK Blocking Buffer include:

- **Nonmammalian**—made from steelhead salmon serum; since protein in the buffer is of nonmammalian origin, the risk of background caused by nonspecific interactions is minimized
- **Convenient**—filtered and stabilized in PBS for compatibility with most assay systems
- **Easy to use**—can be used as supplied or diluted up to 10-fold as needed
- **Flexible**—may be used for many different applications, including as a diluent for antibodies

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Pierce Protein-Free Blocking Buffer

Highlights of Thermo Scientific™ Pierce™ Protein-Free Blocking Buffer include:

- **Protein-free blocker**—minimizes or eliminates cross-reactivity associated with protein-based blocking buffers
- **Application-compatible**—effective in all kinds of protein detection systems, including western blots (membranes), ELISA (microplates), and arrays (coated glass slides)
- **Streptavidin-friendly**—absolutely free of biotin; no possible interference with avidin–biotin detection systems
- **High-performance**—optimized and validated in many protein methods to provide high signal-to-noise ratios (i.e., no quenching of specific binding and signal but eliminating nonspecific binding and background)
- **Flexible**—available in PBS and TBS formulations to suit a variety of applications

## Pierce Fast Blocking Buffer

Highlights of the Thermo Scientific™ Pierce™ Fast Blocking Buffer include:

- **Fast**—use to shorten the typical western blot development by over 2 hours
- **Simple**—optimized protocol makes western blot analysis easier than ever
- **Low background**—provides results comparable to classic western blotting buffers

## Pierce Clear Milk Blocking Buffer (10X)

Highlights of Thermo Scientific™ Pierce™ Clear Milk Blocking Buffer include:

- **Excellent stability**—stable for 1 year stored at 4°C, unlike typical homemade milk buffers
- **Convenient**—concentrated formulation saves storage space and can be diluted easily to obtain optimal blocking results for specific applications
- **Easy to use**—no waiting for powdered milk to dissolve with this ready-to-dilute solution
- **Popular**—nonfat milk has been used for many years in a variety of protein methods, although it is not recommended for avidin-based techniques because it contains some endogenous biotin

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Wash buffers

### BupH dry-blend buffer packs and Pierce concentrate buffers

Thermo Scientific™ BupH™ packs are preblended and premeasured dry mixtures of commonly needed buffers that are easy to prepare; simply empty the contents of a foil envelope pack into a beaker, add ultrapure water, and stir to dissolve. The packs eliminate weighing time and tedious pH adjustments. BupH dry-blend buffer packs and Pierce concentrate buffers are offered for use in a variety of laboratory techniques (Table 19).

#### Highlights:

- **Convenient**—dissolve contents of one envelope in water and the buffer is ready to use
- **Helps save time and trouble**—no weighing, no pH adjustment, no need to stock individual components, and no need to make and store large volumes of stock solution in advance of daily needs
- **Long shelf life**—stocking and storage as dry packs minimizes concerns about long-term stability of stock solutions
- **Eliminate variables**—our quality control helps ensure that every pack will yield consistent buffer

Thermo Scientific™ Pierce™ concentrate buffers are ready to use without having to reconstitute with ultrapure water. The buffers are designed for use in dialysis, crosslinking, enzyme assays, ELISAs, immunohistochemistry, protein plate-coating, biotinylation, and other applications.

#### Highlights:

- **Easy to use**—no packets to open and no powder to dissolve
- **Helps increase accuracy**—minimizes the possibility of powder remaining in a packet
- **Helps save time**—20X concentration minimizes time spent waiting for powder to dissolve
- **Helps save space**—storage as concentrated stock minimizes bench space needed for solutions

Table 19. BupH dry-blend buffer packs and Pierce concentrate buffers.

BupH packs		
Description	Applications	Formulation after reconstitution
Phosphate-buffered saline (PBS)	Crosslinking and biotinylation requiring amine-free buffer	0.1 M sodium phosphate, 0.15 M NaCl, pH 7.2
Modified Dulbecco's PBS	Wash buffers and antibody diluents for ELISA, western blotting, and other immunoassays	8 mM sodium phosphate, 2 mM potassium phosphate, 0.14 M NaCl, 10 mM potassium chloride, pH 7.4
Tris-buffered saline	Wash buffers and antibody diluents for ELISA, western blotting, and other immunoassays	25 mM Tris, 0.15 M NaCl, pH 7.2
Pierce concentrate buffers		
Description	Applications	Formulation after reconstitution
Pierce 20X Modified Dulbecco's PBS Buffer	Used for wash buffers and antibody diluents in applications such as ELISA, western blotting, and other immunoassays	8 mM sodium phosphate, 2 mM potassium phosphate, 0.14 M NaCl, 100 mM KCl, pH 7.4
Pierce 20X Modified Dulbecco's PBS Tween 20 Buffer	A wash buffer for ELISA, western, and other immunoassays, as well as a blocking buffer for plate-based assays	8 mM sodium phosphate, 2 mM potassium phosphate, 0.14 M NaCl, 100 mM KCl, 0.05% Tween 20, pH 7.4
Pierce 20X Phosphate-Buffered Saline	Its ionic strength makes it ideal for crosslinking and biotinylation requiring amine-free buffer	0.01 M sodium phosphate, 0.15 M NaCl, pH 7.5
Pierce 20X PBS Tween 20 Buffer	A wash buffer for ELISA, western, and other immunoassays, as well as a blocking buffer for plate-based assays	0.01 M sodium phosphate, 0.15 M NaCl, 0.05% Tween 20, pH 7.5
Pierce 20X TBS Buffer	Used for wash buffers and antibody diluents in applications such as ELISA, western blotting, and other immunoassays	25 mM Tris, 0.15 M NaCl, pH 7.2
Pierce 20X TBS Tween 20 Buffer	A wash buffer for ELISA, western, and other immunoassays, as well as a blocking buffer for plate-based assays	25 mM Tris, 0.15 M NaCl, 0.05% Tween 20, pH 7.5

# Detergents

## Surfact-Amps Detergent Solutions

Thermo Scientific™ Surfact-Amps™ Detergents are highly purified, precisely diluted (10%) formulations that are ideal for applications or assays that are sensitive to contaminants present in unpurified detergents (Table 20). We test every batch to help ensure that our detergents contain <1.0 µeq/mL peroxides and carbonyls and package them under nitrogen, to prevent oxidization during storage.

### Highlights:

- **Accurate**—precise 10% detergent solution in ultrapure water
- **Easy to use**—solution is simple to dispense and dilute for use
- **Exceptionally pure**—less than 1.0 µeq/mL peroxides and carbonyls
- **Stable**—packaged under inert nitrogen gas in glass ampules or HDPE bottles

Unlike neat detergents, which are extremely viscous, Surfact-Amps 10% Solutions are easy to pipette and accurately dispense. The surfactant solutions are carefully prepared and packaged in nonleaching HDPE bottles under inert nitrogen gas for stability and to minimize the accumulation of peroxides and degradation products. Surfact-Amps Detergent Solutions undergo an extensive multistep purification process to remove deleterious contaminants. This process has been optimized over 20 years and uses purification reagents, phase extraction, extensive washes and other manipulations to significantly reduce the levels of peroxides and carbonyls that can hinder detergent performance and stability and create background effects.

**Table 20. Surfact-Amps Detergents.**

Nonionic detergents	Anionic detergents	Zwitterionic detergents
Triton X-100, 10%	Sodium dodecyl sulfate (lauryl SDS)	CHAPS
Triton X-114, 10%	Sodium dodecyl sulfate (C <sub>12</sub> SDS)	n-Dodecyl-beta-D-maltoside
NP-40, 10%	Sodium dodecyl sulfate (lauryl SDS)	
Brij™-35, 30%	Sodium dodecyl sulfate (lauryl SDS), 20% solution	
Brij-35, 10%	Sodium cholate	
Brij-58, 10%	Sodium deoxycholate	
Tween 20, 10%		
Tween 80, 10%		
Octyl-beta-glucoside		
1-S-Octyl-beta-thioglucoopyranoside (OTG)		

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Protein stabilizers

### Guardian Peroxidase Conjugate Stabilizer/Diluent

Thermo Scientific™ Guardian™ Peroxidase Conjugate Stabilizer/Diluent preserves the functional integrity and activity of HRP-conjugated antibodies and other proteins at very dilute concentrations for long-term storage (Figure 22).

#### Highlights:

- **Preserves HRP activity**—no significant loss of HRP activity over a 6-month period at room temperature (1:1,000 dilution) or 12 months at 4°C
- **Convenient**—store ready-to-use dilutions (1:1,000 to 1:100,000) that maintain enzyme activity in the refrigerator—no aliquoting or freezing necessary
- **Assay-compatible**—simply add your favorite blocking buffer to create the ideal diluent for your HRP-based ELISA system, or store the HRP conjugate as a 1:1,000 stock solution for western blots and dilute in the final assay buffer
- **Helps save money**—less expensive than ordering new HRP conjugates

With Guardian Peroxidase Conjugate Stabilizer/Diluent, typically 1 mg/mL antibody or streptavidin peroxidase conjugates can be diluted as much as 100,000-fold for storage at 4°C. Most ELISA and blotting applications for HRP conjugates require dilution of typical 1 mg/mL stocks at least 1,000-fold (to 1 µg/mL). Sensitive assay systems, such as those using chemiluminescent substrates, often require dilution up to 100,000-fold (10 ng/mL). The Guardian Solution enables these 1X working concentrations of HRP conjugate to be prepared in advance and stored at 4°C for 12 months or at room temperature for 6 months.

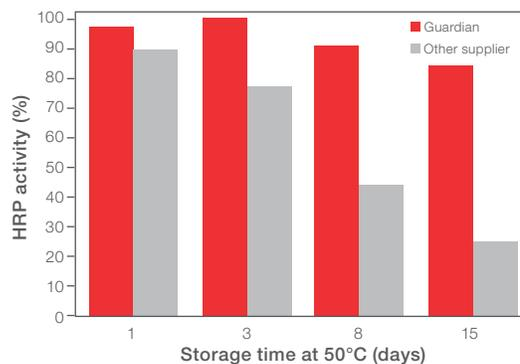
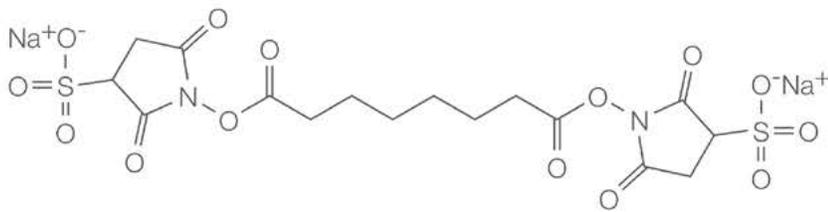


Figure 22. Better HRP conjugate stability with Thermo Scientific™ Guardian™ Peroxidase Conjugate Stabilizer/Diluent. Streptavidin-HRP (1 mg/mL) was diluted 1:1,000 and stored at 50°C in Guardian Stabilizer/Diluent or another supplier's HRP stabilizer. At each time point, the HRP conjugate was diluted 1:5,000,000 with PBS/SuperBlock Blocking Buffer and incubated for 1 hour in a white plate that had been coated with biotinylated BSA. The plate was then washed three times with 200 µL PBST. The Thermo Scientific™ SuperSignal™ ELISA Femto Maximum Sensitivity Substrate (100 µL) was added to the plate and incubated for 1 minute and then measured in a luminometer. Samples at each time point were compared to the control (HRP conjugate that had been stored undiluted). Storing enzymes at 50°C for 2 weeks is equivalent to 12 months at 4°C.



**BS<sup>3</sup>**

Bis(sulfosuccinimidyl)suberate  
MW 572.43  
Spacer arm 11.4 Å



# Detection substrates

Our wide selection of high-quality substrates for use as detection reagents in immunoassays includes chemiluminescent, colorimetric, and chemifluorescent substrates. Substrate choice depends on the platform and sensitivity requirements of the assay.

## Criteria for choosing immunoassay substrates

Optimization of all components is very important for generating the best results with any immunoassay system. Detection substrates differ in their ease-of-use, sensitivity (i.e., lower limit of detection), dynamic range (i.e., logarithmic units of detection), and compatibility with imaging equipment. Table 21 provides some general guidelines for selecting an appropriate substrate. In addition to a wide variety of substrate options, Thermo Fisher can customize your substrate formulations to meet the specific requirements of our immunodiagnostic customers. This includes not only optimizing the performance of a substrate with a customer's platform, but also the price.

Table 21. Comparison of detection substrates.

	Chemiluminescent substrates	Colorimetric substrates	Chemifluorescent substrates
Immunoassay	<ul style="list-style-type: none"> <li>• Highest sensitivity</li> <li>• Rapid signal generation</li> <li>• Largest linear range; enhanced low-end linearity</li> </ul>	<ul style="list-style-type: none"> <li>• Medium/low sensitivity</li> <li>• Slow signal generation</li> <li>• Small linear range; poor low-end linearity</li> <li>• Flexible (stopped, nonstopped, and kinetic assays)</li> </ul>	<ul style="list-style-type: none"> <li>• High sensitivity</li> <li>• Rapid signal generation</li> <li>• Large linear range; enhanced low-end linearity</li> <li>• Flexible (stopped, nonstopped, and kinetic assays)</li> </ul>
Detection equipment	Luminometer	Spectrophotometer	Fluorometer

For more information on all of our ELISA substrates, go to [thermofisher.com/elisasubstrates](https://www.thermofisher.com/elisasubstrates)



## Chemiluminescent substrates

We offer several chemiluminescent substrates for immunoassay development with either HRP, AP, or beta-Gal.

- **Thermo Scientific™ SuperSignal™ ELISA Pico Chemiluminescent Horseradish Peroxidase Substrate** provides excellent performance for a large range of target protein amounts and is easily optimized to detect with greater sensitivity than entry-level colorimetric substrates. Rapid signal generation with 5- to 30-minute signal stability depending on HRP concentration.
- **Thermo Scientific™ SuperSignal™ ELISA Femto Maximum Sensitivity Horseradish Peroxidase Substrate** is one of the most sensitive substrates available for ELISA applications. When properly optimized, the lower detection limit is 1 to 10 orders of magnitude lower than commonly used colorimetric substrates.
- **Invitrogen™ DynaLight™ Alkaline Phosphatase Substrate with RapidGlow™ Enhancer** is our newest ready-to-use chemiluminescent substrate formulation that has been optimized to achieve faster results in solution-based assays. The DynaLight Substrate with RapidGlow Enhancer formulation includes 1,2-dioxetane chemiluminescent substrate and a polymeric enhancer that enables ultrasensitive immunoassay detection by alkaline phosphatase (AP).
- **Applied Biosystems™ AMPPD™, CSPD™, Alkaline Phosphatase, and CDP-Star™ Substrates and Applied Biosystems™ Sapphire™ I, II, & III/Emerald™ I, II, & III generation enhancers** are chemiluminescent AP substrates and enhancers that can also be used in immunoassays. Many ready-to-use formulations of substrate with an enhancer are available, and custom formulations can be developed together to better meet your specifications and price requirements.
- **Invitrogen™ Gal-Star™ Substrate** is a chemiluminescent beta-Gal substrate for use in solution-based assays. Sensitivity can be improved by using an enhancer, such as Invitrogen™ Sapphire-II™ or Invitrogen™ Emerald-II™ Luminescence Enhancer.

## SuperSignal ELISA Pico Chemiluminescent HRP Substrate

SuperSignal ELISA Pico Chemiluminescent Substrate is optimized to generate an intense light signal and provides exceptional performance in luminometer-based assays. SuperSignal ELISA Pico Chemiluminescent Substrate offers greater sensitivity in ELISAs or other solution-based assays for enzyme detection and quantification. The ELISAs, performed in either a test tube or a microplate, are quantified by measuring relative light units (RLU) in a luminometer.

### Highlights:

- **Immediate light generation**—intense signal is produced immediately at room temperature or at 37°C; emits light at 425 nm
- **High signal-to-noise ratios**—minimal background
- **High sensitivity**—detects protein in ELISAs down to picogram levels
- **Convenient**—ambient shipping and room temperature storage conditions
- **High stability**—consistent performance of the working solution over an 8-hour period, with only a 10% decrease in activity at 24 hours
- **Flexible**—signal can be read in black or white opaque plates

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## SuperSignal ELISA Femto Maximum Sensitivity HRP Substrate

SuperSignal ELISA Femto Maximum Sensitivity Substrate is formulated for superior protein detection and low-end linearity in chemiluminescent ELISA applications. The substrate provides greater sensitivity in ELISAs or any other solution-based assay for enzyme detection and quantification. The ELISAs, performed in either a test tube or a microplate, are quantified by measuring relative light units (RLU) in a luminometer.

### Highlights:

- **Immediate light generation**—intense signal generated immediately both at room temperature and 37°C; emits light at 425 nm
- **Improved low-end linearity**—easy detection of low quantities of proteins with high signal-to-noise ratios and low-end linearity of dose response curves
- **High sensitivity**—femtogram-level detection of target proteins in ELISAs (Figure 23)
- **Reduction in assay time**—high sensitivity allows for shorter incubation steps
- **Stability**—storage for 6 months at room temperature or a minimum of 12 months at 4°C with a 6-hour working solution stability

SuperSignal ELISA Femto Maximum Sensitivity Substrate uses an improved enhancer system that meets the needs of high-throughput screening (HTS) applications and diagnostic assay development. SuperSignal ELISA Femto Maximum Sensitivity Substrate generates detectable light within 1 minute of addition to trace amounts of soluble HRP, saving up to 30 minutes per assay. This feature makes SuperSignal ELISA Femto Maximum Sensitivity Substrate ideal for HTS applications with the capability of running as many as 100,000 assays on robotic equipment.

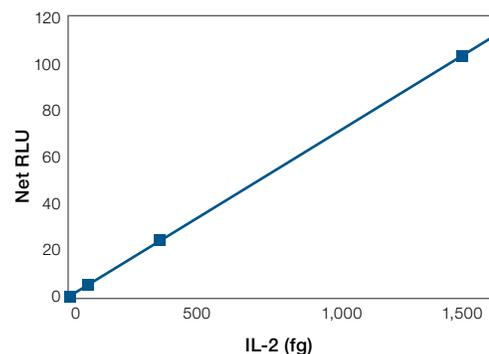


Figure 23. Femtogram-level detection of target protein and superior low-end linearity. The dose response curve generated from an IL-2 ELISA illustrates the exceptional low-end linearity achieved with SuperSignal ELISA Femto Maximum Sensitivity Substrate and the incredible sensitivity attainable. The SuperSignal ELISA Femto Maximum Sensitivity Substrate detected down to 168 fg of IL-2. The  $R^2$  value of the curve was calculated to be 1.00 for signal generated at less than 1,600 fg of IL-2.

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## DynaLight AP Substrate with RapidGlow Enhancer

The DynaLight Substrate with RapidGlow Enhancer is a ready-to-use chemiluminescent substrate formulation that has been optimized to achieve faster results in solution-based assays. The DynaLight Substrate with RapidGlow Enhancer formulation includes 1,2-dioxetane chemiluminescent substrate and a polymeric enhancer that enables ultrasensitive immunoassay detection by AP. DynaLight Substrate with RapidGlow Enhancer is designed to provide superior sensitivity, wider dynamic range, and fast read times that help reduce the overall completion time for an immunoassay.

The substrate provides a robust chemiluminescent reaction catalyzed by AP that achieves maximum light emission within 2 minutes at 37°C and maintains a sustained signal for hours. When combined with DynaLight Trigger Solution, DynaLight Substrate with RapidGlow Enhancer delivers maximum signal within seconds. DynaLight Trigger Solution provides maximum signal within one second of addition and a longer half-life to its ground state than other chemiluminescent flash technologies, allowing ample time to collect data reliably.

### Highlights:

- **Sensitive**—attomole-level detection of purified AP enzyme
- **Fast**—rapid kinetics and higher signal levels than colorimetric, fluorescence, or other chemiluminescence chemistries
- **Optimized**—for immunoassays in either microplate or magnetic bead format
- **Stable**—>1 year of signal stability

### Robust and sensitive assays using multiple formats

DynaLight Substrate with RapidGlow Enhancer can be used both in glow or flash mode, giving attomole-level sensitivity and up to 5 orders of magnitude of dynamic range of detection. As a single reagent addition, the DynaLight Substrate with RapidGlow Enhancer provides a steady chemiluminescence glow signal typically within 2 minutes at 37°C, or within 10 minutes at room temperature (Figure 24). The emission signal lasts several hours, offering maximum flexibility and convenience. The glow option is ideal for microplate immunoassays, where it can take several minutes to read the signal on a luminometer.

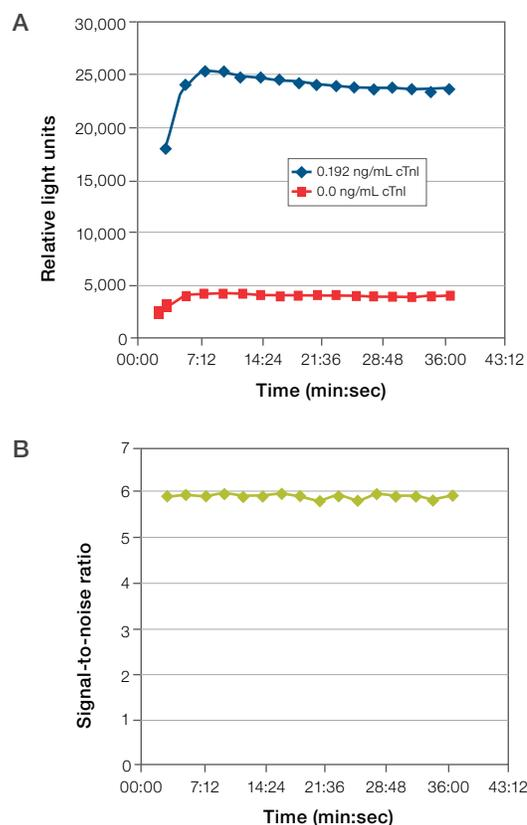


Figure 24. DynaLight Substrate glow light emission kinetics and signal-to-noise ratio. A bead-based sandwich immunoassay was performed at 0 and 0.192 ng/mL human cardiac troponin-I (cTnI) in human serum. DynaLight Substrate with RapidGlow Enhancer (100  $\mu$ L) was added and signal intensity kinetics at 30°C are shown as the assay reaches steady-state light emission. **(A)** Assay signal increases from 3 to 5 minutes after DynaLight Substrate with RapidGlow Enhancer addition until the assay reaches glow emission. **(B)** The signal-to-noise ratio remains stable over time from the initial read at 3 minutes, for a period of over 30 minutes.

For a flash protocol, the DynaLight Trigger Solution may be added as soon as 5 seconds after addition of DynaLight Substrate with RapidGlow Enhancer. The flash chemiluminescence signal has a longer half-life compared to direct-label chemiluminescent labels such as acridinium ester and isoluminol, allowing for further assay detection flexibility. Addition of the optional DynaLight Trigger Solution reduces the read time to seconds and increases assay signal up to 6-fold as quickly as 1 second after its addition (Figure 25).

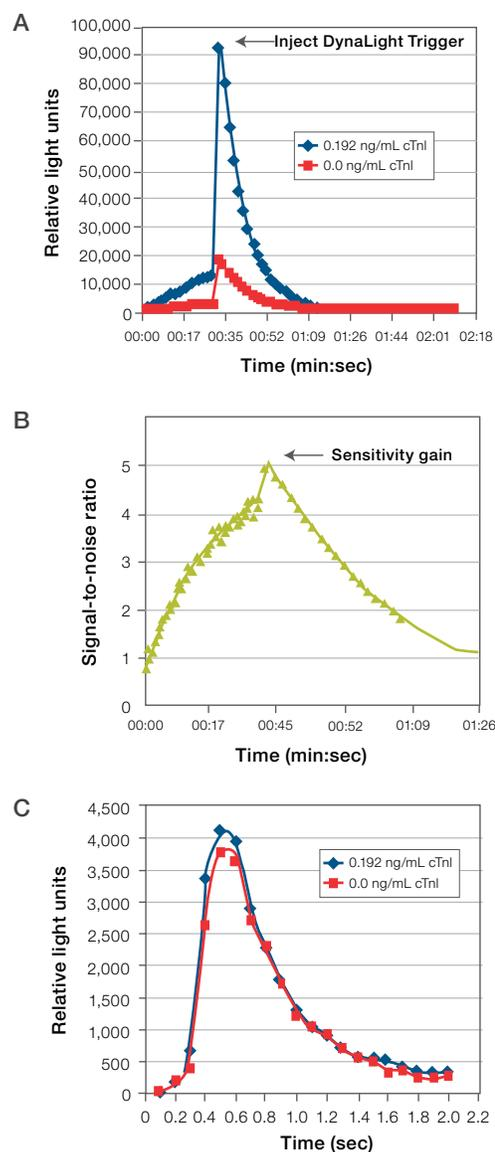


Figure 25. DynaLight Substrate flash light emission kinetics, signal-to-noise ratio with the optional DynaLight Trigger Solution, and acridinium ester flash light emission kinetics. A bead-based sandwich immunoassay was performed at 0 and 0.192 ng/mL human cardiac troponin-I (cTnI) in human serum, using Dynabeads M-280 tosylactivated magnetic beads as the capture surface. DynaLight Trigger Solution (50  $\mu$ L) was added 30 seconds after addition of DynaLight Substrate with RapidGlow Enhancer (50  $\mu$ L). **(A)** Assay signal increases for 30 seconds after DynaLight Substrate with RapidGlow Enhancer addition, with a further 6-fold signal increase within 1 second of DynaLight Trigger Solution addition. **(B)** The signal-to-noise ratio steadily increases over the first 30 seconds, with a further sensitivity improvement after DynaLight Trigger Solution addition. **(C)** Flash light emission kinetics for an acridinium ester-troponin-I conjugate show the more limited sensitivity and smaller read window compared to the DynaLight substrate's kinetic data in (A).

## Convert your colorimetric ELISA to help improve performance, time-to-results, and reduce reagent consumption

Colorimetric, fluorimetric, or other chemiluminescent ELISAs can be converted to a chemiluminescent assay using DynaLight Substrate with RapidGlow Enhancer. Advantages of converting your ELISA include superior sensitivity and dynamic range, combined with the flexibility to help decrease reagent usage and time-to-results, allowing you to tailor your assay to meet your specific needs.

Table 22 shows an example of a colorimetric IL-10 ELISA assay optimized for performance compared to a chemiluminescence assay using DynaLight Substrate with RapidGlow Enhancer optimized for sensitivity, reagent preservation, or minimal incubation times.

**Table 22. Example comparison of colorimetric and DynaLight chemiluminescence assay (IL-10 ELISA) conditions and outcomes.**

Assay conditions	Colorimetric	Chemiluminescence, optimized for:		
		Sensitivity	Reagent preservation	Minimization of incubation times
Capture (µg/mL)	1	1	0.5	1
Detector (µg/mL)	0.16	0.04	0.04	0.08
Streptavidin-enzyme dilution	1/2,500	1/31,250	1/31,250	1/31,250
Total incubation (minutes)	180	70	100	30
Sample and detector Ab	120	30	60	10
Streptavidin-enzyme	30	30	30	10
Substrate	30	10	10	10
Limit of blank (pg/mL)*	2.570	0.090	0.209	0.313
Dynamic range	3 logs	4 logs	4 logs	4 logs

\* Limit of blank = corresponding concentration of mean blank RLU measurements + 2 standard deviations of blank RLU measurements.

Sensitivity: conditions for optimal sensitivity were determined from using reagent titrations. Reagent preservation: the concentration of capture antibody was decreased by 50% to demonstrate reagent preservation. Note that the total assay time is still shorter than that of the colorimetric protocol, despite an increase in the sample and detector antibody incubation times relative to the other chemiluminescence assays. Minimization of incubation times: to minimize total assay time, all incubations were reduced to 10 minutes. The detector antibody concentration is still lower than in the colorimetric protocol, despite an increase relative to the optimal concentration for sensitivity.

For requests or inquiries on our products for immunodiagnostic assay development, please email [idxsupport@thermofisher.com](mailto:idxsupport@thermofisher.com)

## Colorimetric substrates

We offer the following colorimetric (also called chromogenic) substrates for immunoassay development with AP or HRP.

- **PNPP** (*p*-nitrophenyl phosphate, disodium salt) is a widely used substrate for detecting AP in ELISA applications. PNPP produces a yellow water-soluble reaction product that absorbs light at 405 nm. PNPP is available as a crystalline powder, 5 mg tablets, or a ready-to-use formulation.
- **ABTS** (2,2'-azino-bis-[3-ethylbenzothiazoline-6-sulfonic acid]-diammonium salt) is used to detect HRP and yields a water-soluble green end-reaction product. The green product has two major absorbance peaks, at 410 nm and 650 nm. ABTS is less sensitive than the OPD and TMB substrates for HRP detection. Color development is slow (approximately 20 minutes), which may be advantageous if OPD or TMB substrates generate unacceptable background because of their higher sensitivities. ABTS is available in either tablets or a ready-to-use formulation.
- **OPD** (*o*-phenylenediamine dihydrochloride) is used to detect HRP and yields a water-soluble yellow-orange reaction product. The reaction product has an absorbance maximum of 492 nm. OPD is available in either powder or tablet form and easily prepared by dissolving in Thermo Scientific™ Pierce™ Stable Peroxide Substrate Buffer or buffered hydrogen peroxide solution.
- **TMB** (3,3',5,5'-tetramethylbenzidine) soluble substrates yield a blue color when detecting HRP. The major absorbance maxima for the reaction product are 370 nm and 652 nm. The color then changes to yellow with the addition of sulfuric or phosphoric acid, with maximum absorbance at 450 nm. TMB is very sensitive and may produce significant background signal if too much protein or antibody is used. TMB is more quickly oxidized than other HRP substrates, resulting in faster color development.

## Pierce 1-Step Ultra TMB ELISA Substrate

Thermo Scientific™ Pierce™ 1-Step Ultra TMB ELISA Substrate detects HRP activity, yielding a blue color ( $A_{\text{max}} = 370 \text{ nm}$  and  $652 \text{ nm}$ ) that changes to yellow ( $A_{\text{max}} = 450 \text{ nm}$ ) upon addition of a sulfuric or phosphoric acid stop solution. The 1-Step Ultra TMB is the most sensitive of the chromogenic substrates.

### Highlights:

- **Sensitive**—highest sensitivity (2 pg/well) and signal-to-noise ratios of all TMB substrates
- **Ready to use**—single component without DMF or DMSO present in the reagent
- **Stable**—shelf life of 3 years
- **Competitive pricing**—available off-the-shelf in 1 L packaging
- **Safe**—noncarcinogenic
- **Convenient**—no additional reagents or filtering required

TMB is the most popular chromogenic substrate for HRP detection in ELISA and is available in several formats. The 1-Step TMB substrates are single-component substrates that require no preparation before use. Unlike other commercially available substrates, these products do not contain DMF or DMSO.

1-Step Ultra TMB substrate yields the greatest sensitivity among the TMB substrates, followed by Thermo Scientific™ Pierce™ 1-Step Turbo TMB and Thermo Scientific™ Pierce™ 1-Step Slow TMB ELISA substrates. The sensitivity of the 1-Step Turbo TMB substrate is similar to that of OPD-based substrates formulated at approximately 1 mg/mL. The 1-Step Slow TMB substrate has intermediate sensitivity—it is more sensitive than ABTS, but less sensitive than OPD or 1-Step Turbo TMB substrate. The 1-Step Slow TMB is an ideal substrate for kinetic studies.

## Pierce 1-Step NBT/BCIP Substrate Solution

Thermo Scientific™ Pierce™ 1-Step™ NBT/BCIP Substrate Solutions are ready-made solutions of nitro-blue tetrazolium and 5-bromo-4-chloro-3'-indolyphosphate for chromogenic blot and IHC staining with alkaline phosphatase probes. The combination of NBT (nitro-blue tetrazolium chloride) and BCIP (5-bromo-4-chloro-3'-indolyphosphate p-toluidine salt) yields an intense, insoluble black-purple precipitate when reacted with alkaline phosphatase, a popular enzyme conjugate for antibody probes. Pierce NBT and BCIP powders are available individually and in two convenient, ready-to-use solutions (with and without a levamisole suppressor of endogenous phosphatase activity).

### Highlights:

- **AP substrate**—for detection of AP activity on solid media, including nitrocellulose and PVDF membranes and fixed tissue samples
- **Chromogenic**—no special equipment needed for visualization; produces purple-black precipitate that is easy to photograph
- **Package options**—choose individual NBT and BCIP powders, or preformulated, ready-to-use solutions with or without levamisole suppressor

### Highlights of Pierce NBT/BCIP 1-Step Solution:

- Best for immunoblotting applications
- Ready-to-use single component
- Sensitive with low background and minimal assay-to-assay variability

### Highlights of Pierce NBT/BCIP plus Suppressor:

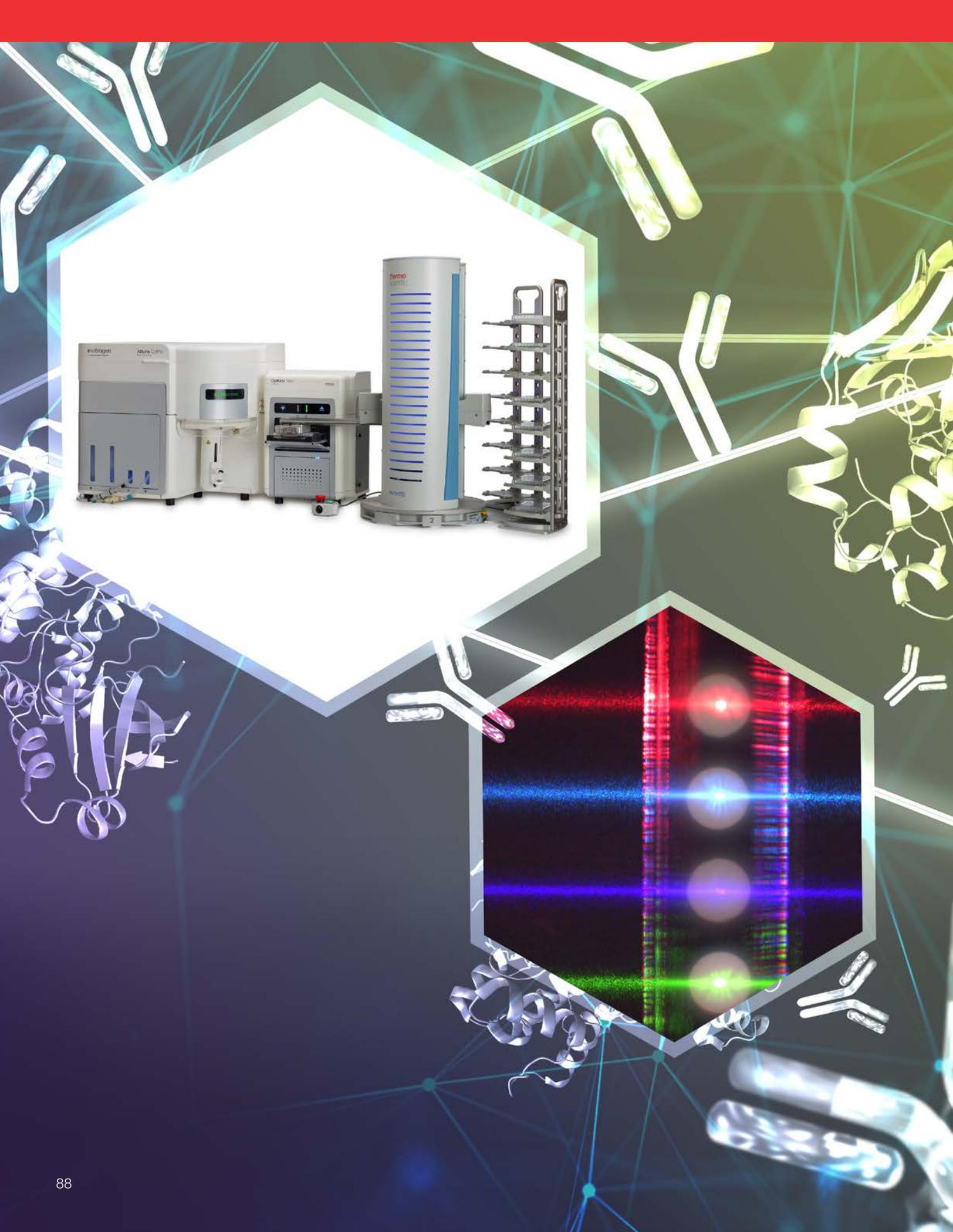
- Best for immunohistochemistry applications
- Ready-to-use single component
- Contains 1 mM levamisole to inhibit endogenous phosphatase activity, while yielding an intense dark purple precipitate in the presence of calf intestinal phosphatase

## Fluorescent substrates

The following chemifluorescent substrates are available for ELISA development with HRP.

- **Thermo Scientific™ QuantaBlu™ Fluorogenic Substrate** has a larger linear detection range with low-end linearity for detection of HRP. The stable fluorescent reaction product has an  $E_{\text{max}}/A_{\text{max}}$  of 420 nm/325 nm, allowing stopped, nonstopped, and kinetic assays to be performed, which is an advantage over the more sensitive chemiluminescent substrates.
- **Thermo Scientific™ QuantaRed™ Enhanced Chemifluorescent Substrate** is the most sensitive fluorescent ELISA substrate available for HRP detection. The fluorescent reaction product (resorufin) is stable for 4 hours, with an  $E_{\text{max}}/A_{\text{max}}$  of 585 nm/570 nm when the reaction is stopped. The red-shifted resorufin reaction product permits detection at a wavelength that has less interference from autofluorescence that can occur in biological samples.
- **Invitrogen™ Amplex™ Red Reagent** a highly sensitive and stable probe for  $\text{H}_2\text{O}_2$ , is one of our best fluorogenic substrates for peroxidase. Because  $\text{H}_2\text{O}_2$  is produced in many different enzymatic reactions, the Amplex Red reagent allows researchers to detect the activity of many different enzymes (Amplex Red Assay Kits).
- **Invitrogen™ Amplex™ UltraRed Reagent** improves upon the performance of our unique Amplex Red reagent, offering brighter fluorescence and enhanced sensitivity on a per-mole basis in peroxidase or peroxidase-coupled enzyme assays.





# Instruments

## Attune Flow Cytometers

### Building a legacy of advancing flow cytometry technology

Technological advancements in flow cytometry are creating an entirely new lineage of reagents and instruments. Innovative flow cytometry solutions enable researchers to create more effective, efficient, and economical solutions. Thermo Fisher Scientific is committed to developing products that fuel groundbreaking research. Our comprehensive suite of flow cytometry solutions, including the Invitrogen™ Attune™ Flow Cytometers and Invitrogen™ CytKick™ Autosamplers and automation options (Figure 26), together with the Invitrogen™ cell health reagents portfolio and Invitrogen™ eBioscience™ antibody conjugates, helps drive discovery of new biological insights for many applications.

Find out more at [thermofisher.com/attune](https://thermofisher.com/attune)

### Designed for efficiency, speed, and accuracy

#### Technology

Acoustic focusing fluidics is the key to the high sensitivity of Attune Flow Cytometers, even at high throughput. The Attune CytPix model adds a high-speed camera that captures brightfield images of cells passing through the flow cell to verify that cell populations consist of single cells and verify morphology.

#### Acoustics-assisted hydrodynamic focusing technology

Attune Flow Cytometers combine ultrasonic waves like those used in medical imaging with hydrodynamic forces to precisely position cells into a single, focused line in the central axis. Enabling cells to be tightly focused at the point of laser interrogation allows the system to collect more photons, helping to ensure data quality regardless of the sample-to-sheath ratio (Figure 27). This allows for a higher degree of data, detail, and throughput that enables processing of a large range of sample types, including large clumpy cells, samples with low concentration of cells, and precious samples, more quickly and accurately than ever before with no loss in data quality.

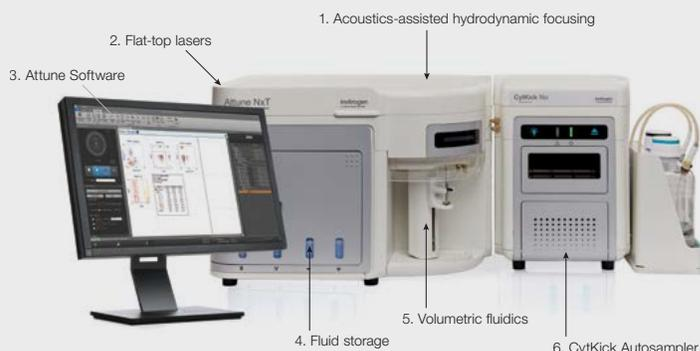


Figure 26. Attune Flow Cytometer components. **(1)** Patented acoustics-assisted hydrodynamic fluidics increase sample input speed while maintaining data integrity. **(2)** Flat-top lasers deliver more even application of light to each cell. **(3)** Invitrogen™ Attune™ Software designed to guide users through complex flow cytometry experiments. **(4)** Fluid storage designed for minimal waste. **(5)** Volumetric fluidics provide cell counting and a resistance to clogging. **(6)** The CytKick Autosampler provides easy one-click transition from tube to plate.

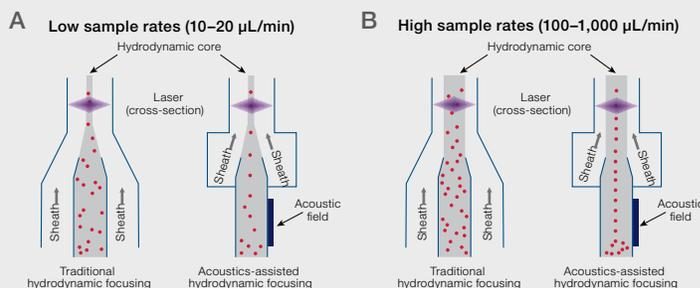


Figure 27. Acoustic focusing vs. traditional hydrodynamic focusing as particles pass through the laser. **(A)** In acoustic focusing, cells remain in tight alignment even at higher sample rates, resulting in less signal variation and improved data quality. **(B)** In traditional hydrodynamic focusing, increasing the sample rate results in widening of the sample core stream, resulting in increased signal variation and compromised data quality.

## Efficient—rapid, accurate acquisition

### Fast, accurate acquisition

Acoustic focusing empowers your lab to rapidly acquire high-quality data (Figure 28). You can achieve sample throughput rates of 12.5  $\mu\text{L}/\text{min}$  or 1,000  $\mu\text{L}/\text{min}$ , up to 10 times faster than traditional hydrodynamic focusing systems and acquisition speeds of up to 35,000 events per second. This means you can process all of your samples—including low-concentration and precious samples—more quickly and accurately with minimal loss in quality.

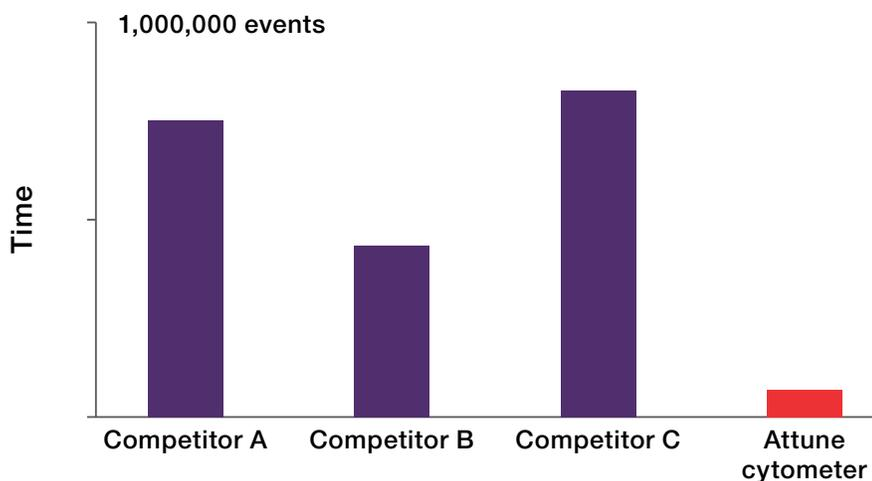


Figure 28. Rapid data acquisition. The time required to acquire 1,000,000 events is compared to three competitor instruments running at maximum sample rates.

“What I like most about the Attune NxT Flow Cytometer is its accuracy and its speed of processing. I could list a range of other things that I like about it because we are building up that list as we discover more aspects of its functions, but the core things are speed and accuracy that are not available with hydrodynamic flow cytometers.”

— Tim Inglis, BM, DM, PhD, FRCPPath, FRCPA, DTM&H  
School of Medicine, University of Western Australia, PathWest Laboratory Medicine

## Rare-event detection

Detection of rare events requires acquisition of high numbers of cells to attain a reliable measure of accuracy (Figures 29 and 30). Attune Flow Cytometers allow dilute samples to be processed quickly at sample input speeds of up to 1 mL/min, significantly faster than conventional cytometers that support maximum sample input rates of 60–100  $\mu\text{L}/\text{min}$ . Acoustic focusing thus offers a unique combination of speed and quality, cutting the time to collect rare events significantly over long acquisition times.

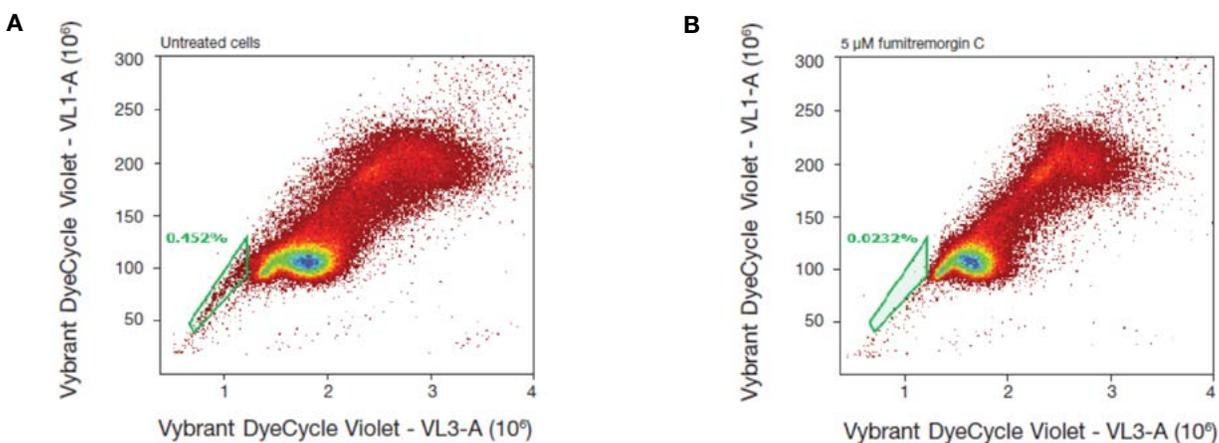


Figure 29. Identifying limbal stem cells (LSC) in a population of differentiated corneal cells. **(A)** Untreated cells have a side population of limbal stem cells (0.452% of total cells) with decreased Invitrogen™ Vybrant™ DyeCycle™ Violet fluorescence due to ABCG2-mediated dye efflux. **(B)** The percentage of cells in the side population is reduced (0.023% of total cells) when the ABCG2 membrane pump is inhibited with 5  $\mu\text{M}$  fumitremorgin C, preventing efflux of DyeCycle Violet dye.

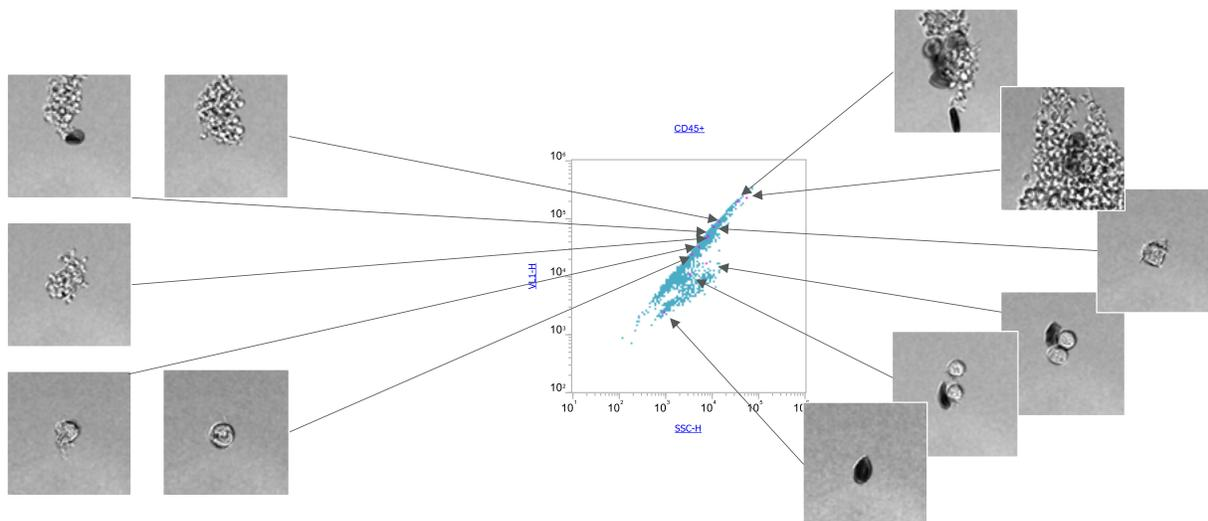


Figure 30. Gain insight with rare-population analysis. Cells were acquired from 24-hour old blood diluted in 1 mM EDTA (<1:4,000). Samples were acquired at 25  $\mu\text{L}/\text{min}$ . The image gate was set to record only CD45<sup>+</sup> events.

# Transformative—break the mold of traditional flow

## Reduce clogging from difficult samples

Your research samples are precious, as they are often difficult to produce. Attune Flow Cytometers are less prone to clogging than traditional flow cytometers, allowing challenging samples such as cardiomyocytes, heterogeneous blood cells, and cancer cells to flow with confidence.

Engineered to actively resist clogging, a syringe-driven system (Figure 31) and larger flow cell help prevent the loss of precious sample such as cancer stem cells from primary pancreatic tumors, and is far less susceptible to clogs. Attune Flow Cytometers employ a nonpressurized system that mechanically decreases the occurrence of clogging.

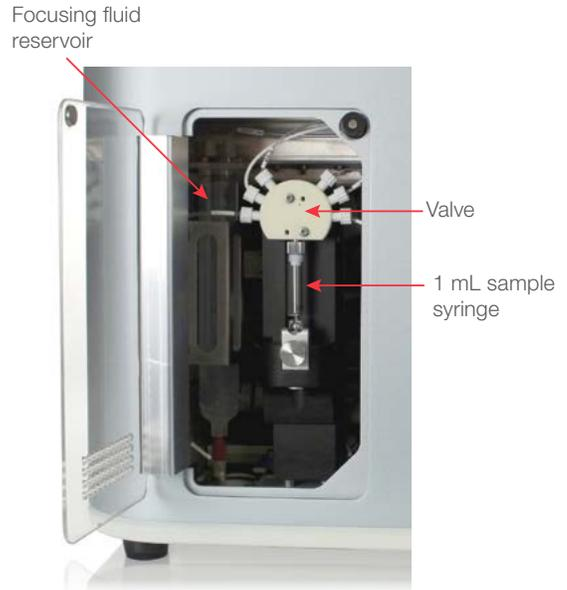


Figure 31. Positive-displacement syringe pump. The syringe is easily removed for cleaning or replacement.

## No-lyse, no-wash, with no compensation

Acoustic focusing allows Attune Flow Cytometers to deliver a no-wash, no-lyse protocol to minimize cell loss, significantly reduce time, and simplify sample preparation (Figure 32).

### Protocol

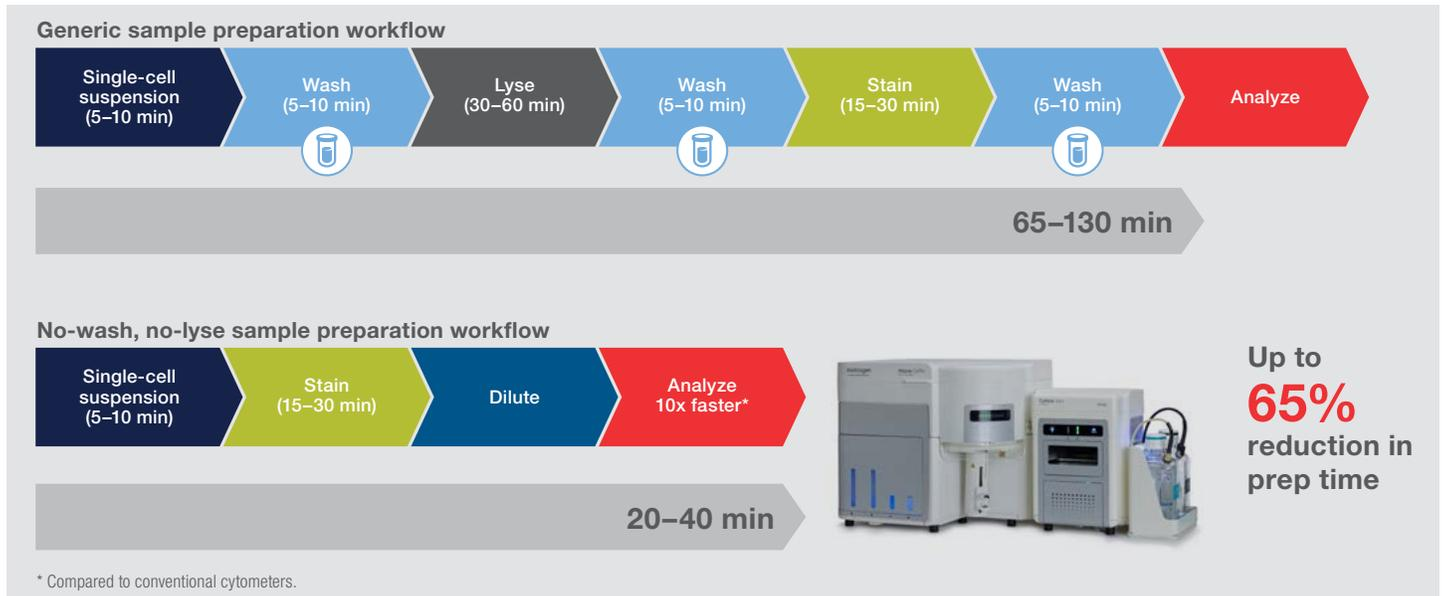


Figure 32. No-wash, no-lyse sample preparation workflow.

## Revealing—brightfield imaging; capture up to 6,000 images per second

With the Attune CytPix Flow Cytometer, you can highlight structural features of large populations in record time (Figure 33). Using image gallery view, you can scan your cell populations rapidly for outliers. This allows you to adjust your gates, if you wish, to include cells of interest while excluding aggregates, unwanted cells, and debris.

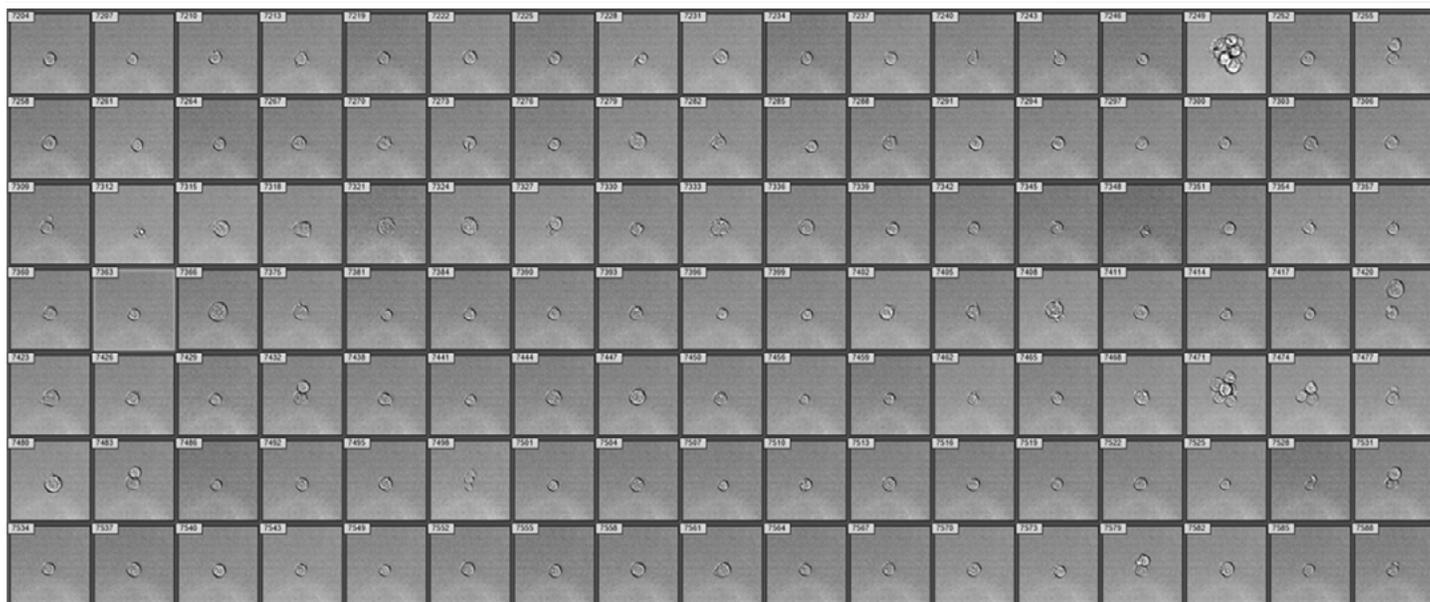


Figure 33. High-level image gallery view. Early log phase Jurkat cells were acquired unfiltered on the Attune CytPix Flow Cytometer at 200  $\mu\text{L}/\text{min}$ ,  $>10^6$  cells/mL. Image gallery view was used to rapidly scan cellular events.

“I can’t believe how quick and easy it is to collect a large body of empirical evidence on the Attune NxT Flow Cytometer to fully support our hypothesis.”

— Jordi Petriz, PhD  
Group Leader, Functional Cytomics Group;  
Josep Carreras Leukaemia Research Institute, Barcelona, Spain

# EVOS M7000 Imaging System

## A powerful, fast, fully automated system

Bring high performance and fast, automated imaging right to your lab bench with the Invitrogen™ EVOS™ M7000 Imaging System. This system has been designed with advanced capabilities to simplify demanding cell-based imaging applications such as live-cell analysis, image tiling, and Z-stacking, so you can focus on acquiring images and data rather than instrument operation.

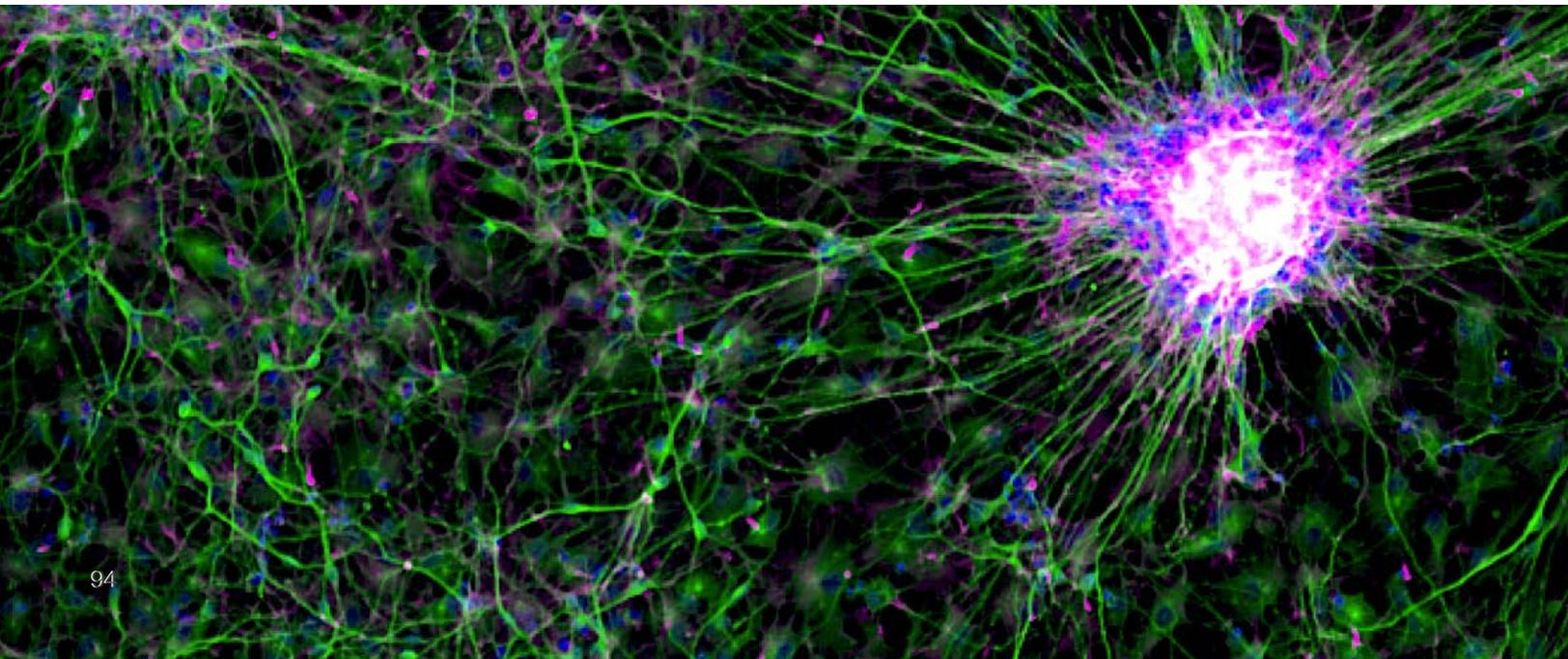
### Features

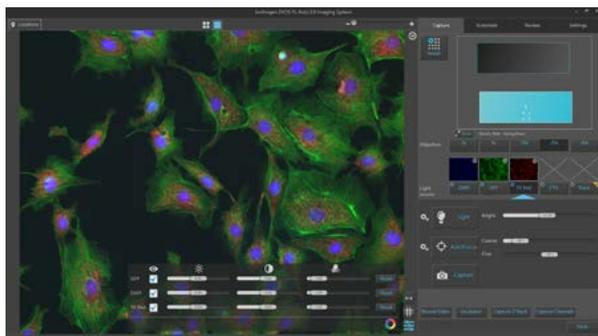
- **Speed**—scan a 96-well plate in three fluorescence channels in less than 5 minutes
- **Flexibility**—customize the system with more than 20 user-changeable LED light cubes, dual cameras (monochrome and color), a variety of objectives ranging from 1.25x to 100x, and multiple vessel holders
- **Time-lapse live-cell imaging**—onstage incubator option for precise control of temperature, humidity, and gases for normoxic or hypoxic conditions allows a wide range of biological studies under physiological conditions
- **Two cameras, no compromises**—all systems come with two cameras: a dedicated high-sensitivity monochrome camera optimized for fluorescence imaging and quantitation, and a dedicated high-resolution color camera optimized for colorimetric imaging
- **Area view**—move rapidly and seamlessly between single-field mode and low- and high-magnification scan modes to easily define and capture the area of interest
- **Automation**—time-saving features such as autofocus, rapid stage movement, and automated routines help reduce time to complete experiments, allowing high throughput, high data quality, and improved experimental reproducibility
- **Data analysis**—extensive quantitative imaging and statistical analysis in combination with Invitrogen™ Celleste™ Image Analysis Software, an optional advanced software package offering powerful tools for image segmentation and classification that can be used for cell counting and for measuring changes in intensity, area, and shape over time

### SmartStart installation and training

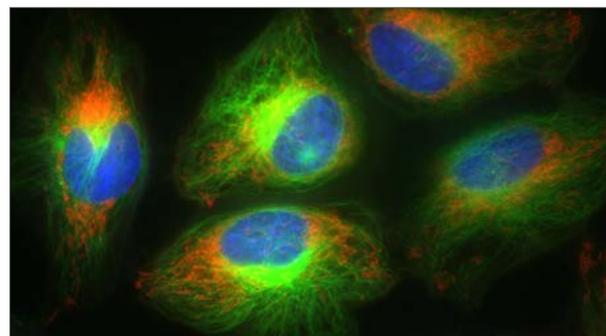
A specialized installation and training program will get you up and running in just one day. One of our dedicated field application scientists comes to your site to provide hands-on workflow training and make sure your lab is quickly enabled to utilize the instrument's powerful features to maximize productivity.

Neural stem cell colony, 10x objective; light cubes: GFP, RFP.





Easy-to-navigate user interface.



Multiplex immunofluorescence staining of HeLa cells with Invitrogen™ NucBlue™ Live ReadyProbes™ Reagent (blue), Tubulin Tracker Green Variety Pack (green), and CellLight™ Mitochondria-RFP BacMam 2.0 (red).

## System highlights

Attribute	Details
Optics	Infinity-corrected optical system; Royal Microscopical Society (RMS) threaded objectives with a 45 mm parfocal distance
Imaging mode	Fluorescence, brightfield, color brightfield, and phase contrast
Illumination	5-position chamber for 4 fluorescence light cubes plus brightfield imaging; light cubes with integrated hard-coated filter set and LED light source with >50,000-hour life; broad selection of standard and specialty light cubes
Imaging methods	Single-color, multicolor, area scan with montage or tile stitch, time lapse, Z-stacking, movie capture
Objective capacity	5-position turret
Objectives (not included)	Wide selection of high-quality long working distance (LWD) and coverslip-corrected objectives
Condenser	60 mm LWD condenser; 4-position turret with a clear aperture and 3 phase annuli
Stage	Motorized X/Y scanning stage; 120 mm x 80 mm travel range with submicron resolution; drop-in inserts to receive vessel holders and lockdown holders to fix sample in place during long scans
Focus mechanism	Automated focus mechanism with submicron resolution
LCD display	23-inch high-resolution touchscreen color monitor (also fully controllable via mouse); 1,920 x 1,080 resolution
Cameras	High-sensitivity 3.2 MP (2,048 x 1,536) monochrome CMOS sensor with 3.45 μm pixel resolution; high-sensitivity 3.2 MP (2,048 x 1,536) color CMOS sensor with 3.45 μm pixel resolution
Computer	External Dell™ PC with an Intel Core™ i7-8700 processor, 32 GB DDR4 RAM, 512 GB PCIe solid-state drive, NVIDIA Quadro™ P1000 graphics card with NVIDIA™ Pascal™ GPU technology and 4 GB memory, and Windows™ 10 software, designed to operate with touchscreen monitor and microscope
Captured images	8-bit TIFF, PNG, JPG; 16-bit RAW monochrome: TIFF, PNG; movies and time-lapse images: AVI, WMV
Output ports	Microscope: USB 3.1 Type B, 4-pin power port Computer: 1 x USB 3.1 Gen 2 Type C; 5 x USB 3.1 Gen 1 Type A; 4 x USB 2.0 Type A; 1 serial; 2 x DisplayPort 1.2; 1 RJ45; 2 PS/2; 1 UAJ; 1 line-out
Networking capability	Connection through Windows/SMB network via an Ethernet cable connection
Power supply	24 V AC adapter with country-specific power cords
Dimensions (L x W x H)	45.7 x 35.6 x 33.0 cm (18 x 14 x 13 in.)
Weight	16 kg (35 lb)

Find out more at [thermofisher.com/evosm7000](https://www.thermofisher.com/evosm7000)

# Plate reader and washers

## Choose the microplate reader that best suits your research needs

Thermo Scientific™ microplate readers provide flexibility, performance, and ease of use for a variety of microplate assays. Whether you need to measure fluorescence, absorbance, luminescence, or time-resolved fluorescence, or perform AlphaScreen™ assays, we offer a microplate reader solution to meet the needs of your specific workflow. With a portfolio of dedicated and modular, upgradeable multimode readers, we also offer solutions that fit your current budget with options to meet your lab's future needs.

Thermo Scientific plate readers have a number of features to help you save time and maximize productivity, including:

- Auto-calibration
- Easy export to Microsoft™ Excel™ format
- Automation readiness with robot compatibility
- No limit to the number of computers on which you can install the intuitive Thermo Scientific™ SkanIt™ Software
- Ready-to-use protocols available in an extensive online protocol library

Considerations	Multiskan FC photometer	Multiskan SkyHigh spectrophotometer	
Applications	Absorbance		
Wavelength range	340–850 nm	200–1,000 nm	
Wavelength selection	Filters	Monochromator	
Plate format	96 wells (384 wells optional)	Thermo Scientific™ $\mu$ Drop™ and $\mu$ Drop™ Duo Plates, 6–48 well plates,* 96- and 384-well plates	
Incubation	Optional	Yes	
Shaking	Yes	Yes	
Dispensers	No	No	
Top/bottom reading	NA	NA	
Cuvettes	No	Optional	
Gas control module	No	No	
			

\* Maximum plate height with lid 19.5 mm.

\*\* Third dispenser requires additional installation.

† Instruments with bottom-read capabilities feature multilocation reads per well.

For more information, go to [thermofisher.com/platereaders](https://www.thermofisher.com/platereaders)

	<b>Fluoroskan fluorometer</b>	<b>Luminoskan luminometer</b>	<b>Fluoroskan FL fluorometer and luminometer</b>	<b>Varioskan LUX scanning multimode reader</b>
	Fluorescence	Luminescence	Fluorescence, luminescence	Absorbance, fluorescence Optional: time-resolved fluorescence (TRF), luminescence, AlphaScreen readout
	Excitation: 320–700 nm Emission: 360–800 nm	270–670 nm	Fluorescence excitation: 320–700 nm emission: 360–670 nm  Luminescence: 270–670 nm	Absorbance and fluorescence excitation: 200–1,000 nm Fluorescence emission: 370–840 nm Luminescence: 360–670 nm TRF excitation: fixed to 334 nm (spectral scanning 200–840 nm) TRF emission: 400–700 nm (spectral scanning 270–840 nm) AlphaScreen excitation: fixed to 680 nm AlphaScreen emission: 400–660 nm
	Filters	Not required for most applications Filters can be used when necessary	Filters	Monochromator for UV-Vis absorbance and fluorescence intensity Filters for luminescence when necessary, TRF, AlphaScreen assays
	6–384 wells	6–384 wells	6–384 wells	6–1,536 wells (fluorometry, TRF, luminometry, AlphaScreen assays) $\mu$ Drop and $\mu$ Drop Duo Plates/6–384 wells (absorbance)
	Yes	Yes	Yes	Yes
	Yes	Yes	Yes	Yes
	Optional (to three)**	Optional (to three)**	Optional (to three)**	Optional (to two)
	Top/bottom†	Top	Top/bottom†	Top (standard) Bottom (optional)†
	No	No	No	With $\mu$ Drop Plate
	No	No	No	Optional
				

# Multiskan SkyHigh Microplate Spectrophotometer

PRODUCT HIGHLIGHT

Spike (trimer) ELISA kit for SARS-CoV-2

## Measure SARS-CoV-2 spike antibodies with this ELISA kit and versatile microplate reader

The Invitrogen™ Human SARS-CoV-2 Spike (Trimer) Ig Total ELISA Kit, along with the Thermo Scientific™ Multiskan™ SkyHigh Microplate Spectrophotometer, provides a sensitive and specific solution for detecting and quantifying the presence of antibodies against SARS-CoV-2 in research samples.

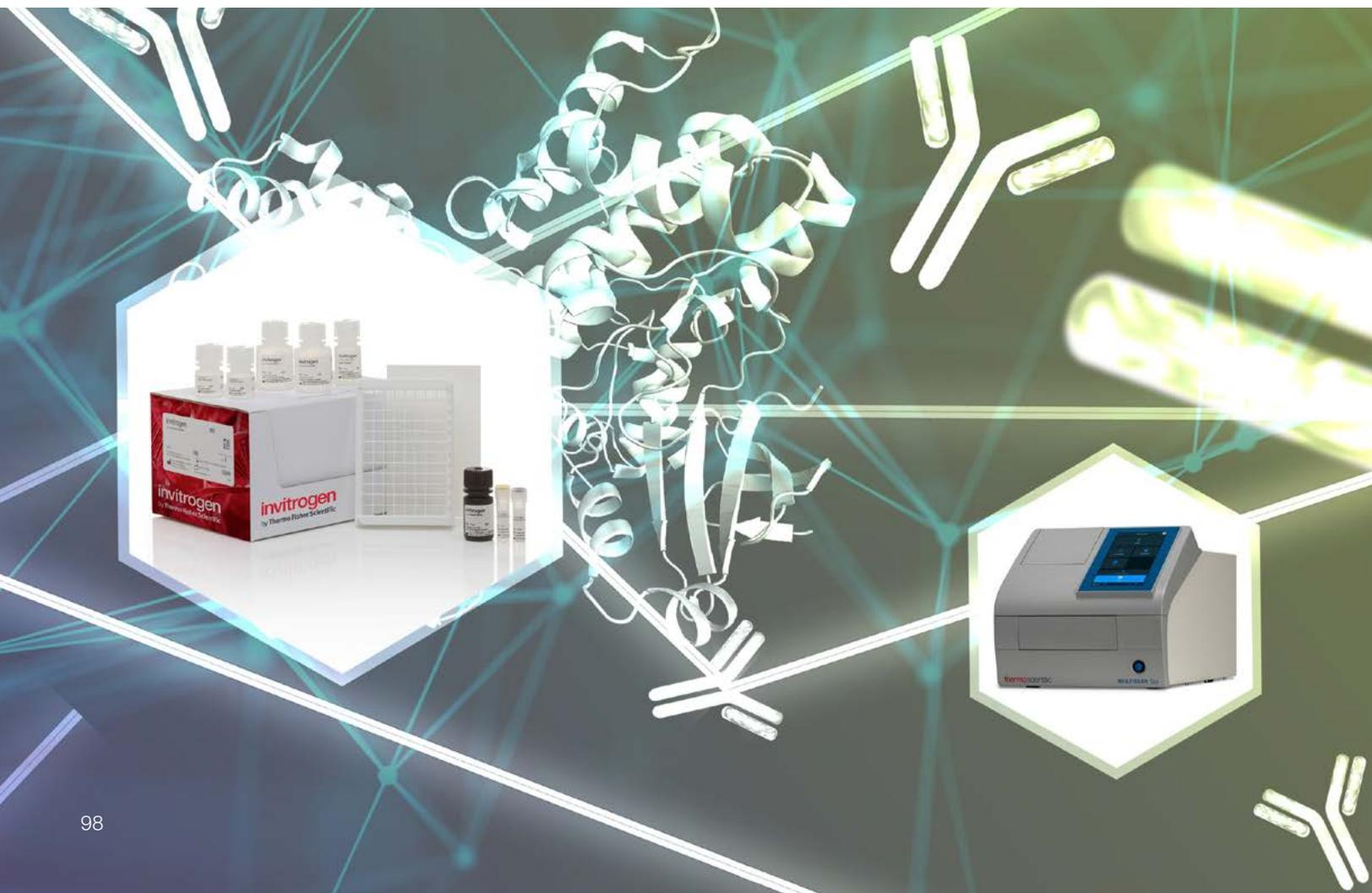
### Ig total ELISA kit highlights:

- **Reliability**—specificity is determined with PCR (+) and PCR (-) samples to minimize false positives, and sensitivity is optimized to detect antibodies to minimize false negatives
- **Flexible results (qualitative or quantitative)**—either positive and negative results, or a 7-point standard curve for quantitation

### The Multiskan SkyHigh plate reader offers:

- **Flexible format**—6, 48, or 96 wells, or 384-well microplates
- **Versatility**—measures absorbance and readouts for a variety of assays
- **Ease of use**—Thermo Scientific™ SkanIt™ software features visual workflows and settings to minimize errors
- **Data security**—remote connectivity and data export available via either Microsoft™ OneDrive™ or Thermo Fisher™ Connect platforms

Order your ELISA kit and/or Multiskan SkyHigh plate reader at [thermofisher.com/sars-serology](https://www.thermofisher.com/sars-serology)



# Qubit Flex Fluorometer

## Accurate and sensitive quantification of DNA, RNA, and protein with flexible throughput

With the same convenience and ease of use as the Qubit 4 Fluorometer, the Invitrogen™ Qubit™ Flex Fluorometer provides improved throughput. The Qubit Flex instrument can selectively and accurately measure the concentrations of up to 8 samples of DNA, RNA, or protein simultaneously.



### Key benefits

- **High sensitivity**—more sensitive than UV absorbance-based quantification
- **Accuracy and speed**—accurately quantifies DNA, RNA, or protein in less than 3 seconds per sample (Figures 34 and 35)
- **Helpful sample prep calculators**—integrated calculators help determine the amount of dye, buffer, and sample needed
- **Specialized calculators for NGS workflows**—easily determine molarity and normalize sample concentrations
- **Ideal for precious samples**—uses as little as 1  $\mu\text{L}$  of sample
- **Plenty of data storage**—stores data for up to 10,000 samples internally
- **Flexible data export**—Wi-Fi, USB drive, or direct connection with a USB cable
- **Flexible, improved throughput**—measure up to 8 samples per run
- **Portable with small footprint**—easy to move and store on shelves in the lab
- **Optimized reagents and tubes**—Qubit reagents and tubes work optimally with Qubit fluorometers
- **Intuitive touchscreen**—quickly get to data generation

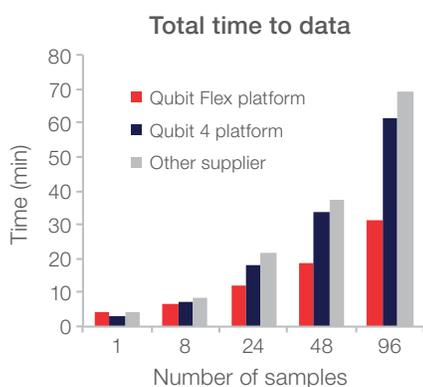


Figure 34. Time trial on 1 to 96 samples shows up to 50% time savings with the Qubit Flex Fluorometer. Time savings began with 8 samples, with more time saved when measuring even more samples.

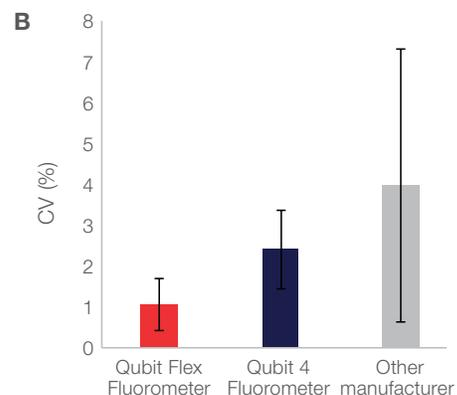
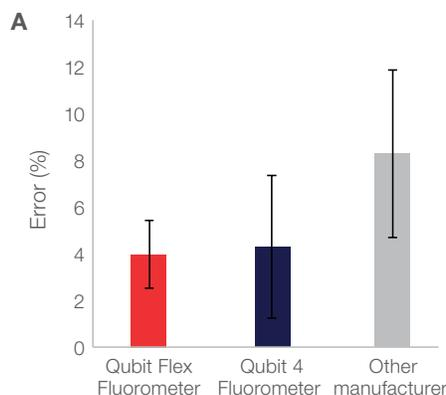


Figure 35. Qubit instruments deliver better accuracy and precision. The Qubit dsDNA BR and 1X dsDNA HS Assay Kits were run with 4 DNA sample concentrations (HS: 0.1, 1, 5, and 10  $\text{ng}/\mu\text{L}$ ; BR: 2, 20, 50, and 100  $\text{ng}/\mu\text{L}$ ) in 8 replicates on the Qubit 4 Fluorometer, Qubit Flex Fluorometer, and another manufacturer's fluorometer. The percent relative error (lower is more accurate) and CV (lower is more precise) were determined for each concentration and averaged across all concentrations for each instrument.

Find out more at [thermofisher.com/qubitflex](https://www.thermofisher.com/qubitflex)

## Summary

Developing and commercializing immunodiagnostic assays are difficult and time-consuming. Optimizing time and assay quality while staying within budget is also challenging, but finding an experienced supplier is the first step in the right direction. From small-scale laboratories to leading immunodiagnostics companies, our team of technical specialists has extensive and relevant experience collaborating with researchers in this area and is familiar with the problems you may face. Therefore, a collaboration with us can aid in overcoming those challenges and help meet your goals.

Partnering with us provides access to our experienced Global OEM and Licensing team that will help develop and commercialize your immunodiagnostic assay. Our breadth and depth of technologies and products for immunodiagnostics development include capture surfaces, antibody conjugation tools and detection probes, linking mechanisms, blocking buffers and detergents, detection substrates, and instrumentation and equipment. Additionally, if our off-the-shelf products do not meet your requirements, we offer customized products according to

your desired specifications. With superior product knowledge and process control, our experienced team of professionals can help select the right products for your workflow, as well as custom products as needed. With over 5,000 patents and licenses in immunodiagnostic products, our business development manager will assist in determining necessary licenses for commercialization.

With world-class manufacturing capabilities, we can help scale up your immunodiagnostic assay. Our ISO 9001– and/or ISO 13485–certified facilities are equipped with large-scale bioprocessing units, HEPA-filtered cleanrooms, chemical synthesis laboratories, custom packaging and labeling capabilities, and thorough QC testing suites to help you in every step of your immunodiagnostic assay workflow. We are the single supplier you need to help accelerate the development and commercialization of your immunodiagnostic assay.

 Find out more at [thermofisher.com/idx](https://thermofisher.com/idx)