

Product Information

Labeled Goat Anti-Human IgG (H+L) Antibodies

Cat. Number	Label	Unit Size
L102A	Andy Fluor 350	100 µg
L102B	Andy Fluor 350	500 µg
L105A	Andy Fluor 405M	100 µg
L105B	Andy Fluor 405M	500 µg
L108A	Andy Fluor 430	100 µg
L108B	Andy Fluor 430	500 µg
L111A	Andy Fluor 488	100 µg
L111B	Andy Fluor 488	500 µg
L115A	Andy Fluor 555	100 µg
L115B	Andy Fluor 555	500 µg
L118A	Andy Fluor 568	100 µg
L118B	Andy Fluor 568	500 µg
L121A	Andy Fluor 594	100 µg
L121B	Andy Fluor 594	500 µg
L127A	Andy Fluor 647	100 µg
L127B	Andy Fluor 647	500 µg
L130A	Andy Fluor 680	100 µg
L130B	Andy Fluor 680	500 µg
L133A	Andy Fluor 750	100 µg
L133B	Andy Fluor 750	500 µg
L136A	Cy3	100 µg
L136B	Cy3	500 µg
L139A	Cy5	100 µg
L139B	Cy5	500 µg
L142A	Cy5.5	100 µg
L142B	Cy5.5	500 µg
L145A	Cy7	100 µg
L145B	Cy7	500 µg
L148A	FITC	100 µg
L148B	FITC	500 µg
L151A	Biotin	100 µg
L151B	Biotin	500 µg

Spectral Properties:

Label	Ex nm	Em nm
Andy Fluor 350	345	440
Andy Fluor 405M	405	450
Andy Fluor 430	430	540
Andy Fluor 488	495	520
Andy Fluor 555	553	565
Andy Fluor 568	578	605
Andy Fluor 594	590	615
Andy Fluor 647	650	665
Andy Fluor 680	680	700
Andy Fluor 750	750	770
Cy3	550	565
Cy5	650	667
Cy5.5	680	700
Cy7	750	770
FITC	490	520

Storage upon receipt:

- 2-8 °C
- Protect from light
- Avoid freeze-thaw cycles

Product Description

Applied BioProbes offers an extensive line of goat anti-human IgG conjugates labeled with a wide selection of fluorescent dyes, biotin, or HRP.

Fluorescent Goat Anti-Human IgG conjugates are ideal for fluorescence microscopy and confocal laser scanning microscopy, flow cytometry, and fluorescent western detection. The breadths of fluorescent markers we offer allow our reagents to be tailored to almost any fluorescent detection system.

The Goat Anti-Human IgG antibody conjugates are prepared from affinity purified antibodies that react with human IgG heavy chains and all classes of human immunoglobulin light chains. To minimize cross-reactivity, the goat anti-human IgG antibodies have been adsorbed against mouse, rabbit and bovine sera prior to conjugation. The degree of labeling for each conjugate is typically 2–8 fluorophore or biotin molecules per IgG molecule. At the time of preparation, the products are certified to be free of unconjugated dyes and are tested in an immunofluorescence experiment to ensure low nonspecific staining.

Product Specification

Physical State: Lyophilized power

Buffer: PBS, pH 7.4

Stabilizer: 0.1% BSA

Preservative: 0.02% Sodium Azide

Reconstitution and Storage: Store lyophilized power at 2-8°C. When ready to use, rehydrate with dH₂O (50 µL for 100 µg antibody or 250 µL for 500 µg antibody) to make 2 mg/mL solution and centrifuge if not clear. Product is stable for about 6 months at 2-8°C as an undiluted liquid. Prepare working dilution fresh each day. For extended storage after rehydration, add an equal volume of glycerol for a final concentration of 50%, and store at -20 °C as a liquid.

Guidelines for Use

Centrifuge the reconstituted protein conjugate solution briefly in a microcentrifuge before use. Add only the supernatant to the experiment. This step eliminates any protein aggregates that may have formed during storage, thereby reducing nonspecific background staining.

Because staining protocols vary with application, the appropriate dilution of antibody should be determined empirically. For the fluorophore- and biotin-labeled antibodies, a final concentration of 1–10 µg/mL should be satisfactory for most immunohistochemical applications. For flow cytometry applications, 0.06–1.0 µg per 1 × 10⁵ cells should yield satisfactory results.

General Protocols

Immunofluorescence Protocol for Microscopy

There are many methods for immunofluorescence staining. The protocol below is a general guideline for staining cells and should be optimized or modified to obtain the best results for each particular application.

1. Coverslip preparation for adherent cells

1.1 Culture cells on slide chambers or sterile glass coverslips. We recommend 18 x 18 mm square coverslips in 6-well plates or 4-well chamber slides.

1.2 Allow cells to adhere and treat as desired.

1.3 Rinse cells gently with PBS.

2. Coverslip preparation for non-adherent cells

2.1 Coat coverslips with 0.01% poly-L-lysine solution for 10 minutes at room temperature.

2.2 Aspirate the poly-L-lysine solution and allow coverslips to dry completely.

2.3 Centrifuge cells in medium and resuspend in PBS. Transfer cells to coverslips.

2.4 Incubate for 30-60 minutes. Check for adherence by microscope.

3. Fixation and Staining

3.1 Fix with 4% paraformaldehyde/PBS, 15 min.

3.2 Rinse twice with PBS to remove traces of fixative.

3.3 Permeabilize with 0.1 - 0.5% TritonX-100/PBS, 5-10 min.

3.4 Block with blocking agent such as with 5% BSA or normal goat serum in PBS, 30 min.

3.5 Dilute primary antibody in dilution buffer as recommended in the specific product's datasheet. Overlay enough diluted antibody to cover cells on coverslip (150-200 μ L is usually sufficient to cover the surface area) or add to each chamber of the chamber slides. Keep slips covered or in a humidified chamber to avoid evaporation.

3.6 Rinse three times with PBS, 5 min each wash.

3.7 Dilute fluorescent secondary antibody in dilution buffer and incubate for 1 hour at room temperature. General range for IgG conjugates is between 1-10 μ g/mL for most applications. Cell samples without primary antibody incubation is recommended for background control. Keep

slips covered or in a humidified chamber to avoid evaporation.

3.8 Rinse three times with PBS, 5 min each wash.

3.9 Additional staining with fluorescent nuclear stains or phalloidins can be done at this step.

3.10 Invert each coverslip onto a pre-cleaned slide with fluorescence anti-fade mounting media. Seal edges with clear polish if desired.

3.11 Store slides in the dark at 4°C.

Staining Protocol for Flow Cytometry

There are many alternative procedures that can be used for specific staining experiments. The protocol below is a general guideline for flow cytometry and should be optimized or modified for each application.

1. Aliquot 1×10^6 cells into 12 X 75 mm polypropylene tubes for flow cytometry.

2. For intracellular staining, cells can be fixed first to ensure stability of soluble antigens or antigens with short half-lives. We recommend a fix and perm kit from reliable manufacturers. Follow manufacturer's instructions.

3. Add the primary antibody or isotype control at the appropriate dilution to the assay tubes. Incubate according to manufacturer's instructions.

4. Rinse cells twice by centrifugation with 2-3 mL incubation buffer.

5. Decant supernatant and re-suspend the pellet in remaining volume of wash.

6. Add fluorescent secondary antibody and incubate for 20-30 minutes. General range for secondary antibodies is between 1-10 μ g/mL for IgG conjugates for most applications.

7. Rinse cells twice by centrifugation with 2-3 mL incubation buffer. Centrifuge to collect cells after each wash. Decant supernatant.

8. Resuspend cells in 0.5 mL of diluent of choice to analyze on flow cytometer. Acquire data using the correct channel.