

APC/iFluor® 700 Goat Anti-human IgG (H+L) Antibody

Catalog Number: 50196

Unit Size: 200 ug

Product Details

| | |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze. |
| Expiration Date | 12 months upon receiving |
| Concentration | 0.2 mg/mL |
| Formulation | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA |

Antibody Properties

| | |
|--------------------|------------|
| Species Reactivity | Human |
| Class | Secondary |
| Clonality | Polyclonal |
| Host | Goat |

Biological Properties

| | |
|-----------------------|--|
| Stabilizer | 0.09% sodium azide, 0.2% (w/v) BSA |
| Appearance | Liquid |
| Preparation | Goat anti-human IgG (H+L) is produced in goat with pooled total human IgG. The antibody is conjugated with APC/iFluor® 700 under optimal conditions. |
| Application | Flow Cytometry (FACS), IF, IHC |
| Recommended Dilutions | Suggested dilutions are only guidelines; users should titrate the product for their specific assay using appropriate controls |

| Application | Recommended dilution |
|-----------------------|----------------------|
| Flow Cytometry (FACS) | 1-5 µg/mL |
| IF | 2 µg/mL |
| IHC | 1-10 µg/mL |

Spectral Properties

| | |
|-----------|-----------------|
| Conjugate | APC/iFluor™ 700 |
|-----------|-----------------|

Excitation Wavelength 651 nm

Emission Wavelength 710 nm

Applications

AAT Bioquest's anti-human secondary antibodies have well-characterized specificity for human immunoglobulins and are useful in the detection, sorting or purification of its specified target. This APC/iFluor® 700-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar APC/iFluor® 700 goat anti-human IgG antibodies from other commercial sources, and thus can significantly increase assay sensitivities. Secondary antibodies offer increased versatility enabling users to use many detection systems (e.g. HRP, AP, fluorescence). They can also provide greater sensitivity through signal amplification as multiple secondary antibodies can bind to a single primary antibody. This antibody was purified through affinity chromatography and conjugated to APC/iFluor® 700 (ex/em = 685/710 nm).