

mFluor™ Violet 510 goat anti-mouse IgG (H+L)

Catalog Number: 49050, 49051

Unit Size: 200 ug, 1 mg

Product Details

| | |
|--------------------|--|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze. |
| Expiration Date | 12 months upon receiving |
| Concentration | 1 mg/mL |
| Formulation | Phosphate-buffered saline (PBS, pH 7.2), 2 mg/mL BSA |

Unit Details

| | | |
|-----------------------|---------------------------|-------------------------|
| Reconstitution Volume | 49050 (200 ug) | 49051 (1 mg) |
| | 200 uL ddH ₂ O | 1 mL ddH ₂ O |

Antibody Properties

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

Chemical Properties

| | |
|------------------|----------|
| Molecular Weight | ~150 kDa |
|------------------|----------|

Biological Properties

| | |
|-----------------------|---|
| Stabilizer | 2 mg/mL BSA |
| Appearance | Solid |
| Preparation | Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG. The antibody is conjugated with mFluor™ Violet 510 under optimal conditions. |
| Application | Flow Cytometry (FACS), IF, IHC, ELISA, WB |
| 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 |
| ELISA | 100 ng/mL |
| WB | 1-10 µg/mL |

Spectral Properties

| | |
|-----------------------|--------------------|
| Conjugate | mFluor™ Violet 510 |
| Excitation Wavelength | 412 nm |
| Emission Wavelength | 505 nm |

Applications

mFluor™ Violet 510 goat anti-mouse conjugates are secondary antibodies designed for optimal performance in immunoassay applications, including flow cytometry, immunofluorescence, and confocal microscopy. These conjugates consist of goat-derived polyclonal antibodies with high affinity and specificity towards mouse IgG, conjugated to the bright and stable mFluor™ Violet 510 fluorochrome. This conjugation is optimized to ensure minimal non-specific binding and enhanced signal clarity, with rigorous purification steps to remove unconjugated components. Provided in a ready-to-use format with a recommended dilution range, the conjugate undergoes stringent quality control tests for performance and specificity. Its compatibility with a wide range of mouse primary antibodies and the exceptional contrast provided by mFluor™ Violet 510 fluorescence makes it a reliable tool for detecting diverse target antigens in multicolor staining protocols. mFluor™ Violet 510 is optimally excited by the violet laser and emits bright green fluorescence maximally at 505 nm. These affinity-purified goat anti-mouse secondary antibodies are valuable for their versatility and sensitivity, enabling efficient detection, sorting, or purification of specific targets through effective signal amplification in research applications.