

mFluor™ Red 780 goat anti-rabbit IgG (H+L)

Catalog Number: 49350, 49351

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 | 49350 (200 ug) | 49351 (1 mg) |
| | 200 uL ddH ₂ O | 1 mL ddH ₂ O |

Antibody Properties

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

Chemical Properties

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

Biological Properties

| | |
|-----------------------|--|
| Stabilizer | 2 mg/mL BSA |
| Appearance | Solid |
| Preparation | Goat anti-rabbit IgG (H+L) is produced in goat with pooled total rabbit IgG. The antibody is conjugated with mFluor™ Red 780 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™ Red 780 |
| Excitation Wavelength | 629 nm |
| Emission Wavelength | 767 nm |

Applications

mFluor™ Red 780 goat anti-rabbit 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 rabbit IgG, conjugated to the bright and stable mFluor™ Red 780 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 rabbit primary antibodies and the contrast provided by mFluor™ Red 780 fluorescence makes it a reliable tool for detecting diverse target antigens in multicolor staining protocols. mFluor™ Red 780 is optimally excited by the red laser and emits NIR fluorescence maximally at 767 nm. These affinity-purified goat anti-rabbit 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.