

## mFluor™ Red 780 goat anti-mouse IgG (H+L) \*Cross-Absorbed\*

Catalog Number: 49550, 49551

Unit Size: 200 ug, 1 mg

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid
Molecular Weight	~150 kDa
Soluble In	Water

### Spectral Properties

---

Excitation Wavelength	629 nm
Emission Wavelength	767 nm

### Applications

---

mFluor™ Red 780 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™ 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 mouse 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-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. To minimize cross-reactivity, these goat anti-mouse IgG whole antibodies have been cross-adsorbed against human, horse, mouse, and bovine IgG.