

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

Catalog Number: 49550, 49551

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	49550 (200 ug)	49551 (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™ 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-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.