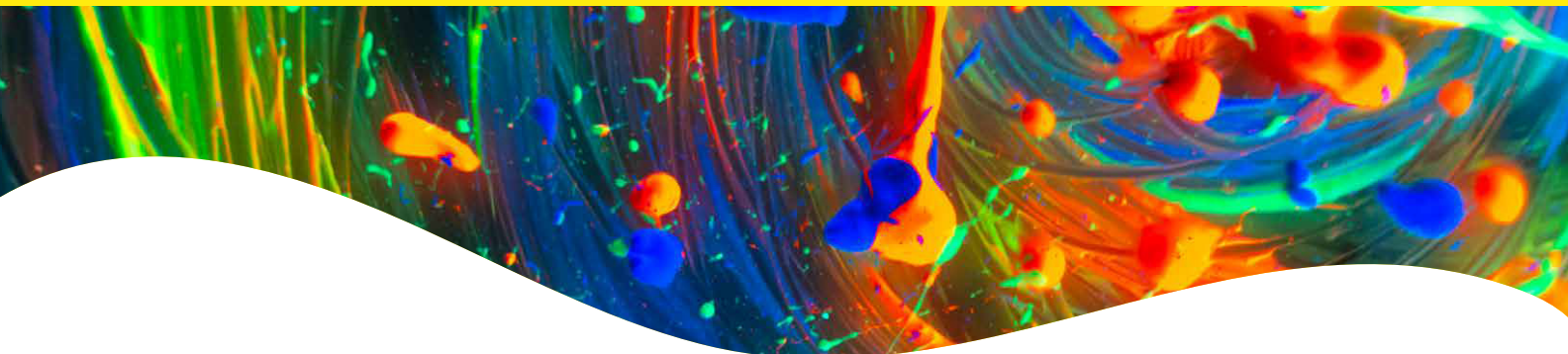


# PhenoVue Fluor - Goat anti-Mouse IgG (H+L) Antibody Conjugates

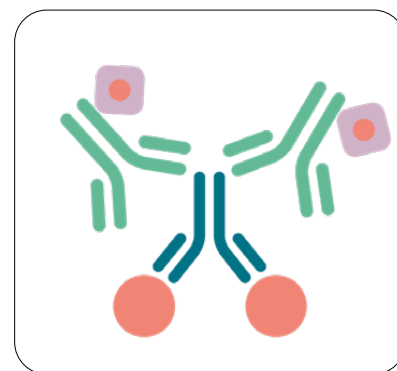


## Overview

Goat anti-mouse IgG (H+L) antibodies are conjugated with our bright PhenoVue™ Fluor dyes.

PhenoVue Fluor dyes Goat anti-mouse IgG (H+L) antibodies, Cross-adsorbed are affinity purified and recognize different mouse IgG isotypes and may display cross-reactivity with other species such as rat or hamster.

PhenoVue Fluor dyes Goat anti-mouse IgG (H+L) antibodies, Highly Cross-adsorbed have been adsorbed against various IgG species such as rat, hamster, bovine, horse, human, rabbit, or chicken to minimize cross-reactivity.



PhenoVue Fluor dyes

## Product information

Product name	Part no.	Numbers of vials per unit	Quantity per vial	Format	Shipping conditions
PhenoVue Fluor 488 - Goat anti-Mouse antibody Cross-adsorbed	2GXM488C1	1	1 mg	Lyophilized	RT
PhenoVue Fluor 555 - Goat anti-Mouse antibody Cross-adsorbed	2GXM555C1				
PhenoVue Fluor 568 - Goat anti-Mouse antibody Cross-adsorbed	2GXM568C1				
PhenoVue Fluor 594 - Goat anti-Mouse antibody Cross-adsorbed	2GXM594C1				
PhenoVue Fluor 647 - Goat anti-Mouse antibody Cross-adsorbed	2GXM647C1				
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed	2GXM488H1	1	1 mg	Lyophilized	RT
PhenoVue Fluor 555 - Goat anti-Mouse antibody Highly Cross-adsorbed	2GXM555H1				
PhenoVue Fluor 568 - Goat anti-Mouse antibody Highly Cross-adsorbed	2GXM568H1				
PhenoVue Fluor 594 - Goat anti-Mouse antibody Highly Cross-adsorbed	2GXM594H1				
PhenoVue Fluor 647 - Goat anti-Mouse antibody Highly Cross-adsorbed	2GXM647H1				

## Storage and stability

- Store lyophilized reagents at 2-8 °C, protected from light.
- The stability of these products is guaranteed until the expiration date provided in the Certificate of Analysis, when stored as recommended and protected from light.
- Allow the powder to warm up to room temperature for 15 min before opening the vials and reconstitution.
- After reconstitution, aliquoted reagents must be stored at -16 °C or below and are stable for 6 months. Avoid repeated freeze / thaw cycles.

## Recommended reconstitution

Product name	Molecular weight	Recommended stock concentration	Working concentration range*
PhenoVue Fluor 488 - Goat anti-Mouse antibody Cross-adsorbed	150000 g/mol	Reconstitution using 1 mL ddH <sub>2</sub> O gives a stock concentration of 1 mg/mL (6.66 µM)	0.1 µg/mL – 10 µg/mL (0.66 nM – 66.6 nM)
PhenoVue Fluor 555 - Goat anti-Mouse antibody Cross-adsorbed			
PhenoVue Fluor 568 - Goat anti-Mouse antibody Cross-adsorbed			
PhenoVue Fluor 594 - Goat anti-Mouse antibody Cross-adsorbed			
PhenoVue Fluor 647 - Goat anti-Mouse antibody Cross-adsorbed			
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed	150000 g/mol	Reconstitution using 1 mL ddH <sub>2</sub> O gives a stock concentration of 1 mg/mL (6.66 µM)	0.1 µg/mL – 10 µg/mL (0.66 nM – 66.6 nM)
PhenoVue Fluor 555 - Goat anti-Mouse antibody Highly Cross-adsorbed			
PhenoVue Fluor 568 - Goat anti-Mouse antibody Highly Cross-adsorbed			
PhenoVue Fluor 594 - Goat anti-Mouse antibody Highly Cross-adsorbed			
PhenoVue Fluor 647 - Goat anti-Mouse antibody Highly Cross-adsorbed			

\* Dilutions can be done in PBS.

## Equivalent number of microplates

Product name	When used at recommended concentration	96-Well microplate (100 µL - 300 µL per well)	384-well microplate (25 µL - 90 µL per well)	1536-well microplate (4 µL - 12 µL per well)
PhenoVue Fluor 488 - Goat anti-Mouse antibody Cross-adsorbed	1.5 µg/mL (10 nM)	Approx. 25-70	Approx. 20-70	Approx. 35-90
PhenoVue Fluor 555 - Goat anti-Mouse antibody Cross-adsorbed				
PhenoVue Fluor 568 - Goat anti-Mouse antibody Cross-adsorbed				
PhenoVue Fluor 594 - Goat anti-Mouse antibody Cross-adsorbed				
PhenoVue Fluor 647 - Goat anti-Mouse antibody Cross-adsorbed				
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed	1.5 µg/mL (10 nM)	Approx. 25-70	Approx. 20-70	Approx. 35-90
PhenoVue Fluor 555 - Goat anti-Mouse antibody Highly Cross-adsorbed				
PhenoVue Fluor 568 - Goat anti-Mouse antibody Highly Cross-adsorbed				
PhenoVue Fluor 594 - Goat anti-Mouse antibody Highly Cross-adsorbed				
PhenoVue Fluor 647 - Goat anti-Mouse antibody Highly Cross-adsorbed				

View our full range of high-quality imaging microplates at [Revvity.com](http://Revvity.com)

## Spectral and photophysical properties

Product name	Maximum excitation wavelength (nm)	Maximum emission wavelength (nm)	Common filter set	Quantum yield ( $\Phi$ )	Epsilon* ( $\epsilon$ in $M^{-1}.cm^{-1}$ at $\lambda$ . max)	Brightness ( $\Phi \times \epsilon$ )
PhenoVue Fluor 488	495	520	FITC	92%	73000	65320
PhenoVue Fluor 555	555	570	Cy3	10%	155000	15500
PhenoVue Fluor 568	578	603	Texas-Red	69%	88000	60720
PhenoVue Fluor 594	590	617	Texas-Red	66%	92000	60720
PhenoVue Fluor 647	650	670	Cy5	30%	240000	72000

\* In methanol

## Cross-reactivity

Product name	Across species	Across IgG isotypes
PhenoVue Fluor 488 - Goat anti-Mouse antibody Cross-adsorbed	Rat: 44% Guinea Pig: 16% Hamster: 50% Horse: 10%	Cross-reactivity with mouse: IgG1 IgG2a IgG2b IgG2c IgG3
PhenoVue Fluor 555 - Goat anti-Mouse antibody Cross-adsorbed		
PhenoVue Fluor 568 - Goat anti-Mouse antibody Cross-adsorbed		
PhenoVue Fluor 594 - Goat anti-Mouse antibody Cross-adsorbed		
PhenoVue Fluor 647 - Goat anti-Mouse antibody Cross-adsorbed		
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed	None	Cross-reactivity with mouse: IgG1 IgG2a IgG2b IgG2c IgG3
PhenoVue Fluor 555 - Goat anti-Mouse antibody Highly Cross-adsorbed		
PhenoVue Fluor 568 - Goat anti-Mouse antibody Highly Cross-adsorbed		
PhenoVue Fluor 594 - Goat anti-Mouse antibody Highly Cross-adsorbed		
PhenoVue Fluor 647 - Goat anti-Mouse antibody Highly Cross-adsorbed		

## Protocols

### Cell Culture

Seed cells in imaging microplates (or any other convenient cell culture vessels). Incubate in the appropriate cell culture conditions, usually 37 °C, 5% CO<sub>2</sub> until 50-70% confluency.

### Fixed-Cell Imaging

Rinse briefly in phosphate-buffered saline (PBS) then proceed with cell fixation.

#### 1. Fixation: 2 options:

- Add ready to use PhenoVue Paraformaldehyde 4% Methanol-Free Solution (PVPFA41) for 10 min at room temperature. Note that paraformaldehyde (PFA) is the most popular fixative reagent.

or

- Add 100% methanol (chilled to -20 °C) at room temperature for 5 min.

#### 2. Washing: Wash three times with PBS.

#### 3. Permeabilization:

- For PFA fixed cells, add ready to use PhenoVue Permeabilization 0.5% Triton X-100 Solution (PVPERM051) for 10 min (for membrane-associated antigens, 100  $\mu$ M digitonin or 0.5% saponin are preferred). Triton X-100 is the most popular detergent for improving the penetration of antibodies. However, it may not be appropriate for some imaging applications since it can destroy membranes.
- Methanol fixed cells do not require permeabilization.

#### 4. Washing: Wash three times with PBS for 5 min.

#### 5. Blocking step: Incubate with PBS + 1% BSA for 60 min at RT.

#### 6. Primary antibody: Incubate with a primary mouse antibody.

#### 7. Washing: Wash three times with PBS for 5 min.

#### 8. Staining: Incubate with 0.1-10 $\mu$ g/mL PhenoVue Fluor Goat anti-Mouse antibody Cross-adsorbed or Highly Cross-adsorbed for 60 min at RT.

9. **Washing:** Wash three times with PBS for 5 min.
10. **Optional:** Incubate with 0.1-2 µg/mL PhenoVue Hoechst 33342 nuclear stain for 10 min.
11. **Washing:** Wash once with PBS for 5 min.
12. Acquire images on an imaging device.

## Tips

- Due to species cross-reactivity of PhenoVue Goat anti-Mouse cross-adsorbed antibodies, especially with rodents, it is preferable to use PhenoVue Goat anti-Mouse highly cross-adsorbed antibodies when performing multiplexing experiments including different primary antibodies (see Figures 4, 5 and 6). Please note that this is not limited to PhenoVue secondary antibodies but rather a general characteristic of antibodies irrespective of the vendor.
- PhenoVue Goat anti-Mouse cross-adsorbed antibodies are well suited for single-plex experiments.

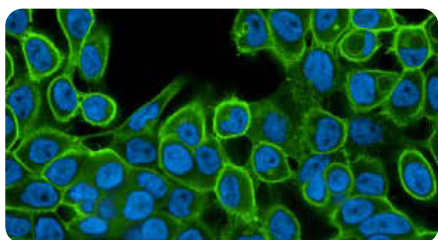
## Safety information

Chemical reagents are potentially harmful, please refer to the Safety Data Sheet (SDS) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Applications

- High-content analysis / high-content screening
- Imaging microscopy
- Flow cytometry

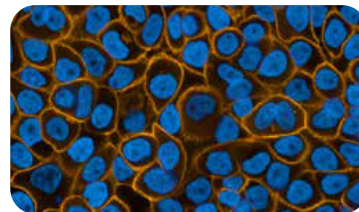
## Validation data



PhenoVue Fluor 488 - Goat anti-Mouse IgG (H+L)  
Cross-adsorbed

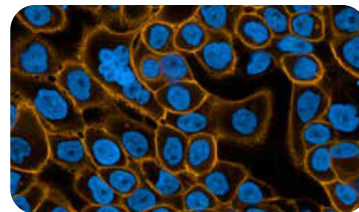
Figure 1: A431 cells were seeded in PhenoPlate™ 96-well microplates (75,000 cells/well) and incubated at 37 °C, 5% CO<sub>2</sub> for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 488 - Goat anti-Mouse IgG (H+L) Cross-adsorbed for 1 hour at RT. Nuclei were stained with 2 µg/mL PhenoVue Hoechst 33342 nuclear stain. Images were acquired on the Operetta CLS™ high-content analysis system.

A



PhenoVue Fluor 555 - Goat anti-Mouse IgG (H+L)  
Cross-adsorbed

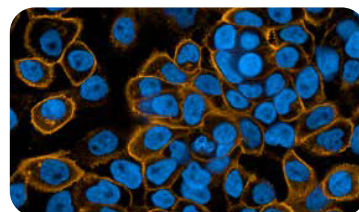
B



PhenoVue Fluor 555 -Goat anti-Mouse IgG (H+L)  
Highly Cross-adsorbed

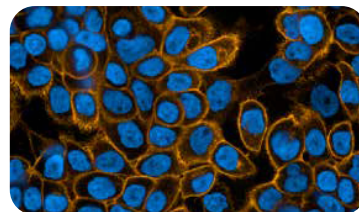
Figure 2: A431 cells were seeded in PhenoPlate 96-well microplates (75,000 cells/well) and incubated at 37 °C, 5% CO<sub>2</sub> for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 555 - Goat anti-Mouse IgG (H+L) Cross-adsorbed (A) or Highly Cross-adsorbed (B) for 1 hour at RT. Nuclei were stained with 2 µg/mL PhenoVue Hoechst 33342 nuclear stain. Images were acquired on the Operetta CLS high-content analysis system.

A



PhenoVue Fluor 568 - Goat anti-Mouse IgG (H+L)  
Cross-adsorbed

B



PhenoVue Fluor 568 Goat anti-Mouse IgG (H+L)  
Cross-adsorbed

Figure 3: A431 cells were seeded in PhenoPlate 96-well microplates (75,000 cells/well) and incubated at 37 °C, 5% CO<sub>2</sub> for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 568 - Goat anti-Mouse IgG (H+L) Cross-adsorbed (A) or Highly Cross-adsorbed (B) for 1 hour at RT. Nuclei were stained with 2 µg/mL PhenoVue Hoechst 33342 nuclear stain. Images were acquired on the Operetta CLS high-content analysis system.

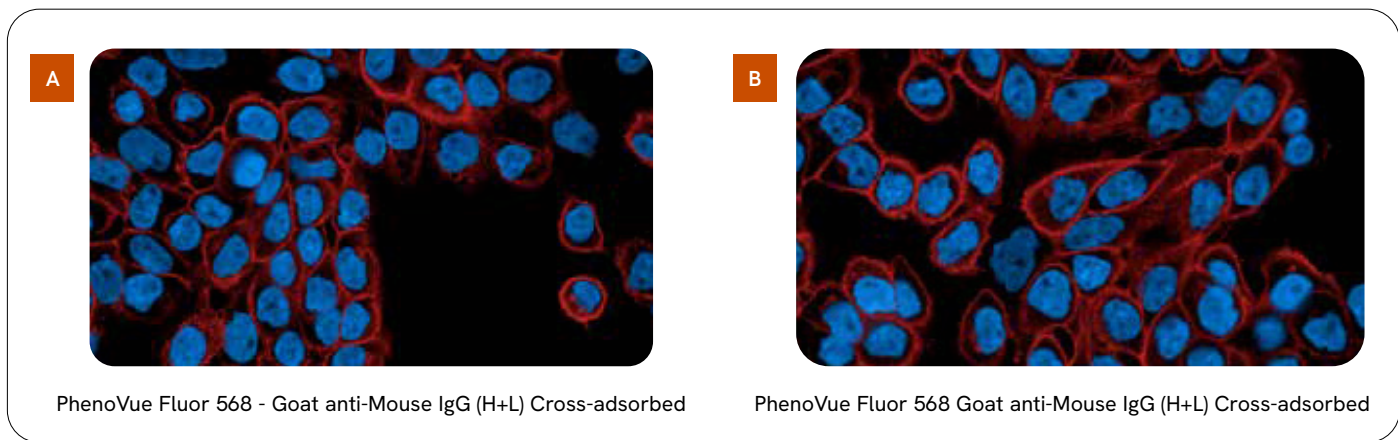


Figure 4: A431 cells were seeded in PhenoPlate 96-well microplates (75,000 cells/well) and incubated at 37°C, 5% CO<sub>2</sub> for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 594 - Goat anti-Mouse IgG (H+L) Cross-adsorbed (A) or Highly Cross-adsorbed (B) for 1 hour at RT. Nuclei were stained with 2 µg/mL PhenoVue Hoechst 33342 nuclear stain. Images were acquired on the Operetta CLS high-content analysis system.

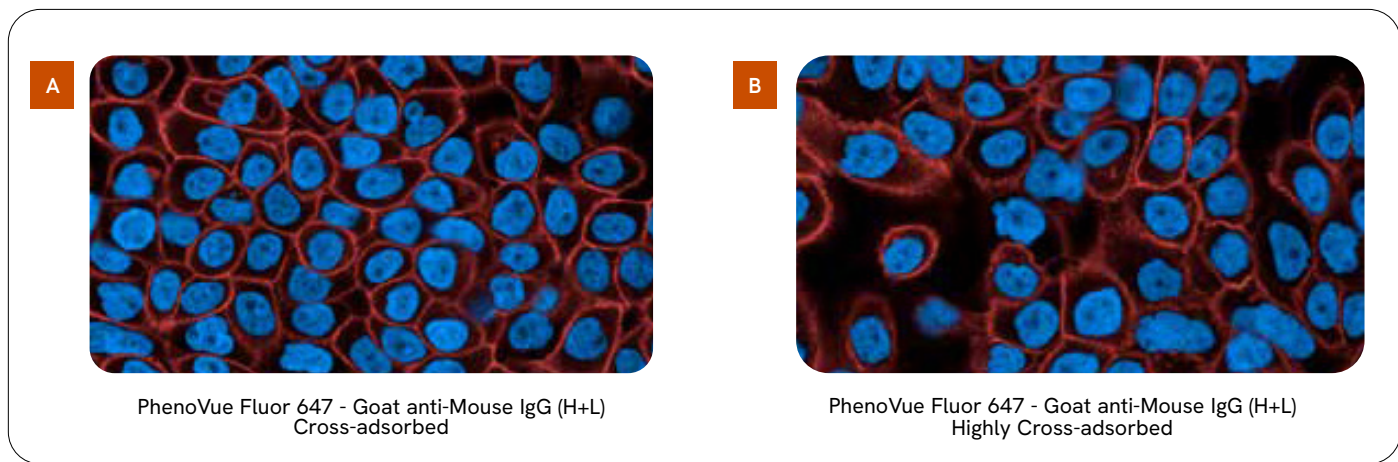
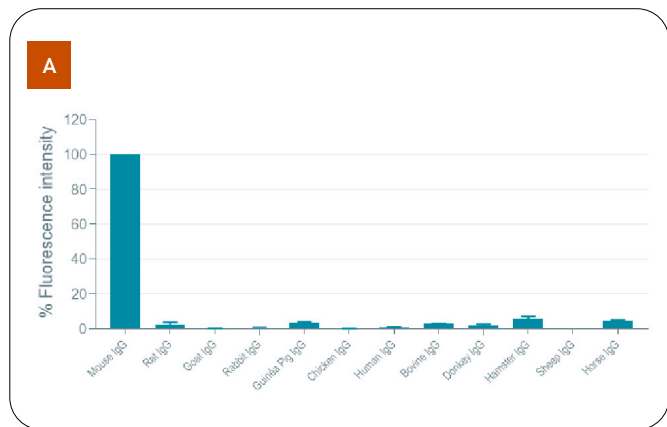
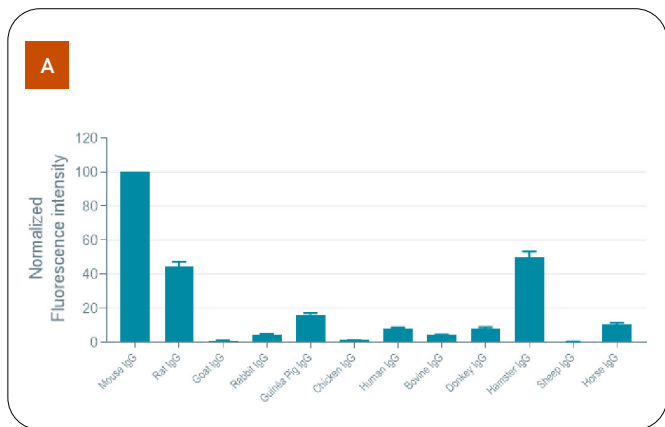


Figure 5: A431 cells were seeded in PhenoPlate 96-well microplates (75,000 cells/well) and incubated at 37°C, 5% CO<sub>2</sub> for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 647 - Goat anti-Mouse IgG (H+L) Cross-adsorbed (A) or Highly Cross-adsorbed (B) for 1 hour at RT. Nuclei were stained with 2 µg/mL PhenoVue Hoechst 33342 nuclear stain. Images were acquired on the Operetta CLS high-content analysis system.



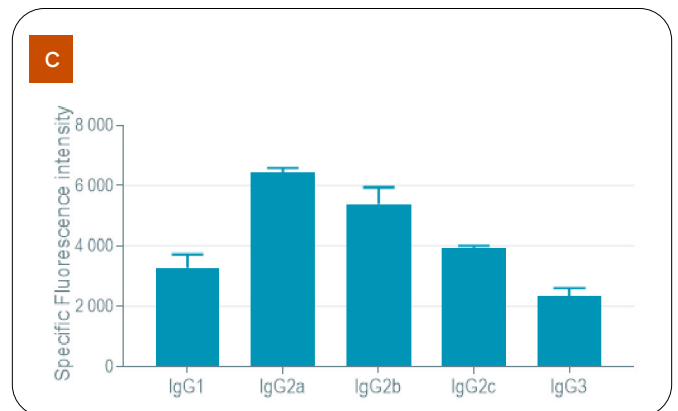
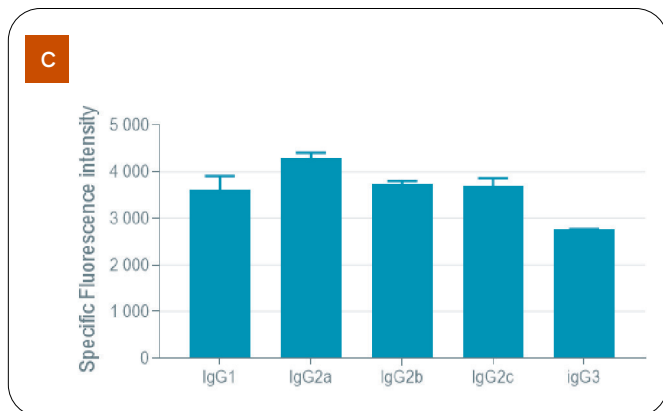
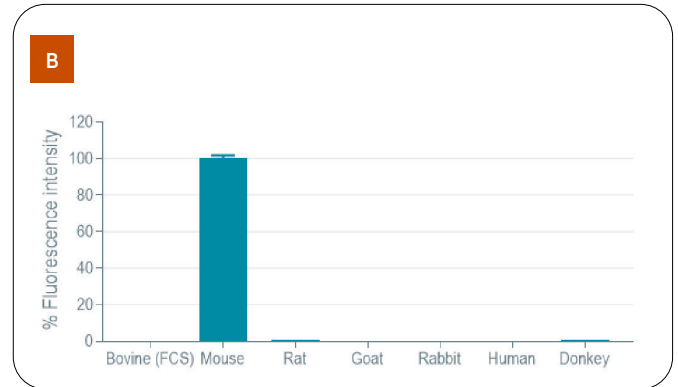
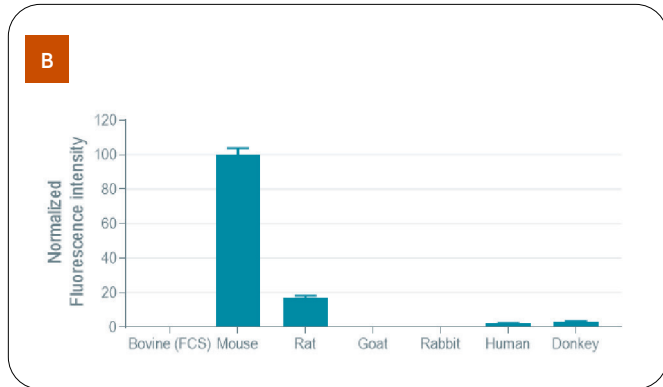
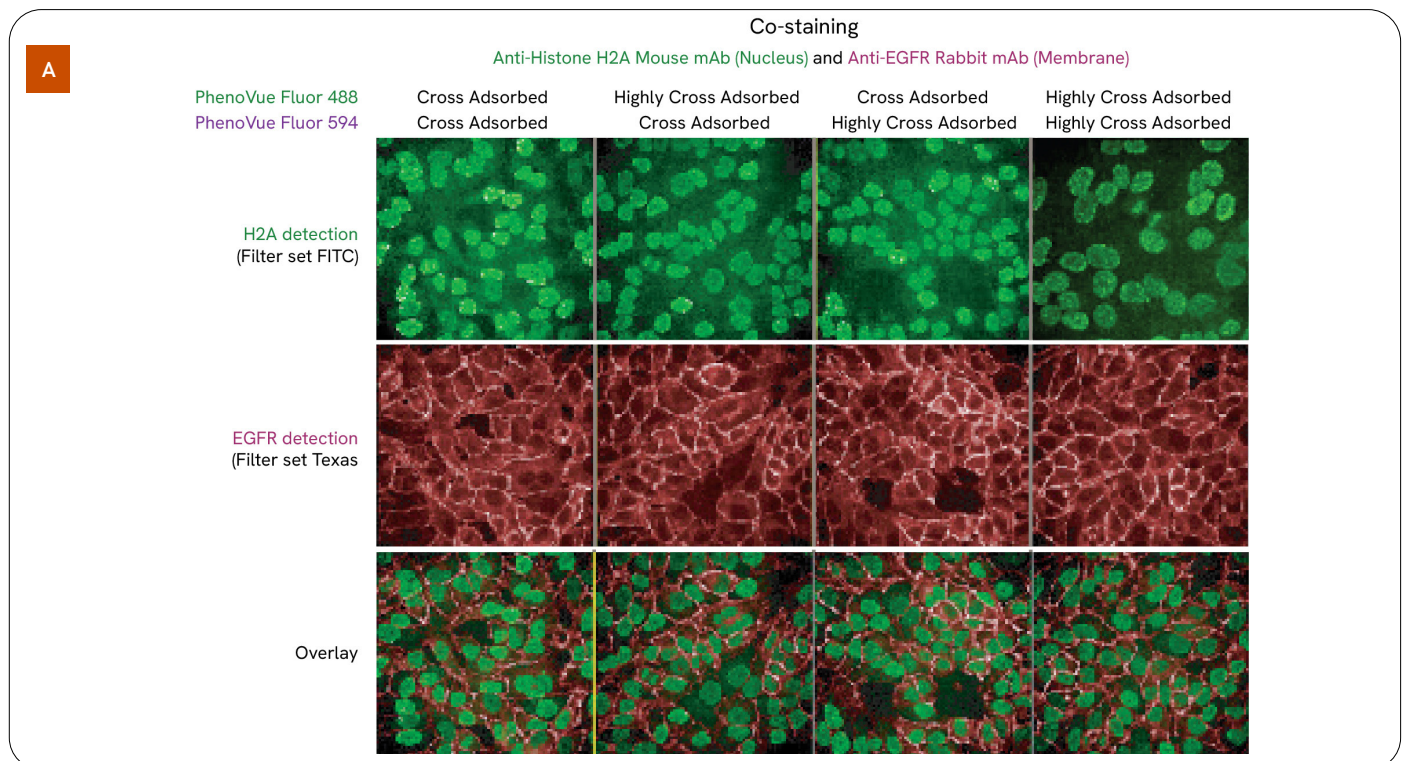


Figure 6.1: F-LISA experiments: different IgG species (A), 10% of serum (B) or IgG isotypes (C) were used to coat a 96-well microplate, then incubated with PhenoVue Fluor 594 - Goat anti-Mouse IgG (H+L) Cross-Adsorbed (1.5 µg/mL). Fluorescence intensity was measured on an EnVision™ multimode plate reader.

Figure 6.2: F-LISA experiments: different IgG species (A), 10% of serum (B) or IgG isotypes (C) were used to coat a 96-well microplate, then incubated with PhenoVue Fluor 594 - Goat anti-Mouse IgG (H+L) Highly Cross-Adsorbed (1.5 µg/mL). Fluorescence Intensity was measured on an EnVision multimode plate reader.

No Cross reactivity of PhenoVue Fluor 488 - Goat anti-Mouse Cross-Adsorbed with primary Rabbit antibody



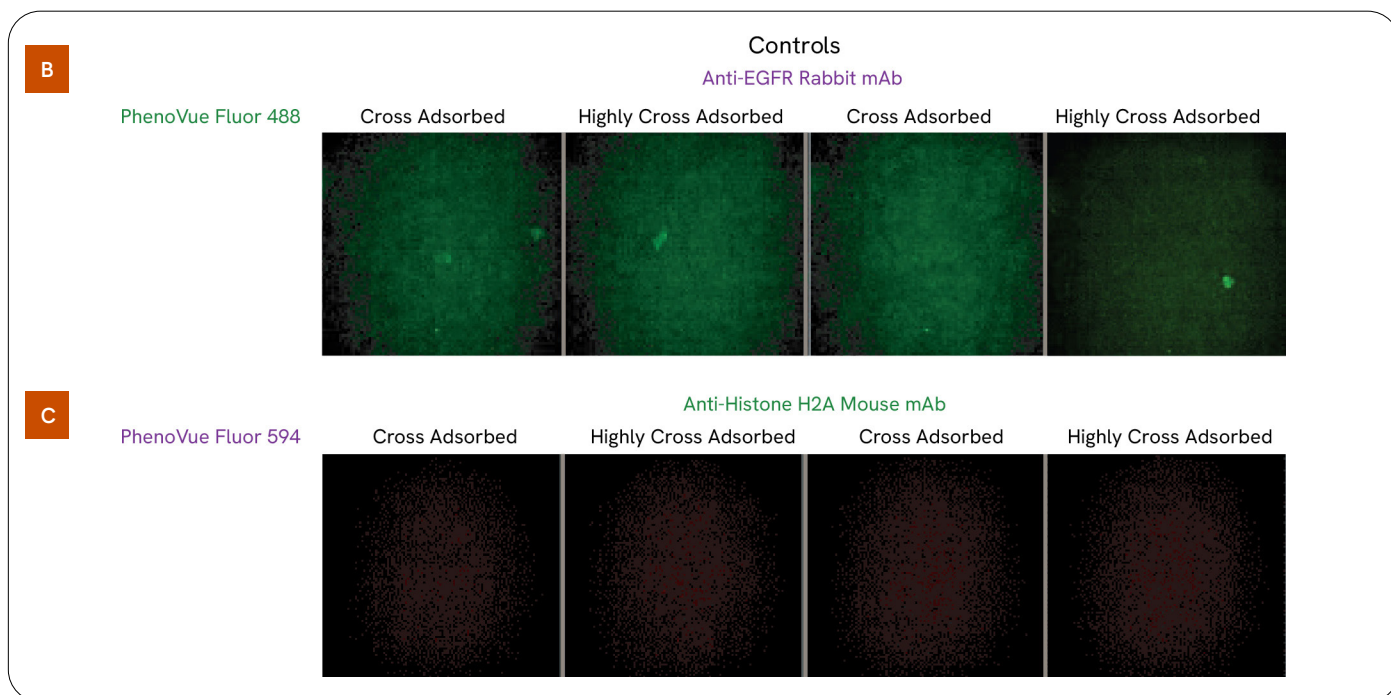
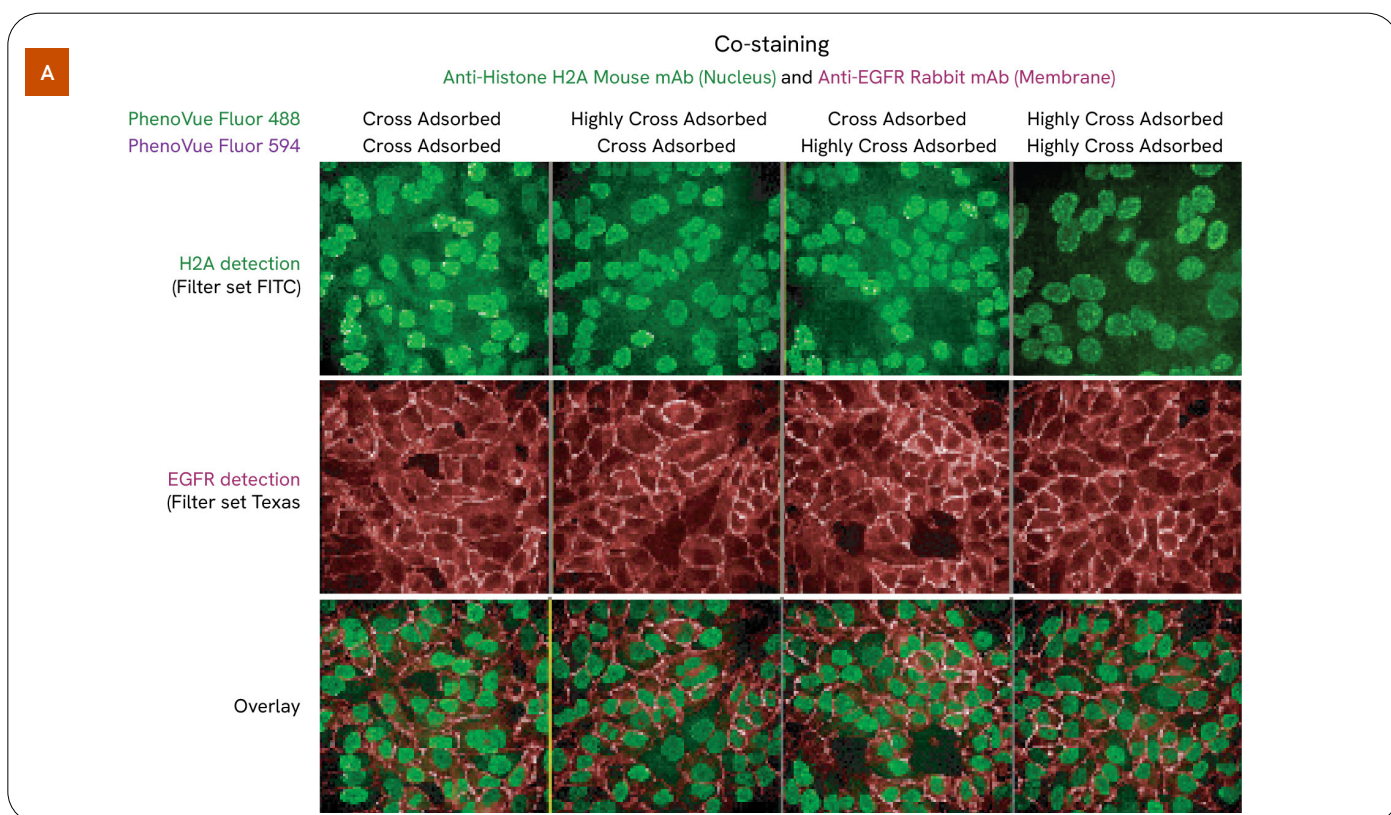


Figure 7: HeLa cells were seeded in PhenoPlate 96-well microplates (50,000 cells/well) and incubated at 37 °C, 5% CO<sub>2</sub> for 24h. Cells were fixed then permeabilized and co-incubated with anti- Histone H2A mouse antibody (5 µg/mL) and anti-EGFR rabbit antibody (0.5 µg/mL). After washing steps, cells were incubated with 1.5 µg/mL PhenoVue Fluor 488 - Goat anti-Mouse IgG (H+L) Cross-Adsorbed or Highly Cross-adsorbed and PhenoVue Fluor 594 - Goat anti-Rabbit IgG (H+L) Cross-Adsorbed or Highly Cross-adsorbed for 1 hour at RT (Panel A). Cross reactivity of PhenoVue Fluor 488 - Goat anti-Mouse antibody either Cross-Adsorbed or Highly Cross-adsorbed was assessed by incubating anti-EGFR rabbit antibody only (Panel B). Cross reactivity of PhenoVue Fluor 594 - Goat anti-Rabbit antibody either Cross-Adsorbed or Highly Cross-adsorbed was assessed by incubating anti-Histone H2A mouse antibody only (Panel C). Images were acquired on the Operetta CLS high-content analysis system.

No Cross reactivity of PhenoVue Fluor 488 - Goat anti-Mouse Highly Cross-Adsorbed with primary Rat antibody



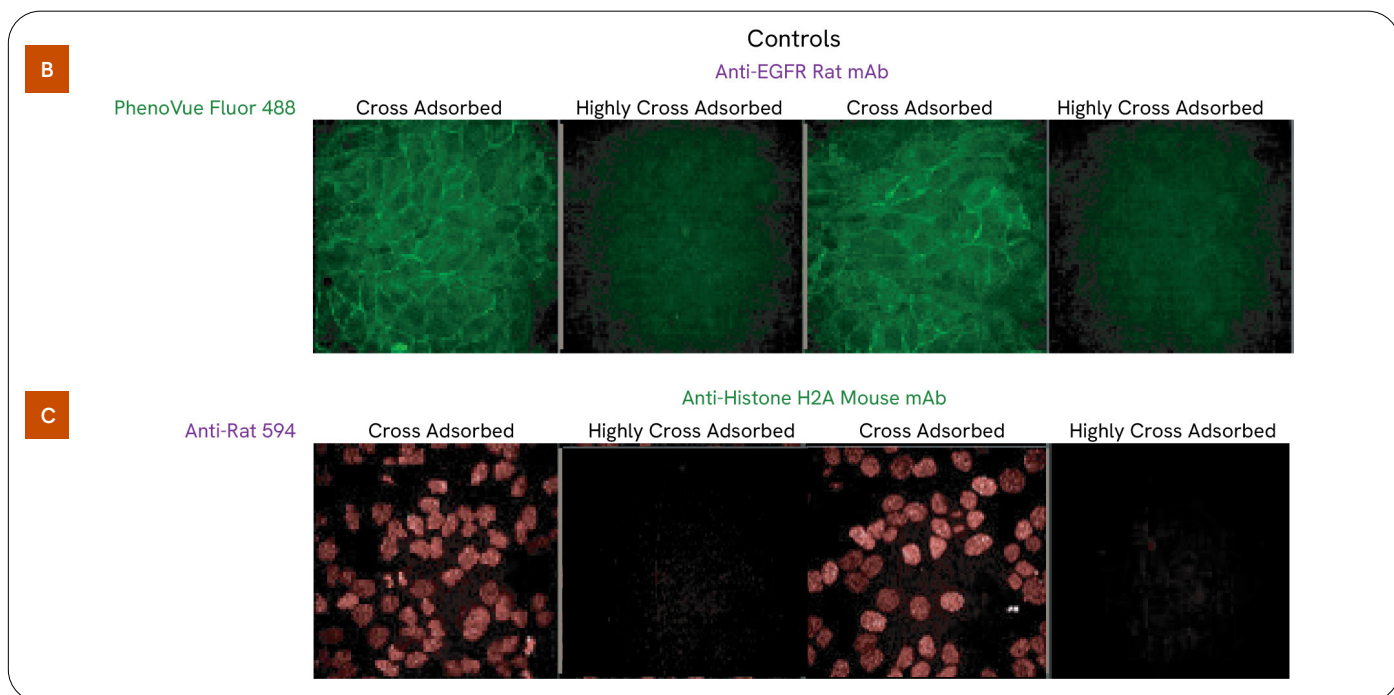


Figure 8: HeLa cells were seeded in PhenoPlate 96-well microplates (50,000 cells/well) and incubated at 37 °C, 5% CO<sub>2</sub> for 24h. Cells were fixed then permeabilized and co incubated with anti- Histone H2A mouse antibody (5 µg/mL) and anti-EGFR Rat antibody (1 µg/mL). After washing steps, cells were incubated with 1.5 µg/mL PhenoVue Fluor 488 - Goat anti-Mouse IgG (H+L) Cross-adsorbed or Highly Cross-adsorbed and Anti-Rat 594 Cross-adsorbed or Highly Cross-adsorbed for 1 hour at RT (Panel A). Cross reactivity of PhenoVue Fluor 488 - Goat anti-Mouse antibody either Cross-adsorbed or Highly Cross-adsorbed was assessed by incubating anti-EGFR rabbit antibody only (Panel B). Cross reactivity of Anti-Rat 594 either Cross-Adsorbed or Highly Cross-adsorbed was assessed by incubating anti-Histone H2A mouse antibody only (Panel C). Anti-Rat 594 is a fluor-labelled goat anti-rat second antibody with a maximum excitation wavelength of 594 nm, and maximum emission wavelength of 617 nm. Images were acquired on the Operetta CLS high-content analysis system.

