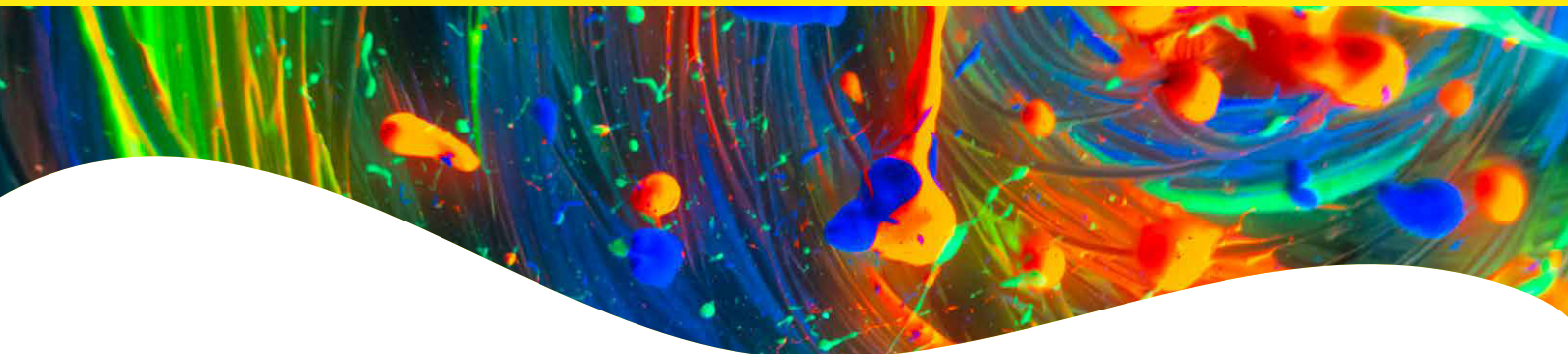


PhenoVue Fluor - Goat anti-Mouse IgG (H+L) Antibody Conjugates

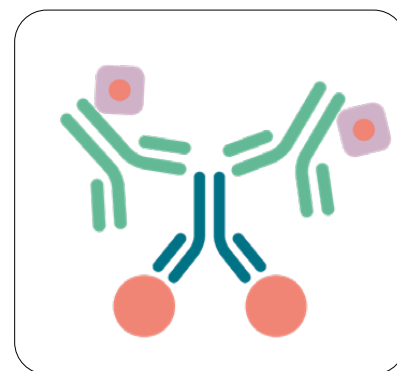


Overview

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 405 - Goat anti-Mouse antibody Highly Cross-adsorbed	2GXM405H1	1	1 mg	Lyophilized	RT
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed	2GXM488H1				
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 ₂ O gives a stock concentration of 1 mg/mL (6.67 µM)	1 µg/mL - 10 µg/mL (6.7 nM - 66.7 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 405 - Goat anti-Mouse antibody Highly Cross-adsorbed	150000 g/mol	Reconstitution using 1 mL ddH ₂ O gives a stock concentration of 1 mg/mL (6.67 µM)	1 µg/mL - 10 µg/mL (6.7 nM - 66.7 nM)
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed			
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	2.5 µg/mL (16.7 nM)	Approx. 14-42	Approx. 14-42	Approx. 22-65
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 405 - Goat anti-Mouse antibody Highly Cross-adsorbed	5 µg/mL (32.3 nM)	Approx. 7-21	Approx. 5.5-20	Approx. 11-33
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed	2.5 µg/mL (16.7 nM)	Approx. 14-42	Approx. 14-42	Approx. 22-65
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

Spectral and photophysical properties

Product name	Maximum excitation wavelength (nm)	Maximum emission wavelength (nm)	Common filter set	Quantum yield (Φ)	Epsilon* (ϵ in $M^{-1}\cdot cm^{-1}$ at λ max)	Brightness ($\Phi \times \epsilon$)
PhenoVue Fluor 405	410	452	Ex: 375-430 nm Em: 440-500 nm	78%	46000	35880
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

Recommended Acquisition settings for Revvity HCS instruments

HCS Instruments		PhenoVue Fluor 405	PhenoVue Fluor 400LS	PhenoVue Fluor 488	PhenoVue Fluor 555	PhenoVue Fluor 568 Or PhenoVue Fluor 594	PhenoVue Fluor 647
Opera Phenix Plus 5 lasers	Excitation laser	425	425	488	561	561	640
	Emission filter	435-480	570-630	500-550	570-630	570-630	650-760
Opera Phenix Plus 4 lasers	Excitation laser	405	405	488	561	561	640
	Emission filter	435-480	570-630	500-550	570-630	570-630	650-760
Operetta CLS 8 LED - 1600	Excitation LED (filter) (nm)	405 (390-420)	405 (390-420)	475 (460-490)	550 (530-560)	550 (530-560)	630 (615-645)
	Emission filter	430-500	570-650	500-550	570-650	570-650	655-760
Operetta CLS 8 LED - 1601	Excitation LED (filter) (nm)	370 (355-385)	440 (435-460)	475 (460-490)	550 (530-560)	580 (560-575)	630 (615-645)
	Emission filter	430-500	600-640 or 570-650	500-550	570-650	585-640	655-760
Operetta CLS 4 LED	Excitation LED (filter) (nm)	370 (355-385)	370 (355-385)	475 (460-490)	550 (530-560)	550 (530-560)	630 (615-645)
	Emission filter	430-500	570-650	500-550	570-650	570-650	655-760

Cross-reactivity

Product name	Across species	Across IgG isotypes
PhenoVue Fluor 488 - Goat anti-Mouse antibody Cross-adsorbed	Mouse:100% 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 405 - Goat anti-Mouse antibody Highly Cross-adsorbed	Mouse:100%	Cross-reactivity with mouse: IgG1 IgG2a IgG2b IgG2c IgG3
PhenoVue Fluor 488 - Goat anti-Mouse antibody Highly Cross-adsorbed		
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% CO₂ until 50-70% confluency.

Fixed-Cell Imaging

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

1. Fixation: 2 options:

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

or

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

2. Washing: Wash three times with PBS.

3. Permeabilization:

1. For PFA fixed cells, add ready to use PhenoVue Permeabilization 0.5% Triton X-100 Solution (PVPERM051) for at least 10 min (for membrane-associated antigens, 100 µM 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.
2. 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 1-10 µg/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 10min or 2-10 µM PhenoVue DRAQ5 Total Cell Nuclear Stain for 15-60 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.

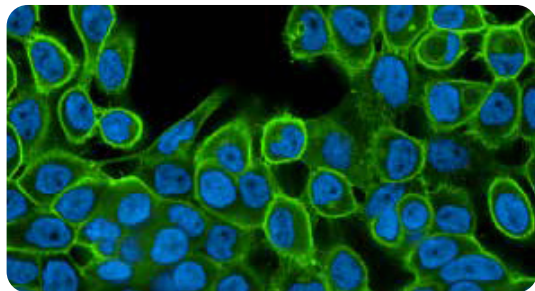
Safety information

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

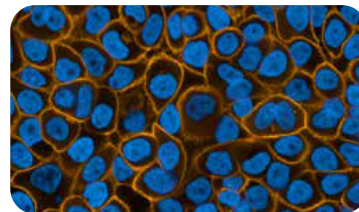
Validation data



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

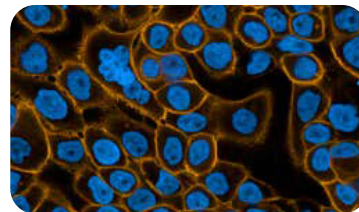
Figure 1: A431 cells were seeded in PhenoPlate™ 96-well microplates (75,000 cells/well) and incubated at 37 °C, 5% CO₂ for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 488 - 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

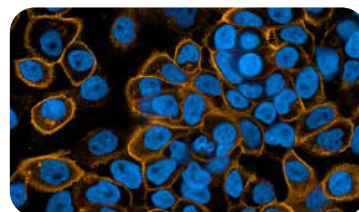
B



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

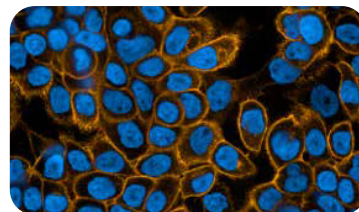
Figure 2: A431 cells were seeded in PhenoPlate 96-well microplates (75,000 cells/well) and incubated at 37 °C, 5% CO₂ for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 555 - 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% CO₂ 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.

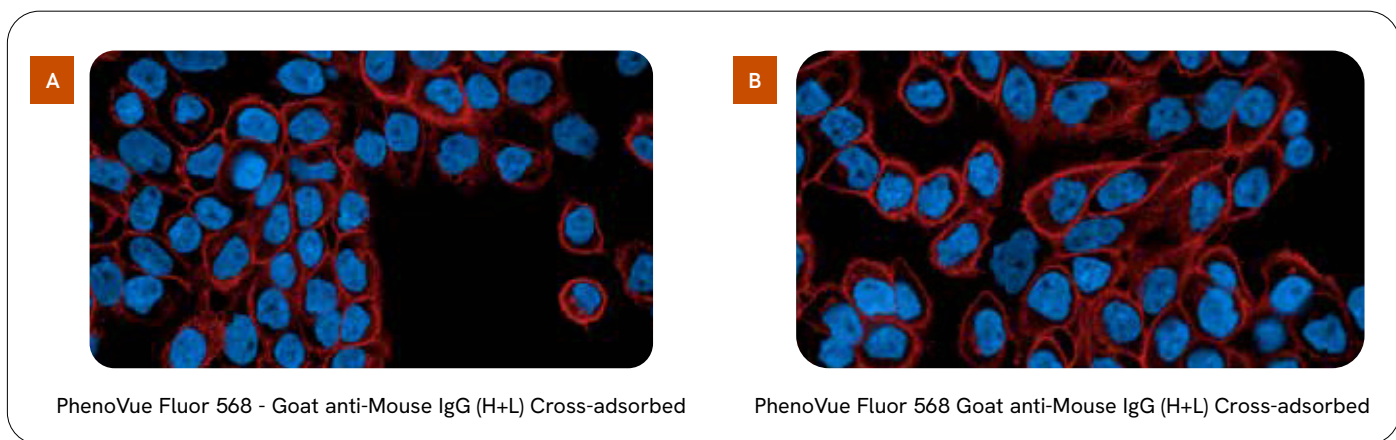


Figure 4: A431 cells were seeded in PhenoPlate 96-well microplates (75,000 cells/well) and incubated at 37°C, 5% CO₂ for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 594 - 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.

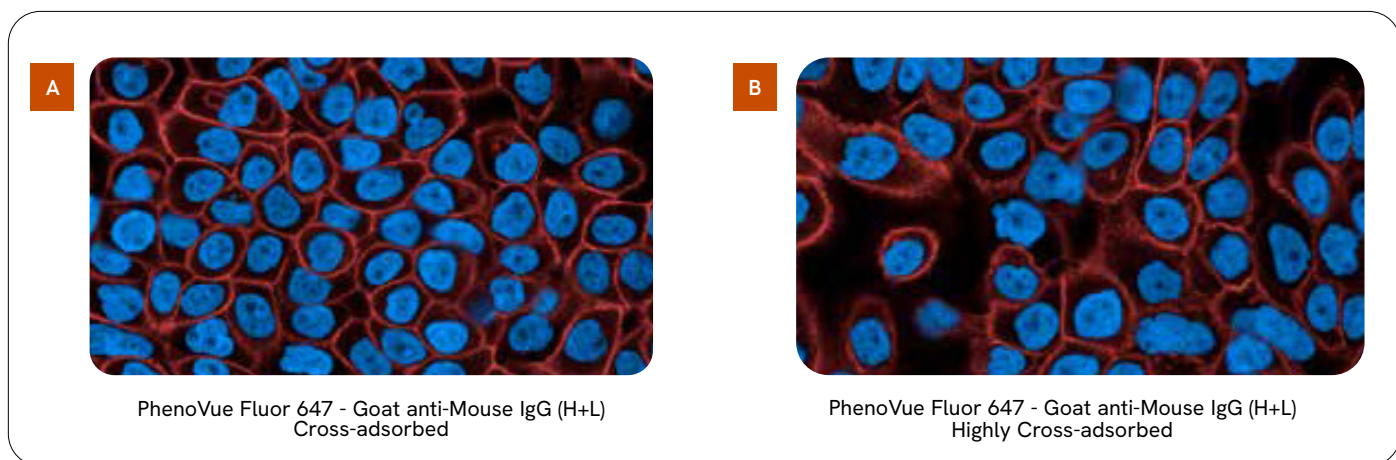


Figure 5: A431 cells were seeded in PhenoPlate 96-well microplates (75,000 cells/well) and incubated at 37°C, 5% CO₂ for 24h. Cells were fixed then permeabilized and incubated with an anti-EGFR Mouse antibody (0.2 µg/mL). After washing steps, cells were incubated with 10 µg/mL of PhenoVue Fluor 647 - 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.

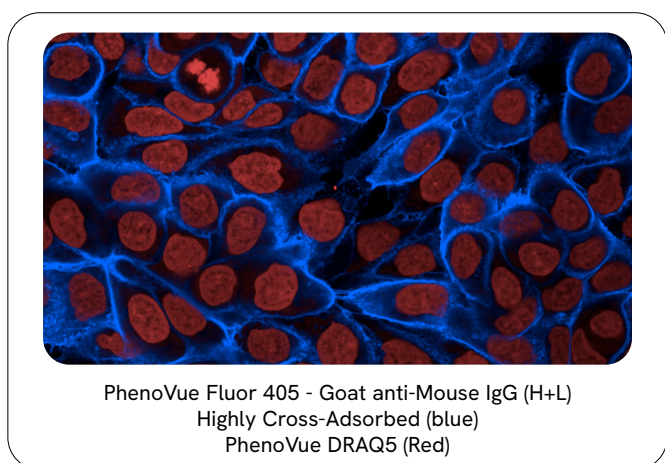


Figure 6: A431 cells were seeded in PhenoPlate™ 96-well microplates at a density of 75,000 cells per well and incubated for 24 hours at 37°C with 5% CO₂. After incubation, the cells were fixed, permeabilized, and treated with a 2 µg/mL concentration of anti-EGFR mouse antibody. Following washing steps, the cells were incubated with 10 µg/mL PhenoVue Fluor 405 - Goat anti-Mouse IgG (H+L), highly cross-adsorbed, for 1 hour at room temperature. The nuclei were stained with 1 µM PhenoVue DRAQ5 Total Cell Nuclear Stain. Images were captured using the Opera Phenix Plus high-content analysis system with 5 lasers, using a 63x water objective in confocal mode.

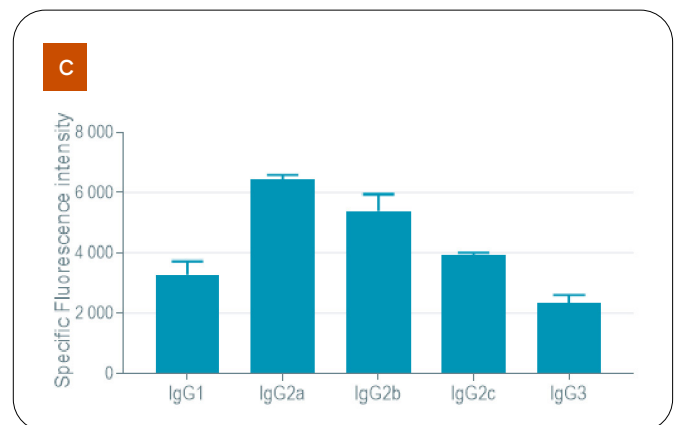
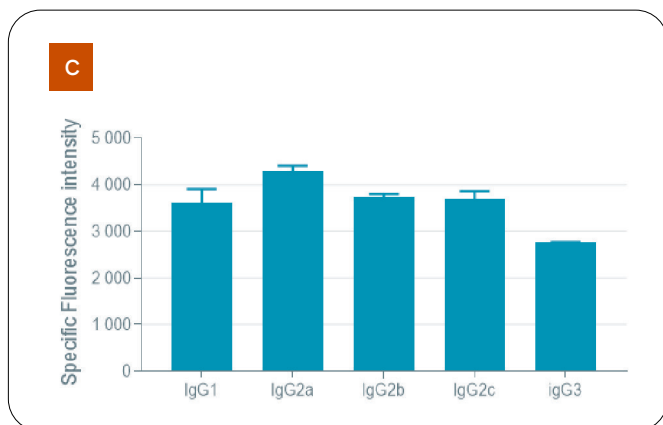
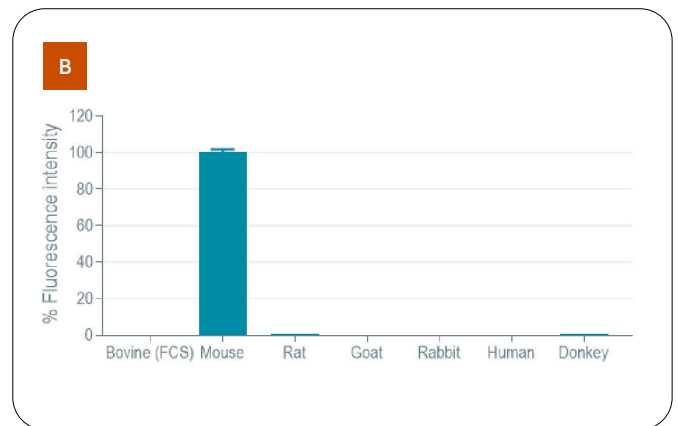
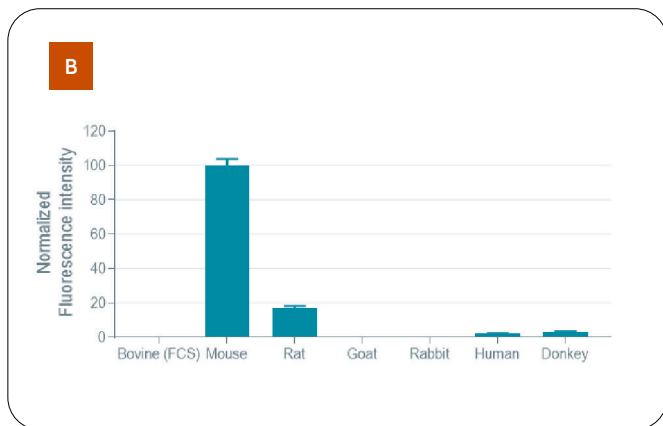
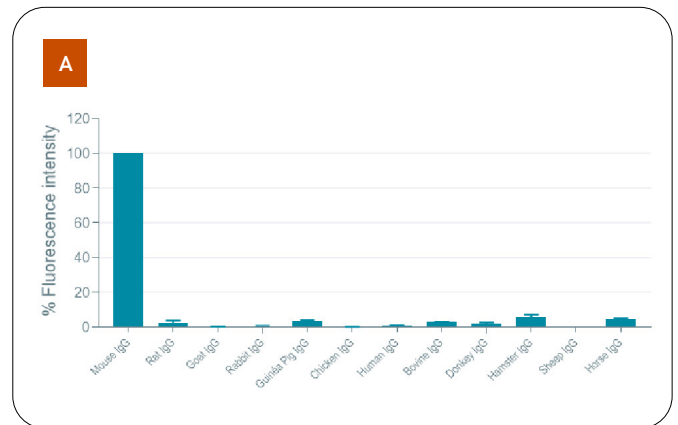
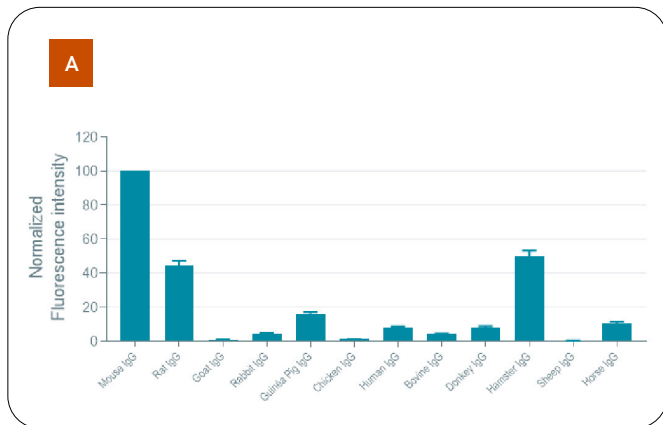


Figure 7.1: F-LISA experiments: different IgG species (A), 10% of serum (B) or IgG isotypes (C) were used to coat a 96-well microplate, then incubated with PhenoVue Fluor 594 - Goat anti-Mouse IgG (H+L) Cross-Adsorbed (1.5 µg/mL). Fluorescence intensity was measured on an EnVision™ multimode plate reader.

Figure 7.2: 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

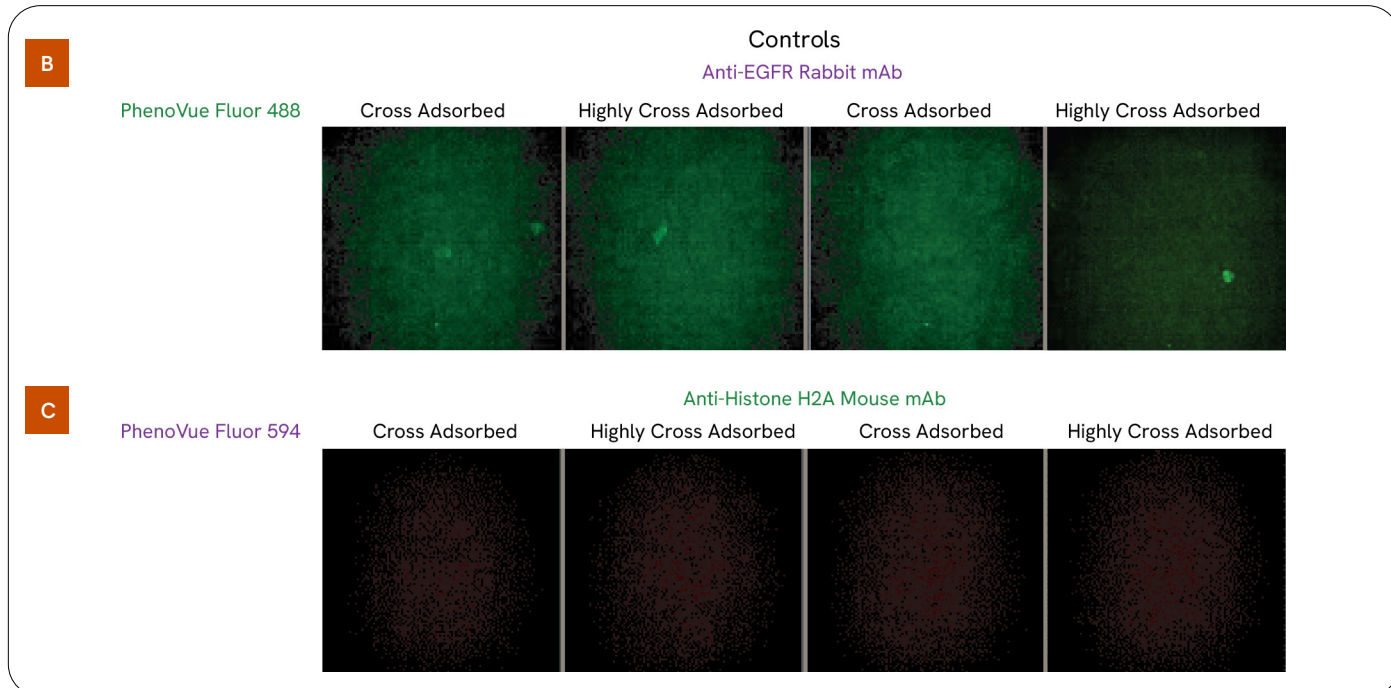
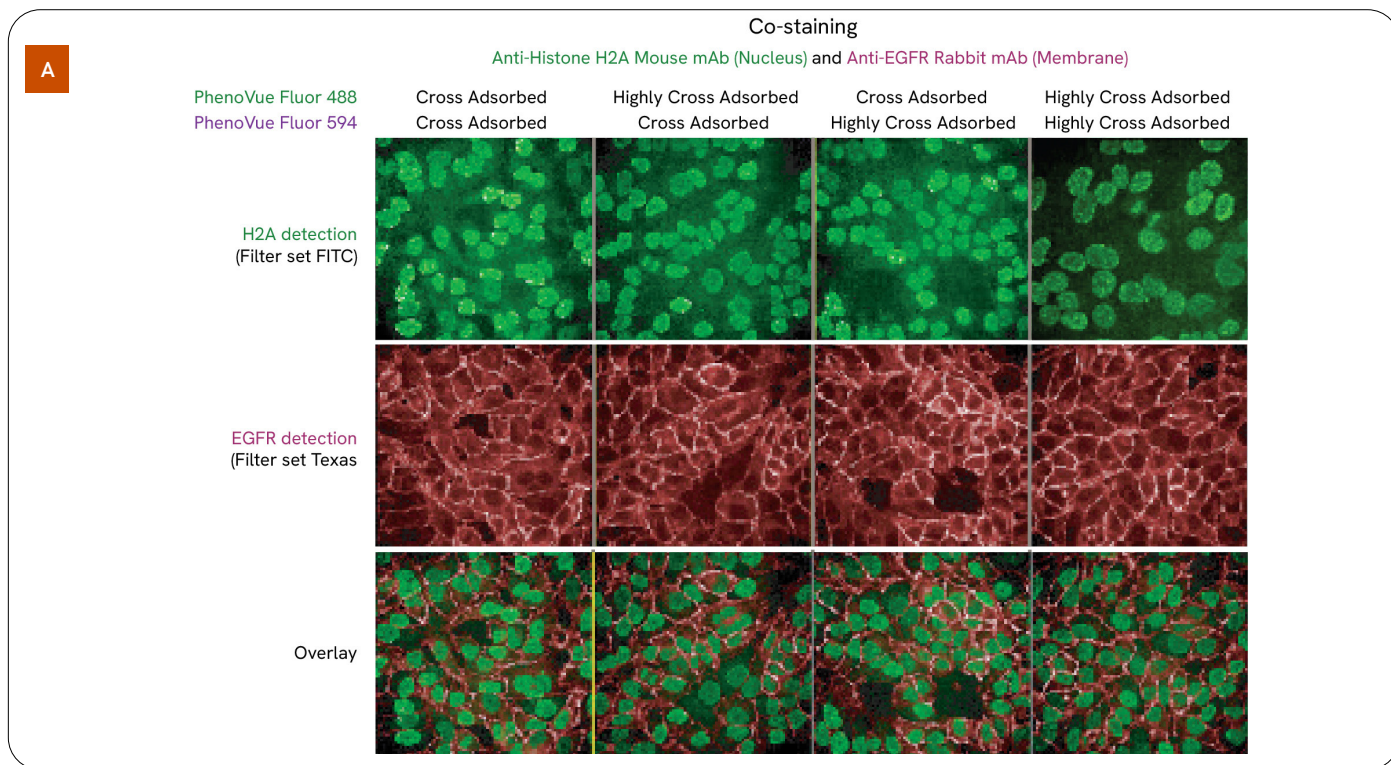


Figure 8: HeLa cells were seeded in PhenoPlate 96-well microplates (50,000 cells/well) and incubated at 37 °C, 5% CO₂ for 24h. Cells were fixed then permeabilized and co-incubated with anti- Histone H2A mouse antibody (5 µg/mL) and anti-EGFR rabbit antibody (0.5 µg/mL). After washing steps, cells were 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 Crossadsorbed 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

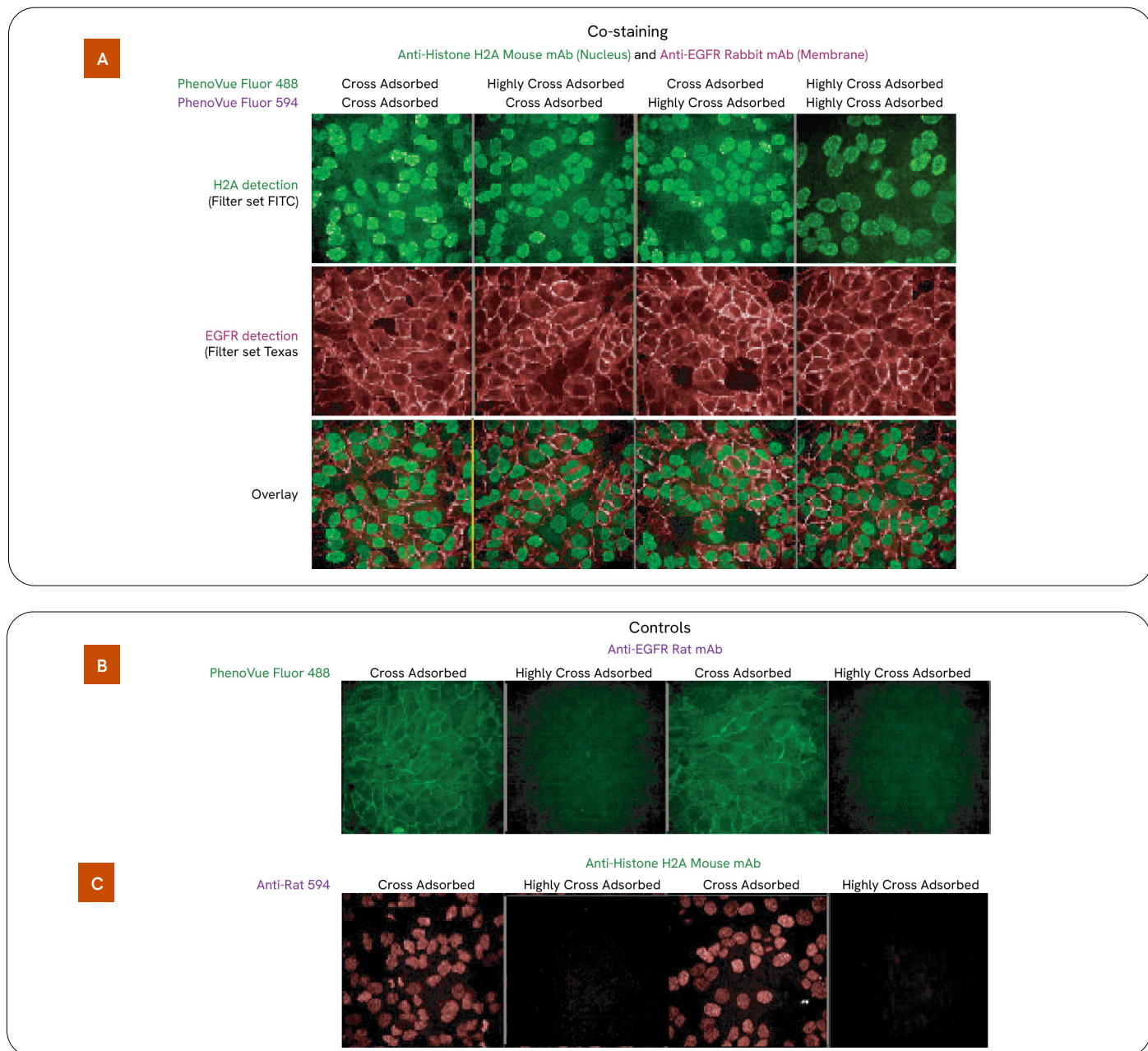


Figure 9: HeLa cells were seeded in PhenoPlate 96-well microplates (50,000 cells/well) and incubated at 37 °C, 5% CO₂ for 24h. Cells were fixed then permeabilized and co incubated with anti- Histone H2A mouse antibody (5 µg/mL) and anti-EGFR Rat antibody (1 µg/mL). After washing steps, cells were incubated 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.

