

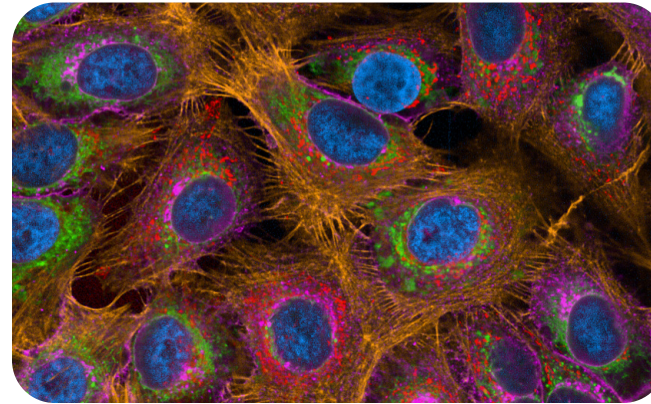


High quality stains for high quality results.

PhenoVue fluorescent reagents and kits for high-content imaging

To get the best results from your high-content assays you need the best cellular imaging reagents.

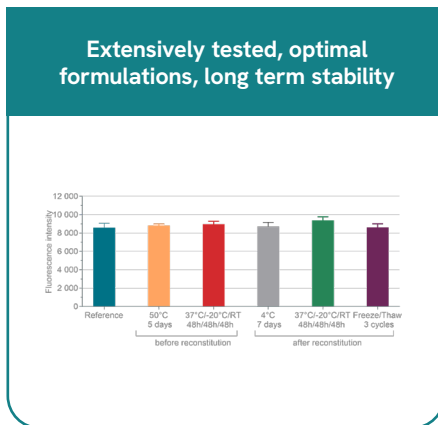
PhenoVue™ kits and reagents are part of our end-to-end solution for high-content workflows, including our market-leading imaging instruments (Opera Phenix™ Plus and Operetta CLS™), automation, and image analysis software, giving you everything you need to streamline your workflow and lead you to breakthroughs faster.



- Organelle and cell compartment stains
- Cell function reagents and kits
- Cell painting kits
- Multi organelle staining kit
- Neuronal differentiation staining kit
- Fluorescent secondary antibodies

Ready to use kits and reagents

Straightforward protocol, easier handling



Performance demonstrated in relevant biological & pharmacological models

iPSC-derived human cortical neurons

Stringent QC High batch to batch reproducibility

Cell painting kit

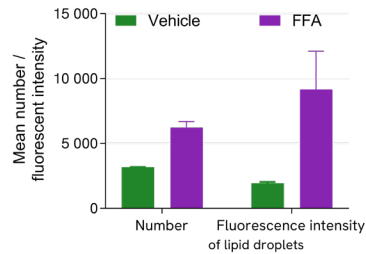
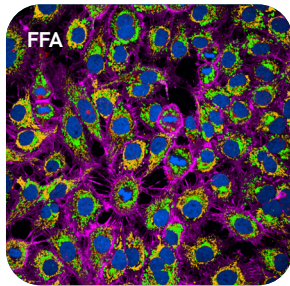
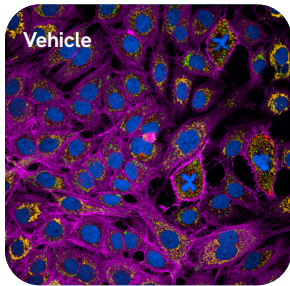
Extensive quality control 87 parameters QC tested

Optimized and verified for phenomenal fluorescent images

Developed by Revvity's highly skilled R&D team with long-standing experience in chemistry, photophysics and assay development, our PhenoVue reagents have been carefully formulated and optimized to deliver **high performance** and **superior image quality** for high-content applications.

With thorough stability testing, and extensive performance verification (including pharmacological studies), our product qualification process is designed to ensure that PhenoVue reagents deliver high performance and reliability, based on our specifications. See our example application data below.

PhenoVue multi organelle staining kit enables to visualize 5 major organelles



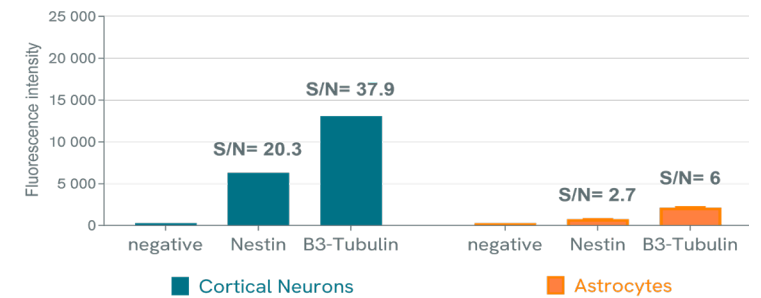
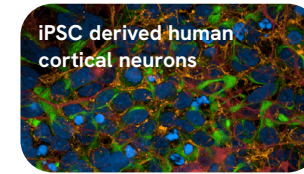
Quantification of fluorescence intensity in lipid droplets

HeLa cells were treated with Free Fatty Acid (FFA) or untreated for 24h and stained using the PhenoVue multi organelle staining kit. Images were acquired on the Opera Phenix HCS system. These results show that exposure to FFA induces an increase in the number of lipid droplets, as expected*.

*Quantification was performed using Signals Image Artist™.

For research use only. Not for use in diagnostic procedures.

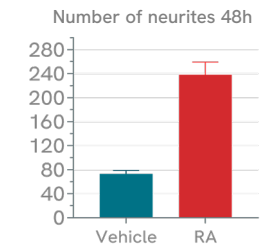
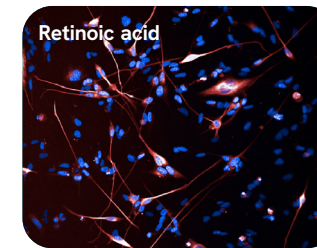
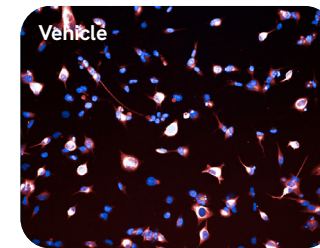
PhenoVue neuronal differentiation staining kit combines markers which are specific to mature neurons



Quantification of fluorescence intensity associated with Nestin (green) and B3 tubulin (red) markers

iPSC-derived human cortical neurons and iPSC-derived human astrocytes were stained using the PhenoVue neuronal differentiation staining kit. Images were acquired on the Operetta CLS HCS system. These results show higher fluorescent signals from the Nestin and B3 tubulin staining in iPSC derived human cortical neurons compared to iPSC-derived human astrocytes, as expected*.

PhenoVue Fluor 647 live cell tubulin enables to visualize microtubules in live cells



Quantification of neurites

SH-SY5Y neuroblastoma cell line treated with retinoic acid or untreated, and incubated with PhenoVue Fluor 647 live cell tubulin stain. Images were acquired on the Opera Phenix Plus HCS system. These results show that exposure to retinoic acid induces an increase in the number of neurites, as expected*.

