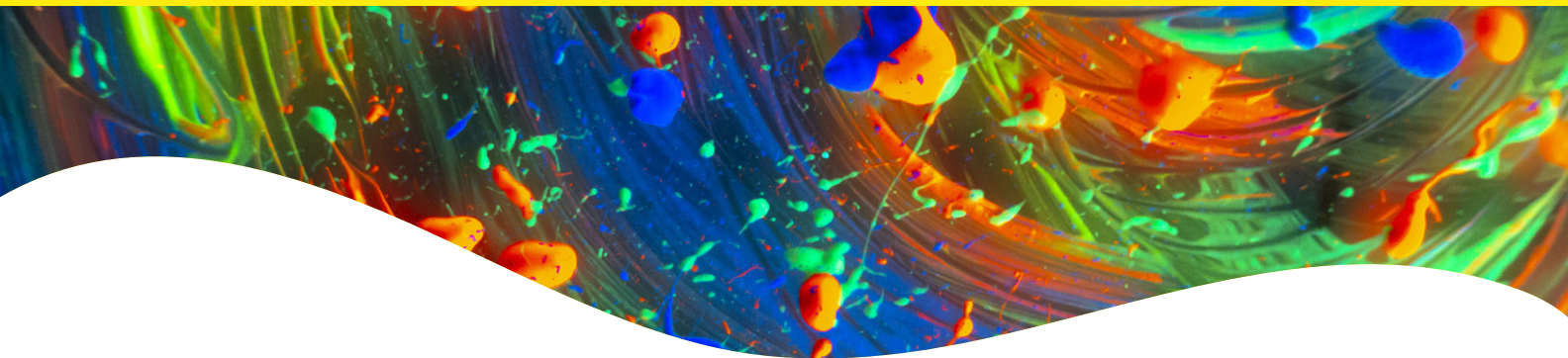


# PhenoVue Paraformaldehyde 4% Methanol-Free Solution



## Overview

Cell fixation is an essential step to preserve cell samples for a wide range of biological assays including cytochemistry and histochemistry. Paraformaldehyde (PFA) has been widely used as a cross-linking fixation agent.

PhenoVue™ paraformaldehyde 4% methanol-free is a ready to use, high purity solution. Together with PhenoVue permeabilization 0.5% Triton X-100 solution, it is ideal for use with PhenoVue Fluor conjugated secondary antibodies, as well as PhenoVue fluorescent probes for organelles & subcellular compartments and PhenoVue cell painting kits.

## Product information

Product name	Part no.	Number of bottles per unit	Quantity per bottle	Format	Shipping conditions
PhenoVue paraformaldehyde, 4% solution, 1 x 25 mL	PVPFA41	1	25 mL	Liquid	RT
PhenoVue paraformaldehyde, 4% solution, 5 x 25 mL	PVPFA42	5	25 mL	Liquid	RT

## Storage and stability

- Store 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.

## Equivalent number of microplates

Product size	96-well microplate (50 $\mu$ L - 100 $\mu$ L per well)	384-well microplate (12 $\mu$ L - 25 $\mu$ L per well)	1536-well microplate (4 $\mu$ L - 12 $\mu$ L per well)
1 x 25 mL	3 to 5	3 to 5	1 to 4
5 x 25 mL	15 to 25	15 to 25	5 to 20

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

## 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

- 1. Rinse** briefly in phosphate-buffered saline (PBS) then proceed with cell fixation.
- 2. Paraformaldehyde fixation:** Add the appropriate volume\* of PhenoVue paraformaldehyde 4% ready to use solution per well for 10 - 15 min at room temperature.  
\*See table above Equivalent number of microplates.
- 3. Washing:** Wash three times with PBS.
- 4. Permeabilization:** Add the appropriate volume of PhenoVue Permeabilization 0.5% Triton X-100 ready to use solution per well for 10 - 15 min at room temperature.
- 5. Washing:** Wash three times with PBS for 5 min.
- 6. Staining:** Perform the staining steps according to your experiment.
- 7. Washing:** Wash three times with PBS for 5 min.
- 8.** Acquire images on an imaging device.

## 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

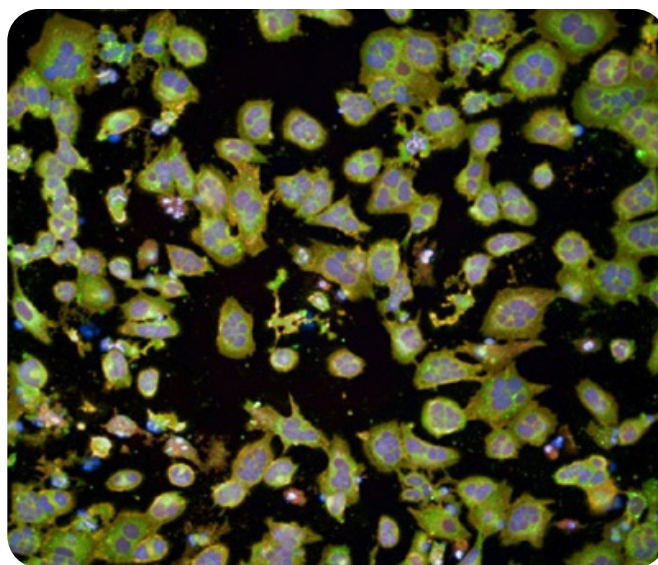


Figure 1: U2OS cells seeded in PhenoPlate™ 384-well, treated with Cytochalasin D and stained according to **PhenoVue Cell Painting JUMP Kit** protocol. Images acquired on the Opera Phenix® high-content screening system.

revvity