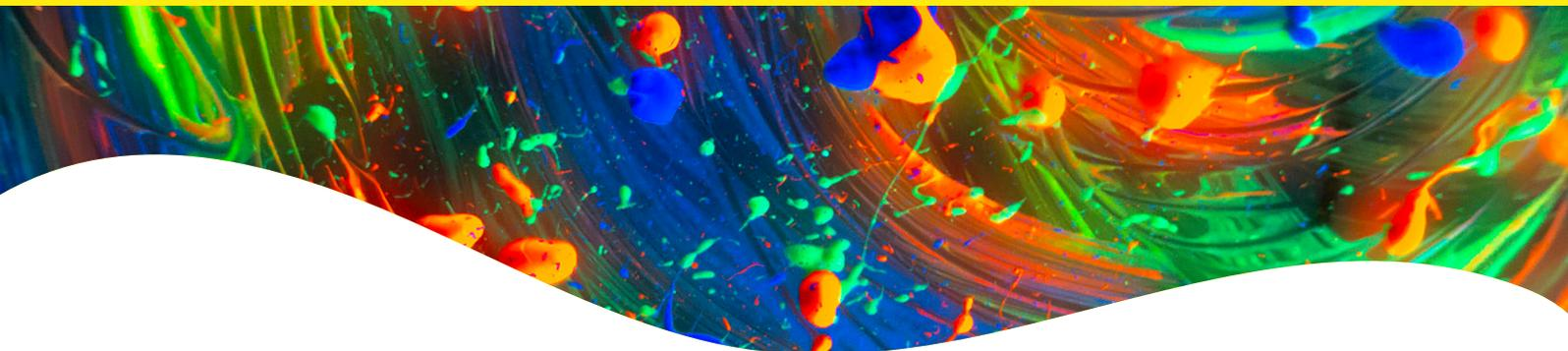


PhenoVue Permeabilization 0.5% Triton X-100 Solution



Overview

After cell fixation, cell permeabilization is an essential step to remove lipids from plasma and nuclear membranes and allow large molecules such as antibodies to penetrate the cell.

PhenoVue™ permeabilization 0.5% Triton X-100 solution is ready to use. Together with PhenoVue paraformaldehyde 4% 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 permeabilization 0.5% Triton X-100 solution, 1 x 25 mL | PVPERM051 | 1 | 25 mL | Liquid | RT |
| PhenoVue permeabilization 0.5% Triton X-100 solution, 5 x 25 mL | PVPERM052 | 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 μ L - 100 μ L per well) | 384-well microplate (12 μ L - 25 μ L per well) | 1536-well microplate (4 μ L - 12 μ 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.revivity.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% CO₂ 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.

REACH European regulations and compliance: This product and/or some of its components include a Triton X-100 concentration of 0.1% or more and as such, it is covered by the REACH European regulations. We recommend researchers using this product to act in compliance with REACH and in particular: to only use the product for *in vitro* research in appropriate and controlled premises by qualified researchers, ii) to ensure the collection and the treatment of subsequent waste, and iii) to make sure that the total amount of Triton X-100 handled does not exceed 1 ton per year.

Applications

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

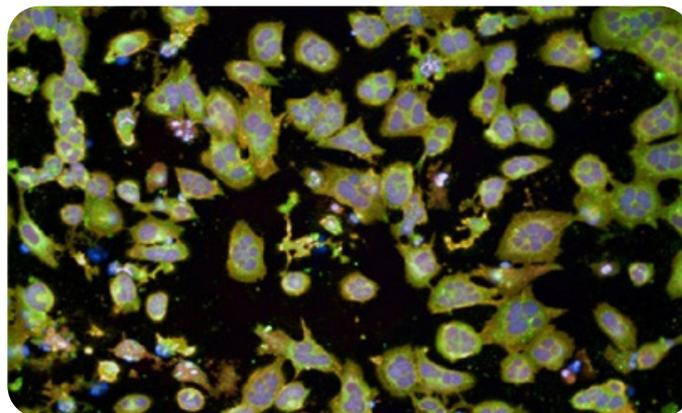


Figure 1: U2OS cells seeded in PhenoPlate™ 384-well, treated with Cytochalasin D and stained according to **PhenoVue Cell Painting JUMP Kit** protocol. Images acquired on the Opera Phenix® high-content screening system.

revvity