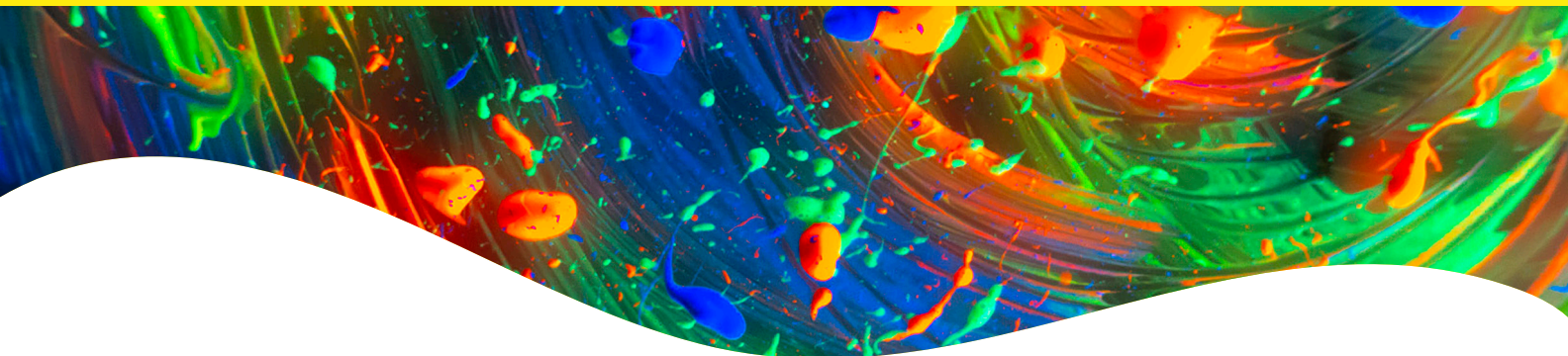


PhenoVue Dye Diluent A (5X)



Overview

PhenoVue™ dye diluent A (5X) is an optimized and concentrated buffer that can be used to dilute PhenoVue fluorescent products.

Product information

Product name	Part no.	Number of bottles per unit	Quantity per bottle	Format	Shipping conditions
PhenoVue dye diluent A (5X)	PVDDA1	1	80 mL	Liquid	Dry ice

Storage and stability

- Store at 2-8 °C, protected from light.
- The stability of this product is guaranteed until the expiration date provided in the Certificate of Analysis, when stored as recommended and protected from light.

Protocol

Dilute PhenoVue dye diluent A 5 times in distilled H₂O to give a 1X ready to use buffer.

Tips

After thawing, the PhenoVue dye diluent A (5X) may display a few aggregates which will not impair product and image quality. If aggregate removal is required, the PhenoVue dye diluent A (5X) can be filtered (0.22 µm filter), prior to dilution. The resulting PhenoVue dye diluent A 1X must be stored at 4 °C and for no more than 2 days.

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

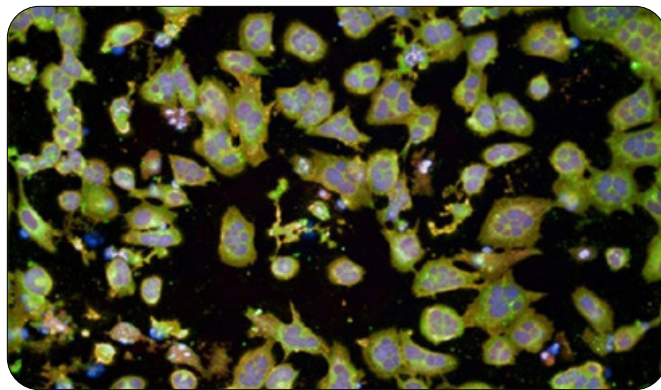
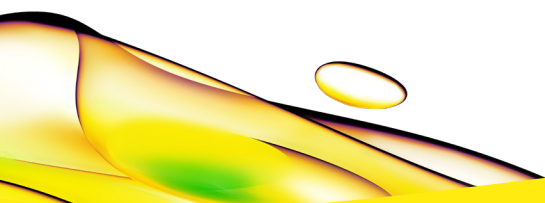


Figure 1: U2OS cells seeded in PhenoPlate™ 384-well, treated with Cytochalasin D and stained according to **PhenoVue cell painting JUMP kit** protocol. Images acquired on the Opera Phenix® high-content screening system.



revvity