

## Screen Quest™ 10X Cell Staining Buffer with Phenol Red Plus™

Ordering Information	Storage Conditions
Product Number: 36300 (10 mL, 10 plates)	Keep in freezer and protect from light

### Typical Assay Protocol (for one 96-well plate)

1. Thaw 10X cell staining buffer at room temperature before use.  
*Note: It is OK to use if the buffer has precipitates.*
2. Prepare 1X Screen Quest™ Cell Staining Buffer: Add **1 mL** of 10X Screen Quest™ cell staining buffer to **9 mL** of HHBS (1X Hank's with 20 mM Hepes buffer, pH 7.0) or buffer of your choice, and mix them well.  
*Note: 10 mL of 1X staining buffer is enough for one plate. The buffer is stable at room temperature. It is recommended to aliquot and store un-used 10X assay buffer at  $\leq -20$  °C. Protect from light and avoid repeated freeze-thaw cycles.*
3. Prepare 2X Assay Solution: Add the cell staining dye stock solution (in general, it is a concentrated DMSO solution) into 1X Screen Quest™ Cell Staining Buffer (from Step 2) to make the final well concentration 2X of the desired concentration.
4. To the microplate well add 2X Assay Solution (from Step 3) which is the same volume as the cell culture medium (e.g., 100  $\mu$ L/well/96-well or 25  $\mu$ L/well/384-well).
5. Incubate the cells in a 37 °C, 5% CO<sub>2</sub> incubator, or as desired.  
*Note: It is possible that the staining dye might interfere with the 1X Screen Quest™ Cell Staining Buffer. In this case, it's recommended to stain the cells by a desired method, and replace the cell staining solution with the cell growth medium or HHBS. Then add 100  $\mu$ L/well/96-well (25  $\mu$ L/well/ 384-well) of 1X Screen Quest™ Cell Staining Buffer into the well.*
6. Observe the cells with a fluorescence microscope or a plate reader as required.

**Disclaimer:** This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.