



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Permeabilization Buffer

Ready-to-use permeabilization buffer for intracellular immunofluorescence staining for microscopy or flow cytometry.



Product attributes

Storage Conditions
Store at room temperature

Product Description

Permeabilization Buffer is a ready to use permeabilization buffer for intracellular immunofluorescence staining for microscopy or flow cytometry.

- Ready-to-use buffered detergent solution
- Permeabilize fixed cells for microscopy and flow cytometry applications
- 100 mL solution

The permeabilization Buffer is also available in the Flow Cytometry Fixation/Permeabilization Kit ([23006](#)).

References

1. DNA Repair (2018) Jun - Jul;66-67:42-49 [doi: 10.1016/j.dnarep.2018.04.007](https://doi.org/10.1016/j.dnarep.2018.04.007)
2. Mol Med Rep (2019) Oct;20(4):3719-3727 [doi:10.3892/mmr.2019.10590](https://doi.org/10.3892/mmr.2019.10590)
3. PLoS One (2019) Oct 9;14(10):e0223725 [doi:10.1371/journal.pone.0223725](https://doi.org/10.1371/journal.pone.0223725)

This datasheet was generated on July 27, 2026 at 10:08:42 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/permeabilization-buffer/>