

10x Glycine buffer pH 3.0 Acidic Glycine Wash Buffer

Cat.-No.	Amount
CPP-A07S	25 ml
CPP-A07L	125 ml

For *in vitro* use only
Quality guaranteed for 12 months
Store at 4°C

Product Information

Form

Concentrated acidic glycine buffer.
Dilute 1:10 with sterile water before use. The diluted
buffer has a concentration of 200 mM.

Description

To remove adsorbed complex from the outside of the membranes a thorough and repeated (3 times) washing of the cells with glycine buffer of pH 3.0 is recommended.

For more information see **general manual**.

Selected References

Kameyama *et al.* (2007) Acid Wash in Determining Cellular Uptake of Fab/Cell-Permeating Peptide Conjugates. *Biopolymers (Peptide Science)* **88**:98.