
Product Information

EasyProbes™ Fixation/Permeabilization Kit

Catalog Number: C065

Table 1. Kit Components and Storage

Material	Amount	Concentration	Storage
Fixative Solution	50 mL	High purity 4% formaldehyde in PBS, pH 7.4	RT
Permeabilization Solution	50 mL	0.5% Triton X-100 in PBS, pH 7.4	RT

Quick Start Guide

1. Remove media from cells.
2. Add 1 ml of Fixative Solution.
3. Incubate 15 min at 20-25°C.
4. Remove Fixative Solution. Wash 3 times with 2 ml of PBS Buffer for 2-5 min.
5. Add 1 ml of Permeabilization Solution.
6. Incubate 15 min at 20-25°C.
7. Remove Permeabilization Solution. Wash 3 times with 2 ml of PBS Buffer for 2-5 min.
8. Add 2 ml of 3% BSA in PBS.
9. Incubate 60 min at 20-25°C.
10. Proceed with desired cell staining and/or antibody labeling.