

PrimeFlow™ Fix/Perm Buffer Set

Catalog Number: 88-17000

GPR: General Purpose Reagents. For Laboratory Use.

Product Information



Contents: PrimeFlow™ Fix/Perm Buffer Set

Catalog Number: 88-17000



Temperature Limitation: Store at 2-8°C.



Batch Code: Refer to vial



Use By: Refer to vial



Contains sodium azide and <10% formaldehyde

Description

The PrimeFlow® Fixation/Permeabilization Buffer Set is a component of the PrimeFlow® RNA Assay (cat. no. 88-18005). These buffers, with or without the PrimeFlow microRNA Pretreatment Buffer (cat. no. 88-18006), can be used to evaluate intracellular staining antibody performance following the fixation and permeabilization steps of the PrimeFlow assay. Note that additional changes to antibody performance may be observed after the hybridization steps of the PrimeFlow RNA assay.

Please refer to the appropriate PrimeFlow RNA Assay protocol on the product data sheet for instructions for use.

Components

PrimeFlow® RNA Fixation Buffer 1A: cat. 00-18100, 50 mL, store at 2-8°C.

PrimeFlow® RNA Fixation Buffer 1B: cat. 00-18200, 50 mL, store at 2-8°C.

PrimeFlow® RNA Permeabilization Buffer (10X): cat. 00-18300, 30 mL, store at 2-8°C. *Note: This buffer has a natural tendency to precipitate, however, its function is not affected by this. To clarify, the solution can be filtered after dilution to a 1X working solution.*

PrimeFlow® RNA Fixation Buffer 2 (8X): cat. 00-18400, 12.5 mL, store at 2-8°C.

PrimeFlow® RNA Wash Buffer: cat. 00-19180, 100 mL, store at 2-8°C.

Applications Reported

This PrimeFlow® Fixation/Permeabilization Buffer Set has been reported for use in flow cytometry.

Applications Tested

This PrimeFlow® Fixation/Permeabilization Buffer Set has been tested by intracellular staining and flow cytometric analysis using the PrimeFlow® RNA Assay and protocol. Note that this buffer set does not include RNase inhibitors, as it is intended only for evaluation of antibody performance.

Related Products

44-17005 PrimeFlow® 96-well plate

88-18005 PrimeFlow® RNA Assay

88-18006 PrimeFlow® microRNA Pretreatment Buffer

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience • info@ebioscience.com