

Catalog No. R37602

Kit contents	Amount
Wash Buffer	500 mL
Fixative Solution	20 mL
Permeabilization Solution	20 mL
Blocking Solution	12 × 10 mL

Required materials not supplied
6-well plate or 35-mm dish (protocol designed for ~10 cm ² growth area)
Dyes
Primary and secondary antibodies



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 Wash 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 Wash Buffer for 2–5 min.



8. Add 2 mL of Blocking Solution.



9. Incubate 60 min at 20–25°C.



10. Proceed with desired cell staining and/or antibody labeling.

For Research Use Only. Not for use in diagnostic procedures.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses.

By use of this product you accept the terms and conditions of all applicable Limited Use Label Licenses.

Manufacturer: Life Technologies Corporation | 29851 Willow Creek | Eugene, OR 97402

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.