



To be used with NaveniFlex – MR Kit | Item No. 39205

GENERAL GUIDELINES:

- Use a PAP pen to draw a sufficiently large border around each cell pellet.
- Reaction volume for control slides is 40 μ l per cell pellet.
- Use best practices when pipetting to minimize reagents consumption.
- Centrifuge vials before pipetting.
- Select the Buffer C vial with the appropriate detection fluorophore for your microscope filter set.
- Completely defrost all buffer mixtures at room temperature, vortex well before use.
- Vortex and spin-down all enzymes (A, B and C) before use.
- Keep enzymes on ice or on a frozen cold block.
- Wait to add enzymes until immediately prior to adding to the sample.
- Buffer C is a light-sensitive reagent. Always keep it protected from light.
- Remove excess washing buffer from samples before adding reagent.
- Do not allow slides/samples to dry.
- Preheat the humidity chamber before each step.
- Incubation times or temperatures other than those specified may give erroneous results.
- NaveniFlex assay might be combined with traditional immunofluorescence techniques, providing that all primary antibodies are from different species and the fluorescence detection systems use different fluorophores.
- As with any product derived from biological sources, proper handling procedures should be used.
- Wear appropriate Personal Protective Equipment to avoid contact with eyes and skin.
- Unused solution should be disposed according to local regulation.

IMPORTANT: Appropriate precautions should be taken to avoid antibodies cross-contamination. Cross-contamination is the main source of unspecific background due to the high sensitivity of the assay.

Avoid bulk washing methods when multiple antibodies are used.



To be used with NaveniFlex – MR Kit | Item No. 39205

CONTROL KIT COMPONENTS:

Material	Cat. No.	Supplier (Cat. No.)	Amount	Storage*
Mouse anti-Her2 antibody (100X)	NF.2.CK.01	Origene (TA503443)	5 µl	From -25 to -15°C
Rabbit anti-Her2 antibody (100X)	NF.2.CK.02	Atlas Antibodies (HPA001383)	5 µl	
Control Slides (BT474)	NF.2.CK.03	Acepix (N/A)	3 slides	

PROTOCOL:

1. Slide Preparation
 - 1.1 Let slides warm up to room temperature.
 - 1.2 Use a PAP pen and draw a border around each cell pellet and let the border dry completely.
 - 1.3 Rehydrate the cells by adding 1x PBS onto each pellet, incubate for 3 min at room temperature.
 - 1.4 Transfer slides into a staining jar with 1x PBS and wash for 2 min with gentle agitation.
2. Blocking
 - 2.1 Add **Blocking Buffer** (1x) to the entire sample area (approximately 40 µl for each 1cm² area).
 - 2.2 Incubate for 60 min at +37 °C in a pre-heated humidity chamber.
3. Primary antibody incubation
 - 3.1 Use the provided **Primary Antibody Diluent** (1x) to dilute the **Mouse and Rabbit Primary Antibodies** according to the table below.

Antibody/ Diluent	Volume
Mouse anti-Her2 antibody (100x)	1 µl
Rabbit anti-Her2 antibody (100x)	1 µl
Naveni Primary Antibody Diluent	98 µl
Total:	100 µl

- 3.2 Decant the Blocking Buffer and add enough of the antibodies to cover one cell pellet – add 40 µl of Primary Antibody Diluent to the second cell pellet as negative control.
- 3.3 Incubate overnight at +4 °C in a humidity chamber.
- 3.4 Aspirate to remove the antibody solution and wash slides for 3x5 min with 1x TBS-T** in a staining jar under gentle agitation.



To be used with NaveniFlex – MR Kit | Item No. 39205

4. Probe incubation
 - 4.1 Prepare the probes by diluting **Probe M1** and **Probe R2** in **Probe Diluent** (1x) (dilute 1:40 each).
 - 4.2 Add enough of the probes to cover the sample area.
 - 4.3 Incubate for 60 min at +37 °C in a pre-heated humidity chamber.
 - 4.4 Decant the solution and wash slides for 3x5 min with 1x TBS-T in a staining jar under gentle agitation.

5. Reaction A
 - 5.1 Start preparing **Reaction A** by diluting **Buffer A** (5x) 1:5 in water. Vortex and spin down.
 - 5.2 Add **Enzyme A** (dilute 1:40) Mix gently by pipetting and spin down.
 - 5.3 Add enough Reaction A to cover the sample area.
 - 5.4 Incubate for 60 min at +37 °C in a pre-heated humidity chamber.
 - 5.5 Decant the solution and wash slides for 2x3 min with 1x TBS-T in a staining jar under gentle agitation.

6. Reaction B
 - 6.1 Start preparing **Reaction B** by diluting **Buffer B** (5x) 1:5 in water. Vortex and spin down.
 - 6.2 Add **Enzyme B** (dilute 1:40). Mix gently by pipetting and spin down.
 - 6.3 Add enough Reaction B to cover the sample area.
 - 6.4 Incubate for 30 min at 37 °C in a pre-heated humidity chamber.
 - 6.5 Wash slides for 2x3 min with 1x TBS-T in a staining jar under gentle agitation.

7. Reaction C
 - 7.1 Select the **Buffer C** vial with the appropriate detection fluorophore for your microscope filter set. Do not use more than one **Buffer C** vial.
 - 7.2 Start preparing **Reaction C** by diluting **Buffer C** (5x) 1:5 in water. Vortex and spin down.
 - 7.3 Add **Enzyme C** (dilute 1:40). Mix gently by pipetting and spin down.
 - 7.4 Add enough Reaction C to cover the sample area.
 - 7.5 Incubate for 90 min at +37 °C in a pre-heated humidity chamber.
 - 7.6 Decant the solution and wash slides for 2 min with 1x TBS in a staining jar under gentle agitation.

Protect from light



To be used with NaveniFlex – MR Kit | Item No. 39205

8. Nuclei staining and mounting
(not provided)

Protect from light

- 8.1 Start preparing a Nuclei staining solution by diluting Dapi in 1xPBS. Vortex and spin down.
- 8.2 Decant wash buffer from the slides.
- 8.3 Add enough Nuclei staining solution to cover the sample area.
- 8.4 Incubate for 5 min at room temperature on the bench.
- 8.5 Decant the solution and wash slides for 2x 10 min with 1x TBS in a staining jar under gentle agitation.
- 8.6 Wash slides for 15 min with 0.1x TBS in a staining jar under gentle agitation.

9. Mounting
(not provided)

Protect from light

- 9.1 Decant excess wash buffer from the slides.
- 9.2 Mount the slides with a coverslip using an anti-fade mounting medium.
- 9.3 Image your slides in fluorescence or confocal microscope, using 20x objective or higher and filter set for DAPI and a corresponding fluorophore (FITC for Atto 488, Texas Red, or Cy5 for Atto 647, respectively).

**TBS-T (Tris-buffered saline supplemented with 0.05% Tween 20)

RECOMMENDED IMAGING SETTINGS:

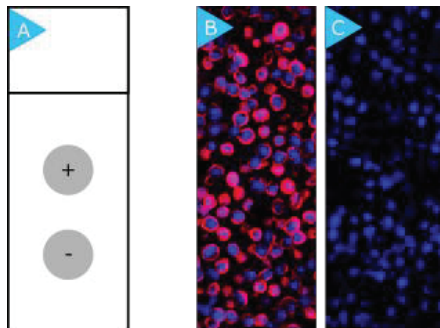


Fig. 1: The recommended slide layout, where “+” indicates both primary antibodies present whilst “-” indicates no primary antibodies present (A). Reference images displaying a typical positive (B) and negative (C) result.