

Let's get started!

This RePlex™ Module provides all the reagents needed to remove primary and secondary antibodies in a RePlex™ assay with two immunoassays or immunoassay with Total Protein assay in a single run. Please note this module can only be used with Jess and Abby.

Components

CLAMSHELL 1, REPLEX™ REAGENT 1 (RP-001-1), STORE AT 18–28 °C

INCLUDES	PART NO
RePlex™ Reagent 1 (12 mL)	043-827

CLAMSHELL 2, REPLEX™ REAGENT 2 (RP-001-2), STORE AT 2–8 °C

INCLUDES	PART NO
RePlex™ Reagent 2 (3 mL)	043-828

Other things you'll need

- Reagents and consumables supplied in the Separation Module and Detection Module.
- Instructions for preparing reagents are provided with the Separation Module and Detection Module inserts, or they can be downloaded from our [website](#).

RePlex™ Module preparation

MIX LUMINOL-S AND PEROXIDE (IF APPLICABLE)

- Combine 450 µL Luminol-S and 450 µL Peroxide in a microcentrifuge tube.
- Vortex to mix and store on ice.

PREPARE REPLEX™ REAGENT

- Combine 1.4 mL of RePlex™ Reagent 1 and 0.35 mL of RePlex™ Reagent 2 in a microcentrifuge tube.
- Close tube and vortex briefly.



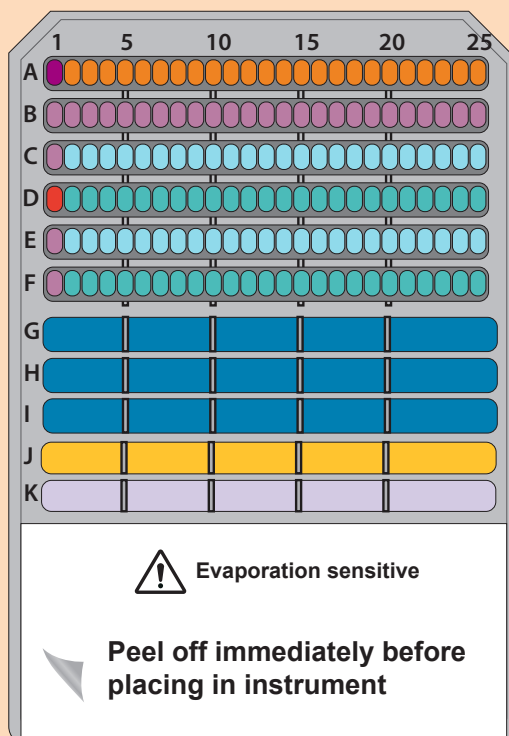
Toll-free: (888) 607-9692
Tel: (408) 510-5500
Fax: (408) 510-5599
info@proteinsimple.com
proteinsimple.com

© 2021 ProteinSimple. ProteinSimple, Simple Western, Jess, Abby and the ProteinSimple logo are trademarks and/or registered trademarks of ProteinSimple.

PL3-0029, Rev B

Pipette your plate

2 IMMUNOASSAYS



● Biotinylated Ladder, 5 μ L; ● Prepared Samples, 3 μ L

● Antibody Diluent (2 or Milk-Free), 10 μ L

● Antibody Diluent (2 or Milk-Free), 10 μ L; ● Primary Antibody for Probe 1, 10 μ L

● Streptavidin-HRP or NIR, 10 μ L; ● Secondary Antibody for Probe 1, 10 μ L

● Antibody Diluent (2 or Milk-Free), 10 μ L; ● Primary Antibody for Probe 2, 10 μ L

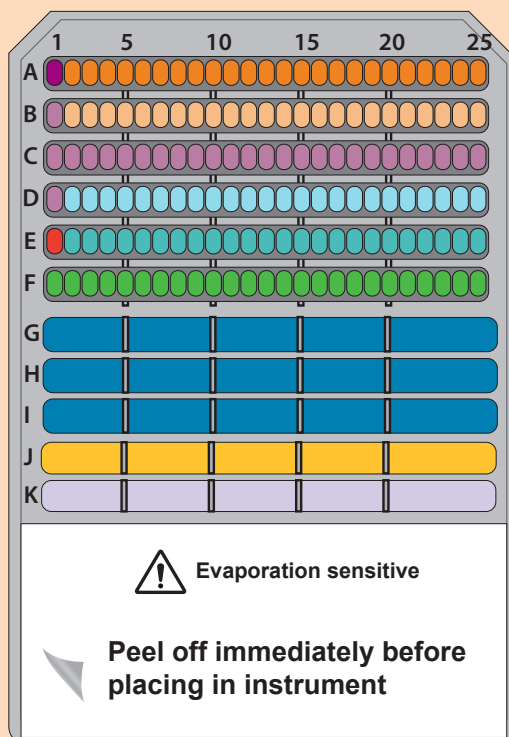
● Antibody Diluent (2 or Milk-Free), 10 μ L; ● Secondary Antibody for Probe 2, 10 μ L

● Wash Buffer
500 μ L/compartments

● Luminol-Peroxide Mix, 170 μ L/compartments (for chemiluminescence only)

● RePlex™ reagent mix, 300 μ L/compartments

IMMUNOASSAY + TOTAL PROTEIN



● Biotinylated Ladder, 5 μ L; ● Prepared Samples, 3 μ L

● Antibody Diluent (2 or Milk-Free), 10 μ L; ● Total Protein labeling reagent, 10 μ L

● Antibody Diluent (2 or Milk-Free), 10 μ L

● Antibody Diluent (2 or Milk-Free), 10 μ L; ● Primary Antibody for Probe 1, 10 μ L

● Streptavidin-HRP or NIR, 10 μ L; ● Secondary Antibody for Probe 1, 10 μ L

● Total Protein Streptavidin-HRP for Probe 2, 8 μ L

● Wash Buffer
500 μ L/compartments

● Luminol-Peroxide Mix, 170 μ L/compartments

● RePlex™ reagent mix, 300 μ L/compartments

For more consistent results, keep the lid on the microplate between reagent additions and minimize bubble formation when adding Wash Buffer to the troughs.

1. Dispense reagents into the assay plate using the volumes shown in the plate diagram.
2. Centrifuge the plate for 5 minutes at 2500 rpm ($\sim 1000 \times g$) at room temperature. Ensure liquid is fully down in all wells.