

ScanLater™ Western Blot Assay Kit

Table 1-1: Available Kits

Assay Kit	Explorer Kit	Bulk Kit
Eu-Labeled Goat Anti-Mouse ScanLater™ Western Blot Assay Kit	R8201	R8205
Eu-Labeled Goat Anti-Rabbit ScanLater™ Western Blot Assay Kit	R8202	R8204
Eu-Labeled Donkey Anti-Goat ScanLater™ Western Blot Assay Kit	R8224	R8227
Eu-Labeled Donkey Anti-Rat ScanLater™ Western Blot Assay Kit	R8225	R8228
Eu-Labeled Streptavidin ScanLater™ Western Blot Assay Kit	R8200	R8203

Quick Start Protocol

1. Dilute the buffer solutions and reconstitute the antibody or streptavidin provided in the assay kit.
2. Transfer the gel to the blotting membrane as required by your application.
3. Soak the membrane in 1X washing buffer for a minimum of 1 minute.
4. Block the membrane in 1X blocking buffer for 1 hour.
5. Add your primary antibody directly into the blocking buffer at the correct dilution for your application and incubate from 90 minutes to 4 hours.
6. Wash the membrane 3 times in 1X wash buffer for 5 minutes each time.
7. Dilute the previously reconstituted antibody or streptavidin in 1X blocking buffer.
 - For the Explorer Kit, the recommended dilution is 1:500.
 - For the Bulk Kit, the recommended dilution is 1:5000.
8. Add the diluted antibody or streptavidin to the membrane and incubate for 60 minutes.
9. Wash the membrane 3 times in 1X wash buffer for 5 minutes each time.
10. Rinse the membrane with water for 15 seconds and let it dry before scanning. The membrane can be scanned wet right after the wash step, but a dry membrane has a higher signal compared to its background.
11. Scan the membrane in a SpectraMax Paradigm or SpectraMax i3 Instrument.

Download Instructions

Complete instructions for using the ScanLater Western Blot Assay Kit can be downloaded from the Knowledge Base on the Molecular Devices support web site: www.moleculardevices.com/support.html

Click **Knowledge Base**, and then click **Manuals** and search for **ScanLater** or **Western Blot**.

Kit Components

Table 1-2: Components of the ScanLater Western Blot Assay Kits

Item	Explorer Kit	Bulk Kit
ScanLater 10X Washing Buffer	100 mL, 1 bottle	300 mL, 3 bottles at 100 mL each
ScanLater 5X Blocking Buffer	40 mL, 1 bottle	120 mL, 1 bottle
ScanLater Eu-Labeled Antibody or Streptavidin	20 µg, 1 vial	60 µg, 1 vial

Before using the components of the assay kit, dilute the buffer solutions and reconstitute and aliquot the Eu-labeled antibody or streptavidin.

If you are using the Explorer Kit, do not aliquot or freeze the reconstituted antibody or streptavidin. Protect the solution from light and store it at 4°C for up to one (1) week.

The Explorer Kit and the Bulk Kit differ in their dilution requirements.

Table 1-3: Dilution Differences in the ScanLater Western Blot Assay Kits

Item	Explorer Kit	Bulk Kit
Reconstitute Eu-labeled antibody or streptavidin	200 µL of 1x blocking buffer	60 µL of reagent-grade water
Recommended dilution of reconstituted antibody or streptavidin for the assay	1:500	1:5000

Table 1-4: Separately Orderable Components

Item	Volume	Part Number
ScanLater 10X Washing Buffer	100 mL bottle	R8206
ScanLater 5X Blocking Buffer	120 mL bottle	R8207
ScanLater Eu-Labeled Goat Anti-Mouse IgG	60 µg vial	R8208
ScanLater Eu-Labeled Goat Anti-Rabbit IgG	60 µg vial	R8209
ScanLater Eu-Labeled Donkey Anti-Goat IgG	60 µg vial	R8226
ScanLater Eu-Labeled Donkey Anti-Rat IgG	60 µg vial	R8229
ScanLater Eu-Labeled Streptavidin	60 µg vial	R8212

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/productpatents>

Product manufactured by Molecular Devices, LLC.
 1311 Orleans Drive, Sunnyvale, California, United States of America 94089.
 Molecular Devices, LLC is ISO 9001 registered.
 © 2014 Molecular Devices, LLC.
 All rights reserved.