

Applications

VRDye 549 Streptavidin (**P/N 926-54030**) can be used as a secondary detection reagent for microscopy, protein arrays, In-Gel Westerns, and Western blotting. Centrifuge briefly before use to eliminate aggregates that may have formed in solution. This will reduce non-specific background staining.

Properties

- Quantity: 0.5 mg
- Storage: 4 °C **protected from light**. Expiration date is listed on the product's label.
- Physical State: Liquid
- Concentration: 1 mg/ml
- Fluorophore: VRDye 549 (1040 g/mole)
- Excitation Wavelength: 555 nm
- Emission Wavelength: 569 nm
- Buffer: 10 mM Phosphate, 183 mM NaCl, and 2.7 mM KCl, pH 7.4 with 0.005% (w/v) sodium azide as a preservative.

Warning: *Contains sodium azide.*

Suggested Dilution Ranges

Suggested working dilutions are given as a guide only. Optimum dilutions will vary and should be determined empirically.

For membrane-based applications and In-Gel Westerns, it is recommended to add SDS (0.02% to 0.1% final concentration) and Tween® 20 (0.1 to 0.2% final concentration) during the detection incubation step to reduce non-specific background staining.

Application	Recommended	Suggested Range
Western blot detection	1:2,000	1:2,000-1:5,000
Other	User Optimized	

Limitation of Liability and Limited Use Label License

For Research Use Only. Not intended for human therapeutic or diagnostic use.

By using this product, you agree to the Limitation of Liability and the Limited Use Label License available online at licor.com/packinserts under the **Pack Inserts** heading (the "License"). The License is incorporated herein by this reference, and the License may be updated from time to time.



© 2021 LI-COR, Inc. LI-COR and VRDye are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com