

**0.5 mg**

## Secondary Antibody Quick Card

**Storage:** Store lyophilized secondary antibody as indicated on the tube. After reconstitution, store stock solution at 4 °C for up to 3 months.

**Reconstitution:** Add 0.5 mL sterile distilled water for a final concentration of 1.0 mg/mL. Mix gently by inverting, and allow to rehydrate for a minimum of 30 minutes before use. Please visit [licor.com/packinserts](https://www.licor.com/packinserts) for detailed product information.

### Recommended Dilutions for Western blotting

Product	Recommended	Suggested Range
IRDye® 800CW	1:20,000	1:5,000 – 1:25,000
IRDye® 680RD	1:20,000	1:5,000 – 1:25,000
IRDye® 680LT	1:20,000	1:10,000 – 1:50,000
VRDye™ 549	1:10,000	1:5,000 – 1:15,000
VRDye™ 490	1:10,000	1:5,000 – 1:15,000

# Secondary Antibody Quick Card

0.5 mg

**Avoid non-specific background staining on PVDF membranes:** If you are making your own antibody diluent, add SDS (final concentration 0.01 – 0.02%) and Tween® 20 (final concentration 0.2%) to the diluent. If using Intercept® T20 Antibody Diluent, only add SDS (final concentration 0.01 – 0.02%).

## Recommended Dilutions for In-Cell Western™ Assay

Product	Recommended	Suggested Range
IRDye® 800CW	1:800	1:200 – 1:1,200
IRDye® 680RD	1:800	1:200 – 1:1,200

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



Doc # 988-19645  
12/21