

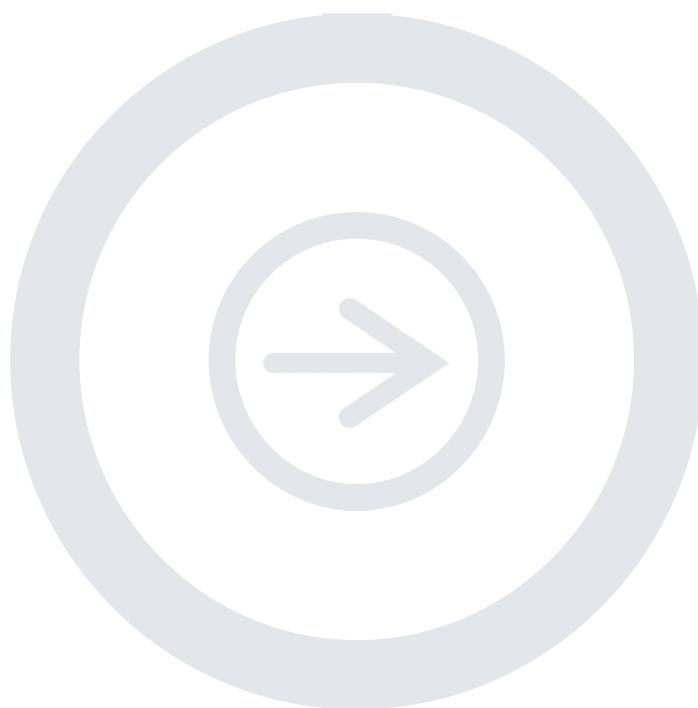
Technical Note

Detection of post-immunoprecipitation proteins by Western blot using the Quick Western Kit–IRDye® 680RD

Developed for:

Aerius,
Odyssey® Classic,
Odyssey CLx, and
Odyssey Sa
Imaging Systems

Please refer to your manual to confirm that this protocol is appropriate for the applications compatible with your Odyssey Imager model.



LI-COR®

Published May 2016. The most recent version of this protocol is posted at <http://www.licor.com/bio/support>

