

Clarity™ Western
ECL Substrate
Transition Guide



Transition Guide. To transition to Clarity, use the following guidelines to adjust exposure times.

Current Substrate	Exposure Using Clarity
GE ECL/Pierce ECL	Decrease 20-fold*
SuperSignal West Pico	Decrease 5-fold*
Bio-Rad® Immun-Star™ HRP	Decrease 5-fold*
Pierce ECL Plus	Equal
Bio-Rad® Immun-Star™ WesternC™	Equal
SuperSignal West Dura	Equal
ECL Prime	Increase 2-fold
SuperSignal West Femto	Increase 5-fold
ECL Select	Increase 5-fold

* Or decrease antibody 2- to 5-fold.

For more information, refer to the full instruction manual available at www.bio-rad.com/clarity.

ECL Select is a trademark of GE Healthcare UK Limited.
SuperSignal is a trademark of Thermo Fisher Scientific Inc.