


In-Cell Western™ Assay – Microplate Setup Guide


Reagents


 Treatment _____
serial dilution from _____ to _____


 Diluent of treatment compound


 Blocking Buffer _____

Primary Antibodies


 Combo #1

 Combo #2

 Combo #3

 Combo #4

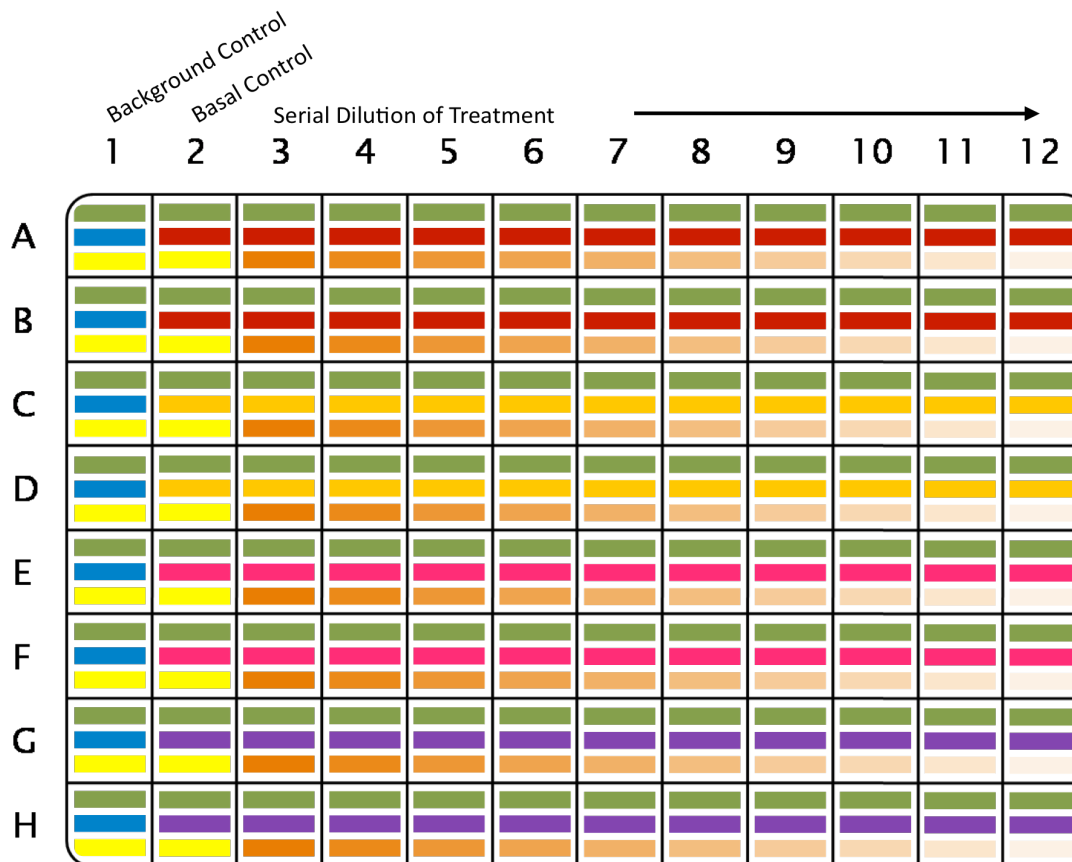
Secondary Antibodies

 _____

LI-COR®

Biosciences

© LI-COR Inc. 2009. LI-COR and In-Cell Western are trademarks or registered trademarks of LI-COR, Inc.



Experimental Conditions:
