



In-Cell Western™ Assay – Microplate Setup Guide

Reagents

 Treatment _____
serial dilution from _____ to _____

 Diluent of treatment compound

 Blocking Buffer

Primary Antibodies

 Combo #1

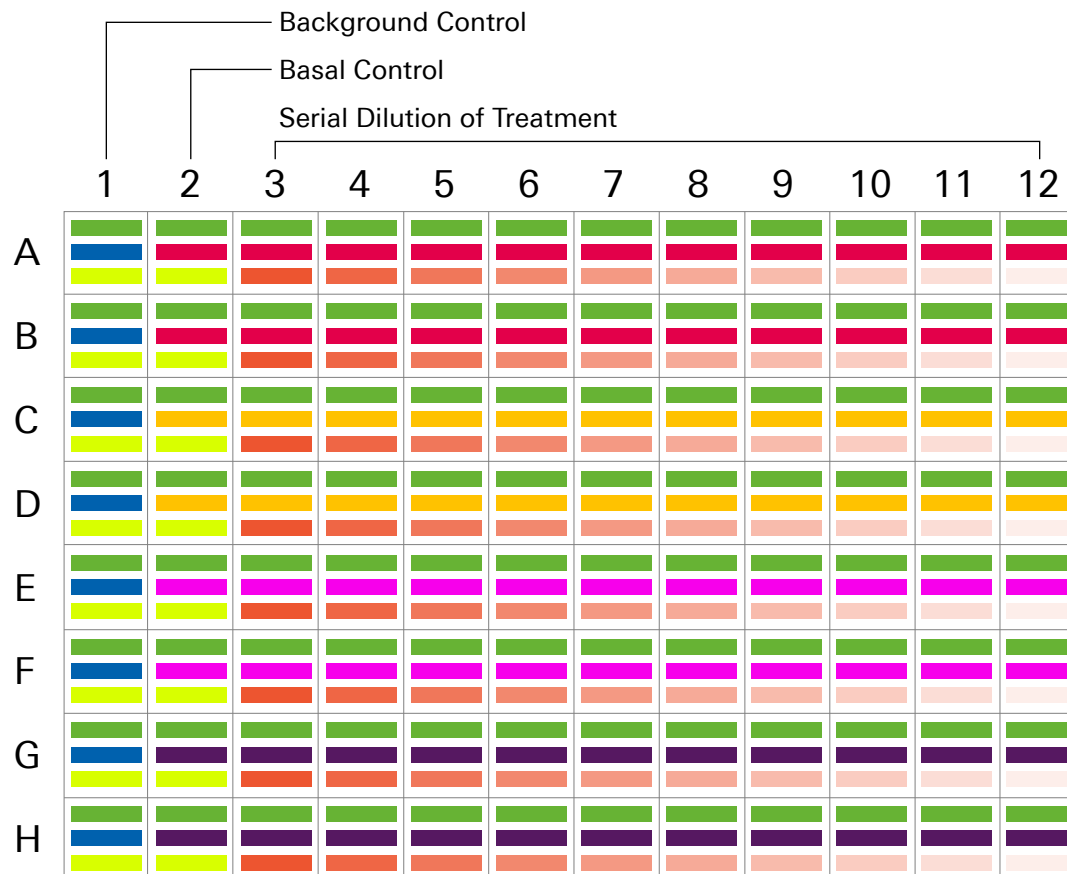
 Combo #2

 Combo #3

 Combo #4

Secondary Antibodies

 _____



Experimental Conditions

