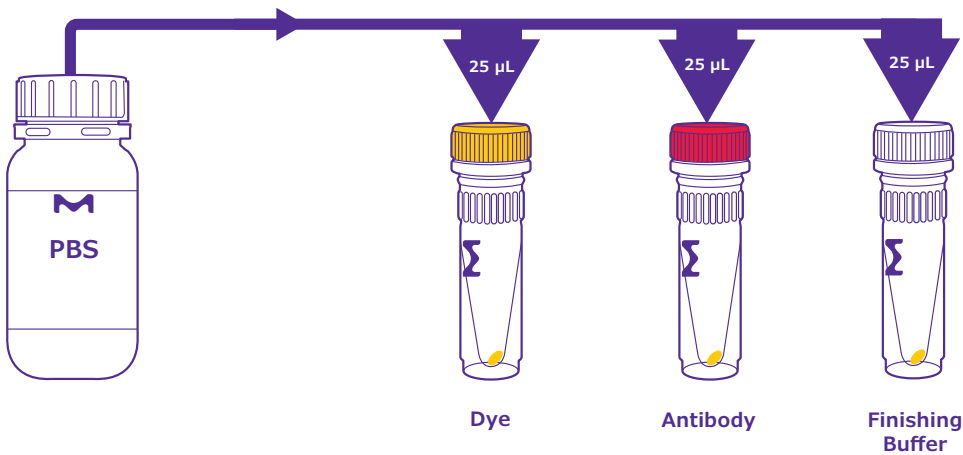
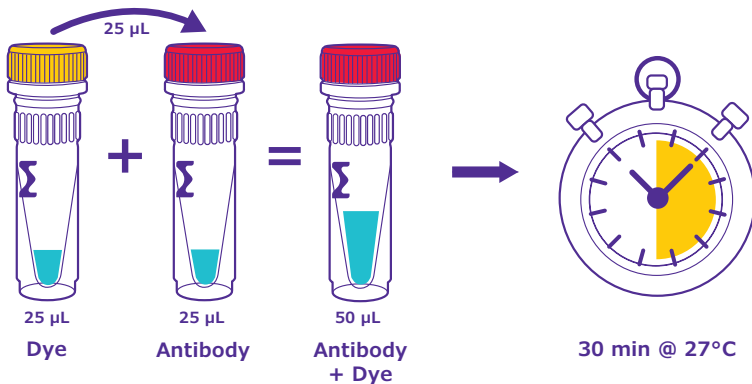


## ColorWheel® 3-Step Protocol

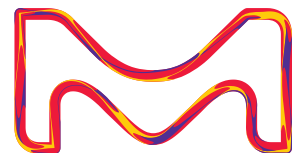
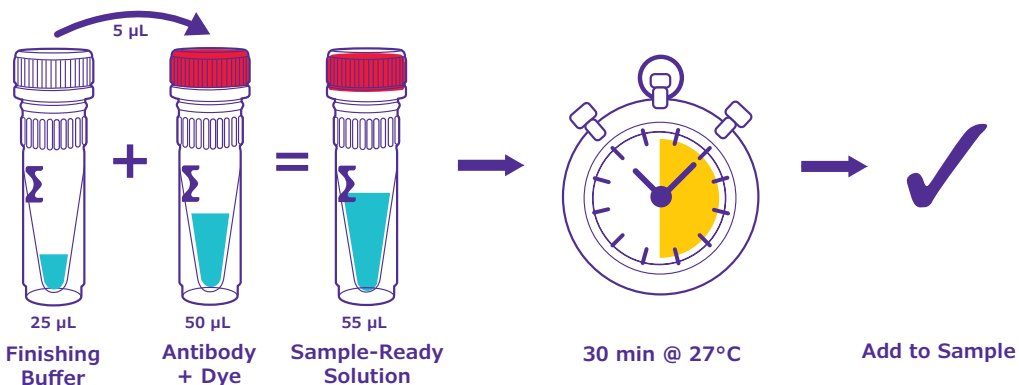
### STEP 1 Spin Down Antibody, Dye, and Finishing Buffer; Resuspend



### STEP 2 Mix Antibody and Dye



### STEP 3 Add Finishing Buffer and Add to Sample



The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

© 2021 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. MilliporeSigma, the vibrant M, Sigma-Aldrich, and CollorWheel are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

MS\_FL8134EN 36633 06/2021