

# Recommended Dilutions and Buffers

The following chart lists many of our reagents, the recommended concentration or dilution, and suggested buffers for various applications. If a specific reagent is not used in one of the listed applications, no buffer will be suggested. Also, although we provide general recommendations, other buffers may often be employed with equivalent results.

Reagent	Working Concentration or Dilution	Tissue Staining	Blots*	ELISAs	Cell Sorting
Unlabeled Antibodies	5-20 µg/ml	A	F	A	
Biotinylated Antibodies	2-10 µg/ml	A	F	A	A
Biotinylated Antibodies (chain specific)	2-5 µg/ml	A	F	A	A
Alkaline Phosphatase Antibodies	1:500-1:1000	A	F	E	
Peroxidase Antibodies	1-5 µg/ml	A	F	F	
Fluorophore Antibodies	5-20 µg/ml	A,B,C	F		B,C
Biotinylated Protein A	2-5 µg/ml	A	F	A	A
Unlabeled Lectins	10-20 µg/ml	B	F	A	B
Biotinylated Lectins	5-20 µg/ml	A,B	F	A	B
Fluorophore Lectins	10-20 µg/ml	A,B	F		B
Anti-Lectins/Antibodies	5-20 µg/ml	A	F	A	A
Biotinylated Anti-Lectins/Antibodies	5-20 µg/ml	A	F	A	A
Biotinylated Alkaline Phosphatase	0.2 units/ml	A	F	E	
Biotinylated Peroxidase	5-10 µg/ml	A	F	F	
Avidin D	10-20 µg/ml	C	F	C	
Avidin DN	10-20 µg/ml	C	F	C	
Alkaline Phosphatase Avidin D	0.2 units/ml	A,C	F	E	
Peroxidase Avidin D	2-5 µg/ml	A,C	F	F	
Fluorescein Avidin DCS	10-20 µg/ml	A,B,C	F		A
Fluorophore Avidin D	10-20 µg/ml	A,B,C	F		A
Anti-Avidin D Antibody	2-5 µg/ml	A	F	A	
Biotinylated Anti-Avidin D	2-5 µg/ml	A	F	A	A
Streptavidin	10-20 µg/ml	A	F	F	
Alkaline Phosphatase Streptavidin	1:500-1:1000	A	F	E	
Peroxidase Streptavidin	2-5 µg/ml	A	F	F	
Fluorophore Streptavidin	10-20 µg/ml	A,C	F		C
Anti-Streptavidin Antibody	2-5 µg/ml	A	F	A	
Biotinylated Anti-Streptavidin Antibody	2-5 µg/ml	A	F	A	A
Anti-Biotin Antibody	2-5 µg/ml	A	F	A	A
Alkaline Phosphatase Anti-Biotin Antibody	1:500-1:1000		F	E	
Peroxidase Anti-Biotin Antibody	1-5 µg/ml	A	F	D	

## Buffers

- |  |                                      |
|--|--------------------------------------|
| A. 10 mM Phosphate, pH 7.5, 0.15 M NaCl          | D. 50 mM TBS, pH 7.5, 0.1% Tween® 20 |
| B. 10 mM HEPES, pH 7.5, 0.15 M NaCl              | E. B + 0.1% Tween 20                 |
| C. 0.1 M Sodium Bicarbonate, pH 8.5, 0.15 M NaCl | F. A + 0.1% Tween 20                 |

\* For chemiluminescence substrates, concentrations are generally lower. Detergents other than Tween 20 may also be used.