

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	Concentration or Dilution	Tissue Staining	Blots*	Elisas	Cell Sorting
Unlabeled Antibodies	5-20 µg/ml	A	G	A	
Biotinylated Antibodies	2-10 µg/ml	A	G	A	A
Biotinylated Antibodies (chain specific)	2-5 µg/ml	A	G	A	A
Alkaline Phosphatase Antibodies	1:500-1:1000	A	G	F	
Peroxidase Antibodies	1-5 µg/ml	A	G	G	
Fluorochrome Antibodies	20 µg/ml	A,B,C	G		B,C
Biotinylated Protein A	2-5 µg/ml	A	G	A	A
Unlabeled Lectins	10-20 µg/ml	B	G	A	B
Biotinylated Lectins	5-20 µg/ml	A,B	G	A	B
Fluorochrome Lectins	10-20 µg/ml	A,B	G		B
Anti-Lectins	5-20 µg/ml	A	G	A	A
Biotinylated Anti-Lectins	5-20 µg/ml	A	G	A	A
Biotinylated IgG	5-20 µg/ml	A	G	A	
Biotinylated Alkaline Phosphatase	0.2 units/ml	A	G	F	
Biotinylated Beta-Galactosidase	0.2 units/ml	A	G	G	
Biotinylated Glucose Oxidase	5 µg/ml	A	G	H	
Biotinylated Peroxidase	5-10 µg/ml	A	G	G	
Avidin D	10-20 µg/ml	C	G	C	
Avidin DN	10-20 µg/ml	C	G	C	
Avidin DX	10-20 µg/ml		G	C	
Alkaline Phosphatase Avidin D	0.2 units/ml	A,C	G	F	
Beta-Galactosidase Avidin D	0.2 units/ml	A,C	G	G	
Glucose Oxidase Avidin D	5 µg/ml	A,C	G	E	
Peroxidase Avidin D	2-5 µg/ml	A,C	G	G	
Phycoerythrin Avidin D	20 µg/ml	A	G		A
Fluorescein Avidin DCS	10-20 µg/ml	A,B,C	G		A
Fluorescein Avidin DN	10-20 µg/ml	A,B,C	G		A
Fluorochrome Avidin D	10-20 µg/ml	A,B,C	G		A
Anti-Avidin D	2-5 µg/ml	A	G	A	
Biotinylated Anti-Avidin D	2-5 µg/ml	A	G	A	A
Peroxidase Anti-Avidin D	1-5 µg/ml	A	G	E	
Alkaline Phosphatase Anti-Avidin D	1:500-1:1000	A	G	F	
Streptavidin	10-20 µg/ml	A	G	G	
Alkaline Phosphatase Streptavidin	1:500-1:1000	A	G	F	
Peroxidase Streptavidin	2-5 µg/ml	A	G	G	
Fluorochrome Streptavidin	10-20 µg/ml	A,C	G		C
Anti-Streptavidin	2-5 µg/ml	A	G	A	
Biotinylated Anti-Streptavidin	2-5 µg/ml	A	G	A	A
Anti-Biotin	2-5 µg/ml	A	G	A	A
Alkaline Phosphatase Anti-Biotin	1:500-1:1000		G	F	
Peroxidase Anti-Biotin	1-5 µg/ml	A	G	E	

Buffers

- A. 10 mM Phosphate, pH 7.8, 0.15M NaCl
- B. 10 mM HEPES, pH 7.5, 0.15M NaCl
- C. 0.1M Sodium Bicarbonate, pH 8.5, 0.15M NaCl
- D. 50 mM Tris, pH 7.5, 0.15M NaCl
- E. D + 0.1% Tween 20
- F. B + 0.1% Tween 20
- G. A + 0.1% Tween 20
- H. 0.1M Sodium Acetate, pH 4.5

* For chemiluminescence, concentrations are generally lower. Blocking reagents other than Tween 20 may also be used.

Vector Laboratories, Inc., 30 Ingold Road, Burlingame, CA 94010 U.S.A., • (650) 697-3600 • Fax: (650) 697-0339

Email: vector@vectorlabs.com • www.vectorlabs.com