

CONSUMABLES

Product	Source	Catalog Number
1.5 mL co-polymer microcentrifuge tubes	USA Scientific	1415-2500
1.5 mL Eppendorf Protein LoBind microcentrifuge tubes	Eppendorf Fisher	022431081 13-698-794
2.0 mL screw-cap microcentrifuge tubes for storage	Fisher	05-669-8
0.2 mL PCR tubes - sterile	Various Vendors	
10 μ L pipette tip refills	Rainin	GPS-10G
250 μ L pipette tip refills	Rainin	GPS-250
1000 μ L pipette tip refills	Rainin	GPS-1000
Zeba Micro Desalt Spin Columns	Pierce	89882 (0.5 mL) 89889 (2 mL)
Slide-A-Lyzer Dialysis Cassettes	Pierce	66330 or 66332
PD-10 column	Amersham Biosciences	17-0851-01
96 well microplate aluminum sealing tape (to reduce light exposure of samples in a plate)	Costar	6570
Aluminum foil	Various Vendors	
Samco transfer pipettes (extra-fine tip, large bulb)	Samco VWR Fisher	233 14670-333 13-711-29
Thermowell 96-well P polycarbonate clear PCR plates - (recommended for XYP heater block)	Costar Fisher	6509 07-200-541
Sealing Mat for 96-well Thermowell P plates	Costar Fisher	6555 07-200-614
Microseal 'A' film	MJ Research	MSA-5001
1.2 μ m PVDF filter microtiter plates (recommended for washed assays)	Millipore	MABVN1250
1 μ m PTFE filter microtiter plates (recommended for washed assays)	Millipore	MAR1N1050
1.2 μ m Supor filter microtiter plates (recommended for washed assays)	Pall Life Sciences VWR	5039 28148-668
96-well black half-area flat bottom plates (recommended for no wash assays)	Costar	3694
96-well half-area flat bottom plates (white, non-treated)	Costar	3693
96-well flat bottom polystyrene solid plates	Costar	3912 (white) 3915 (black)
96-well round bottom polystyrene solid plates	Costar	3789 (white) 3792 (black)