

Each **JBScreen Phosphatase** kit contains 24 unique, sterile filtered reagents, supplied in 10 ml aliquots.

Storage: 4°C

No	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
A1	8 % w/v PEG 4000	None	50 mM HEPES Sodium Salt	7.0	22 mM Lithium Sulfate	0.1 % v/v β -Mercaptoethanol
A2	10 % w/v PEG 4000	None	100 mM HEPES Sodium Salt	7.5	50 mM Ammonium Acetate	None
A3	10 % w/v PEG 4000	10 % v/v Glycerol	50 mM Sodium Succinate	5.5	3 mM Magnesium Chloride	50 mM Tris-Phosphate pH 7.5
A4	12 % w/v PEG 4000	10 % v/v 2-Propanol	100 mM Sodium Citrate	5.6	None	None
A5	12 % w/v PEG 4000	None	200 mM Imidazole Malate	6.0	2 mM Zinc Acetate	None
A6	15 % w/v PEG 4000	None	100 mM HEPES Sodium Salt	7.0	200 mM Magnesium Chloride	None
B1	15 % w/v PEG 4000	15 % v/v PEG 400	50 mM MES Sodium Salt	6.5	200 mM Magnesium Sulfate	50 mM Ammonium Sulfate
B2	15 % w/v PEG 4000	None	85 mM MES Sodium Salt	6.5	170 mM Sodium Acetate	None
B3	20 % w/v PEG 4000	5 % v/v 2-Propanol	100 mM Bicine	9.0	None	None
B4	22 % w/v PEG 4000	10 % v/v 2-Propanol	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate	0.5 % v/v β -Mercaptoethanol
B5	25 % w/v PEG 4000	None	100 mM Bis-Tris Propane	6.5	200 mM di-Ammonium Tartrate	None
B6	30 % w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None	None
C1	30 % w/v PEG 4000	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate	None
C2	30 % w/v PEG 4000	None	100 mM Tris-HCl	7.5	100 mM Cesium Chloride	250 mM Lithium Sulfate
C3	35 % w/v PEG 4000	None	100 mM MES Sodium Salt	6.0	None	None
C4	42 % w/v PEG 4000	None	50 mM Tris-HCl	8.0	100 mM Sodium Perchlorate	None
C5	8 % w/v PEG 6000	10 % v/v 2-Propanol	100 mM Citric Acid	4.0	50 mM Ammonium dihydrogen Phosphate	None
C6	10 % w/v PEG 6000	4 % v/v MPD	100 mM HEPES Sodium Salt	7.5	None	None
D1	10 % w/v PEG 6000	None	100 mM MES Sodium Salt	6.5	None	None
D2	12 % w/v PEG 6000	None	100 mM Sodium Acetate	4.6	None	None
D3	12 % w/v PEG 6000	None	100 mM Tris-Acetate	8.0	None	None
D4	12 % w/v PEG 6000	2% v/v Glycerol	None		25 mM Sodium Phosphate	None
D5	15 % w/v PEG 6000	None	50 mM Sodium Phosphate	7.0	25 mM Potassium Chloride	5 mM Magnesium Sulfate
D6	20 % w/v PEG 6000	None	100 mM Bicine	9.0	None	None