

JBScreen Phosphatase 4

Cat.-No.: CS-174

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	500 mM Ammonium Sulfate	None	100 mM HEPES Sodium Salt	7.5	200 mM Lithium Sulfate	None
A2	1.0 M Ammonium Sulfate	None	100 mM MES Sodium Salt	6.5	100 mM Potassium Chloride	None
A3	1.0 M Ammonium Sulfate	800 mM Potassium Chloride	100 mM HEPES Sodium Salt	7.0	None	None
A4	1.3 M Ammonium Sulfate	None	100 mM CHES	9.5	200 mM Sodium Chloride	None
A5	1.3 M Ammonium Sulfate	None	100 mM Sodium Acetate	5.5	None	None
A6	1.3 M Ammonium Sulfate	5 % v/v Glycerol	100 mM Maleic Acid	6.5	1 mM Magnesium Chloride	5 mM Zinc Chloride
B1	1.5 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate	None
B2	1.7 M Ammonium Sulfate	None	100 mM MES Sodium Salt	6.0	None	None
B3	1.7 M Ammonium Sulfate	4 % v/v MPD	100 mM HEPES Sodium Salt	7.5	None	None
B4	1.7 M Ammonium Sulfate	2 % v/v PEG 400	100 mM Tris-HCl	7.5	None	None
B5	2.0 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	None	None
B6	2.0 M Ammonium Sulfate	10 % v/v PEG 400	50 mM HEPES Sodium Salt	7.0	50 mM Zinc Sulfate	None
C1	2.0 M Ammonium Sulfate	None	100 mM Sodium Citrate	5.6	200 mM Potassium/Sodium Tartrate	None
C2	2.2 M Ammonium Sulfate	6 % v/v PEG 400	100 mM Bis-Tris	6.5	None	None
C3	2.4 M Ammonium Sulfate	None	100 mM HEPES Sodium Salt	7.0	None	None
C4	2.4 M Ammonium Sulfate	None	100 mM Tris-HCl	8.0	10 mM Magnesium Chloride	5 mM Zinc Sulfate
C5	500 mM Lithium Sulfate	10 % v/v PEG 400	100 mM Imidazole-HCl	8.0	None	None
C6	1.5 M Lithium Sulfate	None	100 mM HEPES Sodium Salt	7.5	None	None
D1	2.0 M Lithium Sulfate	2 % v/v PEG 400	100 mM Tris-Acetate	8.0	0.1 % v/v β -Mercaptoethanol	None
D2	20 % v/v MPD	None	100 mM MES Sodium Salt	6.5	None	None
D3	35 % v/v MPD	None	100 mM Sodium Acetate	4.6	None	None
D4	1.4 M Sodium Formate	None	100 mM MES Sodium Salt	6.0	None	None
D5	3.0 M Sodium Formate	None	100 mM Tris-HCl	8.0	None	None
D6	1.0 M Sodium Tartrate	None	100 mM Tris-HCl	8.5	None	None