

JBScreen Phosphatase HTSCat.-No.: **CS-208**

JBScreen Phosphatase comprises 96 unique, sterile filtered reagents, delivered in 96 deep-well blocks, containing either 1.0 ml (S version) or 1.7 ml (L version) aliquots

Storage: 4°C

No	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
A1	15 % v/v PEG 400	None	50 mM Sodium Citrate	5.6	200 mM Sodium Tartrate	200 mM Ammonium Sulfate
A2	22 % v/v PEG 400	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate	None
A3	25 % v/v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Chloride	None
A4	25 % v/v PEG 550 MME	None	100 mM MES Sodium Salt	6.5	10 mM Zinc Sulfate	None
A5	27 % w/v PEG 1500	10 % v/v MPD	10 mM Imidazole-HCl	7.0	220 mM Sodium Nitrate	None
A6	30 % w/v PEG 1500	None	100 mM Bis-Tris	6.5	400 mM Potassium Fluoride	None
A7	8 % w/v PEG 2000 MME	None	10 mM Tris-HCl	7.0	10 mM Nickel (II) Chloride	None
A8	16 % w/v PEG 2000 MME	None	10 mM Sodium Acetate	4.6	400 mM Ammonium Sulfate	None
A9	8 % w/v PEG 3350	None	None		100 mM Sodium Acetate	None
A10	10 % w/v PEG 3350	None	100 mM HEPES Sodium Salt	7.5	200 mM Ammonium Acetate	None
A11	10 % w/v PEG 3350	None	100 mM HEPES Sodium Salt	7.5	50 mM Ammonium Acetate	15 mM Sodium Chloride
A12	10 % w/v PEG 3350	None	100 mM MES Sodium Salt	6.5	20 mM Zinc Acetate	None
B1	15 % w/v PEG 3350	None	None		200 mM Sodium Fluoride	None
B2	15 % w/v PEG 3350	None	100 mM Bis-Tris	6.5	50 mM Glycine	200 mM Sodium Formate
B3	15 % w/v PEG 3350	5 % v/v Glycerol	100 mM Bis-Tris	7.0	200 mM Magnesium Formate	None
B4	20 % w/v PEG 3350	None	None		200 mM Potassium Fluoride	None
B5	20 % w/v PEG 3350	None	100 mM MES Sodium Salt	6.5	100 mM Ammonium Chloride	None
B6	20 % w/v PEG 3350	None	10 mM Tris-HCl	8.0	200 mM Potassium Sulfate	None
B7	20 % w/v PEG 3350	None	None		100 mM Sodium Citrate	None
B8	22 % w/v PEG 3350	None	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Chloride	None
B9	25 % w/v PEG 3350	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride	None
B10	25 % w/v PEG 3350	10 % v/v 2-Propanol	100 mM Sodium Citrate	4.8	100 mM Lithium Chloride	None
B11	30 % w/v PEG 3350	None	None		200 mM Ammonium Nitrate	None
B12	30 % w/v PEG 3350	None	None		200 mM Ammonium dihydrogen Phosphate	None

JBScreen Phosphatase HTS

Cat.-No.: CS-208

No	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
C1	8 % w/v PEG 4000	None	50 mM HEPES Sodium Salt	7.0	22 mM Lithium Sulfate	0.1 % v/v β -Mercaptoethanol
C2	10 % w/v PEG 4000	None	100 mM HEPES Sodium Salt	7.5	50 mM Ammonium Acetate	None
C3	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
C4	12 % w/v PEG 4000	10 % v/v 2-Propanol	100 mM Sodium Citrate	5.6	None	None
C5	12 % w/v PEG 4000	None	200 mM Imidazole Malate	6.0	2 mM Zinc Acetate	None
C6	15 % w/v PEG 4000	None	100 mM HEPES Sodium Salt	7.0	200 mM Magnesium Chloride	None
C7	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
C8	15 % w/v PEG 4000	None	85 mM MES Sodium Salt	6.5	170 mM Sodium Acetate	None
C9	20 % w/v PEG 4000	5 % v/v 2-Propanol	100 mM Bicine	9.0	None	None
C10	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
C11	25 % w/v PEG 4000	None	100 mM Bis-Tris Propane	6.5	200 mM di-Ammonium Tartrate	None
C12	30 % w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None	None
D1	30 % w/v PEG 4000	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate	None
D2	30 % w/v PEG 4000	None	100 mM Tris-HCl	7.5	100 mM Cesium Chloride	250 mM Lithium Sulfate
D3	35 % w/v PEG 4000	None	100 mM MES Sodium Salt	6.0	None	None
D4	42 % w/v PEG 4000	None	50 mM Tris-HCl	8.0	100 mM Sodium Perchlorate	None
D5	8 % w/v PEG 6000	10 % v/v 2-Propanol	100 mM Citric Acid	4.0	50 mM Ammonium dihydrogen Phosphate	None
D6	10 % w/v PEG 6000	4 % v/v MPD	100 mM HEPES Sodium Salt	7.5	None	None
D7	10 % w/v PEG 6000	None	100 mM MES Sodium Salt	6.5	None	None
D8	12 % w/v PEG 6000	None	100 mM Sodium Acetate	4.6	None	None
D9	12 % w/v PEG 6000	None	100 mM Tris-Acetate	8.0	None	None
D10	12 % w/v PEG 6000	2% v/v Glycerol	None		25 mM Sodium Phosphate	None
D11	15 % w/v PEG 6000	None	50 mM Sodium Phosphate	7.0	25 mM Potassium Chloride	5 mM Magnesium Sulfate
D12	20 % w/v PEG 6000	None	100 mM Bicine	9.0	None	None

No	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
E1	8 % w/v PEG 8000	None	None		100 mM Potassium Phosphate	0,2 % v/v β -Mercaptoethanol
E2	10 % w/v PEG 8000	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Acetate	None
E3	10 % w/v PEG 8000	None	100 mM MES Sodium Salt	6.5	200 mM Potassium Chloride	100 mM Magnesium Acetate
E4	10 % w/v PEG 8000	None	100 mM Tris-HCl	8.5	50 mM Sodium Chloride	None
E5	12 % w/v PEG 8000	None	50 mM Sodium Acetate	5.0	300 mM Ammonium Acetate	None
E6	12 % w/v PEG 8000	None	100 mM Tris-HCl	6.3	200 mM Lithium Sulfate	None
E7	12 % w/v PEG 8000	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Acetate	None
E8	13 % w/v PEG 8000	15 % v/v Glycerol	100 mM MOPS	7.0	500 mM Ammonium Sulfate	None
E9	15 % w/v PEG 8000	None	100 mM Potassium Phosphate	4.6	None	None
E10	15 % w/v PEG 8000	None	100 mM Sodium Phosphate	4.6	None	None
E11	15 % w/v PEG 8000	None	100 mM Sodium Succinate	5.5	250 mM Lithium Sulfate	None
E12	18 % w/v PEG 8000	None	100 mM HEPES Sodium Salt	7.0	100 mM Potassium dihydrogen Phosphate	None
F1	18 % w/v PEG 8000	None	None		200 mM Lithium Sulfate	None
F2	18 % w/v PEG 8000	5 % v/v 1,4-Dioxane	100 mM Tris-HCl	8.0	200 mM Sodium Fluoride	None
F3	25 % w/v PEG 8000	5 % v/v Jeffamine M 600	90 mM MES Sodium Salt	6.0	90 mM Magnesium Acetate	None
F4	25 % w/v PEG 8000	None	100 mM Tris-HCl	8.5	200 mM Magnesium Sulfate	None
F5	5 % w/v PEG 10000	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride	None
F6	10 % w/v PEG 10000	None	50 mM Sodium Acetate	5.0	None	None
F7	10 % w/v PEG 10000	None	100 mM HEPES Sodium Salt	7.5	None	None
F8	15 % w/v PEG 10000	None	100 mM Bis-Tris	5.5	100 mM Ammonium Acetate	5 mM LDAO
F9	15 % w/v PEG 10000	None	60 mM Glycine	9.0	20 mM Sodium Citrate	None
F10	20 % w/v PEG 10000	5 % v/v PEG 550 MME	100 mM Sodium Citrate	5.6	None	None
F11	12 % w/v PEG 20000	None	100 mM MES Sodium Salt	6.5	None	None
F12	17 % w/v PEG 20000	None	100 mM Tris-HCl	8.5	100 mM Magnesium Chloride	None

No	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
G1	500 mM Ammonium Sulfate	None	100 mM HEPES Sodium Salt	7.5	200 mM Lithium Sulfate	None
G2	1.0 M Ammonium Sulfate	None	100 mM MES Sodium Salt	6.5	100 mM Potassium Chloride	None
G3	1.0 M Ammonium Sulfate	800 mM Potassium Chloride	100 mM HEPES Sodium Salt	7.0	None	None
G4	1.3 M Ammonium Sulfate	None	100 mM CHES	9.5	200 mM Sodium Chloride	None
G5	1.3 M Ammonium Sulfate	None	100 mM Sodium Acetate	5.5	None	None
G6	1.3 M Ammonium Sulfate	5 % v/v Glycerol	100 mM Maleic Acid	6.5	1 mM Magnesium Chloride	5 mM Zinc Chloride
G7	1.5 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate	None
G8	1.7 M Ammonium Sulfate	None	100 mM MES Sodium Salt	6.0	None	None
G9	1.7 M Ammonium Sulfate	4 % v/v MPD	100 mM HEPES Sodium Salt	7.5	None	None
G10	1.7 M Ammonium Sulfate	2 % v/v PEG 400	100 mM Tris-HCl	7.5	None	None
G11	2.0 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	None	None
G12	2.0 M Ammonium Sulfate	10 % v/v PEG 400	50 mM HEPES Sodium Salt	7.0	50 mM Zinc Sulfate	None
H1	2.0 M Ammonium Sulfate	None	100 mM Sodium Citrate	5.6	200 mM Potassium/Sodium Tartrate	None
H2	2.2 M Ammonium Sulfate	6 % v/v PEG 400	100 mM Bis-Tris	6.5	None	None
H3	2.4 M Ammonium Sulfate	None	100 mM HEPES Sodium Salt	7.0	None	None
H4	2.4 M Ammonium Sulfate	None	100 mM Tris-HCl	8.0	10 mM Magnesium Chloride	5 mM Zinc Sulfate
H5	500 mM Lithium Sulfate	10 % v/v PEG 400	100 mM Imidazole-HCl	8.0	None	None
H6	1.5 M Lithium Sulfate	None	100 mM HEPES Sodium Salt	7.5	None	None
H7	2.0 M Lithium Sulfate	2 % v/v PEG 400	100 mM Tris-Acetate	8.0	0.1 % v/v β- Mercaptoethanol	None
H8	20 % v/v MPD	None	100 mM MES Sodium Salt	6.5	None	None
H9	35 % v/v MPD	None	100 mM Sodium Acetate	4.6	None	None
H10	1.4 M Sodium Formate	None	100 mM MES Sodium Salt	6.0	None	None
H11	3.0 M Sodium Formate	None	100 mM Tris-HCl	8.0	None	None
H12	1.0 M Sodium Tartrate	None	100 mM Tris-HCl	8.5	None	None