

JBScreen PEG/Salt 3

Cat.-No: CS-143

Number	Precipitant	Salt
A 1	20% w/v PEG 5000 MME	200 mM Ammonium Acetate
A 2	20% w/v PEG 5000 MME	200 mM Ammonium Chloride
A 3	20% w/v PEG 5000 MME	200 mM Ammonium Fluoride
A 4	20% w/v PEG 5000 MME	200 mM Ammonium Formate
A 5	20% w/v PEG 5000 MME	200 mM Ammonium Iodide
A 6	20% w/v PEG 5000 MME	200 mM Ammonium Nitrate
B 1	20% w/v PEG 5000 MME	200 mM Ammonium dihydrogen Phosphate
B 2	20% w/v PEG 5000 MME	200 mM di-Ammonium hydrogen Phosphate
B 3	20% w/v PEG 5000 MME	200 mM Ammonium Sulfate
B 4	20% w/v PEG 5000 MME	200 mM Ammonium Sulfite
B 5	20% w/v PEG 5000 MME	200 mM Calcium Acetate
B 6	20% w/v PEG 5000 MME	200 mM Calcium Chloride
C 1	20% w/v PEG 5000 MME	200 mM di-Ammonium Tartrate
C 2	20% w/v PEG 5000 MME	200 mM Potassium Formate
C 3	20% w/v PEG 5000 MME	200 mM Lithium Acetate
C 4	20% w/v PEG 5000 MME	200 mM Lithium Chloride
C 5	20% w/v PEG 5000 MME	200 mM Lithium Citrate
C 6	20% w/v PEG 5000 MME	200 mM Lithium Nitrate
D 1	20% w/v PEG 5000 MME	200 mM Lithium Sulfate
D 2	20% w/v PEG 5000 MME	200 mM Magnesium Acetate
D 3	20% w/v PEG 5000 MME	200 mM Magnesium Chloride
D 4	20% w/v PEG 5000 MME	200 mM Magnesium Formate
D 5	20% w/v PEG 5000 MME	200 mM Magnesium Nitrate
D 6	20% w/v PEG 5000 MME	200 mM Magnesium Sulfate