

JBScreen PEG/Salt 2

Cat.-No: CS-142

Number	Precipitant	Salt
A 1	20% w/v PEG 3350	200 mM Potassium Acetate
A 2	20% w/v PEG 3350	200 mM Potassium Chloride
A 3	20% w/v PEG 3350	200 mM Potassium Fluoride
A 4	20% w/v PEG 3350	200 mM Potassium Iodide
A 5	20% w/v PEG 3350	200 mM Potassium Nitrate
A 6	20% w/v PEG 3350	200 mM Potassium dihydrogen Phosphate
B 1	20% w/v PEG 3350	200 mM di-Potassium hydrogen Phosphate
B 2	20% w/v PEG 3350	200 mM Potassium Sulfate
B 3	20% w/v PEG 3350	200 mM Potassium Thiocyanate
B 4	20% w/v PEG 3350	200 mM Potassium / Sodium Tartrate
B 5	20% w/v PEG 3350	200 mM Sodium Acetate
B 6	20% w/v PEG 3350	200 mM Sodium Chloride
C 1	20% w/v PEG 3350	200 mM Sodium Citrate
C 2	20% w/v PEG 3350	200 mM Sodium Fluoride
C 3	20% w/v PEG 3350	200 mM Sodium Formate
C 4	20% w/v PEG 3350	200 mM Sodium Iodide
C 5	20% w/v PEG 3350	200 mM Sodium Isothiocyanate
C 6	20% w/v PEG 3350	200 mM Sodium Nitrate
D 1	20% w/v PEG 3350	200 mM Sodium dihydrogen Phosphate
D 2	20% w/v PEG 3350	200 mM di-Sodium hydrogen Phosphate
D 3	20% w/v PEG 3350	200 mM Sodium Sulfate
D 4	20% w/v PEG 3350	200 mM di-Sodium Tartrate
D 5	20% w/v PEG 3350	200 mM tri-Potassium Citrate
D 6	20% w/v PEG 3350	200 mM Zinc Acetate