



JBScreen Basic HTS

Cat.-No.: CS-203L

SCREEN FORMULATION



| No. | Precipitant | Precipitant 2 | Buffer | Additive |
|-----|---|------------------------------------|-----------------------------------|---------------------------------------|
| A1 | 25 % v/v Ethylene glycol | none | none | none |
| A2 | 12 % v/v Glycerol | 1.5 M Ammonium sulfate | 100 mM TRIS; pH 8.5 | none |
| A3 | 1 M 1,6-Hexanediol | none | 100 mM Sodium acetate; pH 4.6 | 10 mM Cobalt (II) chloride |
| A4 | 2.5 M 1,6-Hexanediol | none | 100 mM tri-Sodium citrate; pH 5.6 | none |
| A5 | 3.4 M 1,6-Hexanediol | none | 100 mM TRIS; pH 8.5 | 200 mM Magnesium chloride |
| A6 | 30 % v/v 2-Methyl-2,4-pentanediol | none | 100 mM Sodium acetate; pH 4.6 | 200 mM Sodium chloride |
| A7 | 30 % v/v 2-Methyl-2,4-pentanediol | none | 100 mM tri-Sodium citrate; pH 5.6 | 200 mM Ammonium acetate |
| A8 | 30 % v/v 2-Methyl-2,4-pentanediol | none | 100 mM Sodium acetate; pH 4.6 | 20 mM Calcium chloride |
| A9 | 30 % v/v 2-Methyl-2,4-pentanediol | 500 mM Ammonium sulfate | 100 mM HEPES; pH 7.5 | none |
| A10 | 30 % v/v 2-Methyl-2,4-pentanediol | none | 100 mM HEPES; pH 7.5 | 200 mM tri-Sodium citrate |
| A11 | 50 % v/v 2-Methyl-2,4-pentanediol | none | 100 mM TRIS; pH 8.5 | 200 mM Ammonium di-hydrogen phosphate |
| A12 | 70 % v/v 2-Methyl-2,4-pentanediol | none | 100 mM HEPES; pH 7.5 | none |
| B1 | 2 % w/v Ethylene imine polymer | none | 100 mM tri-Sodium citrate; pH 5.6 | 500 mM Sodium chloride |
| B2 | 2 % v/v Polyethylene glycol 400 | 2 M Ammonium sulfate | 100 mM HEPES; pH 7.5 | none |
| B3 | 28 % v/v Polyethylene glycol 400 | none | 100 mM HEPES; pH 7.5 | 200 mM Calcium chloride |
| B4 | 30 % v/v Polyethylene glycol 400 | none | 100 mM TRIS; pH 8.5 | 200 mM tri-Sodium citrate |
| B5 | 30 % v/v Polyethylene glycol 400 | none | 100 mM HEPES; pH 7.5 | 200 mM Magnesium chloride |
| B6 | 30 % v/v Polyethylene glycol 400 | none | 100 mM Sodium acetate; pH 4.6 | 100 mM Calcium chloride |
| B7 | 20 % v/v Polyethylene glycol monomethyl ether 550 | none | 100 mM BICINE; pH 9.5 | 100 mM Sodium chloride |
| B8 | 25 % v/v Polyethylene glycol monomethyl ether 550 | none | 100 mM MES; pH 6.5 | 10 mM Zinc sulfate |
| B9 | 10 % w/v Polyethylene glycol 1,000 | 10 % w/v Polyethylene glycol 8,000 | none | none |
| B10 | 30 % w/v Polyethylene glycol 1,500 | none | none | none |
| B11 | 20 % w/v Polyethylene glycol monomethyl ether 2,000 | none | 100 mM TRIS; pH 8.5 | 10 mM Nickel (II) chloride |
| B12 | 30 % w/v Polyethylene glycol monomethyl ether 2,000 | none | 100 mM Sodium acetate; pH 4.6 | 200 mM Ammonium sulfate |

*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components

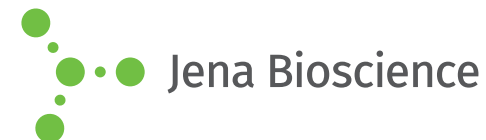




JBScreen Basic HTS

Cat.-No.: CS-203L

SCREEN FORMULATION



| No. | Precipitant | Precipitant 2 | Buffer | Additive |
|-----|---|----------------------------------|-----------------------------------|---------------------------------------|
| C1 | 8 % w/v Polyethylene glycol 4,000 | none | 100 mM Sodium acetate; pH 4.6 | none |
| C2 | 20 % w/v Polyethylene glycol 4,000 | 20 % v/v 2-Propanol | 100 mM tri-Sodium citrate; pH 5.6 | none |
| C3 | 20 % w/v Polyethylene glycol 4,000 | 10 % v/v 2-Propanol | 100 mM HEPES; pH 7.5 | none |
| C4 | 25 % w/v Polyethylene glycol 4,000 | none | 100 mM Sodium acetate; pH 4.6 | 200 mM Ammonium sulfate |
| C5 | 30 % w/v Polyethylene glycol 4,000 | none | none | 200 mM Ammonium sulfate |
| C6 | 30 % w/v Polyethylene glycol 4,000 | none | 100 mM Sodium acetate; pH 4.6 | 200 mM Ammonium acetate |
| C7 | 30 % w/v Polyethylene glycol 4,000 | none | 100 mM tri-Sodium citrate; pH 5.6 | 200 mM Ammonium acetate |
| C8 | 30 % w/v Polyethylene glycol 4,000 | none | 100 mM TRIS; pH 8.5 | 200 mM Sodium acetate |
| C9 | 30 % w/v Polyethylene glycol 4,000 | none | 100 mM TRIS; pH 8.5 | 200 mM Lithium sulfate |
| C10 | 30 % w/v Polyethylene glycol 4,000 | none | 100 mM TRIS; pH 8.5 | 200 mM Magnesium chloride |
| C11 | 30 % w/v Polyethylene glycol monomethyl ether 5,000 | none | 100 mM MES; pH 6.5 | 200 mM Ammonium sulfate |
| C12 | 10 % w/v Polyethylene glycol 6,000 | 2 M Sodium chloride | none | none |
| D1 | 10 % w/v Polyethylene glycol 6,000 | 5 % v/v 2-Methyl-2,4-pentanediol | 100 mM HEPES; pH 7.5 | none |
| D2 | 2 % w/v Polyethylene glycol 8,000 | 1 M Lithium sulfate | none | none |
| D3 | 8 % w/v Polyethylene glycol 8,000 | none | 100 mM TRIS; pH 8.5 | none |
| D4 | 10 % w/v Polyethylene glycol 8,000 | 8 % v/v Ethylene glycol | 100 mM HEPES; pH 7.5 | none |
| D5 | 15 % w/v Polyethylene glycol 8,000 | 500 mM Lithium sulfate | none | none |
| D6 | 18 % w/v Polyethylene glycol 8,000 | none | 100 mM MES; pH 6.5 | 200 mM Calcium acetate |
| D7 | 18 % w/v Polyethylene glycol 8,000 | none | 100 mM MES; pH 6.5 | 200 mM Zinc acetate |
| D8 | 20 % w/v Polyethylene glycol 8,000 | none | none | 50 mM Potassium di-hydrogen phosphate |
| D9 | 20 % w/v Polyethylene glycol 8,000 | none | 100 mM MES; pH 6.5 | 200 mM Magnesium acetate |
| D10 | 30 % w/v Polyethylene glycol 8,000 | none | 100 mM MES; pH 6.5 | 200 mM Sodium acetate |
| D11 | 30 % w/v Polyethylene glycol 8,000 | none | none | 200 mM Ammonium sulfate |
| D12 | 30 % w/v Polyethylene glycol 8,000 | none | 100 mM MES; pH 6.5 | 200 mM Ammonium sulfate |

*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components

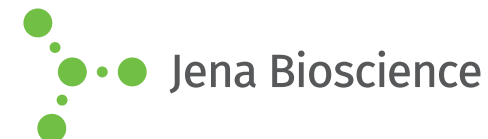




JBScreen Basic HTS

Cat.-No.: CS-203L

SCREEN FORMULATION



| No. | Precipitant | Precipitant 2 | Buffer | Additive |
|-----|-------------------------------------|------------------------|-----------------------------------|----------------------------|
| E1 | 10 % w/v Polyethylene glycol 10,000 | 2 % v/v 1,4-Dioxane | 100 mM BICINE; pH 9.5 | none |
| E2 | 20 % w/v Polyethylene glycol 10,000 | none | 100 mM HEPES; pH 7.5 | none |
| E3 | 12 % w/v Polyethylene glycol 20,000 | none | 100 mM MES; pH 6.5 | none |
| E4 | 5 % v/v 2-Propanol | 2 M Ammonium sulfate | none | none |
| E5 | 20 % v/v 2-Propanol | none | 100 mM HEPES; pH 7.5 | 200 mM tri-Sodium citrate |
| E6 | 20 % v/v 2-Propanol | none | 100 mM Sodium acetate; pH 4.6 | 200 mM Calcium chloride |
| E7 | 30 % v/v 2-Propanol | none | 100 mM HEPES; pH 7.5 | 200 mM Magnesium chloride |
| E8 | 30 % v/v 2-Propanol | none | 100 mM TRIS; pH 8.5 | 200 mM Ammonium acetate |
| E9 | 10 % v/v 1,4-Dioxane | 1.6 M Ammonium sulfate | 100 mM MES; pH 6.5 | none |
| E10 | 35 % v/v 1,4-Dioxane | none | none | none |
| E11 | 10 % v/v Ethanol | 1.5 M Sodium chloride | none | none |
| E12 | 20 % v/v Ethanol | none | 100 mM TRIS; pH 8.5 | none |
| F1 | 25 % v/v 2-Methyl-2-propanol | none | 100 mM TRIS; pH 8.5 | none |
| F2 | 35 % v/v 2-Methyl-2-propanol | none | 100 mM tri-Sodium citrate; pH 5.6 | none |
| F3 | 1 M Imidazole; pH 7.0 | none | none | none |
| F4 | 1 M Lithium sulfate | none | 100 mM TRIS; pH 8.5 | 10 mM Nickel (II) chloride |
| F5 | 1.5 M Lithium sulfate | none | 100 mM HEPES; pH 7.5 | none |
| F6 | 400 mM Potassium Sodium tartrate | none | none | none |
| F7 | 800 mM Potassium Sodium tartrate | none | 100 mM HEPES; pH 7.5 | none |
| F8 | 1.4 M tri-Sodium citrate | none | 100 mM HEPES; pH 7.5 | none |
| F9 | 1.6 M tri-Sodium citrate; pH 6.5 | none | none | none |
| F10 | 10 % v/v Jeffamine® M-600 | none | 100 mM tri-Sodium citrate; pH 5.6 | 10 mM Iron (III) chloride |
| F11 | 20 % v/v Jeffamine® M-600 | none | 100 mM HEPES; pH 7.5 | none |
| F12 | 30 % v/v Jeffamine® M-600 | none | 100 mM MES; pH 6.5 | 50 mM Cesium chloride |

*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components





| No. | Precipitant | Precipitant 2 | Buffer | Additive |
|-----|--|-------------------------------------|-----------------------------------|---|
| G1 | 800 mM Potassium di-hydrogen phosphate | 800 mM Sodium di-hydrogen phosphate | 100 mM HEPES; pH 7.5 | none |
| G2 | 400 mM Ammonium di-hydrogen phosphate | none | none | none |
| G3 | 1 M Ammonium di-hydrogen phosphate | none | 100 mM tri-Sodium citrate; pH 5.6 | none |
| G4 | 2 M Ammonium di-hydrogen phosphate | none | 100 mM TRIS; pH 8.5 | none |
| G5 | 2 M Ammonium formate | none | 100 mM Sodium acetate; pH 4.6 | none |
| G6 | 4 M Ammonium formate | none | 100 mM HEPES; pH 7.5 | none |
| G7 | 2 M Ammonium formate | none | none | none |
| G8 | 500 mM Ammonium sulfate | 1 M Lithium sulfate | 100 mM tri-Sodium citrate; pH 5.6 | none |
| G9 | 1.6 M Ammonium sulfate | none | 100 mM HEPES; pH 7.5 | 100 mM Sodium chloride |
| G10 | 1.8 M Ammonium sulfate | none | 100 mM MES; pH 6.5 | 10 mM Cobalt (II) chloride |
| G11 | 2 M Ammonium sulfate | none | 100 mM TRIS; pH 8.5 | none |
| G12 | 2 M Ammonium sulfate | none | none | none |
| H1 | 2 M Ammonium sulfate | none | 100 mM Sodium acetate; pH 4.6 | none |
| H2 | 2 M Ammonium sulfate | none | 100 mM tri-Sodium citrate; pH 5.6 | 200 mM Potassium Sodium tartrate |
| H3 | 200 mM Magnesium formate | none | none | none |
| H4 | 1.6 M Magnesium sulfate | none | 100 mM MES; pH 6.5 | none |
| H5 | 2 M Magnesium chloride | none | 100 mM BICINE; pH 9.5 | none |
| H6 | 1 M Sodium acetate | none | 100 mM Imidazole; pH 6.5 | none |
| H7 | 1 M Sodium acetate | none | 100 mM HEPES; pH 7.5 | 50 mM Cadmium sulfate |
| H8 | 1.4 M Sodium acetate | none | 100 mM MES; pH 6.5 | none |
| H9 | 500 mM Sodium chloride | 10 mM Magnesium chloride | none | 10 mM Cetyltrimethylammonium bromide |
| H10 | 2 M Sodium chloride | none | 100 mM Sodium acetate; pH 4.6 | none |
| H11 | 2 M Sodium chloride | none | 100 mM MES; pH 6.5 | 100 mM Sodium di-hydrogen phosphate, 100 mM Potassium di-hydrogen phosphate |
| H12 | 4.3 M Sodium chloride | none | 100 mM HEPES; pH 7.5 | none |

*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components