



JBScreen Plus HTS

Cat.-No.: CS-507L

SCREEN FORMULATION



| No. | Additive | Type | CAS Number |
|-----|---------------------------------------|--------------------|------------|
| A1 | 1.00 M Trimethylamine N-oxide | Zwitterion | 62637-93-8 |
| A2 | 1.00 M L-Proline | Zwitterion | 147-85-3 |
| A3 | 100 mM Ectoine | Zwitterion | 96702-03-3 |
| A4 | 1.00 M Glycine | Zwitterion | 56-40-6 |
| A5 | 100 mM Betaine | Zwitterion | 590-47-6 |
| A6 | 100 mM 2-Aminoethanesulfonic acid | Zwitterion | 107-35-7 |
| A7 | 2.00 M Glycerol | Polyalcohol | 56-81-5 |
| A8 | 2.00 M Erythritol | Polyalcohol | 149-32-6 |
| A9 | 2.00 M Xylitol | Polyalcohol | 87-99-0 |
| A10 | 2.00 M Adonitol | Polyalcohol | 488-81-3 |
| A11 | 800 mM Mannitol | Polyalcohol | 69-65-8 |
| A12 | 1.00 M D-Sorbitol | Polyalcohol | 50-70-4 |
| B1 | 200 mM Aluminium chloride | Kosmotropic Cation | 7446-70-0 |
| B2 | 200 mM Magnesium chloride | Kosmotropic Cation | 7791-18-6 |
| B3 | 200 mM Calcium chloride | Kosmotropic Cation | 10035-04-8 |
| B4 | 2.00 M Lithium chloride | Kosmotropic Cation | 7447-41-8 |
| B5 | 1.00 M Manganese (II) chloride | Kosmotropic Cation | 13446-34-9 |
| B6 | 1.00 M Zinc chloride | Kosmotropic Cation | 7646-85-7 |
| B7 | 2.00 M di-Sodium malonate | Kosmotropic Anion | 26522-85-0 |
| B8 | 1.8 M tri-Sodium citrate | Kosmotropic Anion | 6132-04-3 |
| B9 | 1.00 M Sodium fluoride | Kosmotropic Anion | 7681-49-4 |
| B10 | 2.00 M Ammonium sulfate | Kosmotropic Anion | 7783-20-2 |
| B11 | 1.00 M di-Ammonium hydrogen phosphate | Kosmotropic Anion | 7783-28-0 |
| B12 | 1.00 M Ammonium formate | Kosmotropic Anion | 540-69-2 |

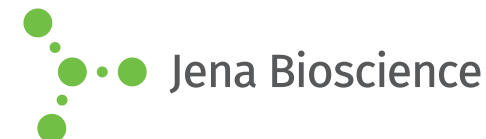




JBScreen Plus HTS

Cat.-No.: CS-507L

SCREEN FORMULATION



| No. | Additive | Type | CAS Number |
|-----|-------------------------------------|-------------------|------------|
| C1 | 1.00 M Urea | Non ionic | 57-13-6 |
| C2 | 4.00 M Dimethyl sulfoxide | Non ionic | 67-68-5 |
| C3 | 4.00 M Dimethylformamide | Non ionic | 68-12-2 |
| C4 | 1.00 M Pyridine | Non ionic | 110-86-1 |
| C5 | 500 mM e-Caprolactam | Non ionic | 105-60-2 |
| C6 | 100 mM Phenol | Non ionic | 108-95-2 |
| C7 | 2.00 M Trimethylamine hydrochloride | Chaotropic Cation | 593-81-7 |
| C8 | 2.00 M Guanidine hydrochloride | Chaotropic Cation | 50-01-1 |
| C9 | 2.00 M Tetramethylammonium Chloride | Chaotropic Cation | 75-57-0 |
| C10 | 1.00 M Cesium chloride | Chaotropic Cation | 7647-17-8 |
| C11 | 1.00 M Rubidium chloride | Chaotropic Cation | 7791-11-9 |
| C12 | 1.00 M Potassium chloride | Chaotropic Cation | 7447-40-7 |
| D1 | 2.00 M Potassium thiocyanate | Chaotropic Anion | 333-20-0 |
| D2 | 2.00 M Potassium Cyanate | Chaotropic Anion | 590-28-3 |
| D3 | 1.00 M Potassium nitrate | Chaotropic Anion | 7757-79-1 |
| D4 | 1.00 M Potassium acetate | Chaotropic Anion | 127-08-2 |
| D5 | 1.00 M Potassium iodide | Chaotropic Anion | 7681-11-0 |
| D6 | 1.00 M Potassium bromide | Chaotropic Anion | 7758-02-3 |
| D7 | 2.00 M Sodium trichloroacetate | Chaotropic Anion | 650-51-1 |
| D8 | 2.00 M Sodium thiocyanate | Chaotropic Anion | 540-72-7 |
| D9 | 2.00 M Sodium perchlorate | Chaotropic Anion | 7601-89-0 |
| D10 | 2.00 M Sodium nitrate | Chaotropic Anion | 7631-99-4 |
| D11 | 1.00 M Sodium iodide | Chaotropic Anion | 7681-82-5 |
| D12 | 2.00 M Sodium bromide | Chaotropic Anion | 7647-15-6 |

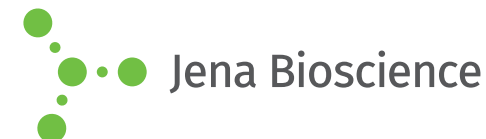




JBScreen Plus HTS

Cat.-No.: CS-507L

SCREEN FORMULATION

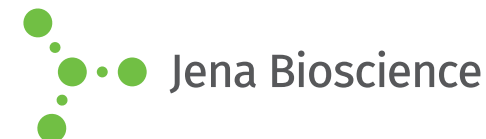


| No. | Additive | Type | CAS Number |
|-----|----------------------------------|--------------------|-------------|
| E1 | 1.00 M tri-Lithium citrate | Lithium Salt | 313222-91-2 |
| E2 | 2.00 M Lithium acetate | Lithium Salt | 6108-17-4 |
| E3 | 2.00 M Lithium bromide | Lithium Salt | 85017-82-9 |
| E4 | 2.00 M Lithium salicylate | Lithium Salt | 552-38-5 |
| E5 | 2.00 M Lithium nitrate | Lithium Salt | 7790-69-4 |
| E6 | 2.00 M Lithium perchlorate | Lithium Salt | 7791-03-9 |
| E7 | 2.00 M Ammonium fluoride | Ammonium Salt | 12125-01-8 |
| E8 | 2.00 M Ammonium chloride | Ammonium Salt | 12125-02-9 |
| E9 | 2.00 M Ammonium bromide | Ammonium Salt | 12124-97-9 |
| E10 | 2.00 M Ammonium nitrate | Ammonium Salt | 6484-52-2 |
| E11 | 2.00 M Ammonium thiocyanate | Ammonium Salt | 1762-95-4 |
| E12 | 2.00 M Ammonium trifluoroacetate | Ammonium Salt | 3336-58-1 |
| F1 | 2.00 M di-Ammonium tartrate | Ammonium Salt | 3164-29-2 |
| F2 | 600 mM Potassium sulfate | Sulfate | 7778-80-5 |
| F3 | 2.00 M Sodium sulfate | Sulfate | 7727-73-3 |
| F4 | 2.00 M Lithium sulfate | Sulfate | 10377-48-7 |
| F5 | 2.00 M Magnesium sulfate | Sulfate | 10034-99-8 |
| F6 | 1.00 M Beryllium sulfate | Sulfate | 13510-49-1 |
| F7 | 100 mM Barium chloride | Multivalent Cation | 10326-27-9 |
| F8 | 100 mM Strontium chloride | Multivalent Cation | 10025-70-4 |
| F9 | 100 mM Cadmium chloride | Multivalent Cation | 10108-64-2 |
| F10 | 100 mM Cobalt (II) chloride | Multivalent Cation | 7791-13-1 |
| F11 | 100 mM Copper (II) chloride | Multivalent Cation | 10125-13-0 |
| F12 | 100 mM Yttrium (III) chloride | Multivalent Cation | 10025-94-2 |



**JBScreen Plus HTS**

Cat.-No.: CS-507L

SCREEN FORMULATION

| No. | Additive | Type | CAS Number |
|-----|--|-----------------------------|---------------------|
| G1 | 2.00 M 6-Aminohexanoic acid | Linker Molecule | 60-32-2 |
| G2 | 2.00 M Ethanolamine | Linker Molecule | 141-43-5 |
| G3 | 2.00 M 1,6-Diaminohexane | Linker Molecule | 124-09-4 |
| G4 | 2.00 M 1,8-Diaminooctane | Linker Molecule | 373-44-4 |
| G5 | 1.00 M 1,2,3-hexanetriol | Polyalcohol | 25323-24-4 |
| G6 | 100 mM Spermidine trihydrochloride | Polyamine | 334-50-9 |
| G7 | 500 mM Polypropylene glycol 400 | Organic Hydrophilic Polymer | 25322-69-4 |
| G8 | 1.00 M Polyethylene glycol 200 | Organic Hydrophilic Polymer | 25322-68-3 |
| G9 | 500 mM Polyethylene glycol 600 | Organic Hydrophilic Polymer | 25322-68-3 |
| G10 | 10.00 % w/v Dextran sulfate sodium salt | Organic Hydrophilic Polymer | 9011-18-1 |
| G11 | 1.00 % w/v Polyvinyl alcohol Type II | Organic Hydrophilic Polymer | 9002-89-5 |
| G12 | 5.00 % w/v Polyvinylpyrrolidone K15 | Organic Hydrophilic Polymer | 9003-39-8 |
| H1 | 1.00 M D-(-)-Fructose | Carbohydrate | 57-48-7 |
| H2 | 1.00 M D-(+)-Glucose | Carbohydrate | 14431-43-7 |
| H3 | 1.00 M D-(+)-Mannose | Carbohydrate | 3458-28-4 |
| H4 | 500 mM Lactose | Carbohydrate | 63-42-3 |
| H5 | 1.00 M D-(+)-Sucrose | Carbohydrate | 57-50-1 |
| H6 | 1.00 M D-(+)-Trehalose | Carbohydrate | 6138-23-4 |
| H7 | 100 mM Tris(2-carboxyethyl)phosphine hydrochloride | Reducing Agent | 51805-45-9 |
| H8 | 10.00 mM L-Glutathione reduced, L-Glutathione oxidized | Reducing Agent | 70-18-8, 27025-41-8 |
| H9 | 100 mM Ethylenediaminetetraacetic acid disodium salt | Chelator | 6381-92-6 |
| H10 | 100 mM Adenosine-5'-triphosphate disodium salt | Co-Factor | 987-65-5 |
| H11 | 1.00 M Benzamidine hydrochloride | Amphiphilic Molecule | 206752-36-5 |
| H12 | 300 mM Glycyl-glycyl-glycine | Amphiphilic Molecule | 556-33-2 |

