



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





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



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	20.00 mM Dextran sulfate sodium salt 5,000	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 mM L-Glutathione reduced, 10 mM 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