



JBScreen Buffers Xtreme

Cat. No.	Amount
CS-215	84 solutions (1,7 ml each)

	1	2	3	4	5	6	7	8	9	10	11	12
A	empty	empty	empty	empty	empty	empty	empty	empty	empty	empty	empty	empty
B	Glycine, pH 3.0	Glycylglycine, pH 3.8	Citrate, pH 3.2	Succinate, pH 3.5	Acetate, pH 4.5	MIB, pH 4.5	BIS-TRIS Propane, pH 8.5	Bicine, pH 8.5	Taurine, pH 8.5	Glycine, pH 10.0	CAPS, pH 9.5	MIB, pH 9.0
C	Formate, pH 3.0	Malate, pH 3.0	Citrate, pH 3.5	Succinate, pH 4.0	Acetate, pH 5.0	MIB, pH 5.0	BIS-TRIS Propane, pH 9.0	Bicine, pH 9.0	Taurine, pH 9.0	Glycine, pH 10.5	CAPS, pH 10.0	MIB, pH 9.5
D	Formate, pH 3.5	Malate, pH 3.5	Citrate, pH 4.0	Succinate, pH 4.5	Acetate, pH 5.5	MIB, pH 5.5	BIS-TRIS Propane, pH 9.5	AMPP, pH 8.5	Taurine, pH 9.5	CHES, pH 8.6	CAPS, pH 9.7	MIB, pH 10.0
E	Formate, pH 4.0	Malate, pH 4.0	Citrate, pH 4.5	Succinate, pH 5.0	Malonate, pH 5.0	SPG, pH 4.0	Glycylglycine, pH 8.9	AMPP, pH 9.0	Troine, pH 8.5	CHES, pH 9.0	CAPS, pH 10.0	SPG, pH 9.5
F	Formate, pH 4.5	Malate, pH 4.5	Citrate, pH 5.0	Succinate, pH 5.5	Malonate, pH 5.5	SPG, pH 4.5	Glycylglycine, pH 8.9	AMPP, pH 9.5	Glycine, pH 8.6	CHES, pH 9.5	CAPS, pH 10.5	SPG, pH 9.9
G	Glycylglycine, pH 5.0	Malate, pH 5.0	Citrate, pH 5.5	Acetate, pH 5.6	MES, pH 5.5	SPG, pH 5.0	TRIS, pH 8.5	TAPS, pH 8.5	Glycine, pH 9.0	CHES, pH 11.0	CAPS, pH 11.0	SPG, pH 9.5
H	Glycylglycine, pH 3.5	Malate, pH 5.5	Succinate, pH 3.2	Acetate, pH 4.0	MIB, pH 4.0	SPG, pH 5.5	TRIS, pH 9.0	TAPS, pH 9.0	Glycine, pH 9.5	CAPS, pH 9.0	MIB, pH 8.5	SPG, pH 10.0
	acidic buffers, pH 3.0 - 5.5						basic buffers, pH 8.5 - 11.0					

Description:

JBScreen Buffers Xtreme contain common buffer stocks (0.5 M concentration) in the extreme pH ranges (3.0 - 5.5 and 8.5 - 11.0) in useful 0.5 pH unit increments.

MIB buffer is produced by mixing Malonic acid:Imidazole:Boric acid in the molar ratios 2:3:3. SPG buffer is produced by mixing Succinic acid:Sodium dihydrogen phosphate:Glycine in the molar ratios 2:7:7. These 'Super Buffers' screen the pH range from 4.0 to 10.0 without changing the chemical composition of the buffer solution.

Related Products:

JBScreen Buffers, #CS-214

JBScreen Thermofluor FUNDAMENT HTS, #CS-332

JBScreen Thermofluor SPECIFIC HTS, #CS-333

Selected References:

Newman (2004) Novel buffer systems for macromolecular crystallization. *Acta Cryst. D* **60**:610.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months