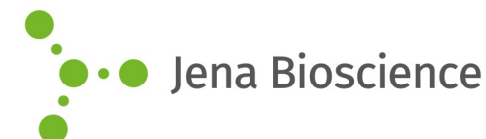


JBScreen Buffers

DATA SHEET



Cat. No.	Amount
CS-214	84 solutions (1.7 ml each)

For *in vitro* use only
Quality guaranteed for 12 months
Store at 8 - 10°C

Buffer concentration

0.5 M

Description

JBScreen Buffers contain common buffer stocks (0.5 M concentration) in the neutral pH range (5.5 - 8.5) in useful 0.5 pH unit increments.

SPG buffer (column 12) is produced by mixing Succinic acid:Sodium dihydrogen phosphate:Glycine in the molar ratios 2:7:7. This 'Super Buffer' screens the pH range from 4.0 to 10.0 without changing the chemical composition of the buffer solution.

Related Products

JBScreen Buffers Xtreme, #CS-215
JBScreen Thermofluor
FUNDAMENT, # CS-332
JBScreen Thermofluor
SPECIFIC, # CS-333

Selected Reference

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

	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	Citrate, pH 5.5	Malonate, pH 5.5	MES, pH 6.0	BIS-TRIS, pH 6.5	Phosphate, pH 7.0	Imidazole, pH 6.2	BIS-TRIS Propane, pH 7.0	MOPS, pH 8.0	TAPSO, pH 7.0	Tricine, pH 7.5	Bicine, pH 8.0	SPG, pH 5.5
C	Citrate, pH 6.0	Malonate, pH 6.0	MES, pH 6.5	BIS-TRIS, pH 7.0	Phosphate, pH 7.5	Imidazole, pH 6.5	BIS-TRIS Propane, pH 7.5	HEPES, pH 7.0	TAPSO, pH 7.5	Tricine, pH 8.0	Bicine, pH 8.5	SPG, pH 6.0
D	Citrate, pH 6.5	Malonate, pH 6.5	MES, pH 6.7	ADA, pH 6.0	Phosphate, pH 8.0	Imidazole, pH 7.0	BIS-TRIS Propane, pH 8.0	HEPES, pH 7.5	TAPSO, pH 8.0	Tricine, pH 8.5	TAPS, pH 7.7	SPG, pH 6.5
E	Citrate, pH 7.0	Malonate, pH 7.0	Maleate, pH 6.0	ADA, pH 6.5	PIPES, pH 6.1	Imidazole, pH 7.5	BIS-TRIS Propane, pH 8.5	HEPES, pH 8.0	TRIS, pH 7.1	Glycylglycine, pH 7.5	TAPS, pH 8.0	SPG, pH 7.0
F	Succinate, pH 5.5	Malonate, pH 7.5	Maleate, pH 6.5	ADA, pH 7.0	PIPES, pH 6.5	Imidazole, pH 7.8	MOPS, pH 6.5	DIPSO, pH 7.0	TRIS, pH 7.5	Glycylglycine, pH 8.0	TAPS, pH 8.5	SPG, pH 7.5
G	Succinate, pH 6.0	Malonate, pH 8.0	Maleate, pH 7.0	Phosphate, pH 6.0	PIPES, pH 7.0	BIS-TRIS Propane, pH 6.3	MOPS, pH 7.0	DIPSO, pH 7.5	TRIS, pH 8.0	Glycylglycine, pH 8.5	AMPD, pH 8.0	SPG, pH 8.0
H	Succinate, pH 6.5	MES, pH 5.5	BIS-TRIS, pH 6.0	Phosphate, pH 6.5	PIPES, pH 7.5	BIS-TRIS Propane, pH 6.5	MOPS, pH 7.5	DIPSO, pH 8.0	TRIS, pH 8.5	Bicine, pH 7.6	AMPD, pH 8.5	SPG, pH 8.5

neutral buffers, pH 5.5 - 8.5

