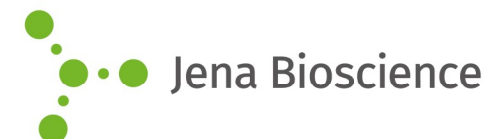


JBScreen Buffers Xtreme

DATA SHEET



Cat. No.	Amount
CS-215	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 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, # 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	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	CAPSO, 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	CAPSO, 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	AMPD, 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 8.5	AMPD, pH 9.0	Tricine, pH 8.5	CHES, pH 9.0	CAPS, pH 10.0	SPG, pH 8.5
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	AMPD, pH 9.5	Glycine, pH 8.6	CHES, pH 9.5	CAPS, pH 10.5	SPG, pH 9.0
G	Glycylglycine, pH 3.0	Malate, pH 5.0	Citrate, pH 5.5	Acetate, pH 3.6	MES, pH 5.5	SPG, pH 5.0	TRIS, pH 8.5	TAPS, pH 8.5	Glycine, pH 9.0	CHES, pH 10.0	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	CAPSO, 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



IFTA AG
Certified QMS and EMS according to
DIN EN ISO 9001 and DIN EN ISO 14001
Reg.-No.: ICV03597 034 and ICV03597 534

Jena Bioscience GmbH | Loebstedter Str. 71 | 07749 Jena, Germany
Phone +49-3641-6285 000 | Fax +49-3641-6285 100
www.jenabioscience.com

Last update: Mar 03, 2021