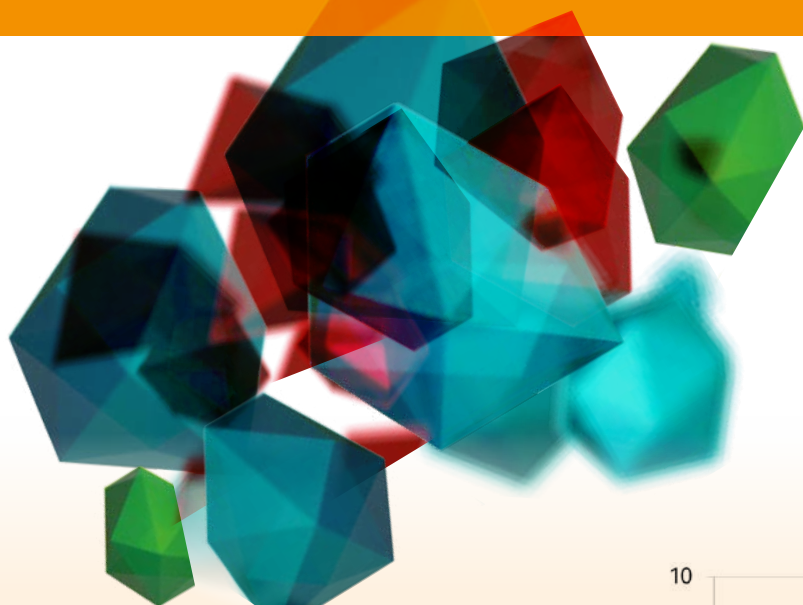




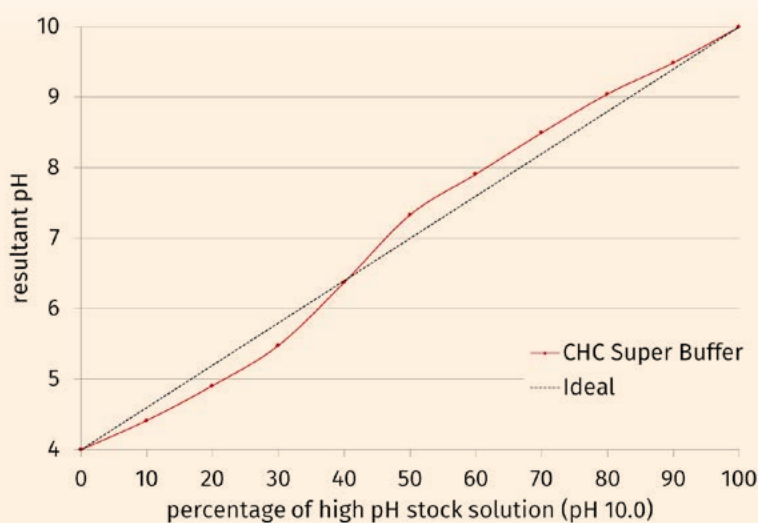
Buffer	pH											
	1	2	3	4	5	6	7	8	9	10	11	
ADA - 1 M CSS-501						pH 6.0 – 7.2						
Bicine - 1 M CSS-502								pH 7.6 – 9.5				
BIS-TRIS - 1 M CSS-503						pH 5.8 – 7.2						
BIS-TRIS Propane - 1 M CSS-504							pH 6.3 – 9.5					
CAPS - 1 M CSS-505										pH 9.7 – 11.1		
CHES - 1 M CSS-506									pH 8.6 – 10.0			
Citrate/Phosphate - 1 M CSS-507			pH 2.6 – 7.0									
Citric Acid - 1 M CSS-508		pH 2.2 – 6.5										
DL-Malic Acid - 1 M CSS-509			pH 2.7 – 6.0									
Glycine - 1 M CSS-510									pH 8.6 – 10.6			
HEPES - 1 M CSS-511							pH 6.8 – 8.2					
Imidazole - 1 M CSS-512							pH 6.2 – 7.8					
Lithium Acetate - 1 M CSS-513								pH 7.5 – 9.5				
MES - 1 M CSS-514						pH 5.5 – 6.7						
MOPS - 1 M CSS-515							pH 6.5 – 7.9					
PIPES - 1 M CSS-516							pH 6.5 – 7.5					
Potassium Phosphate - 1 M CSS-517							pH 5.8 – 8.0					
Sodium Acetate (HCl) - 1 M CSS-518				pH 3.6 – 5.6								
Sodium Acetate - 1 M CSS-519				pH 3.6 – 5.6								
Sodium Phosphate - 1 M CSS-520							pH 6.0 – 7.2					
Sodium Potassium Phosphate - 1 M CSS-521							pH 5.7 – 8.0					
Succinic Acid - 0.5 M CSS-522			pH 3.2 – 6.6									
Tricine - 1 M CSS-523								pH 7.4 – 8.8				
TRIS (TRIS-Acetate) - 1 M CSS-524								pH 7.0 – 9.0				
TRIS - 1 M CSS-525								pH 7.0 – 9.0				
Sodium Citrate - 1 M CSS-526					pH 5.0 – 6.2							



Screen the pH independently from any other parameter.

Super Buffers are composed of a mixture of three individual buffers with distinct pK_a values and cover a broad pH range without changing the chemical environment of the buffer solution.

They are supplied as low and high pH stock solutions, which can be mixed at different ratios to obtain different pH values within the range. Plotting the pH vs. the percentage of high pH stock solution in the mixture results in an almost linear pH function for any JBScreen Super Buffer system.



Super Buffer	Cat. No.	Description
AAB pH 4.0 - 1 M	CSS-404	AAB Buffer System: Sodium Acetate, ADA and Bicine (1:1:1 Molar Ratio)
AAB pH 9.0 - 1 M	CSS-405	
CHC pH 4.0 - 1 M	CSS-402	CHC Buffer System: Citric Acid, HEPES and CHES (2:3:4 Molar Ratio)
CHC pH 10.0 - 1 M	CSS-403	
MIB pH 4.0 - 1 M	CSS-400	MIB Buffer System: Malonic Acid, Imidazole and Boric Acid (2:3:3 Molar Ratio)
MIB pH 10.0 - 1 M	CSS-401	
MMT pH 4.0 - 1 M	CSS-398	MMT Buffer System: Malic Acid, MES and TRIS (1:2:2 Molar Ratio)
MMT pH 9.0 - 1 M	CSS-399	
PCB pH 4.0 - 1 M	CSS-387	PCB Buffer System: Sodium Propionate, Sodium Cacodylate and BIS-TRIS Propane (2:1:2 Molar Ratio)
PCB pH 9.0 - 1 M	CSS-388	
SPG pH 4.0 - 1 M	CSS-389	SPG Buffer System: Succinic Acid, Sodium Dihydrogen Phosphate and Glycine (2:7:7 Molar Ratio)
SPG pH 10.0 - 1 M	CSS-390	
TBG pH 4.0 - 1 M	CSS-384	TBG Buffer System: Sodium Tartrate Dihydrate, BIS-TRIS and Glycylglycine (3:2:2 Molar Ratio)
TBG pH 9.0 - 1 M	CSS-385	
TACS MMF pH 7.0 - 100 %	CSS-420	TACS MMF (aka Tacsimate) is a buffer system containing 1.8305 M Malonic acid, 0.25 M Ammonium citrate tribasic, 0.12 M Succinic acid, 0.3 M DL-Malic acid, 0.4 M Sodium acetate trihydrate, 0.5 M Sodium formate, and 0.16 M Ammonium tartrate dibasic

