



pHat™ Buffer Screen - technical sheet

pH of each formulation shown for each of the 96 well block plate positions

	1M Sodium Lactate / HCl	1M Sodium Acetate / Acetic Acid	1M MES / NaOH	1M Bis-Tris / HCl	1M Imidazole / HCl	1M K ₂ HPO ₄ / NaH ₂ PO ₄	1M HEPES / NaOH	1M Tris / HCl	1M Bicine / NaOH	0.5M CHES / NaOH	1M Glycine / NaOH	0.5M CAPS / NaOH
A1	2.4	3.4	4.6	5.2	5.4	5.8	6.0	6.6	7.0	8.0	8.2	8.8
B1	2.8	3.8	5.0	5.6	5.8	6.2	6.4	7.0	7.4	8.4	8.6	9.2
C1	3.2	4.2	5.4	6.0	6.2	6.6	6.8	7.4	7.8	8.8	9.0	9.6
D1	3.6	4.6	5.8	6.4	6.6	7.0	7.2	7.8	8.2	9.2	9.4	10.0
E1	4.0	5.0	6.2	6.8	7.0	7.4	7.6	8.2	8.6	9.6	9.8	10.4
F1	4.4	5.4	6.6	7.2	7.4	7.8	8.0	8.6	9.0	10.0	10.2	10.8
G1	4.8	5.8	7.0	7.6	7.8	8.2	8.4	9.0	9.4	10.4	10.6	11.2
H1	5.2	6.2	7.4	8.0	8.2	8.6	8.8	9.4	9.8	10.8	11.0	11.6

All formulations are made with ultrapure ASTM Type I water and sterile-filtered stock solutions. Store at 4-25 °C.

© 2004 Emerald BioSystems, Inc., (888) 780-8535, FAX (206) 780-8549, www.emeraldbiosystems.com