

EPIgeneous™ Cell Based Assay Lysis Buffer Guide

Histone Modification Assay	Buffers	Protocols			
	What is included in the kit?	One Plate With Medium (Suspension and Adherent Cells)	One Plate w/o Medium (Adherent cells Only)	Two Plates With Medium (Suspension and Adherent Cells)	Two Plates w/o Medium (Adherent cells Only)
H3K4Me2	Lysis Buffer C part 1 & 2 (4x)	Supplemented Lysis Buffer C (4x)	Supplemented Lysis Buffer C (1x)	Supplemented Lysis Buffer C (4X)	Supplemented Lysis Buffer C (1x)
H3K27Me3	Lysis Buffer A part 1 & 2 (4x)	Lysis Buffer A (4x)	Lysis Buffer A (1x)	Lysis Buffer A (4x)	Lysis Buffer A (1x)
H3K36Me2	Lysis Buffer B (4x) & C part 1 & 2 (4x)	Lysis Buffer B (2.5x)	Supplemented Lysis Buffer C (1x)	Lysis Buffer B (4x)	Supplemented Lysis Buffer C (1x)
H3 Total ▪ used in combination with H3K4Me2, H3K27Me3 or H3K36Me2 for normalization studies	No Buffer included No need to purchase additional buffer	No additional buffer needed Use the lysate from H3K4Me2, H3K27Me3 or H3K36Me2	No additional buffer needed Use the lysate from H3K4Me2, H3K27Me3 or H3K36Me2	No additional buffer needed Use the lysate from H3K4Me2, H3K27Me3 or H3K36Me2	No additional buffer needed Use the lysate from H3K4Me2, H3K27Me3 or H3K36Me2
H3 Total ▪ used “alone”	No Buffer included Need to purchase separately	Lysis Buffer A (4x)	Lysis Buffer A (1x)	Lysis Buffer A (4x)	Lysis Buffer A (1x)