

Essen CellPlayer™ Invasion Assay : Invasion or Migration vs. Time for 3 Different Cell Types

