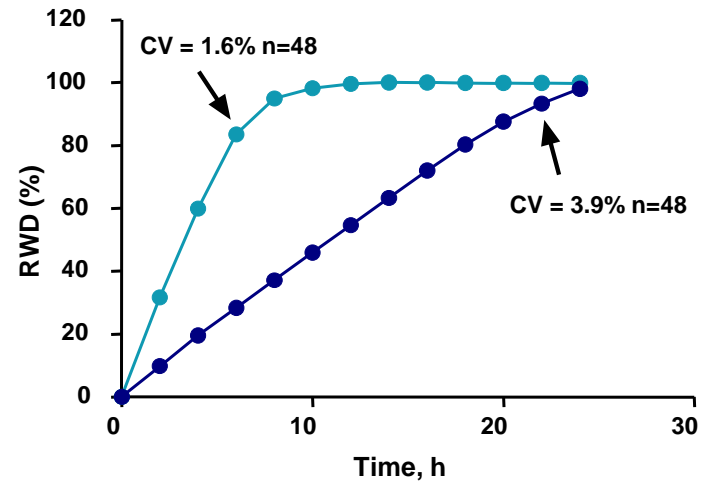


- HT 1080 2D migration on collagen surface
- HT 1080 3D invasion through collagen gel



Summary of statistics from plate on left

Data points represent mean \pm SE, N=48 per condition