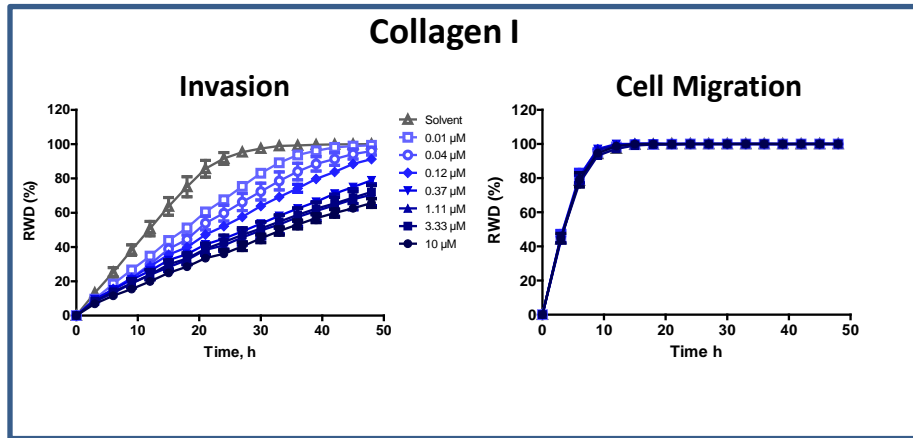
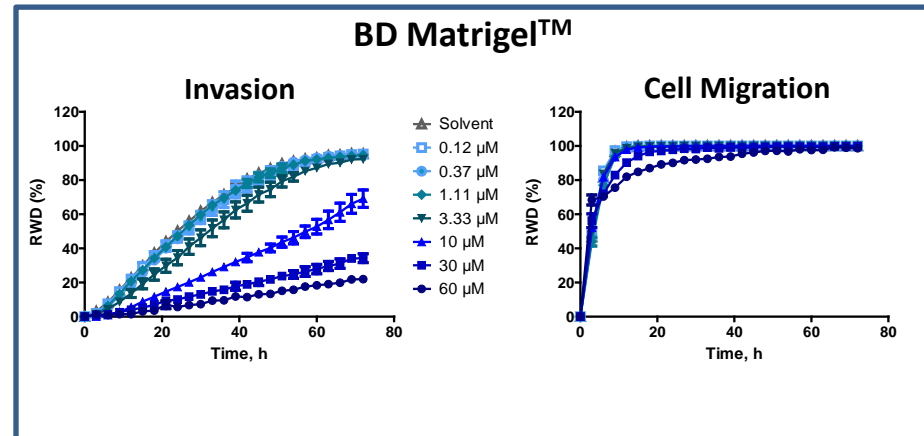
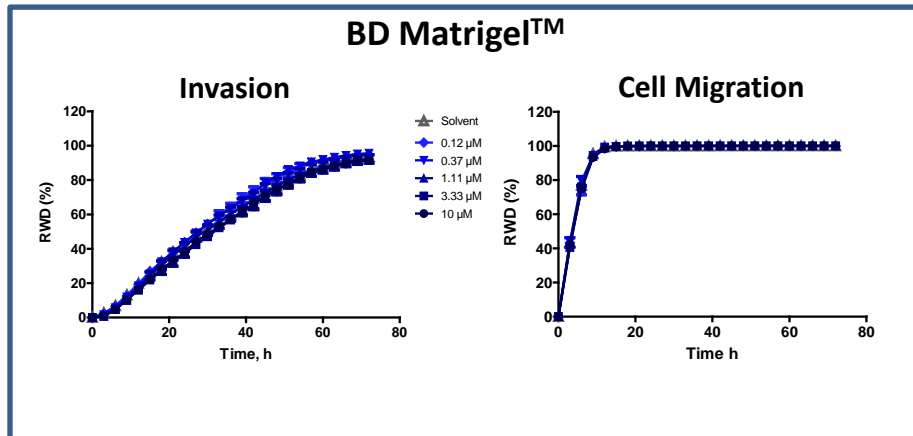
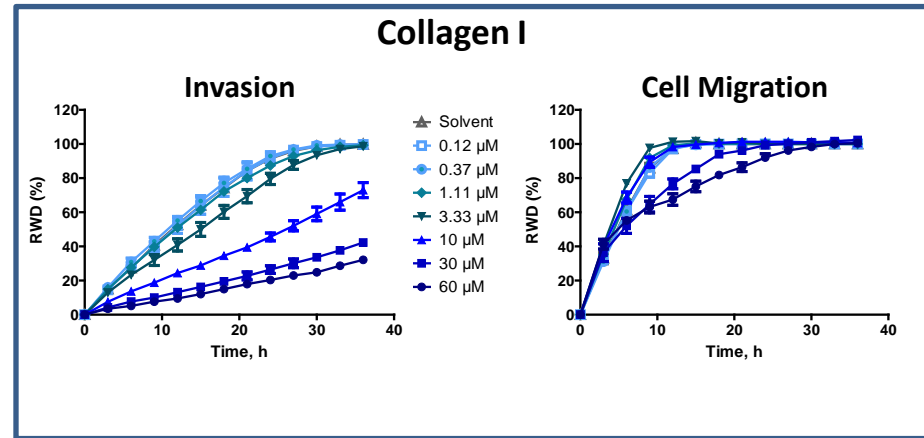


Pharmacological comparison of blebbistatin and GM6001 effect on the invasive phenotype of HT 1080 cells using the Essen BioScience CellPlayer Invasion/Migration Assay

Pharmacology of GM6001: A broad MMP inhibitor



Pharmacology of blebbistatin: A myosin II inhibitor



Conclusion: GM6001 inhibits invasion through Collagen I *but not* Matrigel™

Conclusion: Blebbistatin inhibits invasion through both Collagen I *and* Matrigel™