

Migration and Invasion Assay

Cat #: MIG_HT1080; INV_HT1080

Cells: Human Fibrosarcoma Cancer Epithelial cells, HT-1080

Description:

HT-1080 Cells invade through Matrigel-coated membranes (Invasion) or uncoated membrane (Migration) towards 10% Fetal Bovine Serum in the presence of Doxycycline.

Reference Compound IC50:

Assay	Doxycycline, IC50 (μM)
Migration	29
Invasion	32

