

KINEXUS IN VIVO SERVICES

CHOOSE FROM TWELVE OF THE MOST COMMONLY STUDIED HUMAN TUMOUR CELL LINES. SEND US YOUR DRUG, RNA^{si} OR HORMONE, AND WE WILL PERFORM THE INCUBATIONS AND PREPARE CELL LYSATES FOR KINEX™ ANTIBODY MICROARRAY AND KINETWORKS™ IMMUNOBLOTTING ANALYSIS.

As part of our continuing commitment to ensuring that our clients are able to fully benefit from our proteomics services, we are pleased to offer the convenience and cost savings of our cell incubation and lysate preparation services. We will perform the experimental manipulation of the cells to your specifications, and produce sufficient protein for one Kinex™ and one Kinetworks™ analysis. Our standard charge is just US\$400 for the first lysate and US\$250 for each subsequent lysate from the same cell type.

Cell line: **293**
Origin: Female fetal kidney cells transformed with adenovirus 5



Cell line: **A431**
Origin: Skin epidermoid carcinoma from 85 year old female



Cell line: **A549**
Origin: Lung carcinoma from 58 year old male



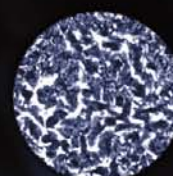
Cell line: **HCT116**
Origin: Colon carcinoma from adult male



Cell line: **HeLa**
Origin: Cervix epithelial adenocarcinoma from 31 year old female



Cell line: **HepG2**
Origin: Liver carcinoma from 15 year old male



Cell line: **HL-60**
Origin: Peripheral blood promyeloblasts from 36 year old female



Cell line: **HUV-EC**
Origin: Umbilical vein endothelial cells from normal adult female



Cell line: **Jurkat**
Origin: T cell leukemia from 14 year old male



Cell line: **MCF-7**
Origin: Breast epithelial adenocarcinoma from 69 year old female



Cell line: **PC-3**
Origin: Prostate adenocarcinoma from bone of 62 year old male



Cell line: **T98G**
Origin: Brain glioblastoma from 61 year old male



Toll free: 1 866 KINEXUS

www.kinexus.ca