

CUSTOM KINASE SUBSTRATE PROFILING (CKSP) SERVICE

List of Available Human Tumour Cell Lysates

Cell Line	Ref. Code	Organ/Tissue	Name	Gender	Comments
A431	CL01	Skin	Human	Female	Skin epidermoid carcinoma from an 85 year old female [ATCC# CRL-1555]
A549	CL02	Lung	Human	Male	Lung carcinoma from a 58 year old male [ATCC# CCL-185]
Daudi	CL03	Blood B Cell	Human	Male	Burkitt's B cell lymphoma from a 16 year old male [ATCC# CCL-213]
HCT116	CL04	Colon	Human	Male	Colon carcinoma from an adult male [ATCC# CCL-247]
HEK 293	CL05	Kidney	Human	Female	Female fetal kidney cells transformed with adenovirus 5 [ATCC# CRL-1573]
HeLa	CL06	Cervix	Human	Female	Cervix epithelial adenocarcinoma from 31 year old female [ATCC# CCL-2]
HL-60	CL07	Blood	Human	Female	Peripheral blood promyeloblasts from a 36 year old female [ATCC# CCL-240]
HepG2	CL08	Liver	Human	Male	Liver carcinoma from a 15 year old male [ATCC# HB-8065]
HU-VEC	CL09	Umbilical Cord	Human	Female	Umbilical vein endothelial cells from a normal adult female [ATCC# CRL-1730]
Jurkat	CL10	Blood T Cell	Human	Male	T cell leukemia from a 14 year old male [ATCC# TIB-152]
MCF7	CL11	Breast	Human	Female	Breast epithelial adenocarcinoma from a 69 year old female [ATCC# HTB-22]
PC3	CL12	Prostate	Human	Male	Prostate adenocarcinoma from bone of 62 year old male [ATCC# CRL-1435]
T98G	CL13	Brain	Human	Male	Brain glioblastoma from a 61 year old male [ATCC# CRL-1690]
THP1	CL14	Blood Monocyte	Human	Male	Acute monocytic leukemia from peripheral blood of a 1 year old male [ATCC# CRL-1435]



KINEXUS

Suite 1, 8755 Ash Street
Vancouver, B.C
Canada V6P 6T3

Phone: 1-866-KINEXUS
Phone: 1-604-323-2547
Facsimile: 1-604-323-2548
E-Mail: info@kinexus.ca
Internet: www.kinexus.ca