



NucView® Validated Cell Lines

[Click on a reference to link to the publication](#)

Immortalized cell lines used with NucView® 405

Cell line	Species	Cell type	References
DLD1	Human	Colon cancer	Grabocka et al. 2016
HeLa	Human	Cervical cancer	Grabocka et al. 2016
NCI-H508	Human	Cecum adenocarcinoma	Grabocka et al. 2016
RH4	Human	Rabdomyosarcoma	Bohm et al. 2016

Immortalized cell lines used with NucView® 530

Cell line	Species	Cell type	References
HCT116	Human	Colorectal carcinoma	Manas et al. 2017
MDA-MB-231	Human	Breast adenocarcinoma	Bueno et al. 2018

Immortalized cell lines used with NucView® 488

Cell line	Species	Cell type	References
293-H	Human	Embryonic kidney	Wu and MacRae 2010
293-T	Human	Embryonic kidney	Ben Salem et al. 2015 ; Boussabbeh et al. 2015 ; Boussabbeh et al. 2015 (2) ; Li et al. 2008 ; Wetzel-Smith et al. 2014
4T1	Mouse	Mammary tumor	Granot et al. 2011
5637	Human	Bladder carcinoma	Brassesco et al. 2013 (A) ; Tang et al. 2013
67NR	Mouse	Mammary carcinoma	Cheng et al. 2009
A172	Human	Glioma	Jeong et al. 2009
A204	Human	Sarcoma	Zhu et al. 2008
A375	Human	Melanoma	Feng et al. 2016 ; Takata et al. 2017 ; Tyciakova et al. 2015
A549	Human	Adenocarcinoma	Yamada et al. 2016
BeWo	Human	Trophoblast	Angeloni et al. 2009
BT549	Human	Breast cancer	Lim et al. 2016
Caco-2	Human	Colorectal	Yoshioka et al. 2014
CCL-134	Human	IPF pulmonary fibroblast	Vuorinen et al. 2007
CCL-190	Human	Pulmonary fibroblast	Vuorinen et al. 2007
CCL-227	Human	Colon	Arsic et al. 2017
CCL-228	Human	Colon	Arsic et al. 2017
CCRF-CEM	Human	Leukemia	Dahlawi et al. 2013 , Pedroza et al. 2014
FU-UR-1	Human	Renal cell carcinoma	Kobos et al. 2013
G-361	Human	Melanoma	Aronchik et al. 2014 (2)
GE11	Mouse	Epithelial	Puigvert et al. 2009
H358	Human	Non-small cell lung cancer	Aronchik et al. 2014
H9c2	Rat	Cardiac myoblast	Ben Salem et al. 2017 ; He et al. 2014 ; Prola et al. 2016 ; Ren et al. 2012
HaCaT	Human	Keratinocyte	Paromov et al. 2011

HaEpi	Insect	Epidermal	Chen et al. 2017 ; Zhao et al. 2016
HCE	Human	Corneal epithelial	Soriano-Romani et al. 2015
HCLE	Human	Corneal epithelial	Mankus et al. 2011
HCT116	Human	Colorectal carcinoma	Aftab et al. 2014 ; Ben Salem et al. 2015 ; Boussabbeh et al. 2015 ; Boussabbeh et al. 2015 (2) ; Johnston et al. 2016 ; Manas et al. 2017
HeLa	Human	Cervical cancer	Angeloni et al. 2009 ; Antczak et al. 2009 ; Balzarini et al. 2013 ; Cen et al. 2008 ; Chiu et al. 2016 ; Gallerne et al. 2013 ; Gao et al. 2014 ; Kasim et al. 2013 ; Moore et al. 2010 ; Sirianant et al. 2016 ; Verma et al. 2013
HepG2	Human	Liver carcinoma	Chiu et al. 2016 ; Liu et al. 2016
HepT1	Human	Hepatoblastoma	Dewerth et al. 2012 ; Eicher et al. 2011
HMEC	Human	Microvascular endothelial	Isherwood et al. 2011
HL-60	Human	Leukemia	Contreras et al. 2018 ; Dahlawi et al. 2013
HOS	Human	Osteosarcoma	Brassesso et al. 2013 ; Takata et al. 2017
HT-29	Human	Colorectal adenocarcinoma	Tyciakova et al. 2015
HT-1080	Human	Breast fibrosarcoma	Wang et al. 2010 ; Zhu et al. 2008
HUH6	Human	Hepatoblastoma	Dewerth et al. 2012 ; Eicher et al. 2011
IMR-90	Human	Lung	Baar et al. 2017
INS-IE	Mouse	Pacreatic islet	Shulga et al. 2015
IOBA-NHC	Human	Conjunctival epithelial	Soriano-Romani et al. 2015
Jurkat	Human	T-lymphocyte	Cen et al. 2008 ; Li and Shively 2013 ; Oldham et al. 2015 ; Stewart et al. 2010
JY	Human	Lymphoblastic leukemia	Berenyi et al. 2011
K562	Human	Myelogenous leukaemia	Sztiller-Sikorska et al. 2009
MCF-7	Human	Breast adenocarcinoma	Balijepalli et al. 2010 ; Hamzeloo-Moghadam et al. 2015 ; Naghibi et al. 2014 ; Renjini et al. 2014
MCF-10A	Human	Breast adenocarcinoma	Hafner et al. 2016 ; Koerner et al. 2013
MDA-MB-231	Human	Breast adenocarcinoma	Aronchik et al. 2011 ; Aronchik et al. 2014 ; Balijepalli et al. 2010 ; Bueno et al. 2018 ; Koerner et al. 2013 ; Palorini et al. 2016 ; Robles-Escajeda et al. 2016 ; Salem et al. 2016 ; Tyciakova et al. 2015
MDA-MB-361	Human	Breast cancer	Gao et al. 2017
MDA-MB-468	Human	Breast cancer	Jangmareddy et al. 2013 (NucView & MitoView Kit) ; Salem et al. 2016
MDCK	Canine	Kidney epithelial	Eisenhoffer et al. 2012 ; Saw et al. 2017
MES-SA	Human	Uterine sarcoma	Zhu et al. 2008
MES-SA/DX	Human	Uterine sarcoma	Zhu et al. 2008
MG-63	Human	Osteosarcoma	Brassesso et al. 2013
Min 6	Mouse	Pancreatic insulinoma	Hofmeister-Brix et al. 2013 ; Pathak et al. 2008
MKN-45	Human	Gastric cancer	Liu et al. 2016
MOLT-3	Human	Leukemia	Dahlawi et al. 2013
MV4-11	Human	Macrophage	Eriksson et al. 2012
Nalm-6	Human	B cell	Robles-Escajeda et al. 2013

N19	Mouse	Oligodendrocyte	Paez et al. 2007 ; Smith et al. 2012
NIH 3T3	Mouse	Fibroblast	Liu et al. 2016
NRK	Rat	Kidney epithelial	Stopper et al. 2009
NRK-52E	Rat	Kidney epithelial	Katsoulieris et al. 2010
OECM-1	Human	Oral cancer	Su et al. 2017
OVK18	Human	Ovarian cancer	De et al. 2018
PC-3	Human	Prostate cancer	Balzarini et al. 2013 ; Bjork et al. 2015 (3D culture) ; Bjorkman et al. 2011 ; Lee et al. 2013 ; Virtanen et al. 2014
PC-9	Human	Non-small cell lung cancer	Dereli-Korkut et al. 2014
PC12	Rat	Pheochromocytoma	Stopper et al. 2009
Ramos	Human	B cell lymphoma	Contreras et al. 2018
R28	Rat	Retinal precursors	Uddin et al. 2015
RD	Human	Rhabdomyosarcoma	Zhu et al. 2008
RINm5F	Rat	Insulinoma	Schmitt et al. 2011
RKO	Human	Colon carcinoma	Cai et al. 2013
RT4	Human	Bladder carcinoma	Brascesco et al. 2013 (A)
RT112	Human	Bladder carcinoma	King et al. 2016
SAS	Human	Oral carcinoma	Su et al. 2017
Saos-2	Human	Osteosarcoma	Lanz et al. 2013
SKBR3	Human	Breast cancer	Jangamreddy et al. 2013 (NucView & MitoView Kit) ; Tyciakova et al. 2015
SKLMS1	Human	Leiomyosarcoma	Wang et al. 2010 ; Zhu et al. 2008
SK-OV-3	Human	Ovarian carcinoma	Tyciakova et al. 2015
SMMC-7721	Human	Hepatocarcinoma	Zhou et al. 2014
SP2/0-AG14	Mouse	Hybridoma	Macsik et al. 2015
STHdh	Mouse	Striatal cells	Lu et al. 2013
SW684	Human	Fibrosarcoma	Zhu et al. 2008
SW872	Human	Liposarcoma	Zhu et al. 2008
T24	Human	Bladder carcinoma	Brascesco et al. 2013 (A)
THP-1	Human	Monocyte	Cestari et al. 2012 ; Dahlawi et al. 2013
TK6	Human	Splenic lymphoblast	Stopper et al. 2009
U2OS	Human	Osteosarcoma	Benetti and Roizman 2007 ; Klotz et al. 2012 ; Lanz et al. 2013
U-251 MG	Human	Glioblastoma	Liu et al. 2015 ; Liu et al. 2016 ; Overmeyer et al. 2008
U-373 MG	Human	Glioblastoma	Jones and Howl 2011
U-87 MG	Human	Glioblastoma	Cribbes et al. 2017 ; Tyciakova et al. 2015
U937	Human	Lymphoma	Wang et al. 2013
VCaP	Human	Prostate cancer	Alinezhad et al. 2014 ; Urbinati et al. 2015
WEHI 7.2	Mouse	Lymphoid	Monaco et al. 2012



Primary cell types used with NucView® 530

Cell type	Species	References
Embryonic fibroblast (MEF)	Mouse	Yao et al. 2018

Primary cell types used with NucView® 488

Cell type	Species	References
Amoeba	Acanthamoeba	Wu et al. 2017
Adipose mesenchymal stem cells	Human	Levy et al. 2014 ; Park et al. 2011
Alveolar epithelial cells	Mouse	Standiford et al. 2012
Astrocyte	Rat	Angelova et al. 2016
B cells	Mouse	Zhang et al. 2013
Brain (whole, ex vivo)	Mouse	Shaw et al. 2015
Cortical neurons	Rat	Kamynina et al. 2013 ; Kovac et al. 2014
Dendritic cells	Mouse	Brodsky and Medzhitov 2008
Embryonic fibroblast (MEF)	Mouse	Chen et al. 2010 ; Handa et al. 2011 ; Yao et al. 2018 ; Zhang et al. 2014
Embryonic stem cells (H9)	Human	Son et al. 2013
Embryo tailbud	Chicken	Olivera-Martinez et al. 2012
Fibroblasts, primary	Human	Dikovskya et al. 2015 ; Kato et al. 2013
Foreskin fibroblasts (HFF)	Human	Cunha et al. 2015 ; Son et al. 2013
Gingival fibroblasts	Human	Tanne et al. 2013
Glia	Rat	Kamynina et al. 2013
Hemocytes	Silkworm (<i>Bombyx mori</i>)	Ishii et al. 2012
Hepatocytes	Rat, Mouse	Kawasaki et al. 2015 ; Kopek et al. 2017 ; Zhang et al. 2011
Hippocampal neurons	Rat	Lefort et al. 2012 ; Kamynina et al. 2013 ; Volosin et al. 2008
Idiopathic pulmonary fibrosis fibroblasts	Human	Vuorinen et al. 2007
Immature B cells	Mouse	Claudio et al. 2009
iPS-derived neurons	Human	Yao et al. 2015
Kidney epithelial cells	Mouse	Leuenroth et al. 2007 ; Schmid et al. 2008
Trypanosome	Leishmania	Awasthi et al. 2016 ; Kathuria et al. 2014
Macrophages	Mouse	Brodsky and Medzhitov 2008 ; Ousingsawat et al. 2015 ; Salpeter et al. 2015
Mammary epithelial cells	Mouse	Jechlinger et al. 2009
Mammary luminal progenitor cells	Human	Knapp et al. 2017
Myoblasts	Pig	Mau et al. 2008
Neutrophils	Human	Amendola et al. 2015 ; Majewska et al. 2012
Neural progenitor cells	Human	Gualda et al. 2014
Neurons, primary	Human	Balez et al. 2016
Neurons, primary	Mouse	Cheli et al. 2015
SVZ neural progenitor cells	Rat	Teng et al. 2008

Oligodendrocytes	Mouse	Clausi et al. 2012 ; Paez et al. 2009
Oligodendrocyte progenitor cells	Mouse	Cheli et al. 2014 ; Guardia et al. 2012
Oocytes	Bovine, mouse	Kujjo et al. 2012
Pancreatic acinar cells	Mouse	Adhikari et al. 2008
Pancreatic beta cells	Rat	Dufer et al. 2011 ; Schmitt et al. 2011
Pancreatic islet cells	Mouse	Gier et al. 2009
Peritoneal macrophages	Mouse	Hanley et al. 2012 ; Yancey et al. 2010
Pollen tubes	Field poppy (<i>Papaver rhoeas</i>)	Bosch and Franklin-Tong 2007
Retinal pigmented epithelial cells	Human, mouse	Yang et al. 2009 ; Yang et al. 2011
Skin fibroblasts	Sand cat (<i>Felis margarita</i>)	Gomez et al. 2008
Skin biopsy cells	Human	De Sica et al. 2016
Smooth muscle cells	Pig	Riches et al. 2013
Stem cells	Human	Choo and Fong 2013 ; Cunha et al. 2015
Thymocytes	Mouse	Tribulatti et al. 2007
T-lymphocytes	Human	Valente et al. 2014
Umbilical vein endothelial cells	Human	Geng et al. 2009
Vascular endothelial cells	Rat	Merlet et al. 2013

3D cell culture used with NucView® 488

Cell system	References
3D PC3 cell culture	Bjork et al. 2015
3D PC9 cell culture	Dereli-Korkut et al. 2014
3D primary mouse epithelium cell culture	Jechlinger et al. 2009
VCaP organoids	Alinezhad et al. 2014
U-87 MG 3D spheroids	Cribbes et al. 2017
HT29 3D spheroids	Santo et al. 2016
Neural 3D aggregates	Gualda et al. 2014
H9c2 cells-on-chip	He et al. 2014
MKN-45 3D cells-on-chip	Liu et al. 2016
Ex vivo mouse brain	Shaw et al. 2015
Ex vivo mouse trachea	Tadokoro et al. 2016