



CYTO-ID[®] Autophagy Detection Kit (ENZ-51031): Cited Samples

Human Cell Lines

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Human	Bladder	Carcinoma	Cell line	5637	D. Lihuan, et al. (2014)
Human	Bladder	Carcinoma	Cell line	T24	D. Lihuan, et al. (2014)
Human	Blood	Leukemia	Cell line	HL60	N. Orfali, et al. (2015) ; N. Sugimori, et al. (2015) ; T. Kozako, et al. (2015) ; T.J. Brown, et al. (2013)
Human	Blood	Leukemia	Cell line	Jurkat	T. Kozako, et al. (2015) ; L.L. Chan, et al. (2012)
Human	Blood	Leukemia	Cell line	K562	S. Jiang, et al. (2016) ; S. Guo, et al. (2015) ; C. Drullion, et al. (2012)
Human	Blood	Leukemia	Cell line	KG-1	S.A.F. Morad, et al. (2019)
Human	Blood	Leukemia	Cell line	MT2	T. Kozako, et al. (2015)
Human	Blood	Leukemia	Cell line	MV4-11	S.A.F. Morad, et al. (2019)
Human	Blood	Leukemia	Cell line	NALM-6	S. Skah, et al. (2018)
Human	Blood	Leukemia	Cell line	NB4	N. Orfali, et al. (2015)
Human	Blood	Leukemia	Cell line	REH	S. Skah, et al. (2018)
Human	Blood	Leukemia	Cell line	S1T	T. Kozako, et al. (2015)
Human	Blood	Leukemia	Cell line	THP-1	S.J. Yu, et al. (2015)
Human	Blood	Multiple myeloma	Cell line	KMS-11	I. Riz, et al. (2015)
Human	Blood	Multiple myeloma	Cell line	KMS-34	I. Riz, et al. (2015)
Human	Blood	Multiple myeloma	Cell line	RPMI 8226	S. Rosebeck, et al. (2016) ; S. Jagannathan, et al. (2015) ; M.A. Abdel Malek, et al. (2014) ; C.M. Thirukkumaran, et al. (2012)
Human	Bone	Osteosarcoma	Cell line	MG-63	N.K. Niu, et al. (2015)
Human	Bone	Osteosarcoma	Cell line	Saos-2	T. Bergeron, et al. (2013)
Human	Bone	Osteosarcoma	Cell line	U-2 OS	N.K. Niu, et al. (2015)
Human	Brain	Glioblastoma	Cell line	LN229	Y. He, et al. (2019) ; A.V. Knizhnik, et al. (2013)
Human	Brain	Glioblastoma	Cell line	LN308	Y. He, et al. (2019)
Human	Brain	Glioblastoma	Cell line	T98G	H. Schmeisser, et al. (2013)
Human	Brain	Glioblastoma	Cell line	U-87	H. Burckel, et al. (2015) ; A.V. Knizhnik, et al. (2013) ; E. Josset, et al. (2013)
Human	Brain	Glioblastoma	Cell line	U-251	H. Wu, et al. (2015)
Human	Brain	Medulloblastoma	Cell line	ONS-76	T.W. Bodnarchuk, et al. (2012)
Human	Breast	Adenocarcinoma	Cell line	BT-474	M.R. Akl, et al. (2015)
Human	Breast	Adenocarcinoma	Cell line	LCC1	J.L. Schwartz-Roberts, et al. (2013)
Human	Breast	Adenocarcinoma	Cell line	LCC9	J.L. Schwartz-Roberts, et al. (2013)
Human	Breast	Adenocarcinoma	Cell line	MCF-7	R. Jakhar, et al. (2016) ; J.P. Li, et al. (2015) ; M. Balan, et al. (2014) ; J.L. Schwartz-Roberts, et al. (2013)
Human	Breast	Adenocarcinoma	Cell line	MDA-MB-231	J.P. Li, et al. (2015) ; T.P. Liu, et al. (2015) ; Q. Tan, et al. (2015) ; J.M. Shi, et al. (2013) ; B. Pandit, et al. (2011)
Human	Breast	Adenocarcinoma	Cell line	MDA-MB-468	Q. Tan, et al. (2015)
Human	Breast	Adenocarcinoma	Cell line	T47D	B.J. Metge, et al. (2015) ; M. Balan, et al. (2014)

Complete citation can be found on product's webpage

Advance Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Human	Breast	Normal epithelial	Cell line	MCF10A	J.P. Li, et al. (2015)
Human	Cervix	Adenocarcinoma	Cell line	HeLa	D. Huang, et al. (2017) ; S. Guo, et al. (2015) ; A. Brüning, et al. (2014) ; A. Brüning, et al. (2013) ; H. Schmeisser, et al. (2013)
Human	Cervix	HeLa contaminant (contains HPV)	Cell line	KB	R. Onodera, et al. (2013)
Human	Cervix	Epithelial	Cell line	VK2/E6E7	R.D. Wagner, et al. (2017)
Human	Colon	Adenocarcinoma	Cell line	Caco-2	J. Keenan, et al. (2018) ; K. Pettersen, et al. (2016) ; L.C. Koay, et al. (2014)
Human	Colon	Adenocarcinoma	Cell line	Caco-2/TC7	A. Brüning, et al. (2014)
Human	Colon	Adenocarcinoma	Cell line	DLD1	I. Lee, et al. (2012)
Human	Colon	Adenocarcinoma	Cell line	HCT-8	I. Lee, et al. (2012)
Human	Colon	Adenocarcinoma	Cell line	HT-29	I. Lee, et al. (2012)
Human	Colon	Adenocarcinoma	Cell line	SW480	H. Kauntz, et al. (2011)
Human	Colon	Adenocarcinoma, metastatic site	Cell line	LoVo	T.C. Chang, et al. (2019) ; J. Keenan, et al. (2018) ; I. Lee, et al. (2012)
Human	Colon	Adenocarcinoma, metastatic site	Cell line	SW620	K. Pettersen, et al. (2016) ; H. Kauntz, et al. (2011)
Human	Colon	Carcinoma	Cell line	HCT-116	T.C. Chang, et al. (2019) ; I. Lee, et al. (2012) ; B. Pandit, et al. (2011)
Human	Esophagus	Carcinoma	Cell line	SNO	E. Wolmarans, et al. (2014)
Human	Esophagus	Squamous cell carcinoma	Cell line	EC109	Y.Y. Wang, et al. (2015)
Human	Esophagus	Squamous cell carcinoma	Cell line	KYSE-450	M.J. Nyhan, et al. (2016) ; T.R. O'Donovan, et al. (2015)
Human	Esophagus	Squamous cell carcinoma	Cell line	KYSE-70	Y.Y. Wang, et al. (2015)
Human	Esophagus	Adenocarcinoma	Cell line	OE19	M.J. Nyhan, et al. (2016) ; T.R. O'Donovan, et al. (2015)
Human	Esophagus	Squamous cell carcinoma	Cell line	OE21	M.J. Nyhan, et al. (2016) ; T.R. O'Donovan, et al. (2015)
Human	Esophagus	Adenocarcinoma	Cell line	OE33	M.J. Nyhan, et al. (2016) ; T.R. O'Donovan, et al. (2015)
Human	Gastric	Adenocarcinoma	Cell line	MKN45	I. Lee, et al. (2012)
Human	Gastric	Adenocarcinoma	Cell line	MKN74	I. Lee, et al. (2012)
Human	Kidney	Adenocarcinoma	Cell line	786-0	P. Banerjee, et al. (2012)
Human	Kidney	Carcinoma	Cell line	A498	R.T. Williams, et al. (2013)
Human	Kidney	Embryonic	Cell line	HEK293	E.D. Laing, et al. (2019) ; N.R. Lee, et al. (2018) ; S. Guo, et al. (2015)
Human	Larynx	Squamous cell carcinoma	Cell line	Tu212	C. Lin, et al. (2015)
Human	Liver	Carcinoma	Cell line	Bel-7402	D.M. Zhang, et al. (2013)
Human	Liver	Carcinoma	Cell line	Hep3B	D.M. Zhang, et al. (2013)
Human	Liver	Carcinoma	Cell line	HepG2	Y.N. Liu, et al. (2015) ; J. Klose, et al. (2014) ; A. Brüning, et al. (2014) ; D.M. Zhang, et al. (2013)
Human	Liver	Carcinoma	Cell line	HepJ5	P.L. Wei, et al. (2019)
Human	Liver	Carcinoma	Cell line	Huh7	J. Klose, et al. (2014)
Human	Liver	Carcinoma	Cell line	Mahlavu	P.L. Wei, et al. (2019)
Human	Lung	Adenocarcinoma	Cell line	H23	Y.C. Li, et al. (2014)
Human	Lung	Carcinoma	Cell line	A549	H.Y. Zhu, et al. (2015) ; Y.C. Yin, et al. (2015) ; I.L. Hsin, et al. (2015) ; Y.C. Li, et al. (2014) ; H. Schmeisser, et al. (2013)
Human	Lung	Carcinoma	Cell line	H1299	Y.C. Yin, et al. (2015)
Human	Lung	Carcinoma	Cell line	H460	H.M. Wahrenius, et al. (2011)
Human	Lymph	Lymphoma	Cell line	BJAB	H. Schmeisser, et al. (2013)
Human	Lymph	Lymphoma	Cell line	Daudi	X. Zeng, et al. (2013) ; H. Schmeisser, et al. (2013)
Human	Lymph	Lymphoma	Cell line	Karpas-422	S. Samuel, et al. (2013)

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Human	Lymph	Lymphoma	Cell line	Raji	J. Fan, et al. (2013) ; X. Zeng, et al. (2013)
Human	Lymph	Lymphoma	Cell line	REC-1	L. Rosich, et al. (2012)
Human	Lymph	Lymphoma	Cell line	U937	H. Schmeisser, et al. (2013)
Human	Lymph	Lymphoma	Cell line	Z138	L. Rosich, et al. (2012)
Human	Ovary	Adenocarcinoma	Cell line	PEO1	F. Simpkins, et al. (2012)
Human	Ovary	Adenocarcinoma	Cell line	SKOV3	Y.H. Ding, et al. (2015)
Human	Ovary	Carcinoma	Cell line	BG-1	F. Simpkins, et al. (2012)
Human	Ovary	Carcinoma	Cell line	OVCAR-3	F. Lang, et al. (2015)
Human	Ovary	Carcinoma	Cell line	OVCAR-4	Y.H. Ding, et al. (2015)
Human	Ovary	Carcinoma	Cell line	OVCAR-8	S. Guo, et al. (2015)
Human	Ovary	Carcinoma	Cell line	OVK18	S.I. Wada, et al. (2015)
Human	Pancreas	Adenocarcinoma	Cell line	BxPC-3	F. Wang, et al. (2015)
Human	Pancreas	Adenocarcinoma	Cell line	Capan-2	J. Gong, et al. (2014)
Human	Pancreas	Carcinoma	Cell line	MIA PaCa-2	D. Uchida, et al. (2014) ; B. Pandit, et al. (2011)
Human	Pancreas	Carcinoma	Cell line	PANC-1	F. Wang, et al. (2015)
Human	Pleural effusion	Myeloma	Cell line	KMS-11	I. Riz, et al. (2015)
Human	Prostate	Adenocarcinoma	Cell line	PC3	Z.W. Zhou, et al. (2015) ; L. Shulak, et al. (2014)
Human	Prostate	Carcinoma	Cell line	DU145	Z.W. Zhou, et al. (2015) ; G. Li, et al. (2013)
Human	Prostate	Carcinoma	Cell line	LNCaP	B. Boutin, et al. (2013) ; G. Li, et al. (2013)
Human	Prostate	Carcinoma (derived from LNCaP)	Cell line	C4-2B	G. Li, et al. (2013)
Human	Prostate	Normal epithelial	Cell line	RWPE-1	G. Li, et al. (2013)
Human	Skin	Keratinocytes	Cell line	HaCaT	Y. Jang, et al. (2016) ; G. Soong, et al. (2015)
Human	Skin	Melanoma	Cell line	A375	Z. Wang, et al. (2014)
Human	Skin	Melanoma	Cell line	Ihara	K. Motoyama, et al. (2015)
Human	Skin	Melanoma, metastatic site	Cell line	Mewo	S.I. Wada, et al. (2015)
Human	Stomach	Adenocarcinoma	Cell line	AGS	C.X. Huan, et al. (2015) ; C.X. Huan, et al. (2015)
Human	Stomach	Carcinoma	Cell line	NCI-N87	C.X. Huan, et al. (2015) ; C.X. Huan, et al. (2015)
Human	Tongue	Squamous cell carcinoma	Cell line	SCC25	S.T. Pan, et al. (2015)

Human Primary Cells

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Human		Alzheimer's disease fibroblasts			D. Ma, et al. (2014)
Human	Blood	Chronic myeloid leukemia	Primary		K. Rothe, et al. (2014)
Human	Blood	Immune cells from patients with RA	Primary		Y.M. Chen, et al. (2018)
Human		Idiopathic pulmonary fibrosis fibroblasts			J. Im, et al. (2015)
Human		Parkinson's disease and control cybrid cells			D.M. Arduino, et al. (2012)
Human	Blood	PBMC	Primary		M.V. Stankov, et al. (2014) ; R. van der Burgh, et al. (2014) ; J. van Loodregt, et al. (2013)
Human	Blood	Sorted CD4+ T cells	Primary		J. van Loodregt, et al. (2013)
Human	Bone marrow	Multiple myeloma	Primary		M.A. Abdel Malek, et al. (2014)
Human	Bone marrow	Leukemic blasts	Primary		S. Skah, et al. (2018)
Human	Brain	Glioma stem-like cells	Primary		R.K. Balvers, et al. (2015)
Human	Umbilical vein	Endothelial cells	Primary	HUVEC	R. Menghini, et al. (2014)

Non-Human Cell Lines

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Canine	Bone	Osteosarcoma	Cell line	D-17	T. Bergeron, et al. (2013)
Hamster	Ovary	Epithelial	Cell line	CHO	J.S. Lee, et al. (2011)
Murine	Blood	Pro B cell line	Cell line	Ba/F3	S. Jiang, et al. (2016)
Murine	Blood	Leukemia	Cell line	C1498	T.J. Brown, et al. (2013)
Murine	Blood	Leukemia	Cell line	RAW 264.7	N. Goldstein, et al. (2019) ; Y. Zhang, et al. (2015)
Murine	Brain	Glioblastoma	Cell line	GL261	K. Suzuki, et al. (2013)
Murine	Brain	Neuroblastoma	Cell line	N2a/APP	L.F. Zhang, et al. (2015)
Murine	Colon	Carcinoma	Cell line	CT26	T.R. O'Donovan, et al. (2015) ; H. Bugaut, et al. (2013)
Murine	Embryo	Blastocyst-derived	Cell line	ES-E14TG2a	N. Ozeki, et al. (2016)
Murine	Liver	Normal hepatocytes	Cell line	AML-12	Y. Zhang, et al. (2015)
Rat	Heart	Cardiomyoblasts	Cell line	H9c2	H.K., Sung, et al. (2019) ; Q. Chen, et al. (2016) ; A. Guilbert, et al. (2014)

Non-Human Primary Cells

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Bovine	Heart	Aortic endothelial cells	Primary	BAEC	T. Shiroto, et al. (2014) ; C.L. Oeste, et al. (2013)
Mouse		Podocytes			C. Li, et al. (2015)
Murine	Blood	B cells (naive, memory and plasmablast)	Primary		A.J. Clarke, et al. (2014)
Murine	Blood	CD4+ T cells	Primary		C.W. Lin, et al. (2014)
Murine	Blood	CD8+ T cells	Primary		C.W. Lin, et al. (2014)
Murine	Blood	iNKT cells	Primary		B. Pei, et al. (2015)
Murine	Embryo	Fibroblasts	Primary		M.M. St. Amad, et al. (2016)
Murine	Epididymis	Adipocytes	Primary		N. Goldstein, et al. (2019)
Murine	Heart	Coronary arterial myocytes	Primary		M. Xu, et al. (2014) ; Y. Zhang, et al. (2014) ; Y.M. Wei, et al. (2013)
Murine	Heart	Vascular smooth muscle cells	Primary		C.F. Michiels, et al. (2015)
Murine	Liver	Hepatocytes	Primary		J. Niemann, et al. (2017)
Murine	Liver	Lymphocytes	Primary		T.E. O'Sullivan, et al. (2016)
Murine	Pancreas	Beta cells	Primary		T. Takatani, et al. (2016)
Murine	Spleen	Splenocytes	Primary		T.E. O'Sullivan, et al. (2016)
Porcine	Ocular	Trabecular meshwork	Primary		K. Porter, et al. (2013)
Rat	Brain	Cortex	Primary		S.N. Weis, et al. (2014)
Rat	Brain	Hippocampus	Primary		S.N. Weis, et al. (2014)
Rat	Heart	Neonatal ventricular myocytes	Primary		A.C. Laurent, et al. (2015)
Rat	Spinal cord	Spinal neurons	Primary		Y. Rong, et al. (2019)

Exotic Cells

SPECIES	CELL	CELL TYPE	NAME	REFERENCES
Beetle	Hemocytes from <i>A. dichotoma</i> larvae			S. Hwang, et al. (2015)
Fish	Grouper brain cells			Y. Huang, et al. (2015)
Fungus	Mycelium from <i>A. parasticus</i>			L.V. Roze, et al. (2015)
Human	Alpha 7 integrin- positive skeletal muscle-derived stems cells		$\alpha 7^+$ hSMSC	N. Ozeki, et al. (2016)
Human	Alzheimer's disease fibroblasts			D. Ma, et al. (2014)
Human	Ataxia telangiectasia fibroblasts	Cell line	FT169A	S.D. Desai, et al. (2013)
Human	Gingival fibroblasts	Cell line	HGF	A.M. Pabst, et al. (2014)

SPECIES	CELL	CELL TYPE	NAME	REFERENCES
Human	iPSC-derived hepatocytes			O. Sirenko, et al. (2014)
Human	Idiopathic pulmonary fibrosis fibroblasts			J. Im, et al. (2015)
Human	Oral keratinocytes	Cell line	HOK	A.M. Pabst, et al. (2014)
Human	Osteoblast-like cells			A.M. Pabst, et al. (2014)
Human	Parkinson's disease and control cybrid cells			D.M. Arduino, et al. (2012)
Human	Tenon tissue fibroblasts			A. Want, et al. (2016)
Mouse	Podocytes			C. Li, et al. (2015)
Oyster	Hemocytes			S. Picot, et al. (2019)
Parasite	<i>T.vaginalis</i>			K.Y. Huang, et al. (2014)
Worm	Hemocytes from <i>T.molitor</i> larvae			H. Tindwa, et al. (2015)