



PROTEOSTAT[®] Aggresome Detection Kit (ENZ-51035): Cited Samples

Cell Lines

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Drosophila	Embryo	Cultured cells	Cell line	Kc167	T.P. Miettinen, et al. (2015)
Human	Blood	Leukemia	Cell line	Jurkat	D. Shen, et al. (2011) ; T.P. Miettinen, et al. (2015)
Human	Blood	Multiple myeloma	Cell line	IM-9	S. Moriya, et al. (2013)
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) ; S. Moriya, et al. (2013)
Human	Blood	Multiple myeloma	Cell line	U266	M.A. Abdel Malek, et al. (2014) ; M.A. Hamouda, et al. (2014) ; S. Moriya, et al. (2013)
Human	Bone	Osteosarcoma	Cell line	U-2 OS	H. Olzscha, et al. (2017) ; T.P. Miettinen, et al. (2015)
Human	Bone marrow	Neuroblastoma	Cell line	SH-SY5Y	Y. Bang, et al. (2016)
Human	Brain	Neuroblastoma	Cell line	SK-N-SH	D. Shen, et al. (2011)
Human	Breast	Adenocarcinoma	Cell line	MCF-7	D.S. Nkandeu, et al. (2013)
Human	Cervix	Adenocarcinoma	Cell line	HeLa	L. Wunderley, et al. (2014) ; A. Kothawala, et al. (2012) ; I. Raju, et al. (2011) ; D. Shen, et al. (2011)
Human	Cervix	Epithelial	Cell line	VK2/E6E7	R.D. Wagner, et al. (2017)
Human	Colon	Carcinoma	Cell line	Caco-2	J. Keenan, et al. (2018)
Human	Colon	Carcinoma	Cell line	HCT-116	P. Magnaghi, et al. (2013)
Human	Colon	Carcinoma	Cell line	HT-29	J. Keenan, et al. (2018)
Human	Embryo	Epithelial stem cells	Cell line	H9	S. Koyuncu, et al. (2018)
Human	Embryo	Mesenchymal stem cells derived from H9 epithelial stem cells	Cell line	H9	S. Wang, et al. (2018)
Human	Esophagus	Carcinoma	Cell line	SNO	E. Wolmarans, et al. (2014)
Human	Kidney	Embryonic	Cell line	HEK293	Y. Nibe, et al. (2018) ; Y. Bang, et al. (2016) ; S. Genier, et al. (2016) ; Y. Kamada, et al. (2015) ; B. Jeric, et al. (2013) ; J. van Hengel, et al. (2012)
Human	Liver	Carcinoma	Cell line	HuH7.5.1	S. Narayanan, et al. (2016)
Human	Lung	Carcinoma	Cell line	A549	I.V. Karagounis, et al. (2016)
Human	Lung	Normal epithelial	Cell line	BEAS-2B	N. Vij, et al. (2016) ; P.C. Shivalingappa, et al. (2016) ; H.W. Chiu, et al. (2015)
Human	Lung	Carcinoma	Cell line	H1299	I.V. Karagounis, et al. (2016)
Human	Pancreas	Carcinoma	Cell line	MIA PaCa-2	M.T. Pacheco, et al. (2014)
Human	Prostate	<i>Adenocarcinoma</i>	Cell line	PC3	M.I. Koukourakis, et al. (2015)
Human	Prostate	Carcinoma	Cell line	DU145	M.I. Koukourakis, et al. (2015)
Human	Skin	Melanoma	Cell line	SK-MEL-28	M.T. Pacheco, et al. (2014)
Human	Stomach	Adenocarcinoma	Cell line	AGS	C.C. Yu, et al. (2013)
Human	Umbilical vein	Endothelial cells	Cell line	HUVEC	T.P. Miettinen, et al. (2015)
Murine	Liver	Hepatocytes	Cell line	Hepa-1c1c7	K. Ohnishi, et al. (2013)
Murine	Pancreas	Insulinoma (beta cells)	Cell line	MIN6	A. Hofmeister-Brix, et al. (2013)
Murine	PNS	Schwann cells	Cell line	TgS1	T. Murakami, et al. (2015)

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

Murine	Blood	Leukemia	Cell line	RAW 264.7	H.W. Chiu, et al. (2015)
Murine	Brain	Neuroblastoma	Cell line	Neuro2a	D. Ishibashi, et al. (2016)
Plant	Tobacco	Cultured cells	Cell line	BY-2	Y. Nakajima, et al. (2013)
Murine	Brain	Neuronally-derived cells	Cell line	ST HDH Q7/7	K. Shen, et al. (2019)
Murine	Brain	Neuronally-derived cells	Cell line	STHDH Q7/111	K. Shen, et al. (2019)
Rat	Heart	Cardiomyoblasts	Cell line	H9c2	J. Liu, et al. (2016) ; J.W. Jahng, et al. (2015)
Rat	Pancreas	Tumor cells	Cell line	AR42J	R.T. Waldron, et al. (2018)
Xenopus	Kidney	Epithelial cells from X.laevis	Cell line	A6	S. Khan, et al. (2014)

Primary Cells

SPECIES	TISSUE	CELL	CELL TYPE	NAME	REFERENCES
Human		Cryopreserved semen	Primary		Y.J. Yi, et al. (2014)
Human	Brain	iPSC neuronal precursor cells	Primary		S.M. Usmani, et al. (2014) ; C.E. Kennedy, et al. (2014)
Murine	Brain	Astrocytes	Primary		M. Polajnar, et al. (2014)
Murine	Brain	Neurons	Primary		B. Schattling, et al. (2019)
Murine	Breast	Epithelial cells	Primary		C. Mukhopadhyay, et al. (2016)
Murine	Heart	Cardiomyocytes	Primary		J. Hu, et al. (2017) ; Y. Hao, et al. (2016)
Murine	Intestine	Peritoneum-derived macrophages	Primary		Y. Nakajima, et al. (2013)
Murine	Liver	Fetal liver hematopoietic stem/progenitor cells	Primary		V. Sigurdsson, et al. (2016)
Murine	Liver	Hepatocytes	Primary		Y. Yamada, et al. (2015) ; U. Tomaru, et al. (2012)
Murine	Skin	Fibroblasts	Primary		A. Hofmeister-Brix, et al. (2013)
Plant	Grapevine	Cultured cells	Primary		Y. Nakajima, et al. (2013)
Plant	Tobacco	Cultured cells	Primary		A. Flierl, et al. (2014)
Porcine		Oocytes	Primary		T. Mocan, et al. (2015)
Rat	Brain	Neurons	Primary		S.S. Sanders, et al. (2019)
Squirrel	Brain	Neurons	Primary		J. Ou, et al. (2018)

Other Samples

SPECIES	TISSUE	SAMPLE TYPE	CELL TYPE	NAME	REFERENCES
Bacteria	<i>E.coli</i>	Cells			S. Navarro, et al. (2016)
Fly	Brain	Tissue section			D.S. Coehlo et al. (2018)
Human	Lung	Tissue section			A.S. Patel, et al. (2012)
Murine	Heart	Tissue section			Y. Maejima, et al. (2013)
Murine	Lung	Tissue section			N. Vij, et al. (2018)
Rat	Brain	Tissue section			H.M. Abdou, et al. (2016)
Rat	Heart	Tissue section			J. Liu, et al. (2016)
Rat	Kidney	Homogenates			M.U. Cheema, et al. (2013)