



Antibodies: Cited Samples

Goat anti-rabbit IgG, polyclonal antibody (HRP conjugate) (ADI-SAB-300) ELISA

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Bacteria				M. Amarasiri, et al. (2016)
Mollusc	Digestive gland	Tissue lysate		F. Gagné, et al. (2013)
Mollusc	Gill	Tissue lysate		F. Gagné, et al. (2013)

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Kidney	Tissue sections		E. Fervenza, et al. (2008)
Rabbit	Blood	Blood cells		I. Guerrero, et al. (2015)

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Coral		Total lysate		D. Seveso, et al. (2017)
Cow	Heart	Bioprosthetic heart valve		S. Lee, et al. (2019)
Cow	Uterus	Endometrial cells		T. Acosta, et al. (2008)
Cuttlefish	Gill	Tissue lysate		J.C. Capaz, et al. (2017)
Cuttlefish	Heart tissue	Tissue lysate		J.C. Capaz, et al. (2017)
Cuttlefish	Mantle tissue	Tissue lysate		J.C. Capaz, et al. (2017)
Fish	Erythrocytes	Tissue lysate		C.D. Ricketts, et al. (2015)
Fish	Heart	Tissue lysate		C.D. Ricketts, et al. (2015)
Fish	Liver	Tissue lysate		C.D. Ricketts, et al. (2015)
Human	Adipose tissue	Adipose-derived stem cells		D. Chae, et al. (2018)
Human	Blood	Mast cell leukemia	HMC-1	E. A. Park, et al. (2019)
Human	Blood	Peripheral Blood Mononuclear Cells	PBMC	W. Wang, et al. (2015)
Human	Blood	Lymphocytes		A. Rothfuss et al. (2001)
Human	Bone marrow	Mesenchymal stem cells		J.Y. Tae, et al. (2018)
Human	Brain	Microglia	HM06	S. K. Jang, et al. (2015)
Human	Breast	Carcinoma	MCF7	S.J. Park, et al. (2015)
Human	Breast	Carcinoma	MCF7-ADR	S.J. Park, et al. (2015)

Complete citation can be found on product's webpage

Global Headquarters
ENZO LIFE SCIENCES, INC.
Ph: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES (ELS) AG
Ph: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg
Ph: +32 3 466 0420
info-be@enzolifesciences.com

France
Ph: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Ph: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Ph: 0845 601 1488 (UK customers)
Ph: +44 1392 825900
info-uk@enzolifesciences.com

For local distributors and detailed product information visit us online: www.enzolifesciences.com

For Research Only. Not for Use in Diagnostic Procedures.

Human	Breast	Carcinoma	MDA-MB-231	S.J. Park, et al. (2015)
Human	Breast	Carcinoma	MDA-MB-231	S.A.E. Ibrahim, et al. (2018)
Human	Breast	Carcinoma	MDA-MB-435s	S.J. Park, et al. (2015)
Human	Breast	Carcinoma	MDA-MB-468	S.A.E. Ibrahim, et al. (2018)
Human	Breast	Carcinoma	SK-BR-3	S.J. Park, et al. (2015)
Human	Cervix	Adenocarcinoma	HeLa	H. Nassour, et al. (2016)
Human	Cervix	Adenocarcinoma; Squamous cell carcinoma	Hela	M.S. Kang, et al. (2018)
Human	Cervix	Adenocarcinoma	HeLa	H.I. Ka, et al. (2020)
Human	Cervix	Squamous cell carcinoma	ME-180	M.S. Kang, et al. (2018)
Human	Colon	Colorectal cancer	HCT-116; SNU-81	S.R. Lee, et al. (2018)
Human	Colon	Colorectal cancer	DLD1	B. Dahal Lamichane, et al. (2019)
Human	Colon	Colorectal cancer	HCT8	B. Dahal Lamichane, et al. (2019)
Human	Colon	Colorectal cancer	HCT116	B. Dahal Lamichane, et al. (2019)
Human	Colon	Colorectal cancer	HT29	B. Dahal Lamichane, et al. (2019)
Human	Colon	Colorectal cancer	LOVO	B. Dahal Lamichane, et al. (2019)
Human	Colon	Colorectal cancer	SNU-81	S.R. Lee, et al. (2018)
Human	Gingiva	Gingiva-derived stem cells		J.Y. Tae, et al. (2018)
Human	Kidney	Embryonic kidney cells	HEK293T	B. Dahal Lamichane, et al. (2019)
Human	Kidney	Embryonic kidney cells	HEK293T	H.I. Ka, et al. (2020)
Human	Kidney	Carcinoma	Ketr-3	Q. Liu, et al. (2018)
Human	Liver	Stellate cells	LX2	S. Wang, et al. (2019)
Human	Liver	Carcinoma	HepG2	H. Nassour, et al. (2016)
Human	Liver	Carcinoma	HepG2	S. Yumnam, et al. (2017)
Human	Liver	Carcinoma	Huh7	H.S. Seol, et al. (2016)
Human	Liver	Carcinoma	SNU449	H.S. Seol, et al. (2016)
Human	Lung	Hembryonic fibroblast	WI-38	S. Raha, et al. (2019)
Human	Muscle			Y.E. Park, et al. (2016)
Human	Muscle	Skeletal Muscle	HSkMC	C.H. Kim, et al. (2016)
Human	Muscle	Skeletal Muscle		J.A. Zoladz, et al. (2014)
Human	Ovary	Adenocarcinoma	A2780	C.H. Choi, et al. (2018)
Human	Ovary	Adenocarcinoma	SKOV-3	C.H. Choi, et al. (2018)
Human	Prostate	Adeonocarcinoma	DU145	U.M. Mazim, et al. (2020)
Human	Prostate	Adeonocarcinoma	DU145	U.M. Mazim, et al. b (2020)
Human	Prostate	Adeonocarcinoma	LNCaP	U.M. Mazim, et al. (2020)
Human	Prostate	Adeonocarcinoma	LNCaP	U.M. Mazim, et al. b (2020)
Human	Prostate	Benign Prostatic Hyperplasia	BPH-1	J.H. Kim, et al. (2018)
Human	Prostate	Stromal cells	WPMY-1	J.H. Kim, et al. (2016)
Human	Saliva	Lysate		F. Ciregia, et al. (2016)
Human	Skin	Keratinocytes	HaCaT	D. Shin, et al. (2018)

Human	Skin	Keratinocytes	HaCaT	D. Shin, et al. b (2018)
Human	Skin	Dermal fibroblasts	CCD986Sk	D. Shin, et al. c (2018)
Human	Skin	Keratinocytes	HaCaT	Y. Oh, et al. (2017)
Human	Skin	Keratinocytes	HaCaT	Y. Oh, et al. (2016)
Human	Skin	Keratinocytes	HaCaT	H.W. Lim, et al. (2016)
Human	Skin	Dermal fibroblasts	CCD986Sk	B. Beomeol, et al. (2016)
Human	Skin	Keratinocytes	HaCaT	S.W. Kim, et al. (2016)
Human	Skin	Keratinocytes	HaCaT	S.J. Oh, et al. (2016)
Human	Skin	Keratinocytes	HaCaT	H.J. Kang, et al. (2016)
Human	Skin	Keratinocytes	HaCaT	S.J. Oh, et al. (2014)
Human	Stomach	Adenocarcinoma	AGS	V. Ventatarama Gowda Saralamma, et al. (2017)
Human	Tongue	Squamous cell carcinoma	CAL 27	S.E. Kim, et al. (2018)
Human	Umbilical vein	Endothelial cells	HUVEC	B.H. Han, et al. (2017)
Mollusc	Gills	Tissue lysate		A. Ravaschiere, et al. (2017)
Mouse	Bone	Femora		Y.J. Moon, et al. (2018)
Mouse	Bone	Fibroblast-like synoviocytes		S.J. Woo, et al. (2015)
Mouse	Bone marrow	Macrophages		D.H. Jeong, et al. (2020)
Mouse	Bone marrow	Total cells		J.Y. Kim, et al. (2019)
Mouse	Brain	Microglia	BV2	S. Raha, et al. (2016)
Mouse	Brain	Microglia	BV2	J.Y. Lee, et al. (2018)
Mouse	Brain	Ischemic region		X. Gui, et al. (2017)
Mouse	Colon			Y. Lee, et al. (2017)
Mouse	Embryo	Fibroblast		D.S. Wheeler, et al. (2003)
Mouse	Embryo	Fibroblasts	MEF	M.N. Kim, et al. (2015)
Mouse	Kidney	Medullary Collecting Duct	mIMCD-3	M.G. Kim, et al. (2017)
Mouse	Pancreas	Adenocarcinoma		J.W. Eng, et al. (2015)
Mouse	Retina			M.C. Trotta, et al. (2019)
Mouse	Spleen			D. Son, et al. (2019)
Mouse	Thymus			D. Son, et al. (2019)
Mouse	Uterus	Tissue lysate and Endothelial cells	HUtMEC	Y.G. Park, et al. (2017)
Pig	Kidney	Proximal Tube	LLC-PK1	K. K. Meldrum, et al. (2001)
Rat	Adrenal gland	Pheochromocytoma	PC-12	Y. Sim, et al. (2016)
Rat	Blood	Erythrocytes		J.S. Kim, et al. (2016)
Rat	Blood	Plasma		J.S. Kim, et al. (2016)
Rat	Bone marrow	Mesenchymal stem cells	MSC	B.W. Song, et al. (2013)
Rat	Brain	Corpus callosum		M.E. Pak, et al. (2018)
Rat	Brain	Hippocampus		H.G. Kim, et al. (2015)
Rat	Heart	Aorta		P. Waghe, et al. (2015)
Rat	Heart	Aortic media		T. Y. Xu, et al. (2015)

Rat	Heart	Cardiomyoblasts	H9c2	M.C. Trotta, et al. (2018)
Rat	Heart	Ventricles		A.A. Fernandes, et al. (2015)
Rat	Heart	Ventricle s		A.A. Fernandes, et al. (2015)
Rat	Intestine			B. Angermayr, et al. (2007)
Rat	Liver	Stellate cells		M.Y. Moon, et al. (2019)
Rat	Skin	Wound (granulation tissue)		M. Ram, et al. (2016)
Shark	Gill	Tissue lysate		A.J. Morash, et al. (2016)
Shark	White muscle	Tissue lysate		A.J. Morash, et al. (2016)
Yeast		Lysate		Y. Shen, et al. (2019)
Yeast		Lysate		Z. Wang, et al. (2014)

For Technical Support:
techserv-usa@enzolifesciences.com
techserv-eu@enzolifesciences.com