

KDEL monoclonal antibody (10C3) (ADI-SPA-827)

ELISA

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Recombinant		KDEL-tagged recombinant protein from plant expression		http://www.ncbi.nlm.nih.gov/pubmed/12151436

EM

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Recombinant		Detection of KDEL-tagged antibody on <i>V. mungo</i> seeds		https://pubmed.ncbi.nlm.nih.gov/10662772/

IF

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Bird		Parasympathetic chick ciliary ganglion cells		https://pubmed.ncbi.nlm.nih.gov/12163544/
Dog		Madin-Darby canine kidney cells	MDCK	https://www.ncbi.nlm.nih.gov/pubmed/26673890
Dog		Madin-Darby canine kidney cells	MDCK	https://www.ncbi.nlm.nih.gov/pubmed/28437467
Hamster	Ovaries	Ovary cells	CHO	http://www.ncbi.nlm.nih.gov/pubmed/26764097
Hamster	Ovaries	Ovary cells	CHO	https://www.ncbi.nlm.nih.gov/pubmed/31662433
Human		iPSC-derived photoreceptors		https://www.ncbi.nlm.nih.gov/pubmed/27889058
Human		Primary fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/31949249
Human	Blood vessel	Primary vascular smooth muscle cells		https://www.ncbi.nlm.nih.gov/pubmed/31242668
Human	Brain	Neuroblastoma cells	SH-SY5Y	http://www.ncbi.nlm.nih.gov/pubmed/26884997
Human	Breast	Mammary gland epithelial cells	T47D	http://www.ncbi.nlm.nih.gov/pubmed/26890638
Human	Cervix	Epithelial Cervix carcinoma cells	HeLa	http://www.ncbi.nlm.nih.gov/pubmed/17387721
Human	Cervix	Carcinoma cells	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/29607936
Human	Embryo	Kidney epithelial cells	HEK293	https://www.pubmed.ncbi.nlm.nih.gov/11938353/
Insect		Drosophila embryonal cells	S2	https://www.pubmed.ncbi.nlm.nih.gov/12581524/

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.

Insect		Drosophila larval wing discs		http://www.ncbi.nlm.nih.gov/pubmed/25942626
Monkey	Kidney	Green monkey fibroblats cells	COS-7	https://pubmed.ncbi.nlm.nih.gov/12368856/
Mouse	Adipose tissue	Fibroblasts	3T3-L1	https://www.ncbi.nlm.nih.gov/pubmed/28111073
Mouse	Blood	Monocytic cells	RAW-D	https://www.ncbi.nlm.nih.gov/pubmed/28791425
Mouse	Blood	Plasmacytoid dendritic cells		https://www.ncbi.nlm.nih.gov/pubmed/25871979
Mouse	Bone marrow	Macrophages		https://www.ncbi.nlm.nih.gov/pubmed/30298952
Mouse	Brain	Cerebellim		https://www.ncbi.nlm.nih.gov/pubmed/30201312
Mouse	Brain	Midbrain		http://www.ncbi.nlm.nih.gov/pubmed/25609609
Mouse	Brain	Neuroblastoma cells	N2a	https://www.ncbi.nlm.nih.gov/pubmed/32376919
Mouse	Brain	Primary cortical neurons		http://www.ncbi.nlm.nih.gov/pubmed/18819914
Mouse	Brain	Primary cortical neurons		http://www.ncbi.nlm.nih.gov/pubmed/25876663
Mouse	Embryo			http://www.ncbi.nlm.nih.gov/pubmed/18547789
Mouse	Embryo			https://www.ncbi.nlm.nih.gov/pubmed/29095923
Mouse	Pancreas	Insulinoma cells	MIN6	http://www.ncbi.nlm.nih.gov/pubmed/23897822
Mouse	Pancreas	Pancreatic islets	MIN6	http://www.ncbi.nlm.nih.gov/pubmed/23897822
Mouse	Pancreas	Pancreatic islets		http://www.ncbi.nlm.nih.gov/pubmed/26046931
Mouse	Spinal cord			http://www.ncbi.nlm.nih.gov/pubmed/17387721
Mouse	Spinal cord			http://www.ncbi.nlm.nih.gov/pubmed/26906090
Rat	Brain			https://www.ncbi.nlm.nih.gov/pubmed/23411409
Rat	Thyroid	Thyrocytes	PCCL3	https://www.ncbi.nlm.nih.gov/pubmed/32241916

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Mouse	Blood vessel	Abdominal aorta		http://www.ncbi.nlm.nih.gov/pubmed/24329494
Human	Blood vessel	Carotid artery		http://www.ncbi.nlm.nih.gov/pubmed/25060744
Human	Blood vessel	Carotid artery		https://www.ncbi.nlm.nih.gov/pubmed/28507169
Mouse	Embryo	Fibroblasts	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/27796797
Mouse	Liver	Liver		http://www.ncbi.nlm.nih.gov/pubmed/26499802
Mouse	Liver			https://www.ncbi.nlm.nih.gov/pubmed/29593095
Mouse	Spinal cord			http://www.ncbi.nlm.nih.gov/pubmed/27030050

IP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Bovine	Kidney	Epithelial cells	MDBK	http://www.ncbi.nlm.nih.gov/pubmed/26939061
Hamster	Kidney	Baby hamster kidney cells	BHK-21	https://pubmed.ncbi.nlm.nih.gov/9557669/
Human	Embryo	Kidney epithelial cells	HEK293	http://www.ncbi.nlm.nih.gov/pubmed/18805786
Recombinant		KDEL sequence as tag for IP in plant expression		http://www.ncbi.nlm.nih.gov/pubmed/12151436

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Hamster	Kidney	Baby hamster kidney cells	BHK-21	https://pubmed.ncbi.nlm.nih.gov/9557669/
Human	Bladder	Carcinoma cells	T24/83	http://www.ncbi.nlm.nih.gov/pubmed/12621026
Human	Blood vessel	Aortic smooth muscle cells	HASMC	http://www.ncbi.nlm.nih.gov/pubmed/12621026
Human	Blood vessel	Human umbilical vein endothelial cells	HUVECs	https://www.ncbi.nlm.nih.gov/pubmed/30396238
Human	Brain	iPSC-derived neurons		https://www.ncbi.nlm.nih.gov/pubmed/29289717
Human	Brain	Neuroblastoma cells	SH-SY5Y	http://www.ncbi.nlm.nih.gov/pubmed/26884997
Human	Cervix	Carcinoma cells	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/22637475
Human	Cervix	Carcinoma cells	HeLa	http://www.ncbi.nlm.nih.gov/pubmed/25880086
Human	Colon	Colorectal tumor biopsy tissue		https://www.ncbi.nlm.nih.gov/pubmed/29374142
Human	Embryo	Kidney epithelial cells	HEK293	http://www.ncbi.nlm.nih.gov/pubmed/18805786
Human	Embryo	Kidney epithelial cells	HEK293	http://www.ncbi.nlm.nih.gov/pubmed/25565208
Human	Embryo	Kidney epithelial cells	HEK293	http://www.ncbi.nlm.nih.gov/pubmed/27079237
Human	Embryo	Kidney epithelial cells	HEK293	https://www.ncbi.nlm.nih.gov/pubmed/30619294
Human	Embryo	Kidney epithelial cells	HEK293	https://www.ncbi.nlm.nih.gov/pubmed/30995221
Human	Liver	Hepatocellular carcinoma cells	HepG2	https://www.ncbi.nlm.nih.gov/pubmed/28827061
Human	Lung	Adenocarcinoma cells	A549	http://www.ncbi.nlm.nih.gov/pubmed/26499802
Human	Lung	Adenocarcinoma cells	A549	https://www.ncbi.nlm.nih.gov/pubmed/26288223
Human	Lung	Non-small cell lung carcinoma cells	H1299	https://www.ncbi.nlm.nih.gov/pubmed/24009855
Human	Pancreas	Pancreatic islets		https://www.ncbi.nlm.nih.gov/pubmed/31184302
Human	Skin	Malignant melanoma cells	A375	https://www.ncbi.nlm.nih.gov/pubmed/27965097
Human	Skin	Malignant melanoma cells	G-361	https://www.ncbi.nlm.nih.gov/pubmed/27965097
Monkey	Kidney	African green monkey kidney cells	37RC	http://www.ncbi.nlm.nih.gov/pubmed/29992955
Mouse	Ascites	Macrophages	J774A.1	https://www.ncbi.nlm.nih.gov/pubmed/28822264
Mouse	Brain	Cerebellum		https://www.ncbi.nlm.nih.gov/pubmed/29289717
Mouse	Brain	Neuroblastoma cells	N2a	http://www.ncbi.nlm.nih.gov/pubmed/17387721
Mouse	Brain	Tissue lysate		https://www.ncbi.nlm.nih.gov/pubmed/27796797
Mouse	Brain	Striatum		https://www.ncbi.nlm.nih.gov/pubmed/29289717
Mouse	Embryo	Fibroblasts	MEFs	http://www.ncbi.nlm.nih.gov/pubmed/26499802
Mouse	Embryo	Fibroblasts	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/27796797

Mouse	Embryo	Fibroblast subcellular fractions	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/31661432
Mouse	Heart	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/12368856/
Mouse	Heart	Tissue lysate		http://www.ncbi.nlm.nih.gov/pubmed/26027784
Mouse	Heart	Tissue lysate		https://www.ncbi.nlm.nih.gov/pubmed/27023784
Mouse	Kidney	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/12368856/
Mouse	Liver	Liver subcellular fractions		https://www.ncbi.nlm.nih.gov/pubmed/23247284
Mouse	Liver	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/12368856/
Mouse	Liver	Tissue lysate		http://www.ncbi.nlm.nih.gov/pubmed/26499802
Mouse	Liver	Tissue lysate		https://www.ncbi.nlm.nih.gov/pubmed/28781602
Mouse	Lung	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/12368856/
Mouse	Muscle	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/12368856/
Mouse	Spleen	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/12368856/
Mouse	Thymus	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/12368856/
Rat	Brain	Adrenal medulla pheochromocytoma cells	PC12	http://www.ncbi.nlm.nih.gov/pubmed/12969271
Rat	Brain	Adrenal medulla pheochromocytoma cells	PC12	https://www.ncbi.nlm.nih.gov/pubmed/26621873
Rat	Brain	Astrocytes		https://www.ncbi.nlm.nih.gov/pubmed/23411409
Rat	Brain	Cerebellum		https://www.ncbi.nlm.nih.gov/pubmed/29289717
Rat	Brain	Cortex		https://www.ncbi.nlm.nih.gov/pubmed/23411409
Rat	Brain	Striatum		https://www.ncbi.nlm.nih.gov/pubmed/29289717
Rat	Brain	Striatum		https://www.ncbi.nlm.nih.gov/pubmed/30366018
Rat	Brain	Substantia nigra		https://www.ncbi.nlm.nih.gov/pubmed/27660269
Rat	Heart	Neonatal ventricular monocytes		https://www.ncbi.nlm.nih.gov/pubmed/30643122
Rat	Heart	Neonatal ventricular cardiomyocytes (NRVCs)		http://www.ncbi.nlm.nih.gov/pubmed/26027784
Rat	Heart	Primary cardiac myocytes		https://www.ncbi.nlm.nih.gov/pubmed/22637475
Rat	Thyroid	Thyocytes	PCCL3	https://www.ncbi.nlm.nih.gov/pubmed/32241916
Yeast		Cell lysates	334	https://pubmed.ncbi.nlm.nih.gov/12363344/
Yeast		Cell lysates	W303a	https://pubmed.ncbi.nlm.nih.gov/12363344/

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

