

Cu/Zn SOD polyclonal antibody (ADI-SOD-100)

Functional

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Embryo	Kidney epithelial cells	HEK293	https://www.ncbi.nlm.nih.gov/pubmed/29991716

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Dog	Spinal Cord			https://www.ncbi.nlm.nih.gov/pubmed/27838005
Dog	Spinal Cord			https://www.ncbi.nlm.nih.gov/pubmed/26162235
Human	Spinal cord			https://www.ncbi.nlm.nih.gov/pubmed/12145308
Mouse	Brain			https://www.ncbi.nlm.nih.gov/pubmed/26643113
Rat	Testicle			https://www.ncbi.nlm.nih.gov/pubmed/27292162

IP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Brain	Neuroglioma cells	H4	https://www.ncbi.nlm.nih.gov/pubmed/26643113
Mouse	Embryo	Fibroblasts	NIH3T3	https://www.ncbi.nlm.nih.gov/pubmed/11278741
Mouse	Spinal cord			https://www.ncbi.nlm.nih.gov/pubmed/29991716

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Fish		Zebrafish whole embryos		https://www.ncbi.nlm.nih.gov/pubmed/30528257
Horse	Bone	Articular chondrocytes		https://www.ncbi.nlm.nih.gov/pubmed/25998424
Human	Bladder	Papilloma/Carcinoma cell nuclear and cytoplasmic extracts	HT1376	https://www.ncbi.nlm.nih.gov/pubmed/27756245
Human	Bladder	Papilloma/Carcinoma cell nuclear and cytoplasmic extracts	RT4	https://www.ncbi.nlm.nih.gov/pubmed/27756245
Human	Bladder	Papilloma/Carcinoma cell nuclear and cytoplasmic extracts	T24	https://www.ncbi.nlm.nih.gov/pubmed/27756245

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	Brain	iPSC-derived Astrocytes		https://www.ncbi.nlm.nih.gov/pubmed/31658977
Human	Brain	Mesencephalic cells	MESC2.10	https://www.ncbi.nlm.nih.gov/pubmed/12145295
Human	Brain	Neuroglioma cells	H4	https://www.ncbi.nlm.nih.gov/pubmed/26643113
Human	Cervix	Carcinoma cells	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/22768276
Human	Embryo	Embryonic kidney cells	HEK293	https://www.ncbi.nlm.nih.gov/pubmed/26297318
Human	Embryo	Embryonic kidney cells	HEK293FT	https://www.ncbi.nlm.nih.gov/pubmed/30341612
Human	Lung	Alveolar type II-like epithelial cells	A549	https://www.ncbi.nlm.nih.gov/pubmed/26073125
Human	Prostate	Adenocarcinoma	DU145	https://www.ncbi.nlm.nih.gov/pubmed/29687779
Human	Prostate	Adenocarcinoma	LNCaP	https://www.ncbi.nlm.nih.gov/pubmed/29687779
Human	Prostate	Adenocarcinoma	PC3	https://www.ncbi.nlm.nih.gov/pubmed/29687779
Human	Skin	Keratinocyte cells	HaCaT	https://www.ncbi.nlm.nih.gov/pubmed/26991844
Human/Dog		Human kidney cells transfected with canine SOD1	HEK293A	https://www.ncbi.nlm.nih.gov/pubmed/26162235
Human/Mouse	Embryo	Mouse fibroblasts overexpressing human SOD1	NIH3T3	https://www.ncbi.nlm.nih.gov/pubmed/11278741
Human/Mouse	Spinal cord	Mouse spinal cords overexpressing human SOD1		https://www.ncbi.nlm.nih.gov/pubmed/11278741
Human/Mouse	Spinal cord	Human-SOD1-transgenic mouse spinal cords		https://www.ncbi.nlm.nih.gov/pubmed/18579581
Mollusc		Clam digestive gland subcellular fraction		https://www.ncbi.nlm.nih.gov/pubmed/17052751
Mouse	Blood vessel	Aorta		https://www.ncbi.nlm.nih.gov/pubmed/22886847
Mouse	Blood vessel	Aorta		https://www.ncbi.nlm.nih.gov/pubmed/30714619
Mouse	Blood vessel	Coronary arteries		https://www.ncbi.nlm.nih.gov/pubmed/27467019
Mouse	Bone	Osteoblast precursor cells	MC3T3 E1	https://www.ncbi.nlm.nih.gov/pubmed/31790903
Mouse	Brain	Brain striatum		https://www.ncbi.nlm.nih.gov/pubmed/27616549
Mouse	Brain	Motor neuronal cells	NSC-34	https://www.ncbi.nlm.nih.gov/pubmed/29415125
Mouse	Brain	Motor neuronal cells	NSC-34	https://www.ncbi.nlm.nih.gov/pubmed/28402699
Mouse	Brain	Neuroblastoma cells	Neuro-2a	https://www.ncbi.nlm.nih.gov/pubmed/22768276
Mouse	Cartilage	Knee joint homogenates		https://www.ncbi.nlm.nih.gov/pubmed/25593283
Mouse	Embryo	Fibroblasts	NIH3T3	https://www.ncbi.nlm.nih.gov/pubmed/26395512
Mouse	Heart			https://www.ncbi.nlm.nih.gov/pubmed/28522596
Mouse	Kidney	Total and nuclear protein extracts		https://www.ncbi.nlm.nih.gov/pubmed/30925687
Mouse	Liver			https://www.ncbi.nlm.nih.gov/pubmed/29237473
Mouse	Liver			https://www.ncbi.nlm.nih.gov/pubmed/30714619
Mouse	Liver	Nuclear extracts		https://www.ncbi.nlm.nih.gov/pubmed/28781602
Mouse	Muscle			https://www.ncbi.nlm.nih.gov/pubmed/29438851
Mouse	Muscle	Gastrocnemius feed		https://www.ncbi.nlm.nih.gov/pubmed/27467019
Mouse	Skin	Dermal fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/23708100
Rat	Blood vessel	Aorta		https://www.ncbi.nlm.nih.gov/pubmed/22796671

Rat	Blood vessel	Cerebral arteries		https://www.ncbi.nlm.nih.gov/pubmed/26945730
Rat	Blood vessel	Feed artery		https://www.ncbi.nlm.nih.gov/pubmed/12381754
Rat	Brain	Nucleus Tractus solitarii (brainstem)		https://www.ncbi.nlm.nih.gov/pubmed/30668894
Rat	Heart			https://www.ncbi.nlm.nih.gov/pubmed/29687779
Rat	Liver			https://www.ncbi.nlm.nih.gov/pubmed/26266409
Rat	Liver	Nuclear extracts		https://www.ncbi.nlm.nih.gov/pubmed/32341232
Rat	Muscle			https://www.ncbi.nlm.nih.gov/pubmed/31306740
Rat	Muscle			https://www.ncbi.nlm.nih.gov/pubmed/30843284
Rat	Ovaries	Whole tissue		https://www.ncbi.nlm.nih.gov/pubmed/26607716
Rat	Spinal cord			https://www.ncbi.nlm.nih.gov/pubmed/31092730

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