

HO-1 polyclonal antibody (ADI-SPA-895)

FC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood	Peripheral Blood Lymphocytes	PBL	https://pubmed.ncbi.nlm.nih.gov/14642388/
Rat	Liver			https://pubmed.ncbi.nlm.nih.gov/9500714/

Immunostaining

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood	Circulating Endothelial Cells	CECs	https://pubmed.ncbi.nlm.nih.gov/11238038/
Human	Blood vessel	Abdominal aortic section		https://pubmed.ncbi.nlm.nih.gov/26296046/
Human	Brain			https://pubmed.ncbi.nlm.nih.gov/8030754/
Human	Kidney			https://pubmed.ncbi.nlm.nih.gov/18805612/
Human	Kidney	Embryonic kidney	HEK-293	https://pubmed.ncbi.nlm.nih.gov/23591596/
Human	Liver	Cholangiocarcinoma		https://pubmed.ncbi.nlm.nih.gov/26726846/
Human	Prostate	Tumor		https://pubmed.ncbi.nlm.nih.gov/27280718/
Mouse	Blood vessel	Aorta		https://pubmed.ncbi.nlm.nih.gov/30108504/
Mouse	Blood vessel	Arteriovenous fistula (AVF)		https://pubmed.ncbi.nlm.nih.gov/21697243/
Mouse	Blood vessel	Infra-renal inferior vena cava		https://pubmed.ncbi.nlm.nih.gov/18988794/
Mouse	Brain			https://pubmed.ncbi.nlm.nih.gov/28276088/
Mouse	Brain	Cerebellum		https://pubmed.ncbi.nlm.nih.gov/28340583/
Mouse	Brain	Hippocampus	HT-22	https://pubmed.ncbi.nlm.nih.gov/12814358/
Mouse	Brain	Striatum		https://pubmed.ncbi.nlm.nih.gov/24952361/
Mouse	Colon			https://pubmed.ncbi.nlm.nih.gov/28257879/
Mouse	Colon	Tissue section and cultured cells	RAW 264	https://pubmed.ncbi.nlm.nih.gov/23188093/
Mouse	Heart	Cardiomyocytes		https://pubmed.ncbi.nlm.nih.gov/28400318/
Mouse	Intestine	Small Intestinal mucosa		https://pubmed.ncbi.nlm.nih.gov/23980531/

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.

Mouse	Kidney	Cortex		https://pubmed.ncbi.nlm.nih.gov/12675844/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/30108504/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/29799837/
Mouse	Liver	Cholangiocarcinoma		https://pubmed.ncbi.nlm.nih.gov/26726846/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/30088830/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/29309760/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/31261663/
Pig	Brain	Hematoma		https://pubmed.ncbi.nlm.nih.gov/27125525/
Rabbit	Eye	Eyeball		https://pubmed.ncbi.nlm.nih.gov/30106111/
Rat	Blood vessel	Femoral vein		https://pubmed.ncbi.nlm.nih.gov/22993073/
Rat	Brain			https://pubmed.ncbi.nlm.nih.gov/22060896/
Rat	Brain			https://pubmed.ncbi.nlm.nih.gov/25054000/
Rat	Brain			https://pubmed.ncbi.nlm.nih.gov/30633838/
Rat	Brain	Mesencephalic cultures		https://pubmed.ncbi.nlm.nih.gov/29154837/
Rat	Kidney			https://pubmed.ncbi.nlm.nih.gov/11799090/
Rat	Kidney			https://pubmed.ncbi.nlm.nih.gov/32235880/
Rat	Liver			https://pubmed.ncbi.nlm.nih.gov/9500714/
Rat	Liver			https://pubmed.ncbi.nlm.nih.gov/27672274/
Rat	Prostate	Tumor		https://pubmed.ncbi.nlm.nih.gov/27550147/
Rat	Prostate	Tumor		https://pubmed.ncbi.nlm.nih.gov/27280718/
Rat	Sciatic nerve	Primary Schwann cells		https://pubmed.ncbi.nlm.nih.gov/31046408/
Rat	Stomach	Gastric mucosa		https://pubmed.ncbi.nlm.nih.gov/18617559/
Sheep	Blood vessels			https://pubmed.ncbi.nlm.nih.gov/9051297/

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Cow	Heart	Aortic Endothelial cells		https://pubmed.ncbi.nlm.nih.gov/10215613/
Cow	Heart	Aortic Endothelial cells		https://pubmed.ncbi.nlm.nih.gov/26825804/
Cow	Heart	Aortic Endothelial cells		https://pubmed.ncbi.nlm.nih.gov/27563876/
Human	Bladder	Carcinoma	T24	https://pubmed.ncbi.nlm.nih.gov/30813460/
Human	Blood	Peripheral blood lymphocytes	PBL	https://pubmed.ncbi.nlm.nih.gov/14642388/
Human	Blood	Peripheral blood mononuclear cells	PBMC	https://pubmed.ncbi.nlm.nih.gov/27694475/
Human	Brain	Glioma	H4	https://pubmed.ncbi.nlm.nih.gov/31586112/
Human	Brain	Microvessels	hCMEC/D3	https://pubmed.ncbi.nlm.nih.gov/26679389/
Human	Brain	Neuroblastoma	SK-N-SH	https://pubmed.ncbi.nlm.nih.gov/30342981/
Human	Breast	Adenocarcinoma	MCF7	https://pubmed.ncbi.nlm.nih.gov/24302665/
Human	Breast	Adenocarcinoma	SKBr3	https://pubmed.ncbi.nlm.nih.gov/24302665/
Human	Colon	Adenocarcinoma	Caco-2	https://pubmed.ncbi.nlm.nih.gov/29791903/

Human	Embryo	Mesenchymal stem cells	hMSCs	https://pubmed.ncbi.nlm.nih.gov/26768768/
Human	Liver	Carcinoma	HepG2	https://pubmed.ncbi.nlm.nih.gov/11707454/
Human	Liver	Cholangiocarcinoma	KKU-100	https://pubmed.ncbi.nlm.nih.gov/22514698/
Human	Liver	Cholangiocarcinoma	KKU-M214	https://pubmed.ncbi.nlm.nih.gov/22514698/
Human	Lung	Alveolar epithelial cells	HPAEpiCs	https://pubmed.ncbi.nlm.nih.gov/32121588/
Human	Prostate	Carcinoma	DU145	https://pubmed.ncbi.nlm.nih.gov/23591596/
Human	Prostate	Carcinoma	PC-3	https://pubmed.ncbi.nlm.nih.gov/23591596/
Human	Skin	Dermal microvascular endothelial cells	HDMEC	https://pubmed.ncbi.nlm.nih.gov/30992483/
Human	Skin	Keratinocytes	HaCaT	https://pubmed.ncbi.nlm.nih.gov/26071936/
Human	Skin	Melanocytes		https://pubmed.ncbi.nlm.nih.gov/26865999/
Human	Trachea	Tracheal smooth muscle	HTSMCs	https://pubmed.ncbi.nlm.nih.gov/31261663/
Human	Umbilical vein	Endothelial cells	HUVEC	https://pubmed.ncbi.nlm.nih.gov/30118148/
Human	Umbilical vein	Endothelial cells	HUVEC	https://pubmed.ncbi.nlm.nih.gov/29610666/
Mouse	Blood	Macrophages	RAW 264.7	https://pubmed.ncbi.nlm.nih.gov/28257879/
Mouse	Blood vessel	Aortic vascular smooth muscle cells	VSMC	https://pubmed.ncbi.nlm.nih.gov/30108504/
Mouse	Blood vessel	Arteriovenous fistula (AVF)		https://pubmed.ncbi.nlm.nih.gov/21697243/
Mouse	Blood vessel	Infra-renal inferior vena cava		https://pubmed.ncbi.nlm.nih.gov/18988794/
Mouse	Blood vessel	Vena cava		https://pubmed.ncbi.nlm.nih.gov/32004074/
Mouse	Bone	Osteoclasts derived from RAW-D cells		https://pubmed.ncbi.nlm.nih.gov/27288918/
Mouse	Bone Marrow	Bone marrow cells		https://pubmed.ncbi.nlm.nih.gov/30018287/
Mouse	Brain			https://pubmed.ncbi.nlm.nih.gov/26724590/
Mouse	Brain	Cerebellum		https://pubmed.ncbi.nlm.nih.gov/28340583/
Mouse	Brain	Cortical neurons		https://pubmed.ncbi.nlm.nih.gov/27112200/
Mouse	Brain	Cortical neurons		https://pubmed.ncbi.nlm.nih.gov/31239473/
Mouse	Brain	Hippocampus	HT-22	https://pubmed.ncbi.nlm.nih.gov/12814358/
Mouse	Brain	Hippocampus		https://pubmed.ncbi.nlm.nih.gov/26698666/
Mouse	Brain	Hippocampus	HT-22	https://pubmed.ncbi.nlm.nih.gov/27939422/
Mouse	Brain	Microglia		https://pubmed.ncbi.nlm.nih.gov/28365213/
Mouse	Brain	Neurons		https://pubmed.ncbi.nlm.nih.gov/27364920/
Mouse	Brain	Striatum		https://pubmed.ncbi.nlm.nih.gov/24952361/
Mouse	Brain	Total brain membrane proteins		https://pubmed.ncbi.nlm.nih.gov/25667042/
Mouse	Colon			https://pubmed.ncbi.nlm.nih.gov/23188093/
Mouse	Colon			https://pubmed.ncbi.nlm.nih.gov/28257879/
Mouse	Heart			https://pubmed.ncbi.nlm.nih.gov/24813593/
Mouse	Heart	Cardiomyoblasts		https://pubmed.ncbi.nlm.nih.gov/31239473/

Mouse	Heart	Cardiomyocytes	HL-1	https://pubmed.ncbi.nlm.nih.gov/23933405/
Mouse	Heart	Cardiomyocytes		https://pubmed.ncbi.nlm.nih.gov/28400318/
Mouse	Heart	Cardiomyocytes		https://pubmed.ncbi.nlm.nih.gov/32376803/
Mouse	Intestine	Peritoneal macrophages		https://pubmed.ncbi.nlm.nih.gov/23980531/
Mouse	Kidney			https://pubmed.ncbi.nlm.nih.gov/12675844/
Mouse	Kidney			https://pubmed.ncbi.nlm.nih.gov/31315267/
Mouse	Kidney	Cortical collecting duct cells	mpkCCD	https://pubmed.ncbi.nlm.nih.gov/31239473/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/26238177/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/27807255/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/30108504/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/29799837/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/32271832/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/26238177/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/30428359/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/29774023/
Mouse	Lung	Alveolar epithelial cells	HPAEpiCs	https://pubmed.ncbi.nlm.nih.gov/30088830/
Mouse	Lung	Alveolar epithelial cells	HPAEpiCs	https://pubmed.ncbi.nlm.nih.gov/29309760/
Mouse	Lung	Alveolar epithelial cells	HPAEpiCs	https://pubmed.ncbi.nlm.nih.gov/30934992/
Mouse	Lung	Microsome		https://pubmed.ncbi.nlm.nih.gov/22490782/
Mouse	Lung	Pleural mesothelium	PMCs	https://pubmed.ncbi.nlm.nih.gov/23399488/
Mouse	Lung	Tracheal smooth muscle cells	HTSMCs	https://pubmed.ncbi.nlm.nih.gov/31261663/
Mouse	Muscle	Quadriceps		https://pubmed.ncbi.nlm.nih.gov/31509445/
Mouse	Spleen			https://pubmed.ncbi.nlm.nih.gov/24813593/
Mouse	Trachea	Tracheal Smooth Muscle; Lysate	HTSMCs	https://pubmed.ncbi.nlm.nih.gov/31261663/
Pig	Brain	Hematoma		https://pubmed.ncbi.nlm.nih.gov/27125525/
Rabbit	Eye	Cornea		https://pubmed.ncbi.nlm.nih.gov/30106111/
Rabbit	Eye	Lens epithelial cells		https://pubmed.ncbi.nlm.nih.gov/10733947/
Rat	Adipose Tissue	Mesenchymal stem cells	ASC	https://pubmed.ncbi.nlm.nih.gov/26836372/
Rat	Adrenal gland	Pheochromocytoma	PC-12	https://pubmed.ncbi.nlm.nih.gov/29154837/
Rat	Blood vessel	Aortic vascular smooth muscle cells	VSMCs	https://pubmed.ncbi.nlm.nih.gov/28542559/
Rat	Blood vessel	Aortic vascular smooth muscle cells	VSMCs	https://pubmed.ncbi.nlm.nih.gov/28274716/
Rat	Blood vessel	Aortic vascular smooth muscle cells	VSMCs	https://pubmed.ncbi.nlm.nih.gov/28257997/
Rat	Brain	Astrocytes		https://pubmed.ncbi.nlm.nih.gov/27005782/
Rat	Brain	Astrocytes	RBA-1	https://pubmed.ncbi.nlm.nih.gov/25148934/
Rat	Brain	Astrocytes	RBA-1	https://pubmed.ncbi.nlm.nih.gov/29209167/
Rat	Brain	Astrocytes	RBA-1	https://pubmed.ncbi.nlm.nih.gov/28497199/
Rat	Brain	Cortical neurons		https://pubmed.ncbi.nlm.nih.gov/27112200/

Rat	Brain	Free-floating sections		https://pubmed.ncbi.nlm.nih.gov/30892744/
Rat	Brain	Neuronal cultures		https://pubmed.ncbi.nlm.nih.gov/31902924/
Rat	Brain	Substantia Nigra		https://pubmed.ncbi.nlm.nih.gov/31832973/
Rat	Heart	Cardiomyoblasts	H9c2	https://pubmed.ncbi.nlm.nih.gov/30986447/
Rat	Heart	Cardiomyocytes		https://pubmed.ncbi.nlm.nih.gov/32376803/
Rat	Kidney		IRPTCs	https://pubmed.ncbi.nlm.nih.gov/11799090/
Rat	Kidney		NRK-52E	https://pubmed.ncbi.nlm.nih.gov/28978536/
Rat	Kidney	Cortex		https://pubmed.ncbi.nlm.nih.gov/26237274/
Rat	Liver			https://pubmed.ncbi.nlm.nih.gov/9500714/
Rat	Liver			https://pubmed.ncbi.nlm.nih.gov/25036135/
Rat	Liver			https://pubmed.ncbi.nlm.nih.gov/27672274/
Rat	Sciatic nerve	Primary Schwann cells		https://pubmed.ncbi.nlm.nih.gov/31046408/
Rat	Stomach	Gastric mucosa		https://pubmed.ncbi.nlm.nih.gov/18617559/
Sheep	Blood vessel	Aorta		https://pubmed.ncbi.nlm.nih.gov/9051297/

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