

HO-1 monoclonal antibody (HO-1-1) (ADI-OSA-110)

ICC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Rat	Brain	Mesencephalic cells		https://pubmed.ncbi.nlm.nih.gov/29154837/

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Kidney	Kidney tissue section		https://pubmed.ncbi.nlm.nih.gov/29636091/
Human	Lung	Lung tissue section		https://pubmed.ncbi.nlm.nih.gov/27526286/
Mouse	Brain	Brain tissue section		https://pubmed.ncbi.nlm.nih.gov/28167258/

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Cow	Kidney	Epithelial cells	MDBK	https://pubmed.ncbi.nlm.nih.gov/31687081/
Human	Blood	Leukemia	THP-1	https://pubmed.ncbi.nlm.nih.gov/30546363/
Human	Blood	Myeloma cell line	INA-6	https://pubmed.ncbi.nlm.nih.gov/27421095/
Human	Blood	PBMC		https://pubmed.ncbi.nlm.nih.gov/14642388/
Human	Blood	PBMC		https://pubmed.ncbi.nlm.nih.gov/28203649/
Human	Blood vessel	Microvascular endothelial cells		https://pubmed.ncbi.nlm.nih.gov/25620054/
Human	Breast	Breast carcinoma	MCF7	https://pubmed.ncbi.nlm.nih.gov/25467193/
Human	Colon	Colorectal adenocarcinoma	SW620	https://pubmed.ncbi.nlm.nih.gov/26585906/
Human	Eye	Retinal pigmented epithelial cells		https://pubmed.ncbi.nlm.nih.gov/26237736/
Human	Kidney	Kidney papilloma	HK-2	https://pubmed.ncbi.nlm.nih.gov/26277392/
Human	Kidney	Kidney papilloma	HK-2	https://pubmed.ncbi.nlm.nih.gov/29039478/
Human	Prostate	Prostate adenocarcinoma	LNCaP	https://pubmed.ncbi.nlm.nih.gov/30077739/
Human	Prostate	Prostate epithelial cells	RWPE-1	https://pubmed.ncbi.nlm.nih.gov/30077739/
Human	Skin	Fibroblast cells		https://pubmed.ncbi.nlm.nih.gov/28947171/
Human	Skin	Keratinocytes	HaCaT	https://pubmed.ncbi.nlm.nih.gov/28142143/
Mouse	Blood	Macrophage cells	RAW264.7	https://pubmed.ncbi.nlm.nih.gov/28820283/

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	Blood vessel	Femoral and mesenteric arteries		https://pubmed.ncbi.nlm.nih.gov/27604527/
Mouse	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/27984168/
Mouse	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/27198537/
Mouse	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/28167258/
Mouse	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/28340583/
Mouse	Brain	Microglial cells	BV-2	https://pubmed.ncbi.nlm.nih.gov/30696869/
Mouse	Embryo	Embryonic stem cells		https://pubmed.ncbi.nlm.nih.gov/29216542/
Mouse	Embryo	Fibroblast cells	3T3-L1	https://pubmed.ncbi.nlm.nih.gov/32540261/
Mouse	Heart	Heart tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/11273999/
Mouse	Kidney	Kidney tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/26232688/
Mouse	Kidney	Kidney tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/29636091/
Mouse	Liver	Liver tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/11493445/
Mouse	Spleen	Spleen tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/29636091/
Rat	Adrenal gland	Pheochromocytoma cell line	PC-12	https://pubmed.ncbi.nlm.nih.gov/26962150/
Rat	Blood vessel	Aorta homogenate		https://pubmed.ncbi.nlm.nih.gov/10904020/
Rat	Blood vessel	Aorta homogenate		https://pubmed.ncbi.nlm.nih.gov/22796671/
Rat	Blood vessel	Aorta homogenate		https://pubmed.ncbi.nlm.nih.gov/25869522/
Rat	Blood vessel	Vascular smooth muscle aortic cells		https://pubmed.ncbi.nlm.nih.gov/30172101/
Rat	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/24723229/
Rat	Brain	Glial cells	OLN-93	https://pubmed.ncbi.nlm.nih.gov/26633426/
Rat	Heart	Cardiac myoblast cells	H9c2	https://pubmed.ncbi.nlm.nih.gov/29689082/
Rat	Heart	Cardiac tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/30428911/
Rat	Intestine	Intestinal epithelial cells	IEC6	https://pubmed.ncbi.nlm.nih.gov/26404472/
Rat	Stomach	Gastric mucosa homogenate		https://pubmed.ncbi.nlm.nih.gov/25263508/
Rat	Testis	Testicular tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/32050098/

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