

Goat anti-mouse IgG F(ab')₂, polyclonal antibody (HRP conjugate) (ADI-SAB-100)

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Kidney	Tissue section		https://pubmed.ncbi.nlm.nih.gov/11238038/

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Coral		Total lysate		https://pubmed.ncbi.nlm.nih.gov/27988888/
Cow	Blood vessel	Aortic endothelial cells		https://pubmed.ncbi.nlm.nih.gov/27800514/
Fish	Erythrocytes	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/25629693/
Fish	Heart	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/25629693/
Fish	Liver	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/25629693/
Fish	Muscle	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/17046286/
Fly	Fat body tissue	Total lysate		https://pubmed.ncbi.nlm.nih.gov/26267495/
Fly	Malpighian tubules	Total lysate		https://pubmed.ncbi.nlm.nih.gov/26267495/
Fly	Oocytes	Total lysate		https://pubmed.ncbi.nlm.nih.gov/26267495/
Fly	Salivary gland	Total lysate		https://pubmed.ncbi.nlm.nih.gov/26267495/
Human	Blood	Peripheral blood mononuclear cells	PBMC	https://pubmed.ncbi.nlm.nih.gov/26021502/
Human	Blood	Serum		https://pubmed.ncbi.nlm.nih.gov/26622380/
Human	Bone marrow	Mesenchymal stem cells		https://pubmed.ncbi.nlm.nih.gov/30186469/
Human	Breast	Carcinoma	MDA-MB-231	https://pubmed.ncbi.nlm.nih.gov/26328244/
Human	Breast	Carcinoma	MDA-MB-435s	https://pubmed.ncbi.nlm.nih.gov/26328244/
Human	Breast	Carcinoma	MCF7	https://pubmed.ncbi.nlm.nih.gov/26328244/
Human	Breast	Carcinoma	MCF7; MCF7-ADR	https://pubmed.ncbi.nlm.nih.gov/26328244/
Human	Breast	Carcinoma	SK-BR-3	https://pubmed.ncbi.nlm.nih.gov/26328244/
Human	Cervix	Carcinoma	HeLa	https://pubmed.ncbi.nlm.nih.gov/27388124/
Human	Cervix; Kidney	Adenocarcinoma; Embryonic Kidney	HeLa; HEK 293T	https://pubmed.ncbi.nlm.nih.gov/32377397/
Human	Gingiva	Gingiva-derived stem cells		https://pubmed.ncbi.nlm.nih.gov/30186469/
Human	Kidney	Embryonic kidney cells	HEK 293	https://pubmed.ncbi.nlm.nih.gov/24291260/

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	Prostate; Kidney	Embryonic kidney cells	HEK 293T	https://pubmed.ncbi.nlm.nih.gov/31178140/
Human	Kidney	Embryonic kidney cells	HEK 293T	https://pubmed.ncbi.nlm.nih.gov/32377397/
Human	Liver	Carcinoma	HepG2	https://pubmed.ncbi.nlm.nih.gov/27388124/
Human	Liver	Carcinoma	HepG2	https://pubmed.ncbi.nlm.nih.gov/28849065/
Human	Liver	Carcinoma	Huh7	https://pubmed.ncbi.nlm.nih.gov/26723877/
Human	Liver	Carcinoma	SNU449	https://pubmed.ncbi.nlm.nih.gov/26723877/
Human	Liver	Stellate cells	LX2	https://pubmed.ncbi.nlm.nih.gov/31374480/
Human	Muscle	Skeletal Muscle	HskMC	https://pubmed.ncbi.nlm.nih.gov/27330287/
Human	Prostate	Adenocarcinoma	DLD1	https://pubmed.ncbi.nlm.nih.gov/31178140/
Human	Prostate	Adeonocarcinoma	DU145	https://pubmed.ncbi.nlm.nih.gov/32319535/
Human	Prostate	Adeonocarcinoma	DU145	https://pubmed.ncbi.nlm.nih.gov/32319589/
Human	Prostate	Adenocarcinoma	HCT-8	https://pubmed.ncbi.nlm.nih.gov/31178140/
Human	Prostate	Adenocarcinoma	HCT-11	https://pubmed.ncbi.nlm.nih.gov/31178140/
Human	Prostate	Adenocarcinoma	HT-29	https://pubmed.ncbi.nlm.nih.gov/31178140/
Human	Prostate	Adeonocarcinoma	LNcaP	https://pubmed.ncbi.nlm.nih.gov/32319535/
Human	Prostate	Adeonocarcinoma	LNcaP	https://pubmed.ncbi.nlm.nih.gov/32319589/
Human	Prostate	Adenocarcinoma	LOVO	https://pubmed.ncbi.nlm.nih.gov/31178140/
Human	Prostate	Stromal cells	WPMY-1	https://pubmed.ncbi.nlm.nih.gov/27325623/
Human	Skin	Keratinocytes	HaCaT	https://pubmed.ncbi.nlm.nih.gov/26648440/
Human	Skin	Dermal endothelium	HMEC	https://pubmed.ncbi.nlm.nih.gov/25620054/
Human	Stomach	Adenocarcinoma	AGS	https://pubmed.ncbi.nlm.nih.gov/28693212/
Human	Umbilical vein	Endothelial cells	HUVEC	https://pubmed.ncbi.nlm.nih.gov/32319555/
Monkey	Kidney	Embryonic kidney cells	COS-7	https://pubmed.ncbi.nlm.nih.gov/24291260/
Monkey	Kidney	Embryonic kidney cells	COS-7	https://pubmed.ncbi.nlm.nih.gov/28322444/
Mouse	Brain	Ischemic region		https://pubmed.ncbi.nlm.nih.gov/27922694/
Mouse	Brain	Microglia	BV2	https://pubmed.ncbi.nlm.nih.gov/30319390/
Mouse	Colon			https://pubmed.ncbi.nlm.nih.gov/28800645/
Mouse	Muscle	Skeletal Myoblast		https://pubmed.ncbi.nlm.nih.gov/25694332/
Mouse	Retina			https://pubmed.ncbi.nlm.nih.gov/30657794/
Mouse	Skin	Melanoma		https://pubmed.ncbi.nlm.nih.gov/31789395/
Mouse	Uterus	Tissue lysate and Endothelial cells		https://pubmed.ncbi.nlm.nih.gov/28849193/
Rabbit	Bone	Articular chondrocytes		https://pubmed.ncbi.nlm.nih.gov/29366902/
Rabbit	Bone	Articular chondrocytes		https://pubmed.ncbi.nlm.nih.gov/30009811/
Rabbit	Bone	Articular chondrocytes		https://pubmed.ncbi.nlm.nih.gov/26851252/
Rabbit	Bone	Articular chondrocytes		https://pubmed.ncbi.nlm.nih.gov/27644154/
Rabbit	Bone	Articular chondrocytes		https://pubmed.ncbi.nlm.nih.gov/27475840/
Rat	Blood vessel	Aortic tissue lysate		https://pubmed.ncbi.nlm.nih.gov/18599595/
Rat	Brain	Corpus callosum		https://pubmed.ncbi.nlm.nih.gov/29199131/
Rat	Heart	Mitochondrial fraction		https://pubmed.ncbi.nlm.nih.gov/11226237/

Rat	Heart	Ventricles		https://pubmed.ncbi.nlm.nih.gov/26176315/
Rat	Liver	Mitochondrial fraction		https://pubmed.ncbi.nlm.nih.gov/11226237/
Rat	Liver	Stellate Cell	HSC-T6	https://pubmed.ncbi.nlm.nih.gov/31173168/
Rat	Muscle	Skeletal Myoblast		https://pubmed.ncbi.nlm.nih.gov/25694332/

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