

α -Gal Epitope (Gal α 1-3Gal β 1-4GlcNAc-R) monoclonal antibody (M86) (ALX-801-090)

ELISA

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Hamster		Cell culture (recombinant protein)	CHO	https://pubmed.ncbi.nlm.nih.gov/16809300/
Chicken	Kidney			https://pubmed.ncbi.nlm.nih.gov/22494361/
Chicken	Muscle			https://pubmed.ncbi.nlm.nih.gov/22494361/
Cow	Eye	Cornea		https://pubmed.ncbi.nlm.nih.gov/30337555/
Cow	Heart			https://pubmed.ncbi.nlm.nih.gov/30337555/
Cow	Kidney			https://pubmed.ncbi.nlm.nih.gov/22494361/
Cow	Kidney			https://pubmed.ncbi.nlm.nih.gov/30337555/
Cow	Muscle			https://pubmed.ncbi.nlm.nih.gov/22494361/
Cow	Skin			https://pubmed.ncbi.nlm.nih.gov/30337555/
Dog	Blood	Serum		https://pubmed.ncbi.nlm.nih.gov/31540167/
Mouse	Heart			https://pubmed.ncbi.nlm.nih.gov/30337628/
Mouse	Kidney			https://pubmed.ncbi.nlm.nih.gov/30337628/
Mouse	Liver			https://pubmed.ncbi.nlm.nih.gov/30337628/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/30337628/
Mouse	Spleen			https://pubmed.ncbi.nlm.nih.gov/30337628/
Pig	Bone			https://pubmed.ncbi.nlm.nih.gov/31342586/
Pig	Eye	Cornea		https://pubmed.ncbi.nlm.nih.gov/30337555/
Pig	Heart			https://pubmed.ncbi.nlm.nih.gov/21118731/
Pig	Heart			https://pubmed.ncbi.nlm.nih.gov/30337555/
Pig	Kidney			https://pubmed.ncbi.nlm.nih.gov/22494361/
Pig	Kidney			https://pubmed.ncbi.nlm.nih.gov/30337555/
Pig	Lung			https://pubmed.ncbi.nlm.nih.gov/21118731/
Pig	Muscle			https://pubmed.ncbi.nlm.nih.gov/22494361/
Pig	Skin			https://pubmed.ncbi.nlm.nih.gov/30337555/

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.

FC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Bacteria				https://pubmed.ncbi.nlm.nih.gov/32344637/
Human	Blood	Leukocytes		https://pubmed.ncbi.nlm.nih.gov/19528506/
Human	Kidney	Renal proximal tubular epithelial cells	RS	https://pubmed.ncbi.nlm.nih.gov/24553072/
Mouse	Skin	Melanoma	BL16	https://pubmed.ncbi.nlm.nih.gov/23740988/
Pig	Embryo	Bone marrow		https://pubmed.ncbi.nlm.nih.gov/12060466/
Pig	Embryo	Kidney		https://pubmed.ncbi.nlm.nih.gov/12060466/
Pig	Heart	Aorta endothelial cells		https://pubmed.ncbi.nlm.nih.gov/22702471/
Pig	Heart	Aorta endothelial cells	PEDSV.15	https://pubmed.ncbi.nlm.nih.gov/15128838/

Functional

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Deer	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/20422448/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/20422448/
Pig	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/20422448/

Immunostaining

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Cow	Heart	Pericardium		https://pubmed.ncbi.nlm.nih.gov/23321301/
Pig	Bone			https://pubmed.ncbi.nlm.nih.gov/31342586/
Pig	Ear	Fibroblasts		https://pubmed.ncbi.nlm.nih.gov/24358349/
Pig	Eye	Retina		https://pubmed.ncbi.nlm.nih.gov/20739475/
Pig	Heart	Aorta endothelial cells		https://pubmed.ncbi.nlm.nih.gov/22702471/
Pig	Heart	Aortic valve		https://pubmed.ncbi.nlm.nih.gov/28763463/
Pig	Heart	Pericardium		https://pubmed.ncbi.nlm.nih.gov/25809726/
Pig	Heart	Pericardium		https://pubmed.ncbi.nlm.nih.gov/29749108/
Pig	Kidney	Tissue Section		https://pubmed.ncbi.nlm.nih.gov/27821126/
Pig	Skin	Derma (ECM and aw reticular dermis)		https://pubmed.ncbi.nlm.nih.gov/20576289/
Rat	Liver	Hepatocytes		https://pubmed.ncbi.nlm.nih.gov/10190554/

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Hamster		Cell culture (recombinant protein)	CHO	https://pubmed.ncbi.nlm.nih.gov/16809300/
Cow	Heart	Pericardium		https://pubmed.ncbi.nlm.nih.gov/27031928/
Cow	Muscle			https://pubmed.ncbi.nlm.nih.gov/24942937/
Cow	Muscle			https://pubmed.ncbi.nlm.nih.gov/24180678/
Horse		Antivenom (MPV, CSL)		https://pubmed.ncbi.nlm.nih.gov/27775867/
Human	Blood	Monocyte-derived dendritic cells	iMDDCs	https://pubmed.ncbi.nlm.nih.gov/30139949/
Pig	Ear	Fibroblasts		https://pubmed.ncbi.nlm.nih.gov/24358349/
Pig	Eye	Cornea		https://pubmed.ncbi.nlm.nih.gov/31284096/