

p62 (human) polyclonal antibody (BML-PW9860)

EM

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Mouse	Brain			https://www.ncbi.nlm.nih.gov/pubmed/29550918

ICC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Breast	Carcinoma	T47D	http://www.ncbi.nlm.nih.gov/pubmed/23227259

IF

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Fish		Zebrafish larvae sections		https://www.ncbi.nlm.nih.gov/pubmed/30718891
Human		Primary fibroblasts		http://www.ncbi.nlm.nih.gov/pubmed/24211199
Human	Brain	Neuroblastoma Cells	SH-SY5Y	https://www.ncbi.nlm.nih.gov/pubmed/32152380
Human	Cervix	Carcinoma	HeLa	http://www.ncbi.nlm.nih.gov/pubmed/23259056
Human	Cervix	Carcinoma	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/23564079
Human	Cervix	Carcinoma	HeLa	http://www.ncbi.nlm.nih.gov/pubmed/26363065
Human	Cervix	Carcinoma	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/28242652
Human	Embryo	Kidney epithelial cells	HEK293	http://www.ncbi.nlm.nih.gov/pubmed/25158237
Human	Embryo	Kidney epithelial cells	HEK293	https://www.ncbi.nlm.nih.gov/pubmed/25311206
Human	Lung	Non-small cell lung carcinoma cells	H1299	https://www.ncbi.nlm.nih.gov/pubmed/25311206
Human	Muscle	Primary myotubes		https://www.ncbi.nlm.nih.gov/pubmed/31341174
Mouse	Brain			https://www.ncbi.nlm.nih.gov/pubmed/29550918
Mouse	Pancreas			https://www.ncbi.nlm.nih.gov/pubmed/32182217
Mouse	Thymus			https://www.ncbi.nlm.nih.gov/pubmed/22719991

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.

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Bone marrow	Bone marrow biopsies		https://www.ncbi.nlm.nih.gov/pubmed/26568842
Human	Brain	Post-mortem autopsies		https://www.ncbi.nlm.nih.gov/pubmed/30591081
Human	Lymph nodes	Post-mortem autopsies		https://www.ncbi.nlm.nih.gov/pubmed/30591081
Human	Stomach	Gastric mucosa biopsies		https://www.ncbi.nlm.nih.gov/pubmed/28983474
Mouse	Brain			https://www.ncbi.nlm.nih.gov/pubmed/29550918
Mouse	Brain			http://www.ncbi.nlm.nih.gov/pubmed/23382946
Mouse	Heart			https://www.ncbi.nlm.nih.gov/pubmed/30771306
Mouse	Liver			http://www.ncbi.nlm.nih.gov/pubmed/23960084
Mouse	Pancreas			http://www.ncbi.nlm.nih.gov/pubmed/22562096
Mouse	Spinal cord			http://www.ncbi.nlm.nih.gov/pubmed/22835604
Rat	Brain			https://www.ncbi.nlm.nih.gov/pubmed/30046966

IP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Cervix	Carcinoma	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/28242652

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Fish		Zebrafish whole larvae homogenates		https://www.ncbi.nlm.nih.gov/pubmed/30718891
Human		Primary fibroblasts	T47D, MCF-7	http://www.ncbi.nlm.nih.gov/pubmed/23227259
Human		Primary fibroblasts		http://www.ncbi.nlm.nih.gov/pubmed/24211199
Human		Primary fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/28378518
Human	Abdomen	Fibroblasts	AG10803	https://www.ncbi.nlm.nih.gov/pubmed/27757122
Human	Blood	B-lymphoblastoid cells	X50-7	http://www.ncbi.nlm.nih.gov/pubmed/27130438
Human	Brain	Neuroblastoma cells	SH-SY5Y	http://www.ncbi.nlm.nih.gov/pubmed/26750654
Human	Brain	Neuroblastoma cells	SH-SY5Y	https://www.ncbi.nlm.nih.gov/pubmed/32152380
Human	Brain	Neuroglioma cells	H4	https://www.ncbi.nlm.nih.gov/pubmed/31586112
Human	Breast	Carcinoma cells	MCF7	http://www.ncbi.nlm.nih.gov/pubmed/23227259
Human	Breast	Carcinoma cells	T47D	http://www.ncbi.nlm.nih.gov/pubmed/23227259
Human	Cervix	Carcinoma cells	HeLa	http://www.ncbi.nlm.nih.gov/pubmed/26363065
Human	Cervix	Carcinoma cells	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/28242652
Human	Colon	Carcinoma cells	Caco-2	http://www.ncbi.nlm.nih.gov/pubmed/26750654
Human	Colon	Carcinoma cells	Caco-2	https://www.ncbi.nlm.nih.gov/pubmed/28475179
Human	Colon	Carcinoma cells	DLD-1	https://www.ncbi.nlm.nih.gov/pubmed/28475179
Human	Colon	Carcinoma cells	HCT-116	https://www.ncbi.nlm.nih.gov/pubmed/28475179
Human	Embryo	Kidney epithelial cells	HEK293	http://www.ncbi.nlm.nih.gov/pubmed/25158237

Human	Embryo	Kidney epithelial cells	HEK293	https://www.ncbi.nlm.nih.gov/pubmed/25311206
Human	Embryo	Kidney epithelial cells	HEK293T	http://www.ncbi.nlm.nih.gov/pubmed/24722279
Human	Lung	Non-small cell lung carcinoma cells	H1299	https://www.ncbi.nlm.nih.gov/pubmed/25311206
Human	Pancreas	Pancreatic cancer cell lines	Panc1	https://www.ncbi.nlm.nih.gov/pubmed/28927939
Human	Pancreas	Pancreatic cancer cell lines	SUIT-2	https://www.ncbi.nlm.nih.gov/pubmed/28927939
Human	Skin	Skin fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/29246495
Human	Stomach	Gastric cancer cell line	AGS	https://www.ncbi.nlm.nih.gov/pubmed/28983474
Mouse		Detection in GST pull-down assays		https://www.ncbi.nlm.nih.gov/pubmed/30771306
Mouse	Bone	Frontal and parietal skull bones		https://www.ncbi.nlm.nih.gov/pubmed/31372669
Mouse	Brain			http://www.ncbi.nlm.nih.gov/pubmed/20947495
Mouse	Brain			https://www.ncbi.nlm.nih.gov/pubmed/29550918
Mouse	Brain	Brain neuron cell line	CATH.a	https://www.ncbi.nlm.nih.gov/pubmed/25902863
Mouse	Brain	Cortical neurons		https://www.ncbi.nlm.nih.gov/pubmed/25216585
Mouse	Brain	Cortical neurons		https://www.ncbi.nlm.nih.gov/pubmed/31076997
Mouse	Brain	Cortical neurons	T4	http://www.ncbi.nlm.nih.gov/pubmed/27112200
Mouse	Brain	Hippocampus		https://www.ncbi.nlm.nih.gov/pubmed/31076997
Mouse	Brain	Primary granule neuron progenitor cells		https://www.ncbi.nlm.nih.gov/pubmed/30301600
Mouse	Brain	Striatum		http://www.ncbi.nlm.nih.gov/pubmed/25583186
Mouse	Breast	Primary Mammary tumor cells		http://www.ncbi.nlm.nih.gov/pubmed/24275666
Mouse	Colon	Colon		https://www.ncbi.nlm.nih.gov/pubmed/29867954
Mouse	Embryo	Fibroblast cells	MEFs	http://www.ncbi.nlm.nih.gov/pubmed/22562096
Mouse	Embryo	Fibroblast cells	MEFs	http://www.ncbi.nlm.nih.gov/pubmed/23508110
Mouse	Embryo	Fibroblast cells	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/25216585
Mouse	Embryo	Fibroblast cells	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/26040720
Mouse	Embryo	Fibroblast cells	MEFs	http://www.ncbi.nlm.nih.gov/pubmed/26750654
Mouse	Embryo	Fibroblast cells	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/28697336
Mouse	Embryo	Fibroblast cells	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/28935812
Mouse	Embryo	Fibroblast cells	MEFs	https://www.ncbi.nlm.nih.gov/pubmed/29514215
Mouse	Embryo	Fibroblast cells	NIH3T3	http://www.ncbi.nlm.nih.gov/pubmed/25880015
Mouse	Eye	Retina		http://www.ncbi.nlm.nih.gov/pubmed/23521856
Mouse	Eye	Retina		http://www.ncbi.nlm.nih.gov/pubmed/27537685
Mouse	Heart			http://www.ncbi.nlm.nih.gov/pubmed/23960084
Mouse	Heart			http://www.ncbi.nlm.nih.gov/pubmed/27246695
Mouse	Heart			https://www.ncbi.nlm.nih.gov/pubmed/30771306
Mouse	Liver			http://www.ncbi.nlm.nih.gov/pubmed/21897364
Mouse	Liver			http://www.ncbi.nlm.nih.gov/pubmed/23960084
Mouse	Liver			https://www.ncbi.nlm.nih.gov/pubmed/28386016
Mouse	Liver	Hepatic stellate cells		https://www.ncbi.nlm.nih.gov/pubmed/22240484
Mouse	Muscle	Skeletal muscle		http://www.ncbi.nlm.nih.gov/pubmed/26244336

Mouse	Skin	Primary keratinocytes		http://www.ncbi.nlm.nih.gov/pubmed/23340736
Mouse	Thymus			https://www.ncbi.nlm.nih.gov/pubmed/22719991
Rat	Brain			http://www.ncbi.nlm.nih.gov/pubmed/23499735
Rat	Brain			http://www.ncbi.nlm.nih.gov/pubmed/23916833
Rat	Brain	Astrocytes		http://www.ncbi.nlm.nih.gov/pubmed/22266730
Rat	Brain	Cortical neurons		http://www.ncbi.nlm.nih.gov/pubmed/27112200
Rat	Liver	Hepatic stellate cells		https://www.ncbi.nlm.nih.gov/pubmed/23485523
Rat	Liver	Hepatic stellate cell line	JS1	https://www.ncbi.nlm.nih.gov/pubmed/23485523

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