

HSP47 monoclonal antibody (M16.10A1) (ADI-SPA-470)

ELISA

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood	Serum		https://www.ncbi.nlm.nih.gov/pubmed/12659832

FC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Breast	Breast cancer cell line	BT474	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MCF7	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MDA-MA-157	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MDA-MB-231	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MDA-MB-468	https://pubmed.ncbi.nlm.nih.gov/32366908/

IF

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Dog	Kidney	Epithelial cells	MDCK	https://www.ncbi.nlm.nih.gov/pubmed/31919357
Dog	Muscle	Primary fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/26004778
Human	Breast	Breast cancer cell line	BT474	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MCF7	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MDA-MA-157	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MDA-MB-231	https://pubmed.ncbi.nlm.nih.gov/32366908/
Human	Breast	Breast cancer cell line	MDA-MB-468	https://pubmed.ncbi.nlm.nih.gov/32366908/
Mouse	Muscle	Primary fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/23234723
Mouse	Muscle	Primary fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/25747978

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Head and neck	FFPE oral squamous cell carcinoma		https://www.ncbi.nlm.nih.gov/pubmed/12710130
Human	Eye	FFPE lacrymal glands		https://www.ncbi.nlm.nih.gov/pubmed/17325149

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	Intestine	FFPE colorectal carcinoma		http://www.ncbi.nlm.nih.gov/pubmed/23265436
Human	Intestine	FFPE colorectal carcinoma		https://www.ncbi.nlm.nih.gov/pubmed/27925182
Human	Lung	FFPE lung tissue section		https://www.ncbi.nlm.nih.gov/pubmed/24583879
Human		FFPE patellar tendons		https://www.ncbi.nlm.nih.gov/pubmed/15314082
Human	Skin	FFPE skin tissue section		https://www.ncbi.nlm.nih.gov/pubmed/28815554
Human	Synovium	FFPE synovial tissue section		https://www.ncbi.nlm.nih.gov/pubmed/27036513
Mouse	Eye	FFPE enucleated eye section		https://www.ncbi.nlm.nih.gov/pubmed/17317786
Mouse	Eye	FFPE enucleated eye section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Intestine	FFPE large intestine section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Intestine	FFPE small intestine section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Lacrimal gland	FFPE lacrimal gland tissue section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Liver	FFPE liver tissue section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Lung	FFPE lung tissue section		https://www.ncbi.nlm.nih.gov/pubmed/24583879
Mouse	Lung	FFPE lung tissue section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Peritoneal tissue	Mesothelial cells		https://www.ncbi.nlm.nih.gov/pubmed/25747978
Mouse	Salivary gland	FFPE salivary gland tissue section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Skin	FFPE skin tissue section		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Rabbit	Eye	FFPE cornea tissue section		https://www.ncbi.nlm.nih.gov/pubmed/27418330
Rat	Blood vessel	FFPE aorta and carotide tissue section		https://www.ncbi.nlm.nih.gov/pubmed/11145931
Rat	Kidney	FFPE kidney tissue section		https://www.ncbi.nlm.nih.gov/pubmed/11141481
Rat	Kidney	FFPE kidney tissue section		https://www.ncbi.nlm.nih.gov/pubmed/19043206
Sheep	Synovium	FFPE joint tissue section		https://www.ncbi.nlm.nih.gov/pubmed/18565987

IP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Connective tissue	HT1080 fibrosarcoma cell line		https://www.ncbi.nlm.nih.gov/pubmed/10329688
Human	Connective tissue	HT1080 fibrosarcoma cell line		https://www.ncbi.nlm.nih.gov/pubmed/12114508
Mouse		Embryonic fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/29351844

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Dog	Muscle	Primary fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/26004778
Human	Adipose tissue			https://www.ncbi.nlm.nih.gov/pubmed/15753122
Human	Blood vessel	Internal thoracic artery	HITB5	https://www.ncbi.nlm.nih.gov/pubmed/12163502
Human	Bone	Osteosarcoma	Saos-2	https://www.ncbi.nlm.nih.gov/pubmed/30840237
Human	Cervical tissue	Cervical cancer cells	HeLa	https://www.ncbi.nlm.nih.gov/pubmed/25742138

Human	Connective tissue	Fibrosarcoma	HT-1080	https://www.ncbi.nlm.nih.gov/pubmed/29276723
Human	Connective tissue	Fibrosarcoma	HT-1080	https://www.ncbi.nlm.nih.gov/pubmed/30840237
Human	Intestine	Colorectal cancer cells		https://www.ncbi.nlm.nih.gov/pubmed/27925182
Human		Rough endoplasmic reticulum fraction		https://www.ncbi.nlm.nih.gov/pubmed/28860186
Human		Recombinant HSP47		https://www.ncbi.nlm.nih.gov/pubmed/16326708
Mouse	Bone	Bone marrow stromal cells		https://www.ncbi.nlm.nih.gov/pubmed/28049723
Mouse	Eye	Enucleated eye lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Intestine	Large intestine lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Intestine	Small intestine lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Lacrimal gland	Lacrimal gland lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Liver	Liver lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Lung	Lung lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Peritoneal tissue	Mesothelial cell lysate		https://www.ncbi.nlm.nih.gov/pubmed/25747978
Mouse	Salivary gland	Salivary gland lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse	Skin	Skin lysate		https://www.ncbi.nlm.nih.gov/pubmed/30307955
Mouse		Embryonic fibroblasts		https://www.ncbi.nlm.nih.gov/pubmed/29351844
Mouse		Fibroblasts	KO-13	https://www.ncbi.nlm.nih.gov/pubmed/26692483
Mouse		Recombinant HSP47		https://www.ncbi.nlm.nih.gov/pubmed/16326708
Rat	Blood vessel	Carotide tissue		https://www.ncbi.nlm.nih.gov/pubmed/11145931

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