

HSP90 monoclonal antibody (AC88) (ADI-SPA-830)

FC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood	Multiple myeloma cell line	OPM-2	https://pubmed.ncbi.nlm.nih.gov/28435516/
Human	Blood	Multiple myeloma cell line	RPML-8226	https://pubmed.ncbi.nlm.nih.gov/28435516/
Human	Skin	Metastatic melanoma	451-LU	https://pubmed.ncbi.nlm.nih.gov/25529535/
Human	Skin	Metastatic melanoma	A375	https://pubmed.ncbi.nlm.nih.gov/25529535/
Human	Skin	Metastatic melanoma	A375	https://pubmed.ncbi.nlm.nih.gov/25762540/
Human	Skin	Metastatic melanoma	M1617	https://pubmed.ncbi.nlm.nih.gov/25529535/
Mouse	Breast	Epithelial carcinoma	EMT6	https://pubmed.ncbi.nlm.nih.gov/30665463/

IF

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Prostate	Prostate adenocarcinoma	PC3	https://pubmed.ncbi.nlm.nih.gov/26921391/
Mouse	Breast	Epithelial carcinoma	EMT6	https://pubmed.ncbi.nlm.nih.gov/30665463/

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Breast	Breast cancer tissue section		https://pubmed.ncbi.nlm.nih.gov/26472975/
Human	Esophagus	Esophageal squamous cell carcinoma section		https://pubmed.ncbi.nlm.nih.gov/28454329/
Rat	Brain	Brain tissue section		https://pubmed.ncbi.nlm.nih.gov/17441507/
Rat	Spinal cord	Spinal cord tissue section		https://pubmed.ncbi.nlm.nih.gov/17441507/

IP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Breast	Breast adenocarcinoma	MDA-MB-231	https://pubmed.ncbi.nlm.nih.gov/23412384/
Human	Cervix	Cervical carcinoma	HeLa	https://pubmed.ncbi.nlm.nih.gov/23412384/
Human	Cervix	Cervical carcinoma	HeLa	https://pubmed.ncbi.nlm.nih.gov/27196783/
Human	Kidney	Embryonic kidney cells	HEK-293	https://pubmed.ncbi.nlm.nih.gov/23412384/
Monkey	Kidney	Fibroblasts	COS-7	https://pubmed.ncbi.nlm.nih.gov/9556607/

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.

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Camel	N/A	Recombinant HSP90 protein		https://pubmed.ncbi.nlm.nih.gov/26234578/
Hamster	Ovary	Ovary cell line	CHO	https://pubmed.ncbi.nlm.nih.gov/31182434/
Hamster	Ovary	Ovary cell line	WCL2	https://pubmed.ncbi.nlm.nih.gov/9195923/
Human	Blood	Leukemia cell supernatant	Jurkat	https://pubmed.ncbi.nlm.nih.gov/30008372/
Human	Blood	Leukemia cell lysate	Jurkat	https://pubmed.ncbi.nlm.nih.gov/30008372/
Human	Blood	Leukemia cell line	K562	https://pubmed.ncbi.nlm.nih.gov/30570667/
Human	Blood	Leukemia cell line	MV4-11	https://pubmed.ncbi.nlm.nih.gov/30570667/
Human	Blood	Leukemia cell line	NB4	https://pubmed.ncbi.nlm.nih.gov/30570667/
Human	Bone marrow	Neuroblastoma	SH-SY5Y	https://pubmed.ncbi.nlm.nih.gov/27163800/
Human	Breast	Breast ductal carcinoma	BT-474	https://pubmed.ncbi.nlm.nih.gov/20111600/
Human	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/26778963/
Human	Brain	Glioblastoma cell line	G55TL	https://pubmed.ncbi.nlm.nih.gov/27488520/
Human	Brain	Glioblastoma cell line	G121	https://pubmed.ncbi.nlm.nih.gov/27488520/
Human	Brain	Glioblastoma cell line	G141	https://pubmed.ncbi.nlm.nih.gov/27488520/
Human	Brain	Glioblastoma cell line	G142	https://pubmed.ncbi.nlm.nih.gov/27488520/
Human	Brain	Primary glioblastoma	NCH644	https://pubmed.ncbi.nlm.nih.gov/27488520/
Human	Breast	Breast carcinoma	MCF7	https://pubmed.ncbi.nlm.nih.gov/18848847/
Human	Breast	Breast adenocarcinoma	MDA-MB-231	https://pubmed.ncbi.nlm.nih.gov/25288732/
Human	Breast	Breast adenocarcinoma	MDA-MB-231	https://pubmed.ncbi.nlm.nih.gov/27788210/
Human	Breast	Breast adenocarcinoma	MDA-MB-231	https://pubmed.ncbi.nlm.nih.gov/27227989/
Human	Cervix	HeLa contaminant	KB	https://pubmed.ncbi.nlm.nih.gov/18848847/
Human	Cervix	Cervical cancer	HeLa	https://pubmed.ncbi.nlm.nih.gov/27788210/
Human	Colon	Healthy colorectal cell line	HCT116	https://pubmed.ncbi.nlm.nih.gov/30487698/
Human	Colon	Healthy colorectal cell line	HCT116	https://pubmed.ncbi.nlm.nih.gov/30570667/
Human	Heart	Cardiomyocytes		https://pubmed.ncbi.nlm.nih.gov/31208058/
Human	Kidney	Embryonic kidney cells	HEK-293	https://pubmed.ncbi.nlm.nih.gov/26778963/
Human	Kidney	Embryonic kidney cells	HEK-293	https://pubmed.ncbi.nlm.nih.gov/27788210/
Human	Kidney	Embryonic kidney cells	HEK-293	https://pubmed.ncbi.nlm.nih.gov/28739485/
Human	Kidney	Embryonic kidney cells	HEK-293	https://pubmed.ncbi.nlm.nih.gov/28986324/
Human	Liver	Hepatocellular carcinoma	HepG2	https://pubmed.ncbi.nlm.nih.gov/32407927/
Human	Prostate	Prostate carcinoma	DU145	https://pubmed.ncbi.nlm.nih.gov/22989880/
Human	Prostate	Prostate adenocarcinoma	LNCaP	https://pubmed.ncbi.nlm.nih.gov/22989880/
Human	Prostate	Prostate adenocarcinoma	PC3	https://pubmed.ncbi.nlm.nih.gov/25288732/
Human	Prostate	Prostate adenocarcinoma	PC3	https://pubmed.ncbi.nlm.nih.gov/26921391/
Human	Prostate	Prostate adenocarcinoma	PC3	https://pubmed.ncbi.nlm.nih.gov/27227989/
Human	Rectum	Rectal cancer cell line	HROC126	https://pubmed.ncbi.nlm.nih.gov/30487698/
Human	Rectum	Rectal cancer cell line	HROC239	https://pubmed.ncbi.nlm.nih.gov/30487698/

Human	Rectum	Rectal cancer cell line	HROC284Met	https://pubmed.ncbi.nlm.nih.gov/30487698/
Human	Skin	Fibroblasts		https://pubmed.ncbi.nlm.nih.gov/32222543/
Human	Umbilical cord	Epithelial cells	HUVECs	https://pubmed.ncbi.nlm.nih.gov/29610666/
Monkey	Kidney	Fibroblasts	COS-7	https://pubmed.ncbi.nlm.nih.gov/9556607/
Mouse	Blood	Plasmacytoid dendritic cells		https://pubmed.ncbi.nlm.nih.gov/25871979/
Mouse	Bone marrow	Bone marrow homogenate		https://pubmed.ncbi.nlm.nih.gov/30022046/
Mouse	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/26778963/
Mouse	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/31596232/
Mouse	Breast	Mammary tumor homogenate		https://pubmed.ncbi.nlm.nih.gov/25659585/
Mouse	Embryo	Embryonic fibroblasts		https://pubmed.ncbi.nlm.nih.gov/18784277/
Mouse	Embryo	Embryonic fibroblasts	NIH-3T3	https://pubmed.ncbi.nlm.nih.gov/26447190/
Mouse	Kidney	Renal adenocarcinoma cell line	Renca	https://pubmed.ncbi.nlm.nih.gov/31932908/
Mouse	Muscle	Myoblasts	C2C12	https://pubmed.ncbi.nlm.nih.gov/28928431/
Mouse	Muscle	Muscle tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/28928431/
Mouse	Pancreas	Islet tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/26397425/
Mouse	Pancreas	Ductal adenocarcinoma cell line	5436	https://pubmed.ncbi.nlm.nih.gov/27838375/
Mouse	Skin	Melanoma	B16-F10	https://pubmed.ncbi.nlm.nih.gov/27683122/
Mouse	Spinal cord	Spinal cord tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/29084947/
Rat	Brain	Brain tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/17441507/
Salmon	Gills	Gills tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/18573527/
Salmon	Liver	Liver tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/18573527/
Salmon	Muscle	Muscle tissue homogenate		https://pubmed.ncbi.nlm.nih.gov/18573527/

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