

Aryl hydrocarbon receptor polyclonal antibody (BML-SA210)

ChIP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Bird	Embryo	Chichen embryo hepatocytes		http://www.ncbi.nlm.nih.gov/pubmed/20876576
Human	Colon	Epithelial cell line	Caco-2	https://www.ncbi.nlm.nih.gov/pubmed/27837168
Mouse	Blood	B-cells		https://www.ncbi.nlm.nih.gov/pubmed/27875245
Mouse	Bone marrow	Macrophages		https://www.ncbi.nlm.nih.gov/pubmed/30962630
Mouse	Brain	primary astrocytes		https://www.ncbi.nlm.nih.gov/pubmed/27158906
Mouse	Colon			https://www.ncbi.nlm.nih.gov/pubmed/30119997
Mouse	Colon	Epithelial cell line	YAMC	https://www.ncbi.nlm.nih.gov/pubmed/27837168
Mouse	Liver			https://www.ncbi.nlm.nih.gov/pubmed/29752288
Mouse	Liver	Hepatoma cells	Hepa-1c1c7	http://www.ncbi.nlm.nih.gov/pubmed/27101761
Mouse	Skin	Skin fibroblast cells	JB6	https://www.ncbi.nlm.nih.gov/pubmed/30578876

ICC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Mouse	Heart	Cardiomyocytes		http://www.ncbi.nlm.nih.gov/pubmed/27163630

IF

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Bladder	Normal urothelial cells		https://www.ncbi.nlm.nih.gov/pubmed/29323757
Human	Brain	Brain with multiple sclerosis lesions		https://www.ncbi.nlm.nih.gov/pubmed/27158906
Human	Brain	Ex vivo glioblastoma samples		https://www.ncbi.nlm.nih.gov/pubmed/30962630
Human	Prostate	Benign prostatic hyperplasia cells	BPH-1	http://www.ncbi.nlm.nih.gov/pubmed/22560882

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.

IP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human		Recombinant protein expressed in yeast		http://www.ncbi.nlm.nih.gov/pubmed/26923060
Mouse	Liver			http://www.ncbi.nlm.nih.gov/pubmed/24150244
Mouse	Liver			http://www.ncbi.nlm.nih.gov/pubmed/25224401

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human		Recombinant protein expressed in yeast		http://www.ncbi.nlm.nih.gov/pubmed/26923060
Human	Bladder	Normal urothelial cells		https://www.ncbi.nlm.nih.gov/pubmed/29323757
Human	Blood	Myeloid leukaemia cells	KBM7	https://www.ncbi.nlm.nih.gov/pubmed/32248720
Human	Cervix	Carcinoma cells	HeLa	http://www.ncbi.nlm.nih.gov/pubmed/24265133
Human	Colon	Epithelial cell line	Caco-2	https://www.ncbi.nlm.nih.gov/pubmed/27837168
Human	Lung	Adenocarcinoma cells	CL1-5	http://www.ncbi.nlm.nih.gov/pubmed/22273977
Human	Lung	Adenocarcinoma cells	NCI-H1355	http://www.ncbi.nlm.nih.gov/pubmed/23229511
Human	Prostate	Benign prostatic hyperplasia cells	BPH-1	http://www.ncbi.nlm.nih.gov/pubmed/22560882
Mouse	Blood	B lymphocytes		https://www.ncbi.nlm.nih.gov/pubmed/27875245
Mouse	Blood	Macrophage cells	RAW264.7	https://www.ncbi.nlm.nih.gov/pubmed/30962630
Mouse	Colon	Colonic crypts		https://www.ncbi.nlm.nih.gov/pubmed/31905023
Mouse	Colon	Epithelial cell line	YAMC	https://www.ncbi.nlm.nih.gov/pubmed/27837168
Mouse	Liver			http://www.ncbi.nlm.nih.gov/pubmed/24150244
Mouse	Liver	Hepatoma cells	Hepa-1c1c7	http://www.ncbi.nlm.nih.gov/pubmed/24265133
Mouse	Liver	Hepatoma cells	Hepa-1c1c7	https://www.ncbi.nlm.nih.gov/pubmed/27105554
Mouse	Ovaries	Primary granulosa cells		http://www.ncbi.nlm.nih.gov/pubmed/25500125
Mouse	Ovaries	Ovary antral follicles		http://www.ncbi.nlm.nih.gov/pubmed/25500125
Rat		Uterus		http://www.ncbi.nlm.nih.gov/pubmed/22127542

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