

Ceramide monoclonal antibody (MID 15B4) (ALX-804-196)

Dot Blot

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Skin	Stratum corneum		https://pubmed.ncbi.nlm.nih.gov/28295641/
Mouse	Brain	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/28706199/

FC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/15130896/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/21923134/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/26584281/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/26021265/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27442249/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27960155/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27978526/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27544243/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27889774/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27050423/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27889758/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27855368/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27197532/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27394133/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/26824457/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/27395049/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/26937949/
Human	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/29035889/
Human	Blood	Lymphocytes		https://pubmed.ncbi.nlm.nih.gov/11279185/
Human	Blood	Lymphocytes		https://pubmed.ncbi.nlm.nih.gov/11751974/

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	Blood	Lymphocytes		https://pubmed.ncbi.nlm.nih.gov/12011072/
Human	Blood	Monocytes		https://pubmed.ncbi.nlm.nih.gov/27870891/
Human	Blood	T Cells		https://pubmed.ncbi.nlm.nih.gov/19834551/
Human	Cervix	Adenocarcinoma	HeLa	https://pubmed.ncbi.nlm.nih.gov/21406729/
Human	Cervix	Adenocarcinoma	HeLa	https://pubmed.ncbi.nlm.nih.gov/30389666/
Human	Prostate	Adenocarcinoma	DU-145	https://pubmed.ncbi.nlm.nih.gov/21406729/
Human	Prostate	Adenocarcinoma	PPC-1	https://pubmed.ncbi.nlm.nih.gov/21406729/
Mouse	Blood	Erythrocytes		https://pubmed.ncbi.nlm.nih.gov/14615798/
Mouse	Blood	Platelet		https://pubmed.ncbi.nlm.nih.gov/30187671/
Mouse	Neck	Squamous cell carcinoma		https://pubmed.ncbi.nlm.nih.gov/22354109/

Functional

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
		Lipids		https://pubmed.ncbi.nlm.nih.gov/12454264/
Human	Blood	Lymphocytes		https://pubmed.ncbi.nlm.nih.gov/11279185/
Human	Colon			https://pubmed.ncbi.nlm.nih.gov/27702691/
Mouse	Lung	Edema		https://pubmed.ncbi.nlm.nih.gov/14704790/

Immunostaining

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Cow	Heart	Aortic Endothelial Cells	BAEC	https://pubmed.ncbi.nlm.nih.gov/27702691/
Guinea pig	Colon			https://pubmed.ncbi.nlm.nih.gov/30389666/
Human	Brain	Glioma	U-373 MG	https://pubmed.ncbi.nlm.nih.gov/31788096/
Human	Blood	Leukemia	Jurkat	https://pubmed.ncbi.nlm.nih.gov/12767914/
Human	Blood	Leukemia	Jurkat	https://pubmed.ncbi.nlm.nih.gov/15849201/
Human	Blood	Lymphocytes		https://pubmed.ncbi.nlm.nih.gov/11279185/
Human	Blood	Lymphocytes		https://pubmed.ncbi.nlm.nih.gov/11751974/
Human	Blood	Lymphocytes		https://pubmed.ncbi.nlm.nih.gov/12011072/
Human	Blood	Monocytes		https://pubmed.ncbi.nlm.nih.gov/27870891/
Human	Blood	T cells		https://pubmed.ncbi.nlm.nih.gov/19834551/
Human	Breast	Adenocarcinoma	MCF7	https://pubmed.ncbi.nlm.nih.gov/21538359/
Human	Cervix	Adenocarcinoma	HeLa	https://pubmed.ncbi.nlm.nih.gov/29507116/
Human	Cervix	Adenocarcinoma	HeLa	https://pubmed.ncbi.nlm.nih.gov/30389666/
Human	Larynx	Carcinoma		https://pubmed.ncbi.nlm.nih.gov/15023837/
Human	Lung			https://pubmed.ncbi.nlm.nih.gov/30500419/
Human	Lung	Fibroblasts		https://pubmed.ncbi.nlm.nih.gov/12563314/
Human	Lung	Carcinoma	H157	https://pubmed.ncbi.nlm.nih.gov/22764099/
Human	Neck	Squamous cell carcinoma	SCC-17B	https://pubmed.ncbi.nlm.nih.gov/25635908/
Human	Neck	Squamous cell carcinoma	SCC-17B	https://pubmed.ncbi.nlm.nih.gov/25739041/

Human	Trachea	Epithelial cells		https://pubmed.ncbi.nlm.nih.gov/12563314/
Human	Umbilical Vein	Endothelial cells	HUVEC	https://pubmed.ncbi.nlm.nih.gov/26909363/
Mouse	Blood vessel	Aorta		https://pubmed.ncbi.nlm.nih.gov/31743567/
Mouse	Bone Marrow	Mesenchymal stromal cells	MSCs	https://pubmed.ncbi.nlm.nih.gov/26857229/
Mouse	Brain			https://pubmed.ncbi.nlm.nih.gov/21615590/
Mouse	Brain			https://pubmed.ncbi.nlm.nih.gov/25270989/
Mouse	Brain			https://pubmed.ncbi.nlm.nih.gov/28706199/
Mouse	Embryo	Inner cell mass	J1-ES	https://pubmed.ncbi.nlm.nih.gov/12885759/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/26487756/
Mouse	Lung			https://pubmed.ncbi.nlm.nih.gov/30500419/
Mouse	Lung	Fibroblasts		https://pubmed.ncbi.nlm.nih.gov/12563314/
Mouse	Prostate	Tumor (Xenograft)	PC-3	https://pubmed.ncbi.nlm.nih.gov/23643061/
Mouse	Prostate	Tumor (Xenograft)	PC-3	https://pubmed.ncbi.nlm.nih.gov/25051356/
Mouse	Skin	Epidermis		https://pubmed.ncbi.nlm.nih.gov/31564077/
Mouse	Trachea	Epithelial cells		https://pubmed.ncbi.nlm.nih.gov/12563314/
Pig	Ovary	Oocyte		https://pubmed.ncbi.nlm.nih.gov/29385411/
Pig	Ovary	Oocyte-granulosa cell complexes	OGCs	https://pubmed.ncbi.nlm.nih.gov/31518729/
Rat	Heart	Cardiomyoblasts	H9C2	https://pubmed.ncbi.nlm.nih.gov/25529920/
Rat	Heart	Left ventricle cells		https://pubmed.ncbi.nlm.nih.gov/29309922/
Rat	Skin			https://pubmed.ncbi.nlm.nih.gov/30368923/

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Mouse	Brain	Tissue lysate		https://pubmed.ncbi.nlm.nih.gov/25270989/

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