

IL-33 monoclonal antibody (Nessy-1) (ALX-804-840)

ELISA/ELISPOT

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Mouse	Peritoneum	Macrophages	F4/80	https://pubmed.ncbi.nlm.nih.gov/19933859/

ICC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Mouse	Bone marrow	Mast cells		https://pubmed.ncbi.nlm.nih.gov/19933859/

IF

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood vessel	Carotid tissue sections		https://pubmed.ncbi.nlm.nih.gov/21737781/
Human	Blood vessel	Endothelial cells	HUVECs	https://pubmed.ncbi.nlm.nih.gov/18787100/
Human	Blood vessel	Endothelial cells	HUVECs	https://pubmed.ncbi.nlm.nih.gov/27694941/
Human	Blood vessel	Temporal artery		https://pubmed.ncbi.nlm.nih.gov/32286393/
Human	Breast	Ductal mammary carcinoma sections		https://pubmed.ncbi.nlm.nih.gov/18787100/
Human	Colon	Adenocarcinoma sections		https://pubmed.ncbi.nlm.nih.gov/18787100/
Human	Colon	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/27748438/
Human	Colon	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/28249897/
Human	Kidney	Renal clear cell carcinoma sections		https://pubmed.ncbi.nlm.nih.gov/18787100/
Human	Kidney	Renal papillary carcinoma sections		https://pubmed.ncbi.nlm.nih.gov/18787100/
Human	Lung	Bronchoalveolar lavage cells		https://pubmed.ncbi.nlm.nih.gov/30973786/
Human	Lung	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/18787100/
Human	Skin	Keratinocytes		https://pubmed.ncbi.nlm.nih.gov/25425229/
Human	Skin	Keratinocytes		https://pubmed.ncbi.nlm.nih.gov/33011245/
Human	Skin	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/18787100/

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	Small intestine	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/18787100/
Mouse		Fibroblasts		https://pubmed.ncbi.nlm.nih.gov/22215666/
Mouse	Eye	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/27701734/
Mouse	Kidney	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/21949094/
Rat	Brain	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/27022724/
Rat	Skin	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/18787100/

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Colon	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/27748438/
Human	Colon	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/26679377/
Human	Kidney	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/29763892/
Human	Lung	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/19801525/
Human	Lung	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/19841166/
Human	Lung	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/23219168/
Human	Lung	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/25003325/
Human	Ovary	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/31322193/
Human	Tendon	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/25857925/
Mouse	Lung	Tissue sections		https://pubmed.ncbi.nlm.nih.gov/21422152/

IP

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood vessel	Lung arterial endothelial cells		https://pubmed.ncbi.nlm.nih.gov/25003325/
Human	Blood vessel	Lung microvascular endothelial cells		https://pubmed.ncbi.nlm.nih.gov/25003325/
Mouse	Embryo	Fibroblasts	MEFs	https://pubmed.ncbi.nlm.nih.gov/21734074/
Mouse	Peritoneum	Macrophages	F4/80	https://pubmed.ncbi.nlm.nih.gov/19933859/

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human		Recombinant IL-33 protein		https://pubmed.ncbi.nlm.nih.gov/19439663/
Human	Blood vessel	Lung arterial endothelial cells		https://pubmed.ncbi.nlm.nih.gov/25003325/
Human	Blood vessel	Lung microvascular endothelial cells		https://pubmed.ncbi.nlm.nih.gov/25003325/
Human	Blood vessel	Umbilical vein endothelial cells	HUVECs	https://pubmed.ncbi.nlm.nih.gov/18787100/

Human	Blood vessel	Umbilical vein endothelial cells	HUVECs	https://pubmed.ncbi.nlm.nih.gov/27694941/
Human	Blood vessel	Umbilical vein endothelial cells	HUVECs	https://pubmed.ncbi.nlm.nih.gov/28715472/
Human	Colon	Carcinoma	HCT116	https://pubmed.ncbi.nlm.nih.gov/28715472/
Human	Colon	Carcinoma	LS513	https://pubmed.ncbi.nlm.nih.gov/28715472/
Human	Embryo	Kidney cells	293T	https://pubmed.ncbi.nlm.nih.gov/19465481/
Human	Embryo	Kidney cells	293T	https://pubmed.ncbi.nlm.nih.gov/28715472/
Human	Muscle	Bronchial smooth muscle cells	BSMCs	https://pubmed.ncbi.nlm.nih.gov/33043044/
Mouse		Recombinant IL-33 protein		https://pubmed.ncbi.nlm.nih.gov/19933859/
Mouse	Brain	Oligodendrocytes	OPC	https://pubmed.ncbi.nlm.nih.gov/31291887/
Mouse	Colon	Carcinoma	CT26	https://pubmed.ncbi.nlm.nih.gov/26679377/
Mouse	Eye	Tissue extract		https://pubmed.ncbi.nlm.nih.gov/27701734/
Mouse	Kidney	Tissue extract		https://pubmed.ncbi.nlm.nih.gov/21949094/
Mouse	Lung	Tissue extract		https://pubmed.ncbi.nlm.nih.gov/19933859/
Mouse	Lung	Tissue extract		https://pubmed.ncbi.nlm.nih.gov/22331917/
Mouse	Lung	Tissue extract		https://pubmed.ncbi.nlm.nih.gov/28666799/
Mouse	Ovary	Tissue extract		https://pubmed.ncbi.nlm.nih.gov/24860190/
Mouse	Ovary	Cancer cells	ID8	https://pubmed.ncbi.nlm.nih.gov/31322193/
Mouse	Spinal cord	Tissue extract		https://pubmed.ncbi.nlm.nih.gov/26363151/
Rat	Brain	Oligodendrocytes	OPC	https://pubmed.ncbi.nlm.nih.gov/27022724/
Rat	Brain	Oligodendrocytes	OPC	https://pubmed.ncbi.nlm.nih.gov/31291887/

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