

Human IgG1 Isotype Control (ALX-804-133)

FC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Hamster	Ovary	Epithelial cell	CHO-K1	https://www.ncbi.nlm.nih.gov/pubmed/26914241
Human	Blood	CAR T cells		https://www.ncbi.nlm.nih.gov/pubmed/25358763
Human	Blood	Leukemia	K562	https://www.ncbi.nlm.nih.gov/pubmed/23530123
Human	Kidney	Embryonic kidney cell	293F	https://www.ncbi.nlm.nih.gov/pubmed/19264769
Human	Kidney	Embryonic kidney cell	293T	https://www.ncbi.nlm.nih.gov/pubmed/19264769
Mouse	Colon	Colon carcinoma cell	CT26	https://www.ncbi.nlm.nih.gov/pubmed/27653549

Function

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood	PBMCs		https://www.ncbi.nlm.nih.gov/pubmed/23894722
Human	Blood	Leukemia	THP-1	https://www.ncbi.nlm.nih.gov/pubmed/22937039
Human	Colon	Fibroblasts	CCD-18Co	https://www.ncbi.nlm.nih.gov/pubmed/22937039
Human		Fibroblasts	Ko77	https://www.ncbi.nlm.nih.gov/pubmed/22937039
Mouse	Skin	T cells		https://www.ncbi.nlm.nih.gov/pubmed/26048148

IHC

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Colon	FFPE colorectal cancer tissue section		https://www.ncbi.nlm.nih.gov/pubmed/21197415
Human	Lung	FFPE lung cancer tissue section		https://www.ncbi.nlm.nih.gov/pubmed/21197415
Human	Pancreas	FFPE pancreatic cancer tissue section		https://www.ncbi.nlm.nih.gov/pubmed/21197415

In vivo

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Mouse			SCID	https://www.ncbi.nlm.nih.gov/pubmed/17431103

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

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.