

## FLIP (human) monoclonal antibody (7F10) (ALX-804-961)

WB

SPECIES	TISSUE	SAMPLE TYPE	NAME	REFERENCES
Human	Blood	B cell lymphoma cell line	Bjab	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26467384">https://www.ncbi.nlm.nih.gov/pubmed/26467384</a>
Human	Blood	B cell lymphoma cell line	OCI-Ly7	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26467384">https://www.ncbi.nlm.nih.gov/pubmed/26467384</a>
Human	Bone	Osteosarcoma cell line	U2OS	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27210546">https://www.ncbi.nlm.nih.gov/pubmed/27210546</a>
Human	Brain	Glioblastoma cell line	M059J	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30696705">https://www.ncbi.nlm.nih.gov/pubmed/30696705</a>
Human	Breast	Breast carcinoma cell line	BT-20	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29434473">https://www.ncbi.nlm.nih.gov/pubmed/29434473</a>
Human	Breast	Breast carcinoma cell line	MCF-7	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29434473">https://www.ncbi.nlm.nih.gov/pubmed/29434473</a>
Human	Breast	Breast carcinoma cell line	MCF-7	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29363524">https://www.ncbi.nlm.nih.gov/pubmed/29363524</a>
Human	Breast	Breast carcinoma cell line	MDA-MB-231	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30696705">https://www.ncbi.nlm.nih.gov/pubmed/30696705</a>
Human	Breast	Breast carcinoma cell line	T47D	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29363524">https://www.ncbi.nlm.nih.gov/pubmed/29363524</a>
Human	Cervix	Cervical cancer cell line	HeLa	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27062399">https://www.ncbi.nlm.nih.gov/pubmed/27062399</a>
Human	Cervix	Cervical cancer cell line	HeLa	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28212752">https://www.ncbi.nlm.nih.gov/pubmed/28212752</a>
Human	Cervix	Cervical cancer cell line	HeLa	<a href="https://www.ncbi.nlm.nih.gov/pubmed/32027418">https://www.ncbi.nlm.nih.gov/pubmed/32027418</a>
Human	Colon	Colorectal adenocarcinoma cell line	HT29	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28106882">https://www.ncbi.nlm.nih.gov/pubmed/28106882</a>
Human	Colon	Colon carcinoma cell line	SW620	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29254146">https://www.ncbi.nlm.nih.gov/pubmed/29254146</a>
Human	Kidney	Kidney carcinoma cell line	293FT	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31046799">https://www.ncbi.nlm.nih.gov/pubmed/31046799</a>
Human	Kidney	Kidney carcinoma cell line	A498	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30008897">https://www.ncbi.nlm.nih.gov/pubmed/30008897</a>
Human	Kidney	Renal cell carcinoma cell line	Caki	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27528031">https://www.ncbi.nlm.nih.gov/pubmed/27528031</a>
Human	Kidney	Renal cell carcinoma cell line	Caki	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26725939">https://www.ncbi.nlm.nih.gov/pubmed/26725939</a>
Human	Kidney	Renal cell carcinoma cell line	Caki	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29565452">https://www.ncbi.nlm.nih.gov/pubmed/29565452</a>
Human	Kidney	Renal cell carcinoma cell line	Caki	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31624721">https://www.ncbi.nlm.nih.gov/pubmed/31624721</a>
Human	Lung	Lung adenocarcinoma cell line	A549	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25682199">https://www.ncbi.nlm.nih.gov/pubmed/25682199</a>
Human	Lung	Lung adenocarcinoma cell line	A549	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26448608">https://www.ncbi.nlm.nih.gov/pubmed/26448608</a>
Human	Lung	Lung adenocarcinoma cell line	A549	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31046799">https://www.ncbi.nlm.nih.gov/pubmed/31046799</a>
Human	Lung	Lung carcinoma cell line	Calu-1	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31046799">https://www.ncbi.nlm.nih.gov/pubmed/31046799</a>

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](http://www.enzolifesciences.com)

For Research Only. Not for Use in Diagnostic Procedures.

Human	Lung	Lung carcinoma cell line	H1299	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30736843">https://www.ncbi.nlm.nih.gov/pubmed/30736843</a>
Human	Lung	Lung carcinoma cell line	H1299	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31046799">https://www.ncbi.nlm.nih.gov/pubmed/31046799</a>
Human	Lung	Lung carcinoma cell line	H157	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30736843">https://www.ncbi.nlm.nih.gov/pubmed/30736843</a>
Human	Lung	Lung carcinoma cell line	H157	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31046799">https://www.ncbi.nlm.nih.gov/pubmed/31046799</a>
Human	Lung	Lung carcinoma cell line	A1792	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31046799">https://www.ncbi.nlm.nih.gov/pubmed/31046799</a>
Human	Lung	Lung carcinoma cell line	H226	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29345617">https://www.ncbi.nlm.nih.gov/pubmed/29345617</a>
Human	Lung	Lung carcinoma cell line	H460	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30736843">https://www.ncbi.nlm.nih.gov/pubmed/30736843</a>
Human	Lung	Lung carcinoma cell line	H460	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31046799">https://www.ncbi.nlm.nih.gov/pubmed/31046799</a>
Human	Stomach	Gastric adenocarcinoma cell line	AGS	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24603337">https://www.ncbi.nlm.nih.gov/pubmed/24603337</a>

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