

MultiView[®] Universal Tissue Microarray

Catalog #: ADI-950-020

5 slides, 25 slides and 50 slides

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.



Please read
entire booklet
before
proceeding.



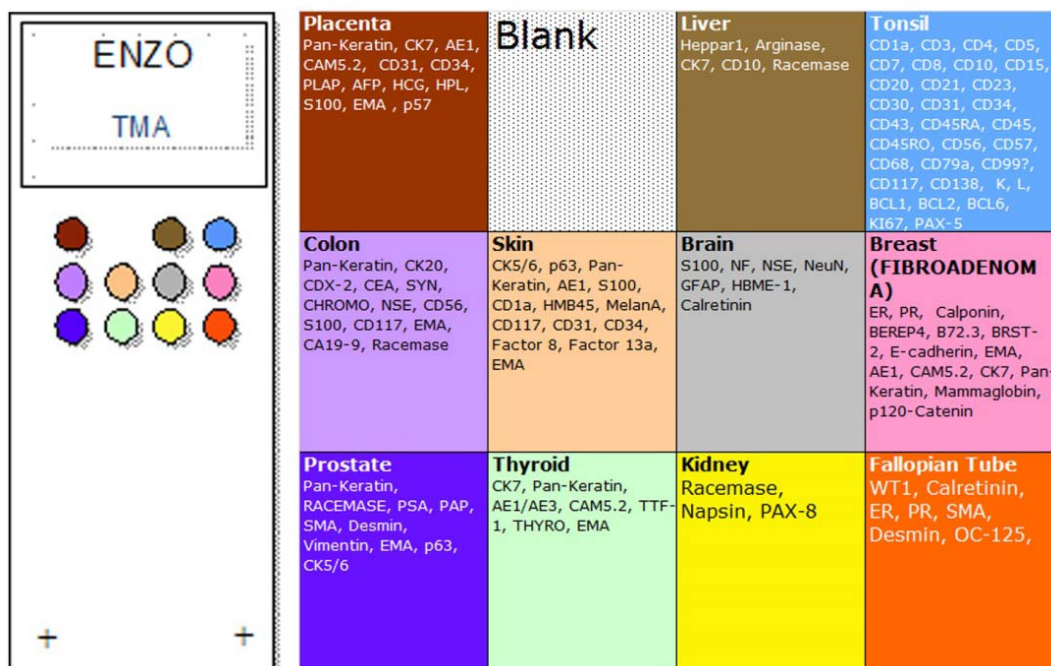
Please contact
Enzo Life
Sciences
Technical
Support if
necessary.

TABLE OF CONTENTS

Description	2
Uses and Advantages	2
Tissue Reactivity Chart	3
Quality Control	8
Handling	8
Storage and Stability	8
Shipping	8
References	9
Contact Information	10

DESCRIPTION

The MultiView[®] Universal Tissue Microarray consists of 11 tissue specimens, including placenta, liver, tonsil, colon, skin, brain, breast (fibroadenoma), prostate, thyroid, kidney and fallopian tube. The samples are fixed in formalin and paraffin embedded, and the cores are approximately 2.0mm in diameter. The sections are cut at 3µm and placed on a SuperFrost Plus slides.



Please refer to the tissue reactivity chart below for details.

USES AND ADVANTAGES

This MultiView[®] Universal Tissue Microarray slide is intended to provide broad-spectrum reactivity in manual and fully automated procedures for:

- Histochemical stains** *example: hematoxylin and eosin (H&E)*
- Special stains** *example: trichrome, PAS*
- Immunohistochemical stains** *example: CK7, CK20, etc.*
- In situ hybridization** *example: 17q*

The use of tissue microarrays for controls has many advantages, including uniformity of slides and improved operational efficiency (with fewer control types to maintain). The selected tissues are designed to cover the majority of commonly used immunohistochemical stains. When used for the optimization or validation of stains, tissue microarrays provide a rapid and relatively inexpensive screening tool. Usual protocols for FFPE (formalin fixed paraffin embedded) tissue on charged slides may be followed.

TISSUE REACTIVITY CHART

	PLACENTA	FALLOPIAN TUBE	BREAST	THYROID	LIVER	KIDNEY	COLON	BRAIN	SKIN	TONSIL	PROSTATE	Count of reactive tissue cores
Androgen Receptor									y		y	2
a-1-A												0
ACTH												0
AE1	y	y	y	y	y	y	y		y	y	y	10
AE1/AE3	y	y	y	y	y	y	y		y	y	y	10
AFP	y											1
ALK-1												0
Annexin A1										y		1
B72.3	y						y					2
Bcl-1			y									1
Bcl-2			y									1
Bcl-6										y		1
BER-EP4		y					y					2
Beta-Catenin							y		y			2
BOB-1										y		1
Breast Mix			y									1
Breast Triple			y									1
BRST-2			y									1
CA19-9					y		y					2
Calcitonin												0
Calponin	y	y	y		y	y	y	y	y	y	y	10
Calretinin		y						y				2
CAM5.2	y	y										2
CD10	y					y						2
CD117	y								y			2
CD138	y									y		2
CD15	y									y		2
CD163	y	y								y		3
CD1a									y	y		2
CD2			y							y		2
CD20										y		1
CD20/CD3										y		1

	PLACENTA	FALLOPIAN TUBE	BREAST	THYROID	LIVER	KIDNEY	COLON	BRAIN	SKIN	TONSIL	PROSTATE	Count of reactive tissue cores
CD21										y		1
CD23										y		1
CD25										y		1
CD3										y		1
CD30										y		1
CD31										y		1
CD34	y	Y	y	y	y	y	y	y	y	y	y	11
CD4										y		1
CD43										y		1
CD45 (LCA)										y		1
CD5										y		1
CD56								y		y		2
CD68										y		1
CD7										y		1
CD79a										y		1
CD8										y		1
CD99/12E7										y		1
CD99/HBA-71										y		1
CDX-2							y					1
CEAm	Y	Y	y	y	y	y	y		y	y	y	10
Chromogranin							y	y			y	3
CK19		Y										1
CK20							y					1
CK5/6		y	y						y	y	y	5
CK7	Y	Y	y	y	y	y			y	y		8
CMV												0
CXCL13										y		1
D2-40	Y	Y	y			y	y		y	y	y	8
DBA44										y		1
Desmin	Y	Y							y	y	y	5
DOG-1			y				y					2
E-Cadherin			y									1
EBV (EBER)												0

	PLACENTA	FALLOPIAN TUBE	BREAST	THYROID	LIVER	KIDNEY	COLON	BRAIN	SKIN	TONSIL	PROSTATE	Count of reactive tissue cores
EGFR	Y	Y										2
EMA	Y	Y	y						Y	Y		5
ER		Y	y									2
Factor VIII	Y	Y	y									3
Factor XIIIa	Y	Y	y									3
Galectin-3			y				y		y	y	y	5
GFAP							y	y				2
Glycophorin	y									y		2
Glypican-3												0
H. pylori												0
HBME-1								y				1
HCG	Y											1
Hepatocyte				y								1
HepBCore												0
HepBsAg												0
HER2												0
HgH												0
HHV-8												0
HMB45									Y			1
HPL	Y											1
HPV HR												0
HPV LR												0
HSV												0
Inhibin												0
INI-1	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
Insulin												0
Kappa										y		1
Ki-67	Y	Y	y							y		4
Lambda										y		1
Lysozyme	Y	Y	y									3
Mammaglobin			y									1
Melan-A									y			1
MITF									y			1

	PLACENTA	FALLOPIAN TUBE	BREAST	THYROID	LIVER	KIDNEY	COLON	BRAIN	SKIN	TONSIL	PROSTATE	Count of reactive tissue cores
MLH1	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
MOC-31		Y	Y				Y					3
MPO	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
MSA												0
MSH2	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
MSH6	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
Myogenin												0
NeuN								y				1
Neurofilament		Y						y				2
NSE							y	y				2
OC-125		Y										1
OCT-2										y		1
OCT-4												0
p120 Catenin			y									1
p16			y				y			y		3
p40								y	y			2
p53			y									1
p57	Y											1
p63		y	y					y	y	y		5
Pan Cytokeratin	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	10
PAPm											y	1
Parvovirus B19												0
PAX-5										y		1
PD1										y		1
PLAP	Y											1
PMS2	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
PR		Y	y									2
Prolactin												0
PSA											y	1
PSAP											y	1
PTH												0
Racemase					Y	y						2
S100	Y		y									2

	PLACENTA	FALLOPIAN TUBE	BREAST	THYROID	LIVER	KIDNEY	COLON	BRAIN	SKIN	TONSIL	PROSTATE	Count of reactive tissue cores
SMA	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
Synaptophysin							Y	Y				2
TdT												0
Thyroglobulin				Y								1
TSH												0
TTF-1				Y								1
Tyrosinase									Y			1
Vimentin	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	Y	11
WT-1		Y										1
Total covered	107											
%	0.78102											

NOTE: Individual results may vary and validation is the responsibility of the end-user.

QUALITY CONTROL

Initial and final sections are stained by H&E and the presence of appropriate tissue is confirmed. The ultimate quality control and validation of the control tissue array will be performed by the end user.

STORAGE AND STABILITY

This product may be stored and maintained at room temperature (20-25 °C). Excessive heating for more than one hour at 60 °C may cause a significant decrease or loss of immunoreactivity. To preserve reactivity with some stains, slides should be stained within 12 weeks of sectioning when stored at room temperature (20-25 °C), although stability studies show excellent reproducibility after 6 months at room temperature (4 stains tested: TTF-1, Pankeratin, ER and CD34).

HANDLING

Standard universal precautions should be taken when handling formalin-fixed paraffin- embedded tissue, including both tissue blocks and slide-mounted tissue. Standard procedures for disposal of glass products should be observed with the disposition of this product after use.

REFERENCES

1. Lundgaard HB, Winther H, Moller K. Excessive section drying of breast cancer tissue prior to deparaffinisation and antigen retrieval causes a loss in HER2 immunoreactivity, *Immunocytochemistry* 2008; 6,119-22.
2. Henwood AF. Effect Of Slide Drying At 80°C On Immunohistochemistry. *J Histotechnology* 2005; 28:45- 46.
3. Karlsson C, Karlsson MG. Effects of Long-Term Storage on the Detection of Proteins, DNA, and mRNA in Tissue Microarray Slides *J Histochem Cytochem* 2011 59: 1113
4. Hewitt, SM, et al, Validation of proteomic-based discovery with tissue microarrays. *Proteomics Clin. Appl.* 2008, 2, 1460–1466.
5. Fedor, HI and DeMarzo AM, Practical Methods for Tissue Microarray Construction, *Methods in Molecular Medicine, Vol. 103: Pancreatic Cancer: Methods and Protocols* Edited by: G. Su © Humana Press Inc., Totowa, NJ, 89-101.

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com

