

MULTIVIEW®

Universal Tissue Microarray

MULTIVIEW® Universal Tissue Microarray with 11 different human tissue specimens for a broad spectrum of immunohistochemical staining procedures.

The MULTIVIEW® Universal Tissue Microarray consists of 11 human 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 slide.

Ordering Information

[Order Online »](#)

ADI-950-020-0025	25 slides
ADI-950-020-0050	50 slides

Manuals, SDS & CofA

[View Online »](#)

- Diverse set of 11 tissues designed to cover the majority of commonly used immunohistochemical stains
- Provides a rapid, inexpensive screening tool for optimization or validation of stains
- Reactivity with over 130 different certified markers provided for each tissue
- High-consistency specimen & slide preparation reduce control slide maintenance
- Suitable for use with common FFPE tissue protocols



ENZO TMA		Placenta Pan-Keratin, CK7, AE1, CAM5.2, CD31, CD34, PLAP, AFP, HCG, HPL, S100, EMA, p57	Blank	Liver HepPar1, Arginase, CK7, CD10, Racemase	Tonsil CD1a, CD3, CD4, CD5, CD7, CD8, CD10, CD11, CD20, CD21, CD23, CD30, CD31, CD34, CD43, CD45RA, CD45, CD45RO, CD56, CD57, CD68, CD79a, CD99, CD117, CD138, X, L, BCL1, BCL2, BCL6, Ki67, PAX-5
		Colon Pan-Keratin, CK20, CDK-2, CEA, SYN, CHROMO, NSE, CD56, S100, CD117, EMA, CA19-9, Racemase	Skin CK5/6, p63, Pan-Keratin, AE1, S100, CD1a, HMB45, MelanA, CD117, CD31, CD34, Factor 8, Factor 13a, EMA	Brain S100, NF, NSE, NeuN, GFAP, HBME-1, Calretinin	Breast (FIBROADENOMA) ER, PR, Calponin, BEREPA, 872.3, BREST-2, E-cadherin, EMA, AE1, CAM5.2, CK7, Pan-Keratin, Mammaglobin, p120-Catenin
Prostate Pan-Keratin, RACEMASE, PSA, PAP, SMA, Desmin, Vimentin, EMA, p63, CK5/6		Thyroid CK7, Pan-Keratin, AE1/AE3, CAM5.2, TTF-1, THYRO, EMA	Kidney Racemase, Napsin, PAX-8	Fallopian Tube WT1, Calretinin, ER, PR, SMA, Desmin, OC-125	

Example H&E stained slide to highlight the different morphologies of the tissues. The slides sold are unstained.

The MULTIVIEW® Universal Tissue Microarray array consists of 11 tissue specimens, including placenta, liver, tonsil, colon, skin, brain, breast (fibroadenoma), prostate, thyroid, kidney and fallopian tube. Please refer to the product manual for details of the tissue reactivity.

Handling & Storage

Handling	To preserve reactivity with some stains, slides should be stained within 12 weeks of sectioning when stored at room temperature (20-25 °C).
Short Term Storage	Ambient
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes

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

- (1) histochemical stains (example: hematoxylin and eosin (H&E));
- (2) special stains (example: trichrome, PAS);
- (3) immunohistochemical stains (example: CK7, CK20) and
- (4) *in situ* hybridization (example: 17q).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)