

FEATURED IMAGE

MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C MaTek's Cells and Media, NHEK-CRY-AD Echo Microscopes, Revolve Normal Human Epidermal Keratinocytes (NHEK-CRY-AD) were cultured on MatTek P35G-1.5-14-C Glass Bottom dishes. Cells were fixed and stained via immunocytochemistry for cytokeratin 19 (red), phalloidin (green), and dapi (blue) and imaged at 10x on the Revolve ECHO microscope.



January

2023

SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
01	02	03	04	05	06	07
New Year's Day						
08	0.9	10	11	12	13	14
		10		12	15	
15	16	17	18	19	20	21
	Martin Luther King Day					
22	23	24	25	26	27	28
29	30	31	01	02	03	04

FEATURED IMAGE

Elly Liao

University of Minnesota MatTek's Glass Bottom 12-well Plate, P12G-1.5-14-F Our lab developed this SHSY5Y neurosphere model to study the molecular mechanisms that are at work in neurite outgrowth. Introduction of mutant proteins, like tau and alpha synuclein, change the cellular microenvironment and influence neurite extension. Using this assay, we can screen and evaluate compounds that promote or rescue neurite growth. For evaluation, the neurospheres were plated on MatTek 12-well plates coated with poly-L-ornithine/laminin and stained with β III-tubulin (red) and DAPI (blue).





SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
29	30	31	01	02	03	04
05	06	07	08	09	10	11 International Day of Women and Girls in Science
12	13	14 Valentine's Day	15	16	17	18
19	20 Presidents' Day	21	22	23	24	25
26	27	28	01	02	03	04

FEATURED IMAGE

Julien Courchet

Inserm / University of Lyon MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C Our ability to interact with the world comes from the delicate interaction between neurons and muscle cells. This picture features one of the first neuron (green)/muscle (red) cocultures I performed. Wandering around with the microscope, I found in the glass bottom dish that spinal cord neurons self-organized into a heart-shaped spheroid. Nuclei were marked with DAPI (blue).





SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
26	27	28	01	02	03	04
05	06	07	08	09	10	11
12	13	14	15	16	17 St. Patrick's Day	18
19	20	21	22	23	24	25
26	27	28	29	30	31	01

FEATURED IMAGE

Nicole Heiberger

Genemarkers MatTek's EpiDermFT™ Model, EFT-400 Skin tissues exposed to HEV light for 2hr each day for up to 5-days, resulted in a significant increase in IL-6 production. IL-6 expression is visible in nuclear as well as in the cytoplasmic regions. The EFT-400 model was used as it has a well-defined stratum corneum, epidermal, and dermal layer separations.





SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
26	27	28	29	30	31	01
02	03	04	05	06	07	08
09	10	11	12	13	14	15
16	17	18	19	20	21	22 Earth Day
23	24	25 National DNA Day	26	27	28	29
30	01	02	03	04	05	06

FEATURED IMAGE

MatTek, Advanced BioMatrix, and Discover Echo Kaitlyn Coen

ECM Select Array Kit Cat # 5170 & MatTek NHBF-CRY

MatTek's Normal Human Bronchial Fibroblasts (NHBF-CRY) stained with Vimentin (red), Phalloidin (green), and DAPI (blue) grown on Advanced Biomatrix ECM Select Array Kit Collagen I/Collagen IV demonstrating successful cell growth after 96hrs.





SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
30	01	02	03	04	05	06
					Cinco De Mayo	
07	08	09	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30	31	01	02	03
	Memorial Day					

FEATURED IMAGE

Geoffrey Williams

Brown University MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C Lily of the valley Convallaria majalis, whole flower fixed, dehydrated into 100% ethanol, then stained with Acridine orange.

Trichome found on a Russian Olive (Elaeagnus angustifolia) flower. Stained with Acridine Orange.



June

2023

SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
28	29	30	31	01	02	03
04	05	06	07	08	09	10
11	12	13	14	15	16	17
18	19	20	21	22	23	24
25	26	27	28	29	30	01

FEATURED IMAGE

Sarah Guest

UC Davis MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C This is a Rhizarian amoeba called Filoreta, which grows and moves in a mechanism similar to how neurons develop their branches. They are multinucleate cells (syncytia) that make a complex branched network as they grow to macroscopic sizes. In this image, a small syncutium has recently fused with its neighbor. These delicate syncytia were grown and immunostained on glass-bottom Mattek dishes. Nuclei = dapi, microtubules = alexa-594 antibody (red), actin = phalloidin 488.





SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
25	26	27	28	29	30	01
02	03	04 Independence Day	05	06	07	08
09	10	11	12	13	14	15
16	17	18	19	20	21	22
23	24	25	26	27	28	29
30	31	01	02	03	04	05

FEATURED IMAGE

Laurens Kimps

KU Leuven MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C Mouse embryonic fibroblast(s) with fluorescently labeled Vinculin cultured on glass substrate. The stained Vinculin reveals the beautiful focal adhesions of the cells, through which they adhere to the glass substrate.





SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
30	31	01	02	03	04	05
06	07	08	09	10	11	12
13	14	15	16	17	18	19
20	21	22	23	24	25	26
27	28	29	30	31	01	02

FEATURED IMAGE

Priya Gohel

Stockholm University MatTek's 35 mm Glass Bottom Dish, P35GC-1.5-14-C Live imaging of sfGFP-tagged protein along with His2Av-RFP during syncytial mitotic divisions in Drosophila embryo resulting in "TulipOsomes" - dividing chromosomes (red) and a DNA-binding transcription factor (green) that look like tulips.



September

2023

SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
27	28	29	30	31	01	02
03	04 Labor Day	05	06	07	08	09
10	11 Patriot Day	12	13	14	15	16
17	18	19	20	21	22	23
24	25	26	27	28	29	30

FEATURED IMAGE

Sergey Pryshchep

Rensselaer Polytechnic institute MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C Three-dimensional confocal image of THP-1 cells migration on Human umbilical vein endothelial cells (HUVEC) monolayer. THP-1 cells marked with 488ViaFluor® Microtubule Stains (green), HUVEC cells marked with DAPI dye (blue) for nucleic acid and 647CF® Dyes-phalloidin(red) for actin filamets.





SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
01	02	03	04	05	06	07
08	09	10	11	12	13	14
45		17	10	10	20	21
15	10	17	18	19	20	21
22	23	24	25	26	27	28
		24	0.1			
29	30	31	01	02	03	04
		Halloween				

FEATURED IMAGE

Discover Echo - A BICO COMPANY

Courtesy of Benedicte Gobert, Dynacure, Illkirch, France. Their mission is Changing the Lives of Patients with Myotubular and Centronuclear Myopathies Human immortalized myoblasts were plated on coated coverslips and differentiated for 4 days in order to obtain multinucleated fibers (green : myosin heavy chain staining, blue: nuclei staining) A Mosaic 4x4 Stich with the Revolution Microscope.



Novemeber

2023

SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
29	30	31	01	02	03	04
05	06	07	08 International STEM Day	09	10	11 Veterans' Day
12	13	14	15	16	17	18
19	20	21	22	23 Thanksgiving	24	25
26	27	28	29	30	01	02

FEATURED IMAGE

Noa Martin

Instituto de Investigacion Sanitaria Hospital Universitario de la Princesa

MatTek's 35 mm Glass Bottom Dish, P35G-1.5-10-C

T lymphocytes are contacting antigen-presenting cells through the recognition of specific antigens. Cytoskeltal components such as microtubules (green) reorganize to bring the centrosome near the contact area, forming an immunological synapse. There, the T cell receptor and CD3 complex (red) accumulate to sustain T cell activation. Intracellular compartment, such as endosomes, contain signaling molecules delivered to the synaptic contact such as LAT (blue).



DEC December

2023

SUNDAY	MONDAY	TUESDAY	WEDNESDAY	THURSDAY	FRIDAY	SATURDAY
26	27	28	29	30	01	02
03	04	05	06	07	08	09
				First Night		
10	11	10	12	of Hanukkah	45	
10	11	12	13	14	15	16
17	18	19	20	21	22	23
24	25	26	27	28	29	30
	Christmas Day	First Night				
		of Kwanzaa				
31	01	02	03	04	05	06
New Year's Eve	New Year's Day					
	rear a bay					

FEATURED IMAGE

Caterina Ciacci

University of Urbino MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C

Representative images of HaCaT Keratinocytes. The image was obtained using a Leica TCS SP5 confocal setup mounted on a Leica DMI6000 CS inverted microscope (Leica Microsystems, Heidelberg, Germany) using a 63 x 1.4 oil objective. HaCaT cells were stained with Alexa Fluor 647 phalloidin for F-actin and with SYBR Green I for nuclei. Cells were grown on MatTek Glass bottom chambers, fixed with 3% paraformaldehyde for 10min and treated with 0.1% Nonidet (NP-40) for cell permeabilization before staining.



CALENDAR

FEATURED IMAGE

MatTek's 35 mm Glass Bottom Dish, P35G-1.5-14-C MaTek's Cells and Media, NHDF-CRY-AD Normal Human Dermal Fibroblasts (NHDF-CRY-AD) were cultured on MatTek P35G-1.5-14-C Glass Bottom dishes. Cells were fixed and stained via immunocytochemistry for vimentin (red), phalloidin (green), and dapi (blue) and imaged at 100x on the Olympus fv100 confocal microscope.