



H I G H C O N T E N T I M A G I N G G A L L E R Y

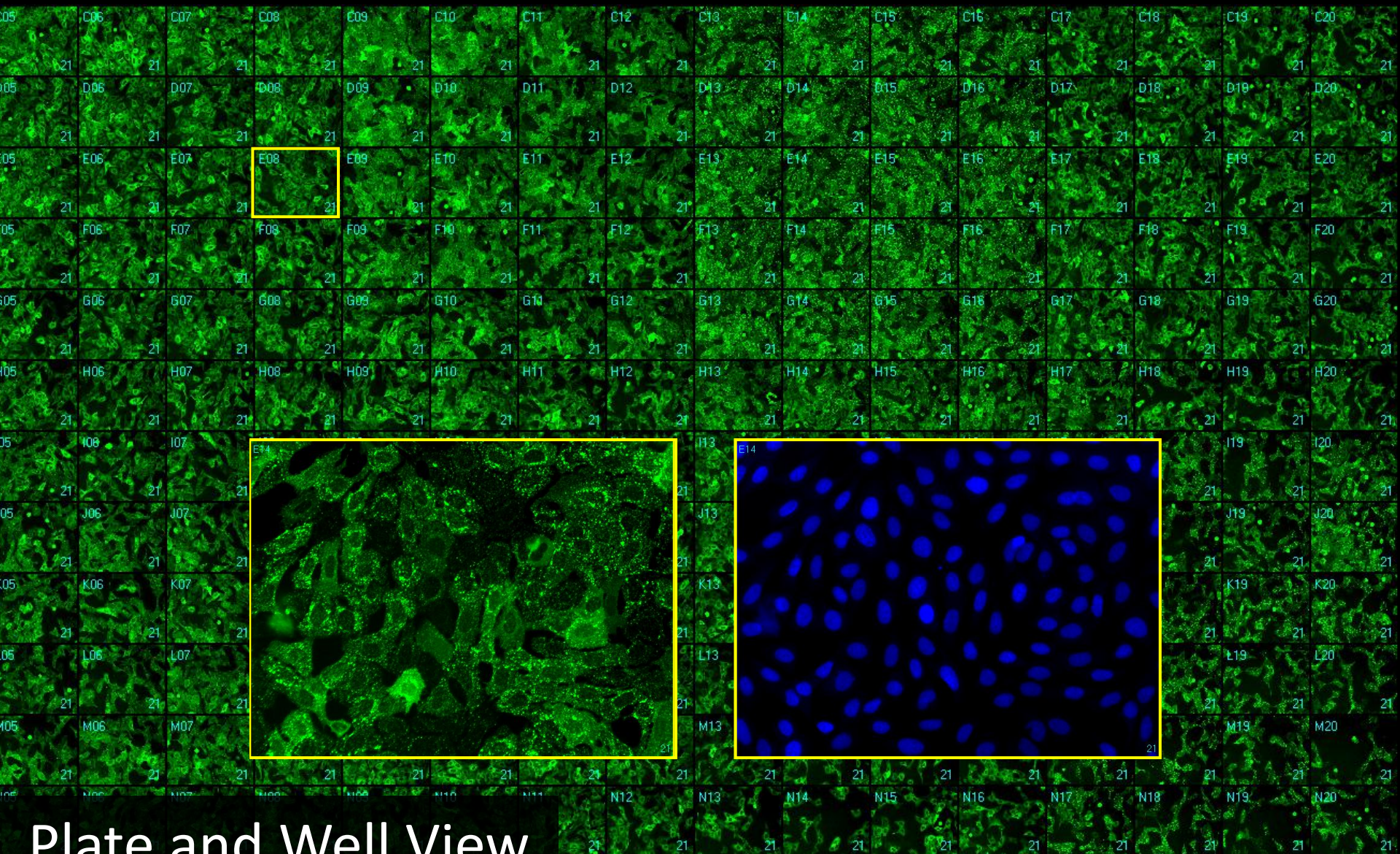
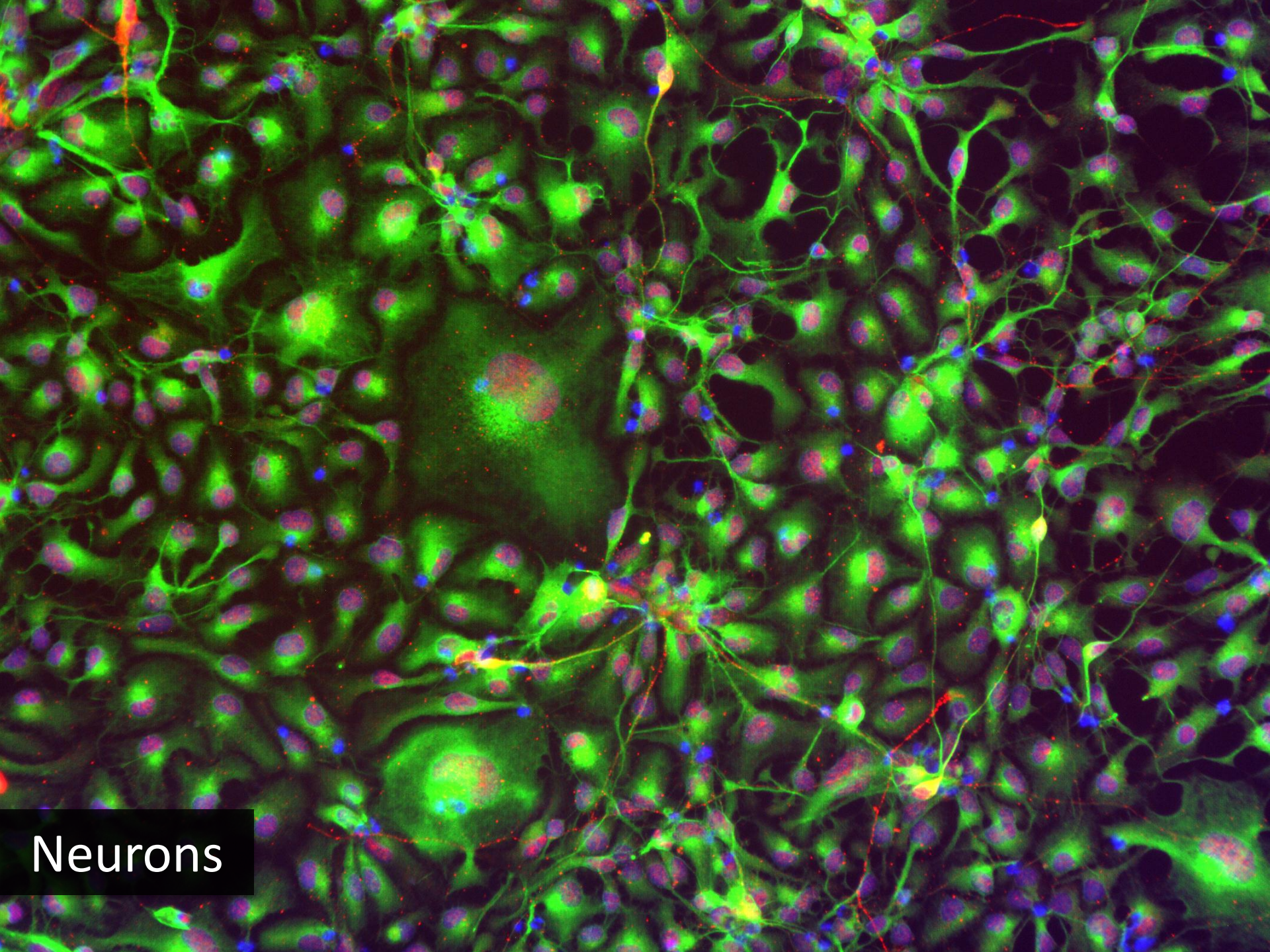
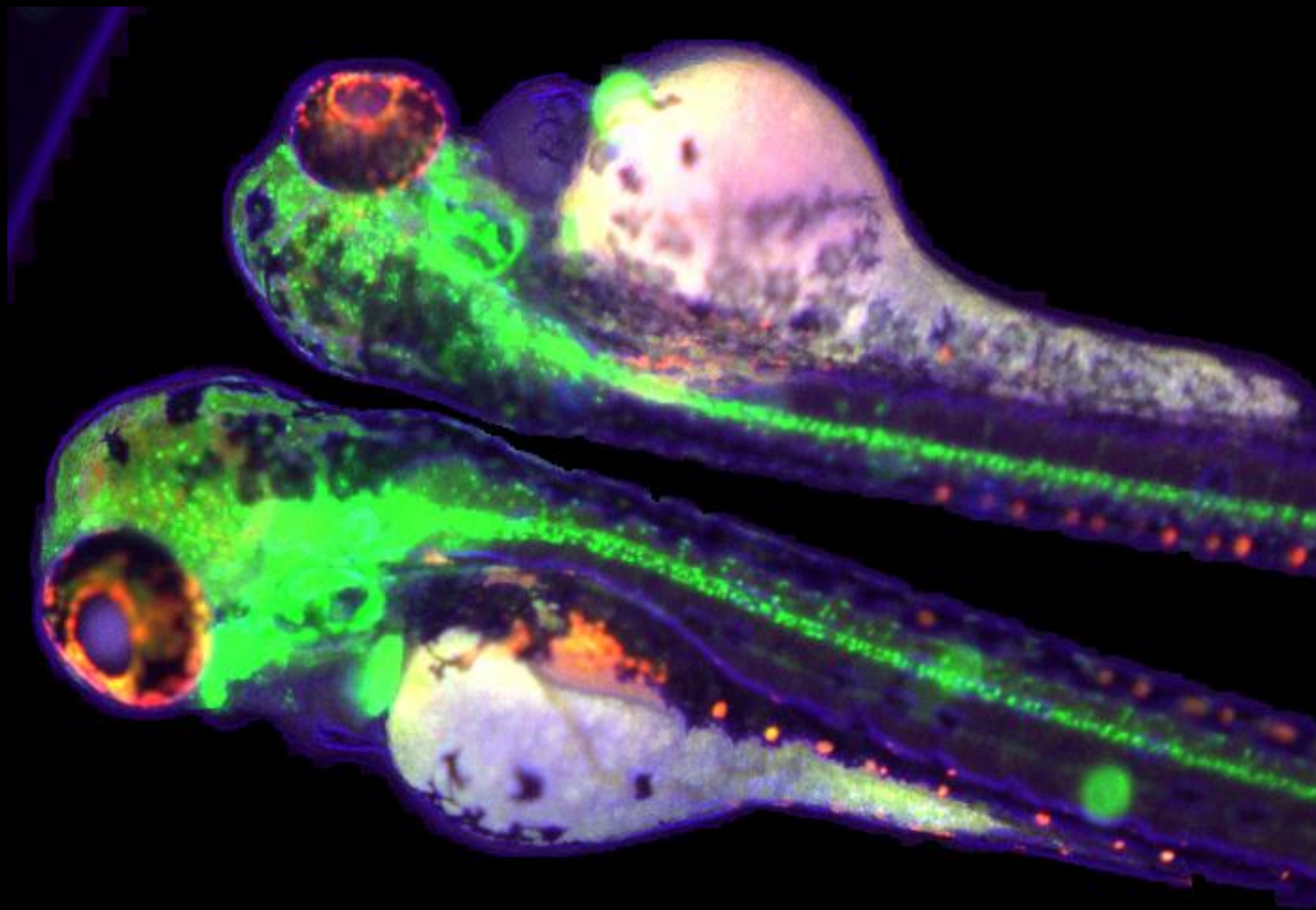


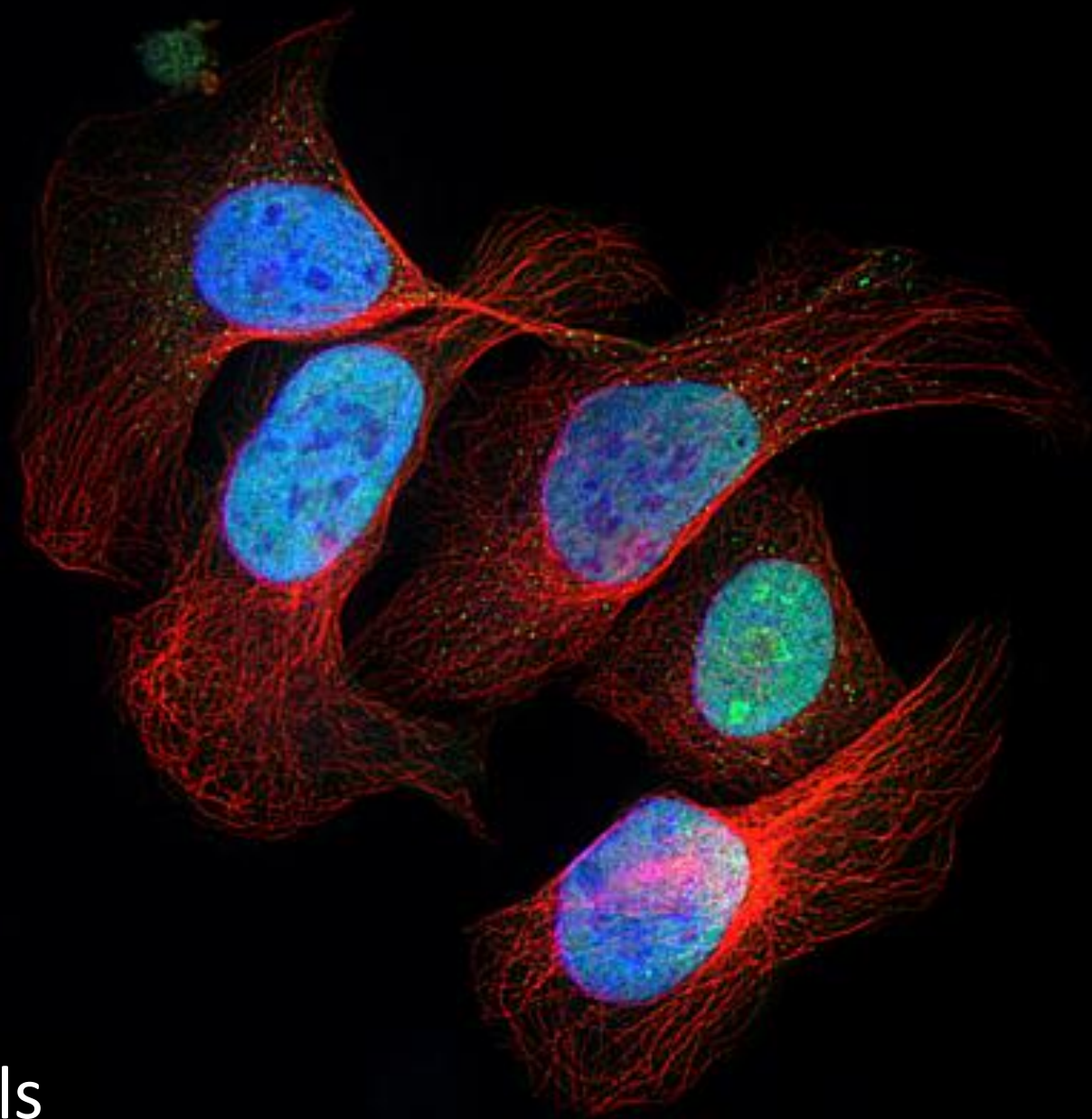
Plate and Well View



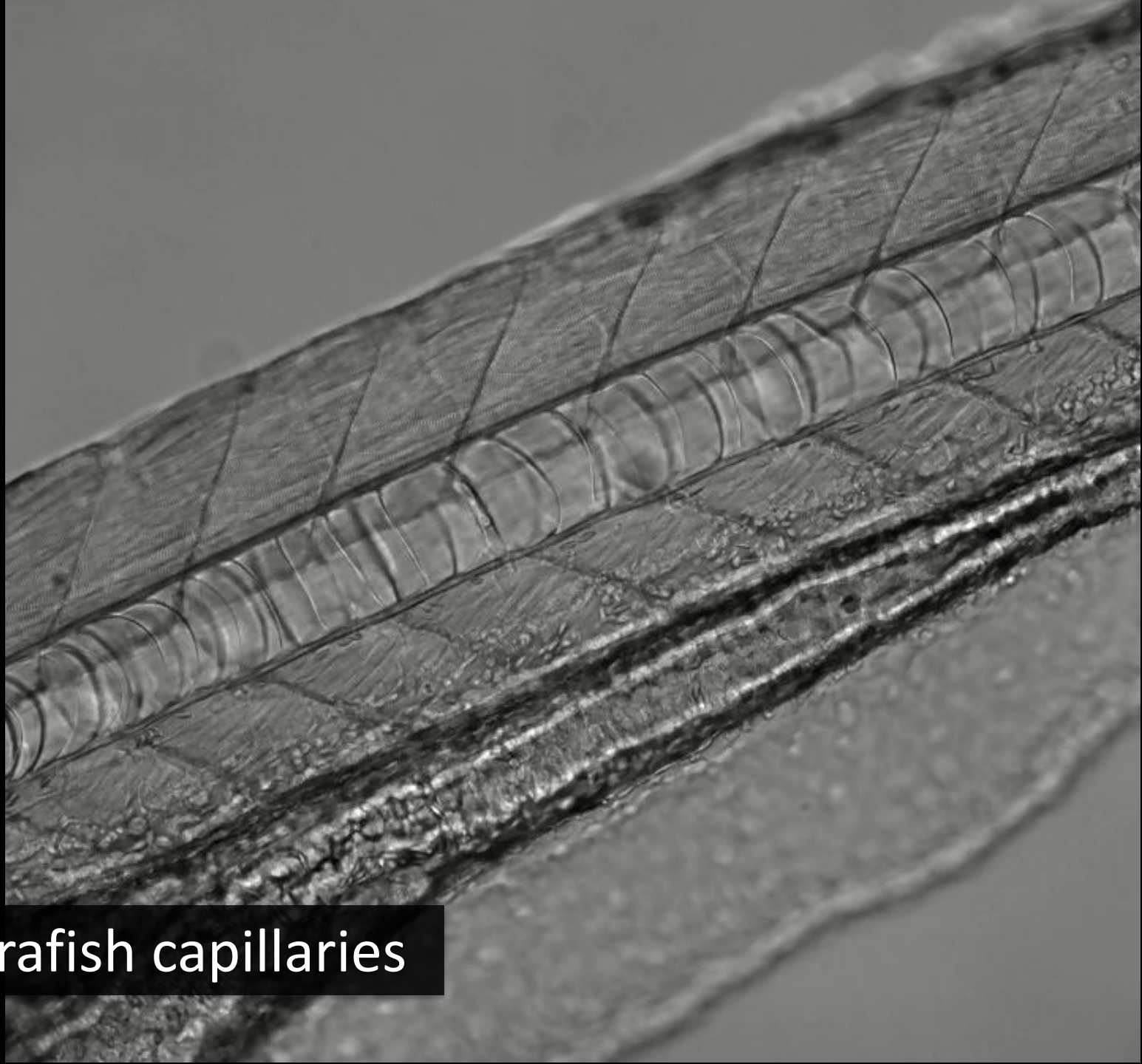
Neurons



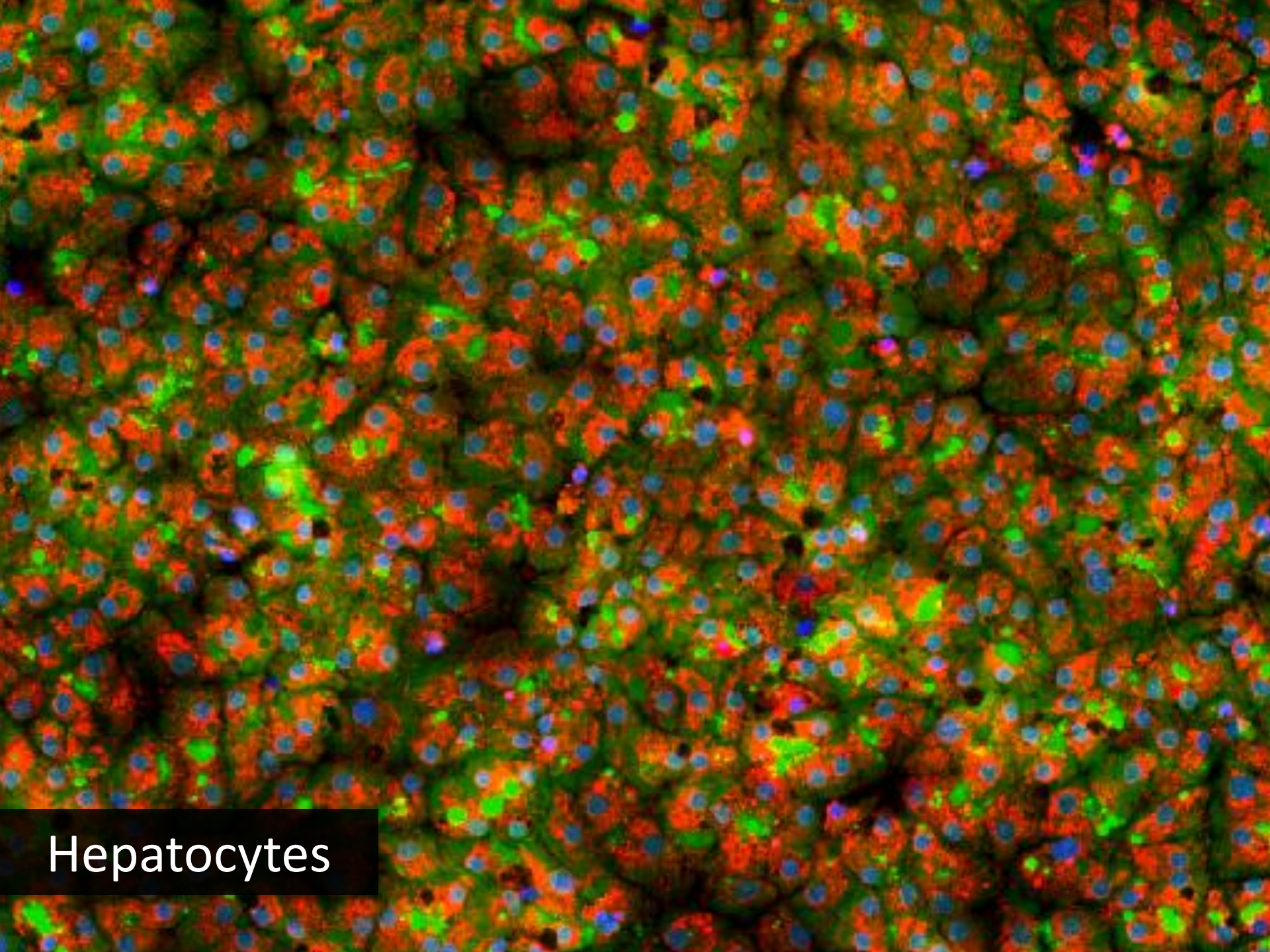
Zebrafish



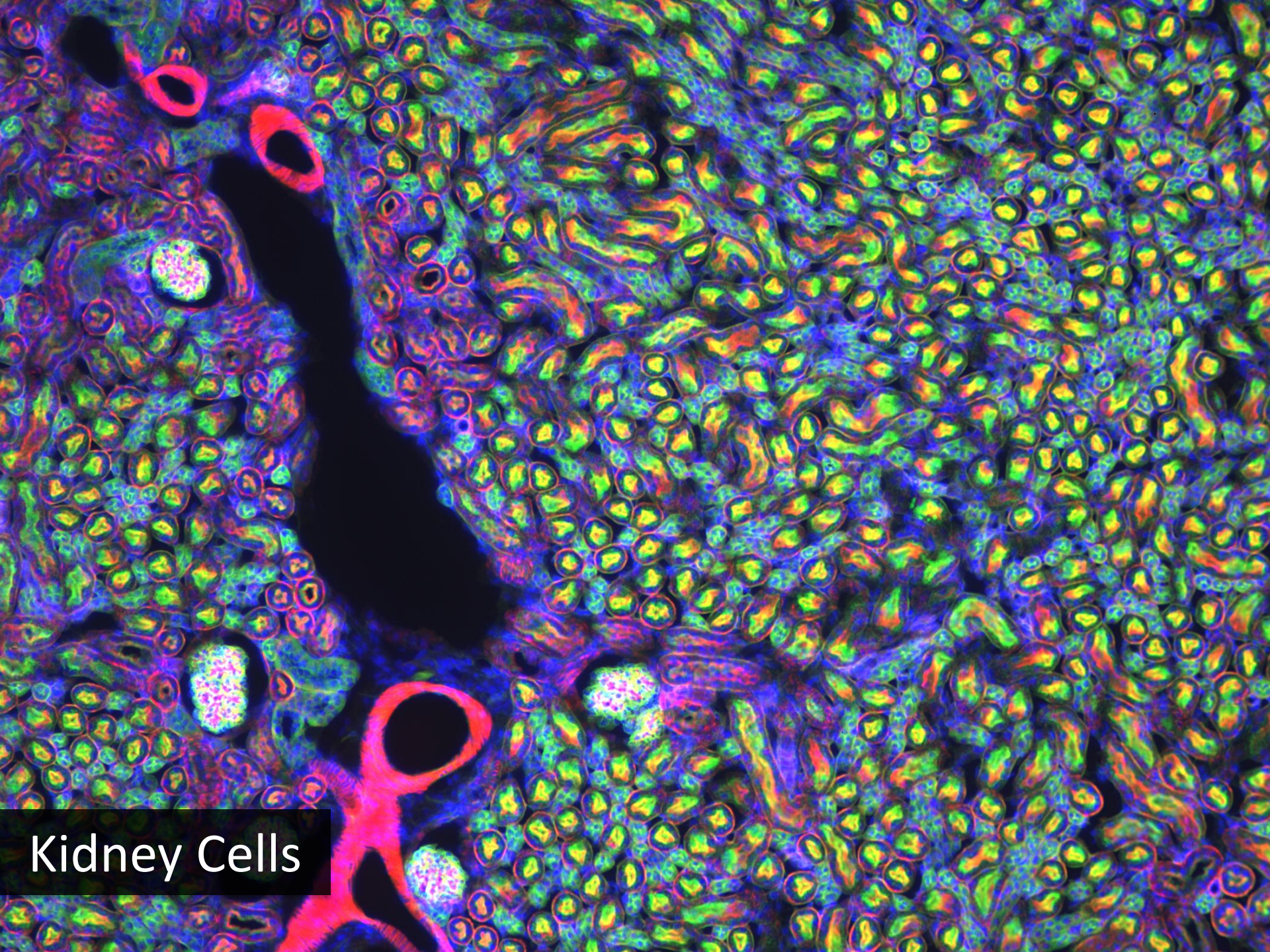
U2OS cells



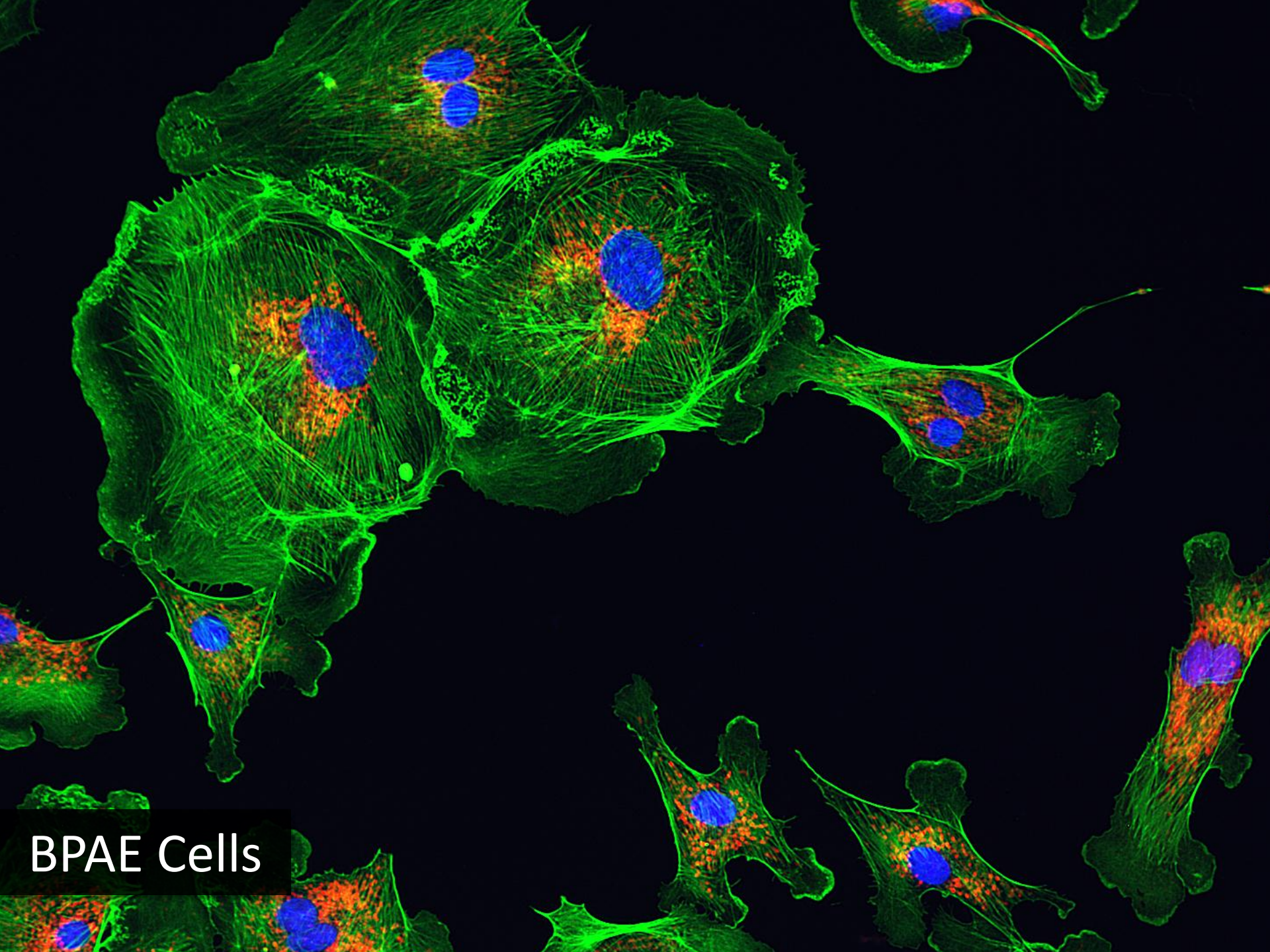
Zebrafish capillaries



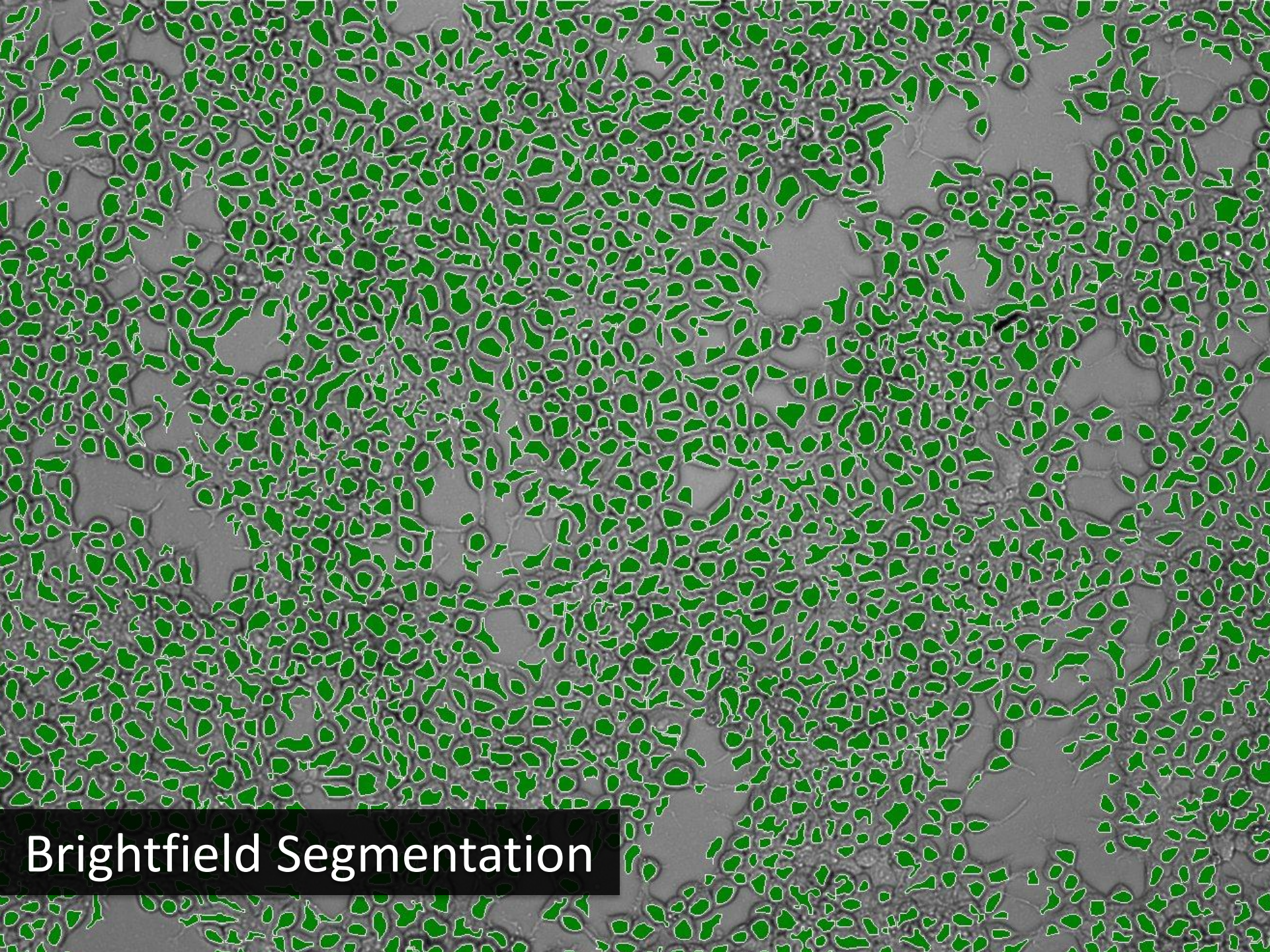
Hepatocytes



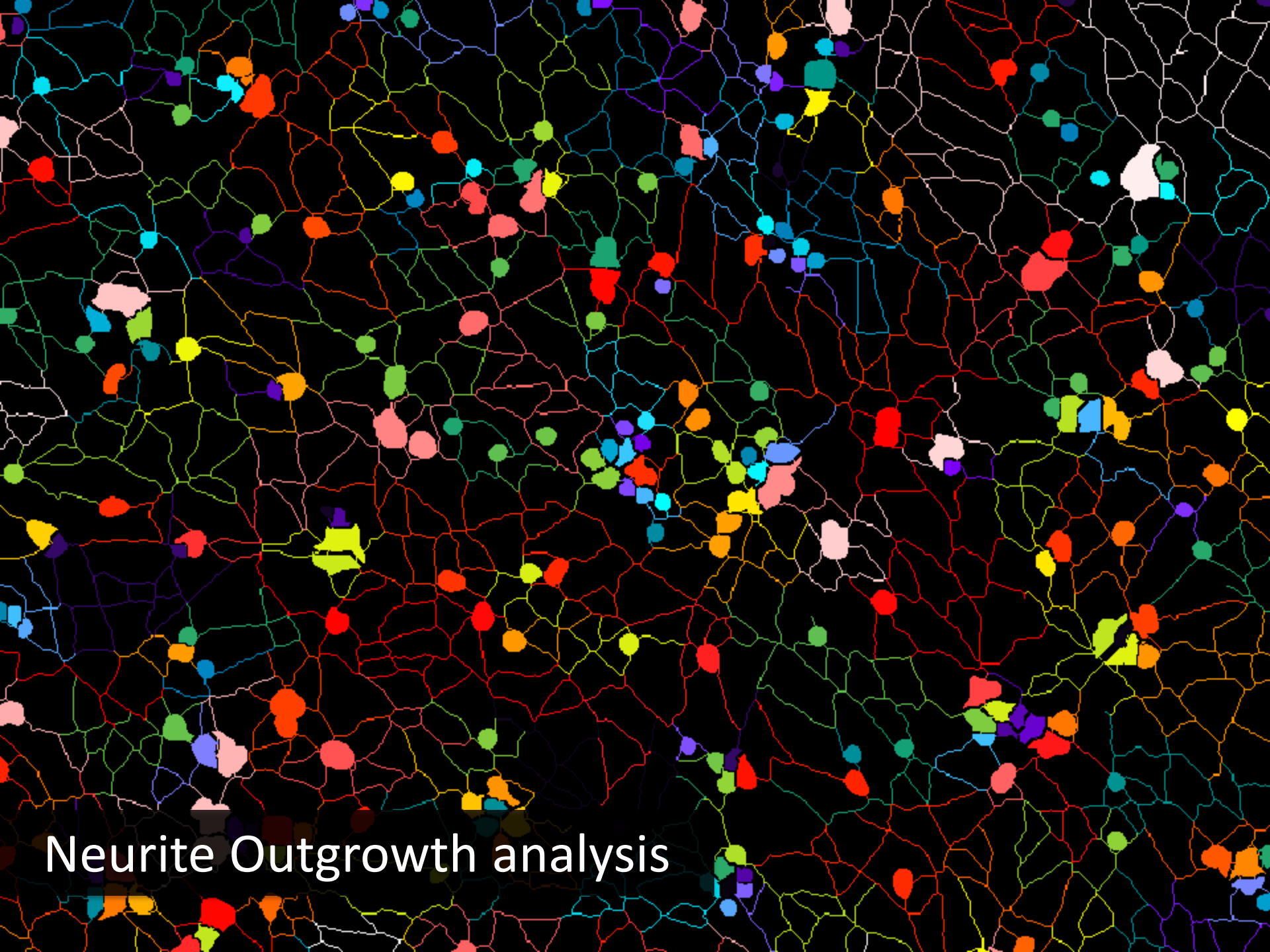
Kidney Cells



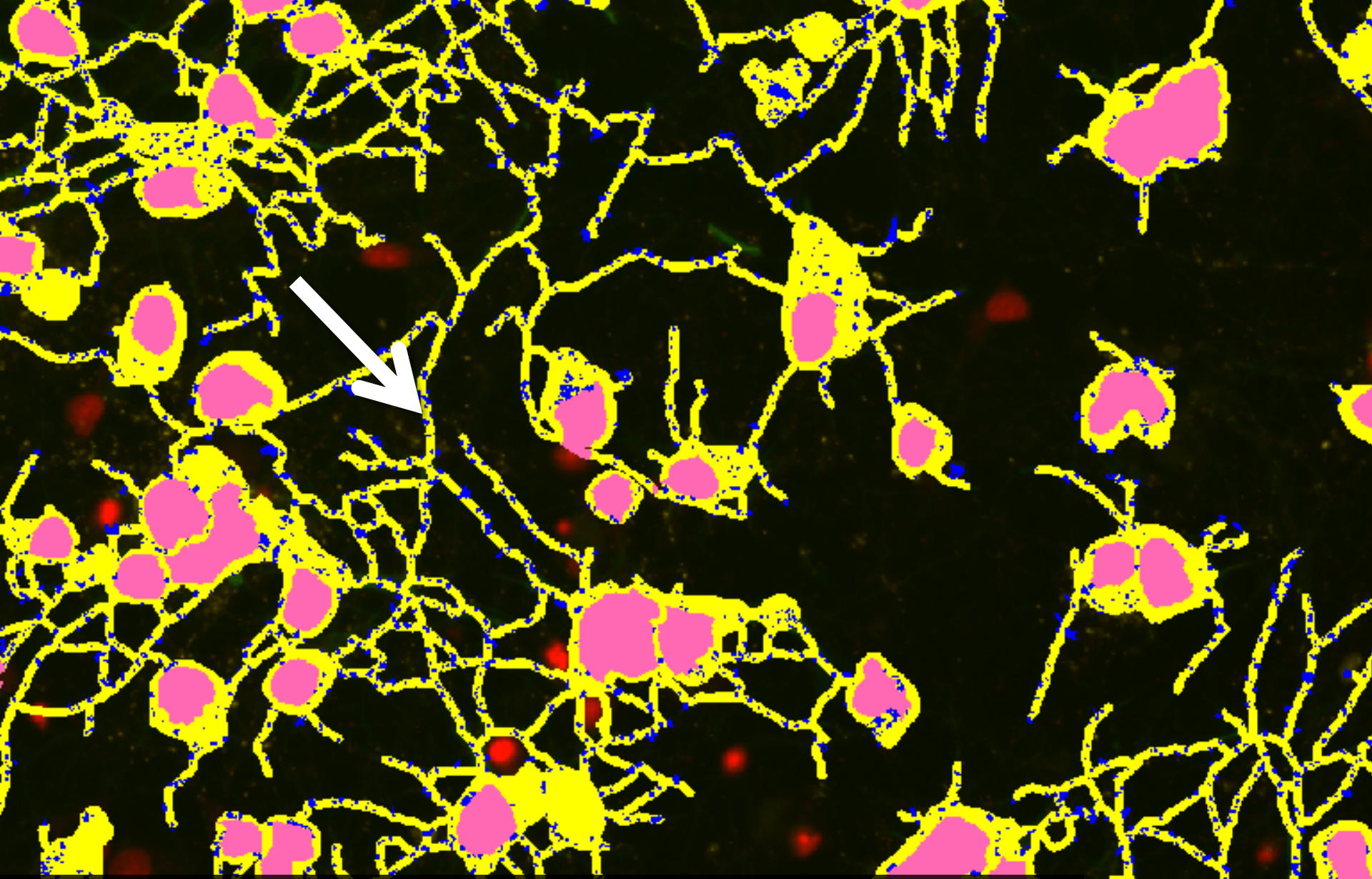
BPAE Cells



Brightfield Segmentation



Neurite Outgrowth analysis

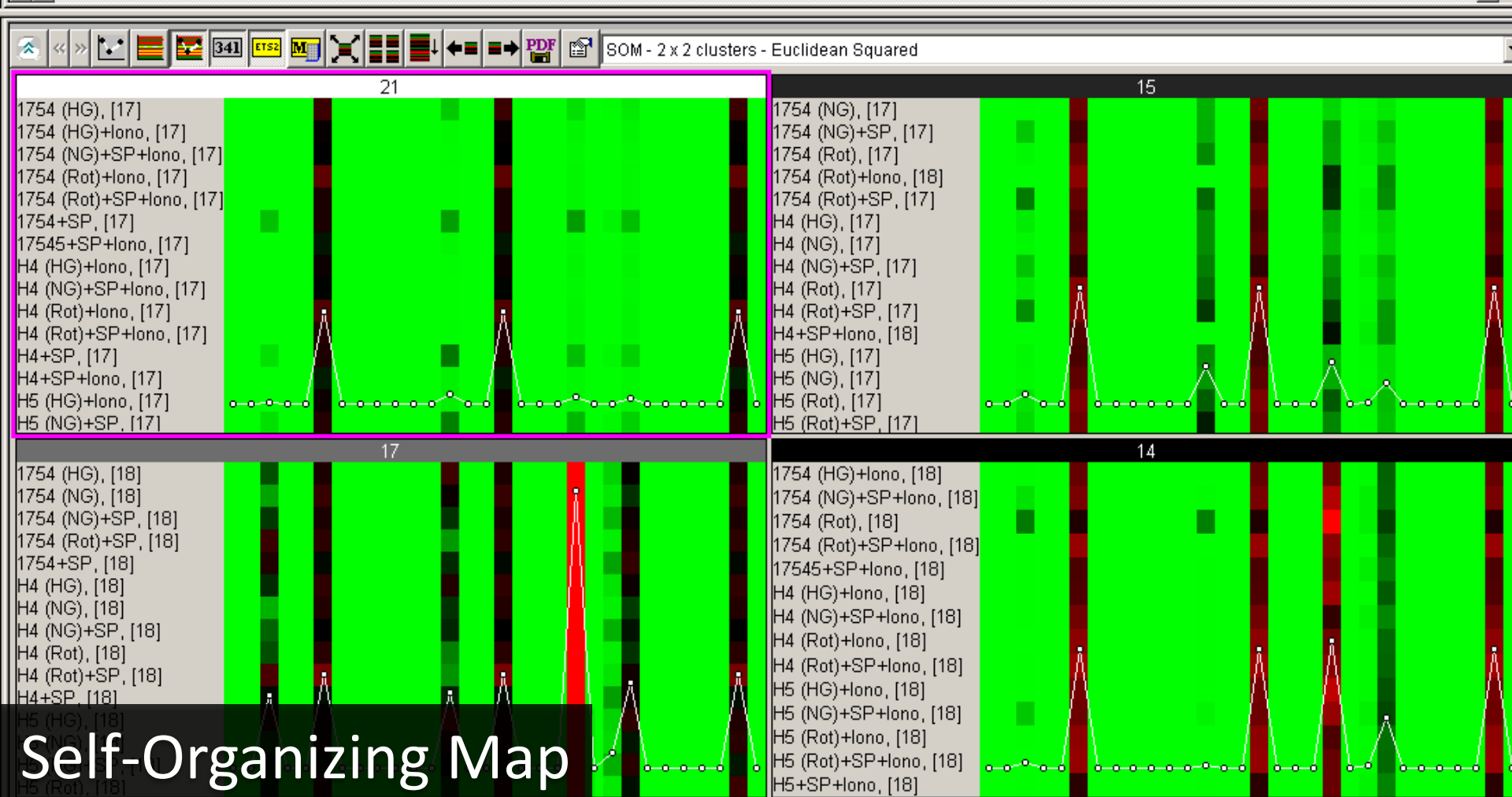


Punctae in Neurites with Custom Modules

	01	02	03	04	05	06	07	08	09	10	11	12	13	14	15	16	17	18	19	20	21	22	23
A																							
B																							
C			27	29	37	36	39	39	14	26	2	30	63	68	58	59	25	26	18	27	24	24	
D			35	31	34	31	41	35	33	30	7	32	63	62	64	62	25	27	24	29	24	27	
E			35	27	32	29	30	31	29	34	21	32	59	68	62	61	20	24	21	25	23	24	
F			37	25	37	36	34	42	34	31	27	32	61	63	58	59	23	26	28	27	28	24	
G			31	33	33	35	31	25	32	30	32	25	53	50	52	53	49	51	45	51	27	21	
H			38	31	22	35	33	33	35	33	34	28	33	36	33	26	65	65	62	62	22	25	
I			34	30	36	30	33	35	39	28	28	29	29	26	23	30	64	65	62	63	25	27	
J			42	35	36	39	34	35	36	31	33	34	27	25	21	16	67	60	54	65	22	22	
K			43	35	36	38	36	34	43	35	32	33	28	24	15	0	63	65	64	62	24	24	
L			37	30	36	37	34	42	29	32	32	34	21	24	25	27	64	60	66	64	22	25	
M			43	35	37	42	34	34	25	36	30	24	21	20	26	17	23	65	66	68	23	26	
N			46	35	37	31	42	35	40	38	34	33	23	24	4	19	10	3	68	66	20	21	
O																							
P																							

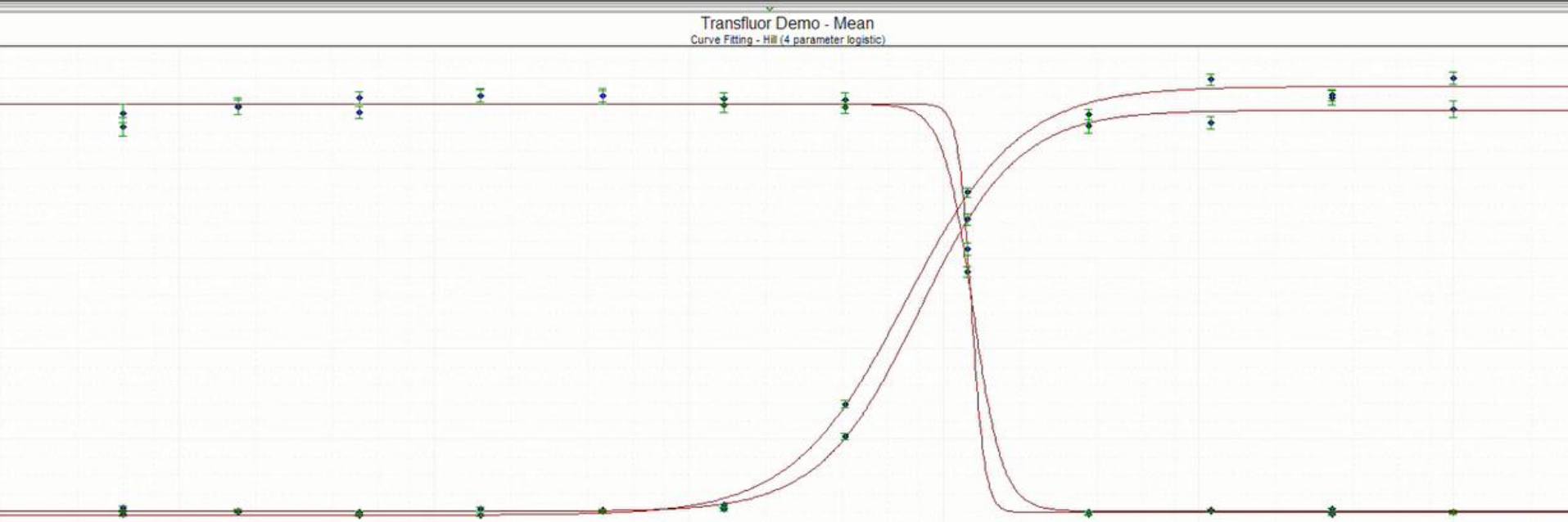
Heat Map

Data		Row Descriptors		Web Links		Statistics [Mean]		Advanced		Curve Fit				
Row: Compound [Well]		Column: Concentration		Measurement:		Cell: Vesicle Count		Calculation: Mean						
Compound [Well Annotation], Plate ID [Cell Measurement]		0 - [Cell: Vesicle Count ...]	0 - [Cell: Vesicle Total ...]	0 - [Cell: Vesicle Integrated Intensity ...]	0 - [Cell: Vesicle Average ...]	0 - [Cell: Nuclear Total Area ...]	0 - [Cell: Nuclear Integrated ...]	0 - [Cell: Nuclear Average ...]	0 - [Cell: Texture Index ...]	0 - [Cell: Gradient Index ...]	0 - [Cell: Laplacian Index ...]	0 - [Cell: Granule Count ...]	0 - [Cell: Granule Total ...]	0 - [Cell: Granule Integrated ...]
1754 (HG), [17]		0.00193	0.087	7692.954	16.538	136.342	7345196.500	5815.083	407.710	99.847	73.613	10.818	22.204	1231949
1754 (HG), [18]		0.586	35.385	4105755.250	6107.340	140.862	8649820	5160.035	2585.222	479.038	340.797	13.901	47.364	7464454
1754 (HG)+Iono, [17]		0	0	0	0	121.932	6149482	5460.573	16.115	30.916	22.496	1.899	2.433	118672.453
1754 (HG)+Iono, [18]		0.00803	0.419	46311.961	92.368	137.306	8358854	6703.698	627.422	184.328	132.903	0.317	0.335	17999.520



Self-Organizing Map

ound [Well Annotation], Plate ID [Cell Measurement]	6e-006 - [Cell: Vesicle Total Area ...]	1.7e-005 - [Cell: Vesicle Total Area ...]	5.1e-005 - [Cell: Vesicle Total Area ...]	0.000152 - [Cell: Vesicle Total Area ...]	0.000457 - [Cell: Vesicle Total Area ...]	0.001372 - [Cell: Vesicle Total Area ...]	0.004115 - [Cell: Vesicle Total Area ...]	0.012346 - [Cell: Vesicle Total Area ...]	0.037037 - [Cell: Vesicle Total Area ...]	0.111111 - [Cell: Vesicle Total Area ...]	0.333333 - [Cell: Vesicle Total Area ...]	1 - [Cell: Vesicle Total Area ...]	C50	Correlation Coefficient
#14 Propranolol DIC, [201]	47.524	45.498	48.248	47.535	47.099	48.602	47.287	37.072	4.725	2.675	2.316	2.472	0.016988	0.981341
#14 Propranolol DIC, [202]	50.164	47.507	49.536	47.899	47.759	46.471	49.311	28.717	2.661	2.045	2.042	1.975	0.012597	0.975066
#14 Propranolol DIC, [203]	38.720	43.672	47.783	47.822	46.348	45.994	44.886	22.733	1.819	2.159	2.514	2.578	0.012294	0.936259
#14 Propranolol DIC, [204]	44.639	46.842	47.821	48.137	48.043	47.057	47.639	31.079	1.684	2.067	1.629	1.903	0.012803	0.976700
#14 Propranolol DIC, [205]	46.120	46.958	46.258	48.113	48.072	47.755	46.818	28.562	1.840	2.005	2.186	1.921	0.012952	0.983972
#14 Isoproterenol DRC, [200]	2.206	2.544	1.952	2.286	1.867	4.565	14.104	37.242	45.964	48.840	47.535	48.898	0.006968	0.983968
#14 Isoproterenol DRC, [201]	2.115	2.395	1.738	1.734	2.310	3.002	15.683	40.880	46.437	48.921	48.477	49.751	0.006100	0.981103
#14 Isoproterenol DRC, [202]	2.696	1.821	2.119	2.357	2.016	4.039	13.267	37.033	47.563	50.443	50.781	50.557	0.007645	0.989125
#14 Isoproterenol DRC, [203]	2.320	2.019	1.821	2.224	2.065	2.168	10.259	34.354	44.749	45.141	47.888	46.585	0.007990	0.979175
#14 Isoproterenol DRC, [204]	1.626	1.923	1.589	1.555	1.999	2.626	13.886	37.351	46.004	49.903	48.190	50.061	0.007092	0.978112
#14 Isoproterenol DRC, [205]	1.506	1.398	1.909	2.734	1.832	2.569	11.265	33.984	45.177	48.680	48.062	48.403	0.008242	0.984203



Curve Fit Statistics