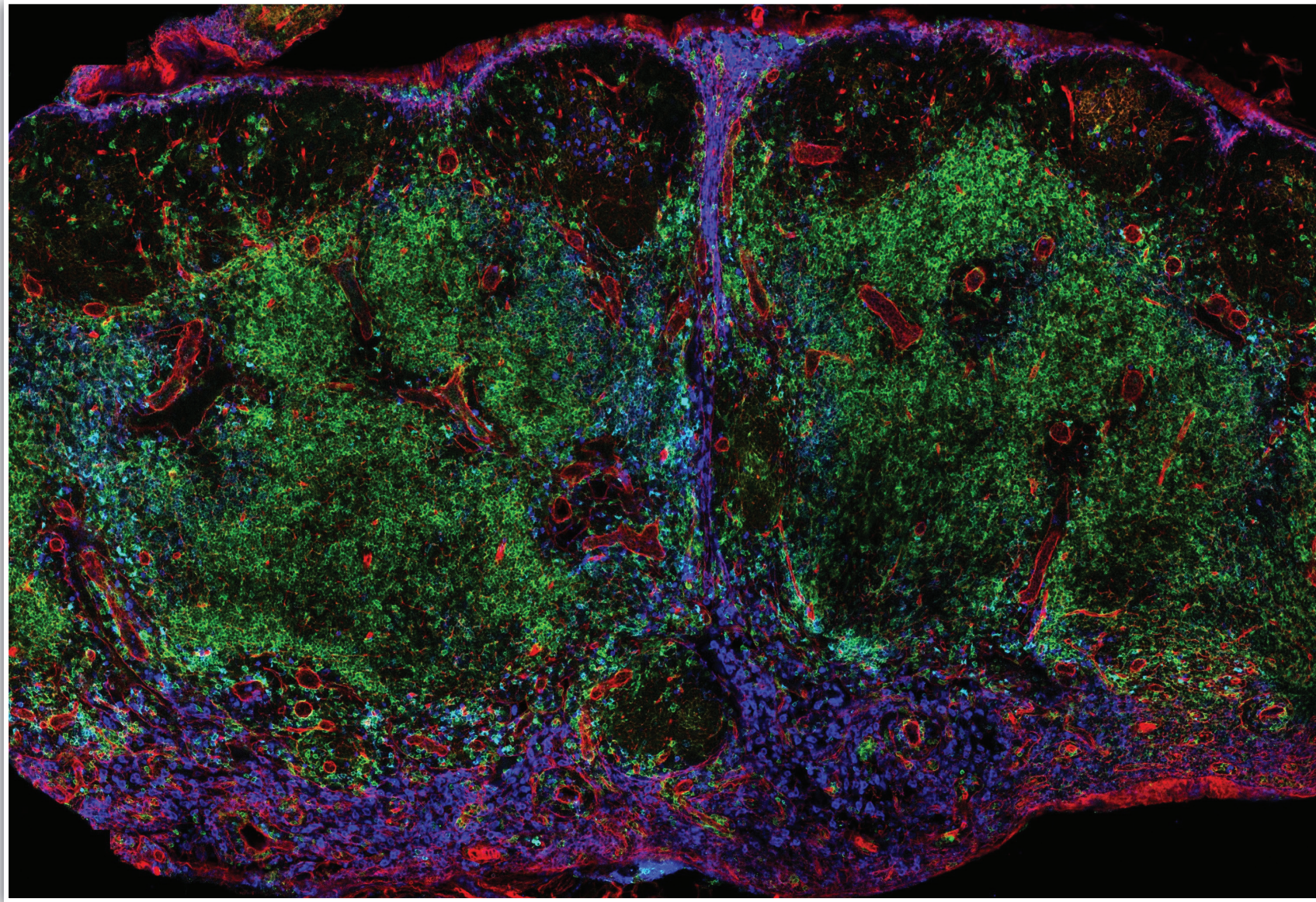
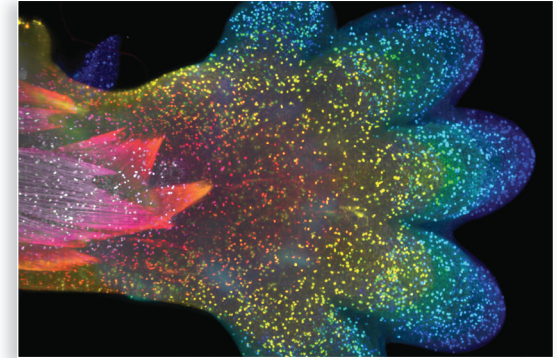
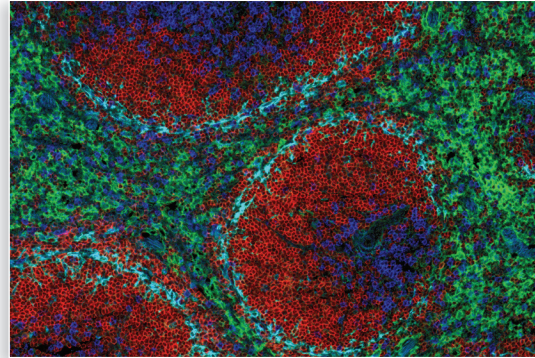
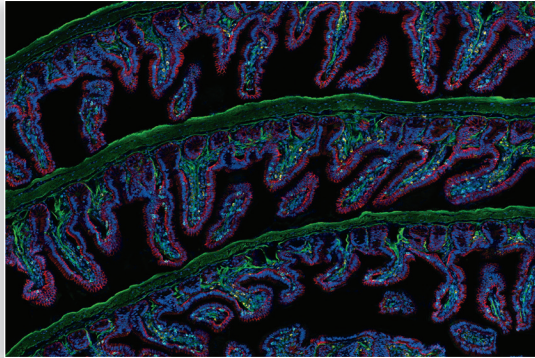


Legendary Microscopy 2024





For over 20 years, BioLegend has partnered with scientists around the world to develop novel reagents. Our beginnings were rooted in flow cytometry solutions for immunology, but we have since expanded our portfolio into a variety of applications and research areas. From cell culture, multiomics, and immunoassay applications to cancer, neuroscience, and stem cell research, our growth would not have been possible without your input and collaboration. For our 2024 calendar, we are focusing on how our tools are advancing spatial biology with a showcase of customer-generated microscopy images. We hope you enjoy the spectacular visuals in this calendar as much as we enjoyed seeing how our reagents support your research. As we continue to pave our way in the world of spatial biology, we are excited to continue our partnership with you.

2024 CALENDAR

2025 CALENDAR

JANUARY						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30	31			

FEBRUARY						
S	M	T	W	T	F	S
				1	2	3
4	5	6	7	8	9	10
11	12	13	14	15	16	17
18	19	20	21	22	23	24
25	26	27	28	29		

MARCH						
S	M	T	W	T	F	S
					1	2
3	4	5	6	7	8	9
10	11	12	13	14	15	16
17	18	19	20	21	22	23
24	25	26	27	28	29	30
31						

JANUARY						
S	M	T	W	T	F	S
			1	2	3	4
5	6	7	8	9	10	11
12	13	14	15	16	17	18
19	20	21	22	23	24	25
26	27	28	29	30	31	

FEBRUARY						
S	M	T	W	T	F	S
						1
2	3	4	5	6	7	8
9	10	11	12	13	14	15
16	17	18	19	20	21	22
23	24	25	26	27	28	

MARCH						
S	M	T	W	T	F	S
						1
2	3	4	5	6	7	8
9	10	11	12	13	14	15
16	17	18	19	20	21	22
23	24	25	26	27	28	29
30	31					

APRIL						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30				

MAY						
S	M	T	W	T	F	S
		1	2	3	4	
5	6	7	8	9	10	11
12	13	14	15	16	17	18
19	20	21	22	23	24	25
26	27	28	29	30	31	

JUNE						
S	M	T	W	T	F	S
						1
2	3	4	5	6	7	8
9	10	11	12	13	14	15
16	17	18	19	20	21	22
23	24	25	26	27	28	29
30						

APRIL						
S	M	T	W	T	F	S
	1	2	3	4	5	
6	7	8	9	10	11	12
13	14	15	16	17	18	19
20	21	22	23	24	25	26
27	28	29	30			

MAY						
S	M	T	W	T	F	S
		1	2	3		
4	5	6	7	8	9	10
11	12	13	14	15	16	17
18	19	20	21	22	23	24
25	26	27	28	29	30	31

JUNE						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30				

JULY						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30	31			

AUGUST						
S	M	T	W	T	F	S
				1	2	3
4	5	6	7	8	9	10
11	12	13	14	15	16	17
18	19	20	21	22	23	24
25	26	27	28	29	30	31

SEPTEMBER						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30				

JULY						
S	M	T	W	T	F	S
	1	2	3	4	5	
6	7	8	9	10	11	12
13	14	15	16	17	18	19
20	21	22	23	24	25	26
27	28	29	30	31		

AUGUST						
S	M	T	W	T	F	S
					1	2
3	4	5	6	7	8	9
10	11	12	13	14	15	16
17	18	19	20	21	22	23
24	25	26	27	28	29	30
31						

SEPTEMBER						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30				

OCTOBER						
S	M	T	W	T	F	S
	1	2	3	4	5	
6	7	8	9	10	11	12
13	14	15	16	17	18	19
20	21	22	23	24	25	26
27	28	29	30	31		

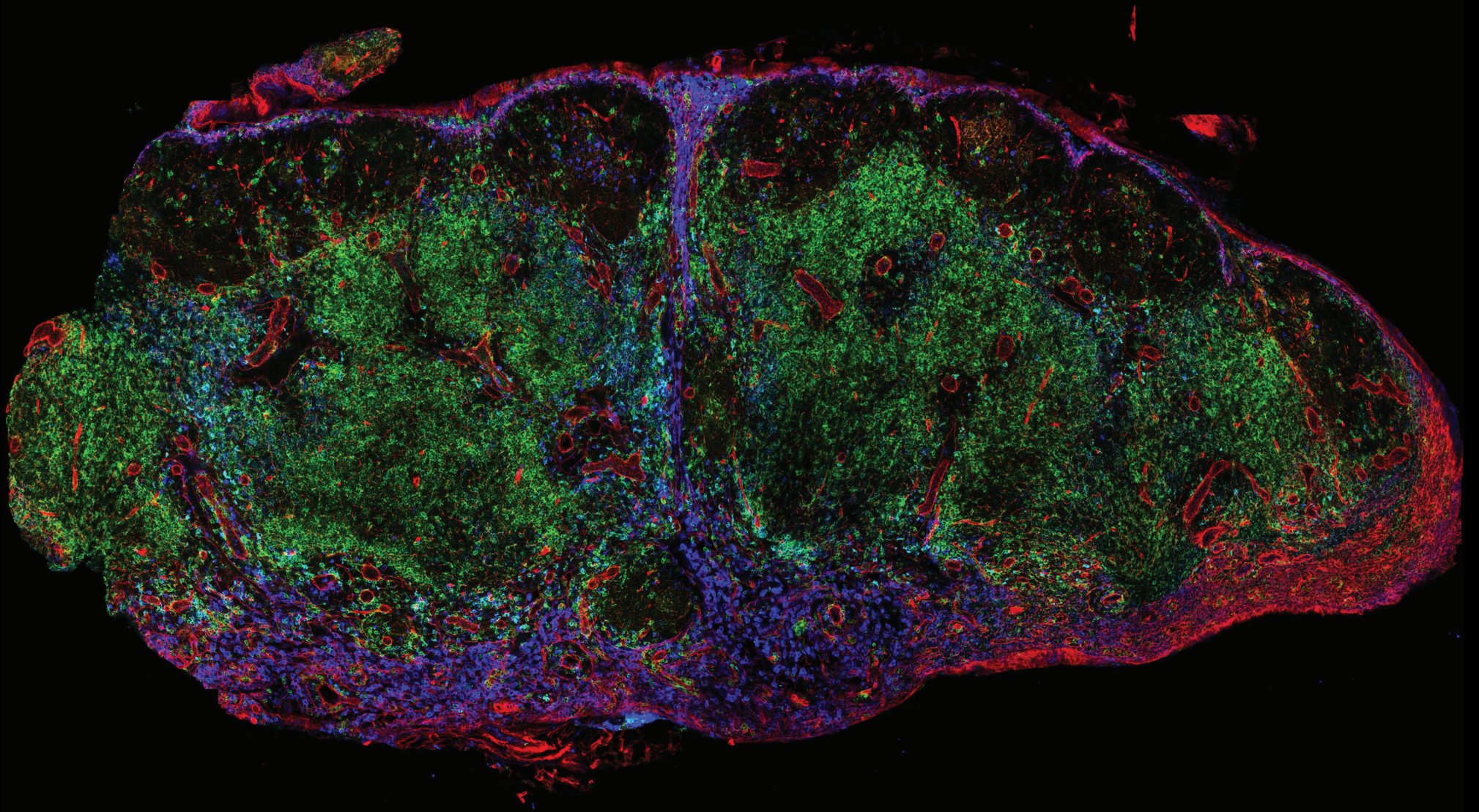
NOVEMBER						
S	M	T	W	T	F	S
					1	2
3	4	5	6	7	8	9
10	11	12	13	14	15	16
17	18	19	20	21	22	23
24	25	26	27	28	29	30

DECEMBER						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30	31			

OCTOBER						
S	M	T	W	T	F	S
			1	2	3	4
5	6	7	8	9	10	11
12	13	14	15	16	17	18
19	20	21	22	23	24	25
26	27	28	29	30	31	

NOVEMBER						
S	M	T	W	T	F	S
						1
2	3	4	5	6	7	8
9	10	11	12	13	14	15
16	17	18	19	20	21	22
23	24	25	26	27	28	29
30						

DECEMBER						
S	M	T	W	T	F	S
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30	31			

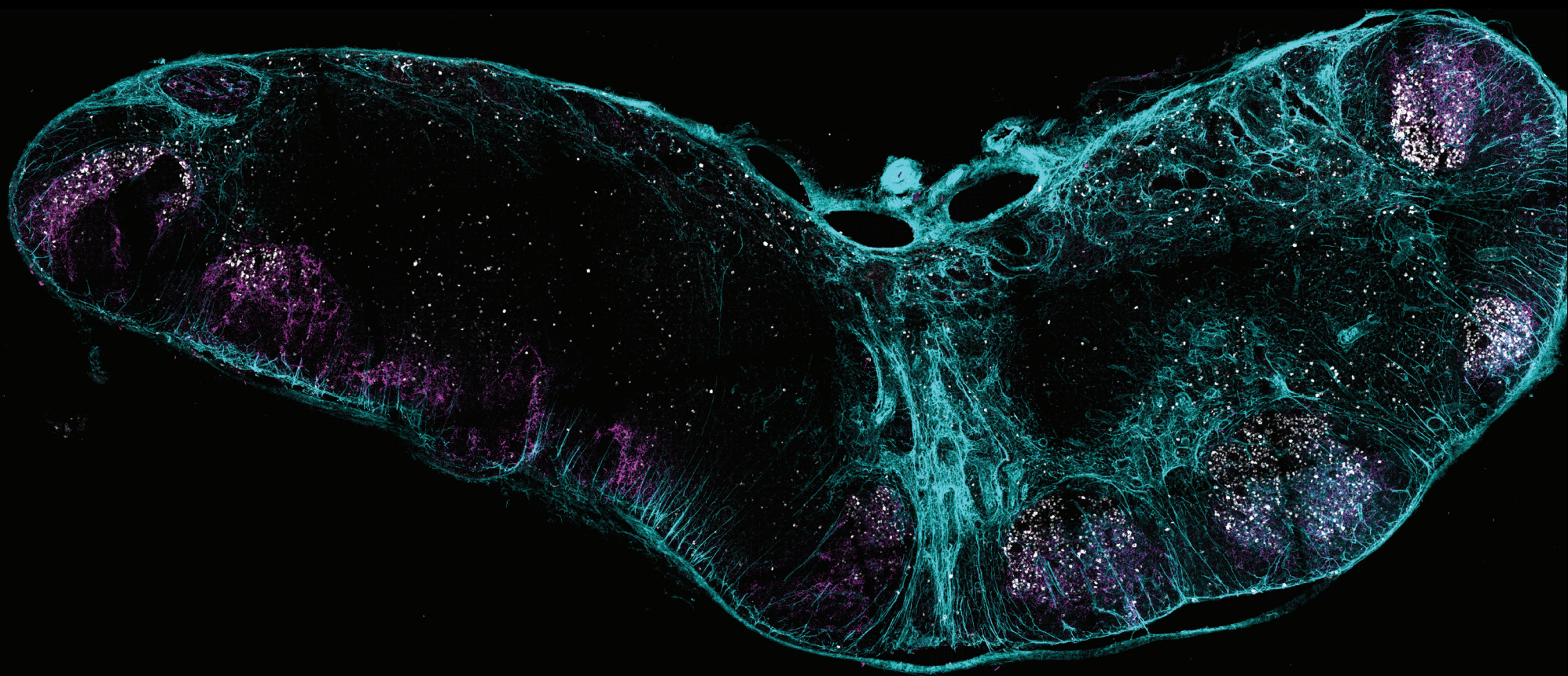


Jake Colautti
McMaster University

Mouse lymph node tissue stained with Alexa Fluor[®] 488 anti-mouse CD3 (green; Cat. No. 100212), PE anti-mouse/human CD11b (blue; Cat. No. 101208), and anti-CD29 (red).

JANUARY 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																																	
<p>DECEMBER</p> <table border="1"> <thead> <tr> <th>S</th> <th>M</th> <th>T</th> <th>W</th> <th>T</th> <th>F</th> <th>S</th> </tr> </thead> <tbody> <tr> <td></td> <td></td> <td></td> <td></td> <td>1</td> <td>2</td> <td></td> </tr> <tr> <td>3</td> <td>4</td> <td>5</td> <td>6</td> <td>7</td> <td>8</td> <td>9</td> </tr> <tr> <td>10</td> <td>11</td> <td>12</td> <td>13</td> <td>14</td> <td>15</td> <td>16</td> </tr> <tr> <td>17</td> <td>18</td> <td>19</td> <td>20</td> <td>21</td> <td>22</td> <td>23</td> </tr> <tr> <td>24</td> <td>25</td> <td>26</td> <td>27</td> <td>28</td> <td>29</td> <td>30</td> </tr> <tr> <td>31</td> <td></td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S					1	2		3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31							1	2	3	4	5	6
S	M	T	W	T	F	S																																																	
				1	2																																																		
3	4	5	6	7	8	9																																																	
10	11	12	13	14	15	16																																																	
17	18	19	20	21	22	23																																																	
24	25	26	27	28	29	30																																																	
31																																																							
7	New Year's Day	9	10	11	12	13																																																	
14	15	16	17	18	19	20																																																	
21	22	23	24	25	26	27																																																	
28	Martin Luther King Jr. Day	30	31	<p>FEBRUARY</p> <table border="1"> <thead> <tr> <th>S</th> <th>M</th> <th>T</th> <th>W</th> <th>T</th> <th>F</th> <th>S</th> </tr> </thead> <tbody> <tr> <td></td> <td></td> <td></td> <td></td> <td>1</td> <td>2</td> <td>3</td> </tr> <tr> <td>4</td> <td>5</td> <td>6</td> <td>7</td> <td>8</td> <td>9</td> <td>10</td> </tr> <tr> <td>11</td> <td>12</td> <td>13</td> <td>14</td> <td>15</td> <td>16</td> <td>17</td> </tr> <tr> <td>18</td> <td>19</td> <td>20</td> <td>21</td> <td>22</td> <td>23</td> <td>24</td> </tr> <tr> <td>25</td> <td>26</td> <td>27</td> <td>28</td> <td>29</td> <td></td> <td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S					1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29											
S	M	T	W	T	F	S																																																	
				1	2	3																																																	
4	5	6	7	8	9	10																																																	
11	12	13	14	15	16	17																																																	
18	19	20	21	22	23	24																																																	
25	26	27	28	29																																																			

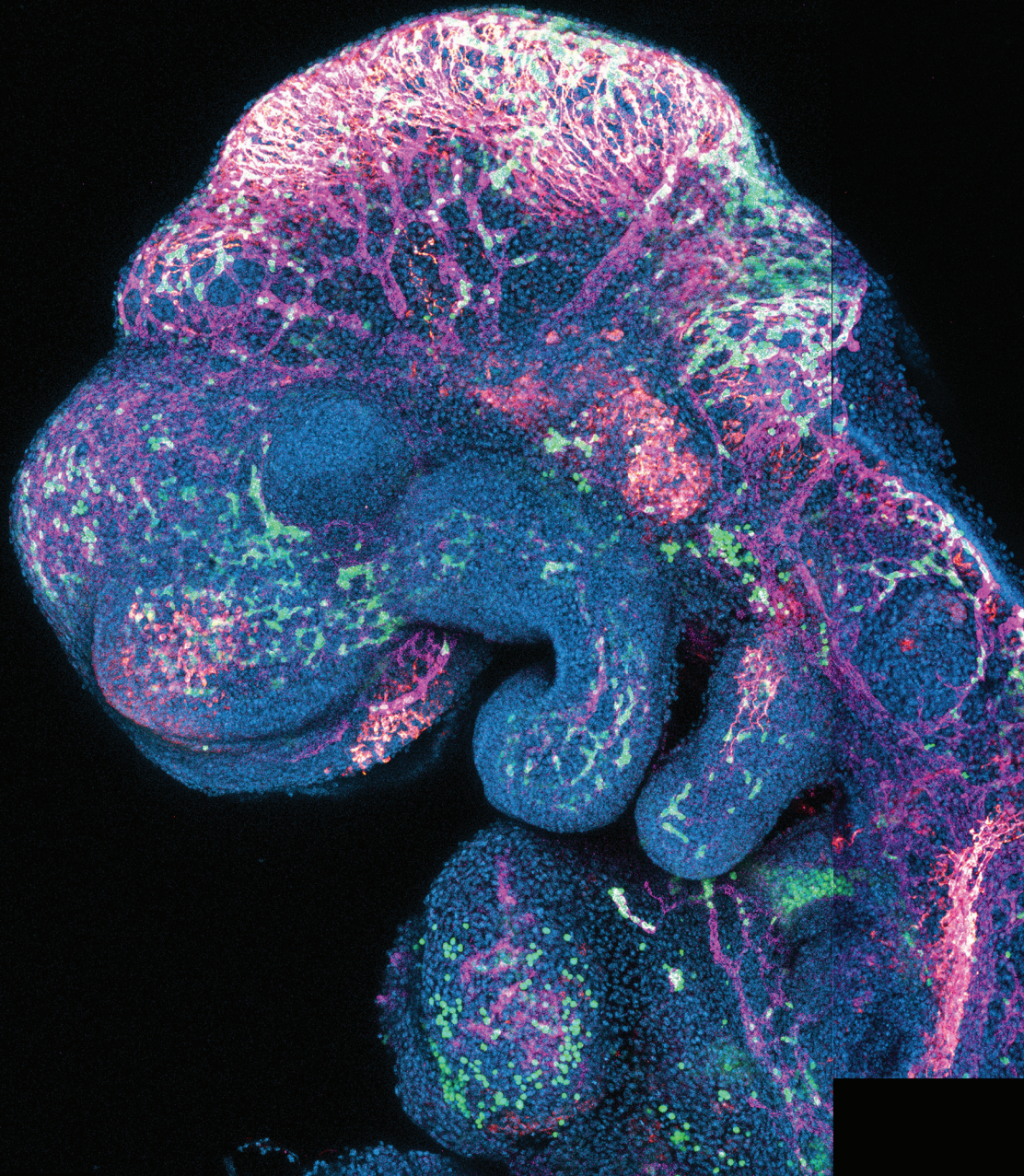


Darelynn Oo
University of Calgary

Mouse inguinal lymph node tissue stained with Ultra-LEAF™ Purified anti-mouse CD21/CD35 (CR2/CR1)(magenta; Cat. No. 123430), anti-Ki-67 (grey), and FITC tracer (cyan).

FEBRUARY 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																																	
		JANUARY <table border="1"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr><td></td><td>1</td><td>2</td><td>3</td><td>4</td><td>5</td><td>6</td></tr> <tr><td>7</td><td>8</td><td>9</td><td>10</td><td>11</td><td>12</td><td>13</td></tr> <tr><td>14</td><td>15</td><td>16</td><td>17</td><td>18</td><td>19</td><td>20</td></tr> <tr><td>21</td><td>22</td><td>23</td><td>24</td><td>25</td><td>26</td><td>27</td></tr> <tr><td>28</td><td>29</td><td>30</td><td>31</td><td></td><td></td><td></td></tr> </tbody> </table>	S	M	T	W	T	F	S		1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31						1	2	3						
S	M	T	W	T	F	S																																																	
	1	2	3	4	5	6																																																	
7	8	9	10	11	12	13																																																	
14	15	16	17	18	19	20																																																	
21	22	23	24	25	26	27																																																	
28	29	30	31																																																				
4	5	6	7	8	9	10 Lunar New Year																																																	
11	12	13	14 Valentine's Day	15	16	17																																																	
18	19 Presidents' Day	20	21	22	23	24																																																	
25	26	27	28	29	MARCH <table border="1"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr><td></td><td></td><td></td><td></td><td></td><td>1</td><td>2</td></tr> <tr><td>3</td><td>4</td><td>5</td><td>6</td><td>7</td><td>8</td><td>9</td></tr> <tr><td>10</td><td>11</td><td>12</td><td>13</td><td>14</td><td>15</td><td>16</td></tr> <tr><td>17</td><td>18</td><td>19</td><td>20</td><td>21</td><td>22</td><td>23</td></tr> <tr><td>24</td><td>25</td><td>26</td><td>27</td><td>28</td><td>29</td><td>30</td></tr> <tr><td>31</td><td></td><td></td><td></td><td></td><td></td><td></td></tr> </tbody> </table>	S	M	T	W	T	F	S						1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31							
S	M	T	W	T	F	S																																																	
					1	2																																																	
3	4	5	6	7	8	9																																																	
10	11	12	13	14	15	16																																																	
17	18	19	20	21	22	23																																																	
24	25	26	27	28	29	30																																																	
31																																																							



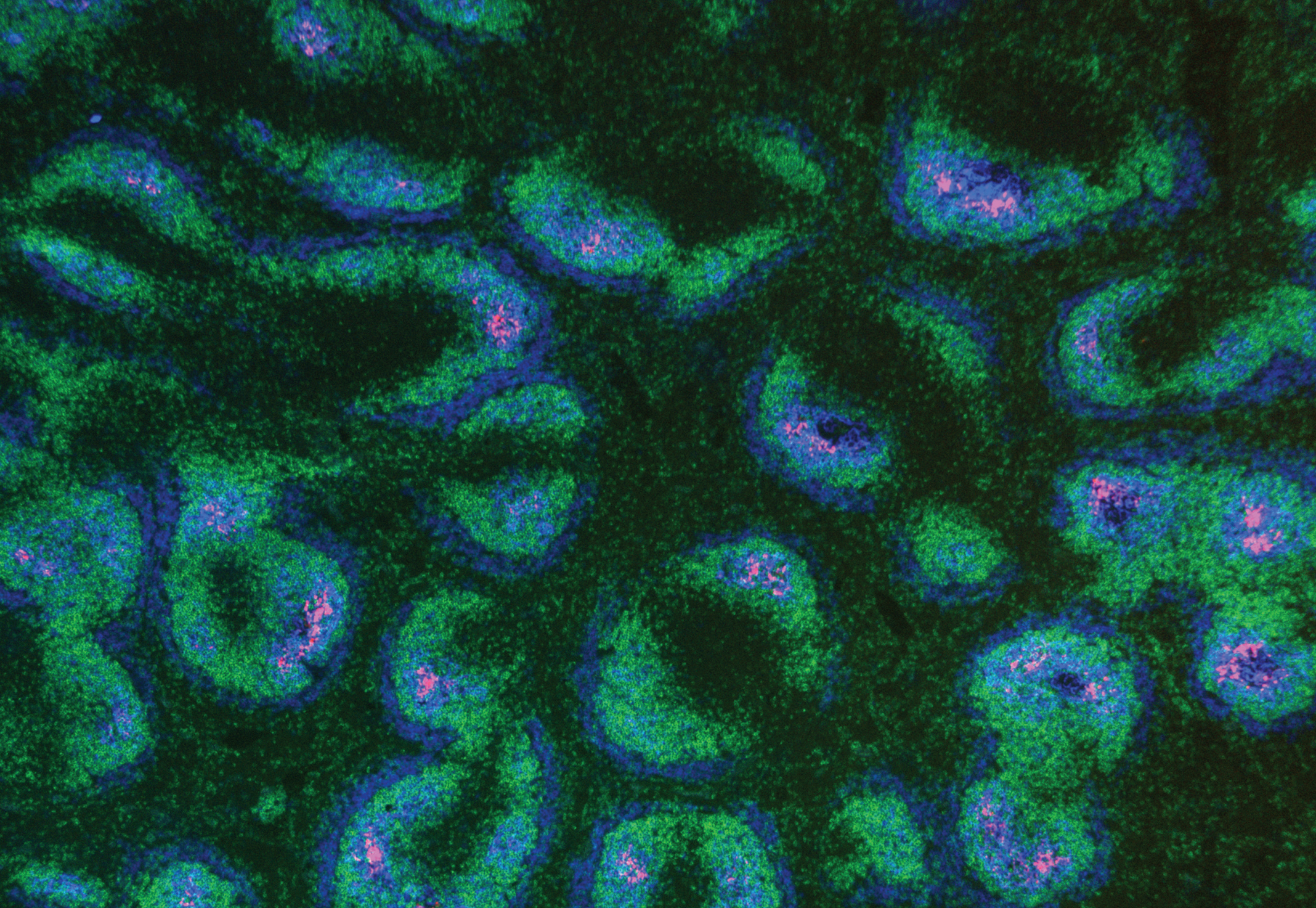
Rachel Keuls

Baylor College of Medicine

Mouse embryo tissue stained with Purified anti-Tubulin β 3 (TUBB3)(orange; Cat. No. 801213), anti-erythroblasts (green), anti-CD31 (magenta), and DAPI (blue).

MARCH 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																										
				FEBRUARY <table border="1"> <thead> <tr> <th>S</th> <th>M</th> <th>T</th> <th>W</th> <th>T</th> <th>F</th> <th>S</th> </tr> </thead> <tbody> <tr> <td></td> <td></td> <td></td> <td></td> <td>1</td> <td>2</td> <td>3</td> </tr> <tr> <td>4</td> <td>5</td> <td>6</td> <td>7</td> <td>8</td> <td>9</td> <td>10</td> </tr> <tr> <td>11</td> <td>12</td> <td>13</td> <td>14</td> <td>15</td> <td>16</td> <td>17</td> </tr> <tr> <td>18</td> <td>19</td> <td>20</td> <td>21</td> <td>22</td> <td>23</td> <td>24</td> </tr> <tr> <td>25</td> <td>26</td> <td>27</td> <td>28</td> <td>29</td> <td></td> <td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S					1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29			1	2
S	M	T	W	T	F	S																																										
				1	2	3																																										
4	5	6	7	8	9	10																																										
11	12	13	14	15	16	17																																										
18	19	20	21	22	23	24																																										
25	26	27	28	29																																												
3	4	5	6	7	8	9																																										
		AD/PD Lisbon, Portugal (until 3/9)																																														
10	11	12	13	14	15	16																																										
Daylight Saving Time Begins			WIRM XVIII Davos, Switzerland (until 3/16)																																													
17	18	19	20	21	22	23																																										
St. Patrick's Day		First Day of Spring																																														
24	25	26	27	28	29	30																																										
31	APRIL <table border="1"> <thead> <tr> <th>S</th> <th>M</th> <th>T</th> <th>W</th> <th>T</th> <th>F</th> <th>S</th> </tr> </thead> <tbody> <tr> <td></td> <td>1</td> <td>2</td> <td>3</td> <td>4</td> <td>5</td> <td>6</td> </tr> <tr> <td>7</td> <td>8</td> <td>9</td> <td>10</td> <td>11</td> <td>12</td> <td>13</td> </tr> <tr> <td>14</td> <td>15</td> <td>16</td> <td>17</td> <td>18</td> <td>19</td> <td>20</td> </tr> <tr> <td>21</td> <td>22</td> <td>23</td> <td>24</td> <td>25</td> <td>26</td> <td>27</td> </tr> <tr> <td>28</td> <td>29</td> <td>30</td> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S		1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30									
S	M	T	W	T	F	S																																										
	1	2	3	4	5	6																																										
7	8	9	10	11	12	13																																										
14	15	16	17	18	19	20																																										
21	22	23	24	25	26	27																																										
28	29	30																																														
Easter																																																



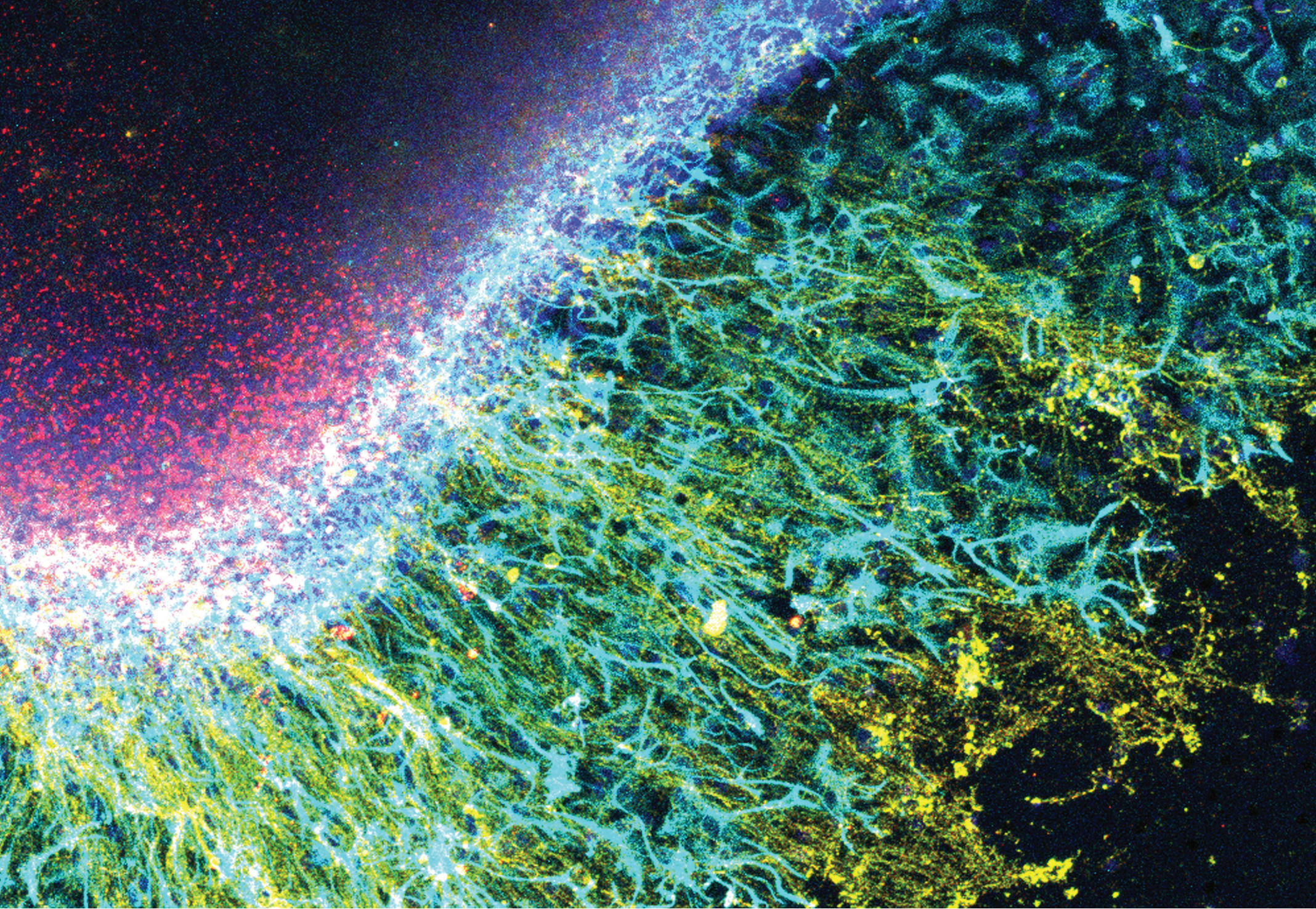
Hanover Matz, PhD

Washington University in St. Louis

Mouse spleen tissue stained with FITC anti-mouse IgD (green; Cat. No. 405704), Brilliant Violet 421™ anti-mouse CD21/CD35 (CR2/CR1)(magenta; Cat. No. 123422), and DAPI (blue).

APRIL 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																																	
MARCH <table border="1"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr> <td></td><td></td><td></td><td></td><td>1</td><td>2</td><td></td> </tr> <tr> <td>3</td><td>4</td><td>5</td><td>6</td><td>7</td><td>8</td><td>9</td> </tr> <tr> <td>10</td><td>11</td><td>12</td><td>13</td><td>14</td><td>15</td><td>16</td> </tr> <tr> <td>17</td><td>18</td><td>19</td><td>20</td><td>21</td><td>22</td><td>23</td> </tr> <tr> <td>24</td><td>25</td><td>26</td><td>27</td><td>28</td><td>29</td><td>30</td> </tr> <tr> <td>31</td><td></td><td></td><td></td><td></td><td></td><td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S					1	2		3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31							1	2	3	4	5	6
S	M	T	W	T	F	S																																																	
				1	2																																																		
3	4	5	6	7	8	9																																																	
10	11	12	13	14	15	16																																																	
17	18	19	20	21	22	23																																																	
24	25	26	27	28	29	30																																																	
31																																																							
7	8	9	10	11	12 AACR San Diego, USA (until 4/10)	13																																																	
14	15	16	17	18	19	20																																																	
21	22	23	24	25	26	27																																																	
28	29	30	MAY <table border="1"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr> <td></td><td></td><td></td><td>1</td><td>2</td><td>3</td><td>4</td> </tr> <tr> <td>5</td><td>6</td><td>7</td><td>8</td><td>9</td><td>10</td><td>11</td> </tr> <tr> <td>12</td><td>13</td><td>14</td><td>15</td><td>16</td><td>17</td><td>18</td> </tr> <tr> <td>19</td><td>20</td><td>21</td><td>22</td><td>23</td><td>24</td><td>25</td> </tr> <tr> <td>26</td><td>27</td><td>28</td><td>29</td><td>30</td><td>31</td><td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S				1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31											
S	M	T	W	T	F	S																																																	
			1	2	3	4																																																	
5	6	7	8	9	10	11																																																	
12	13	14	15	16	17	18																																																	
19	20	21	22	23	24	25																																																	
26	27	28	29	30	31																																																		



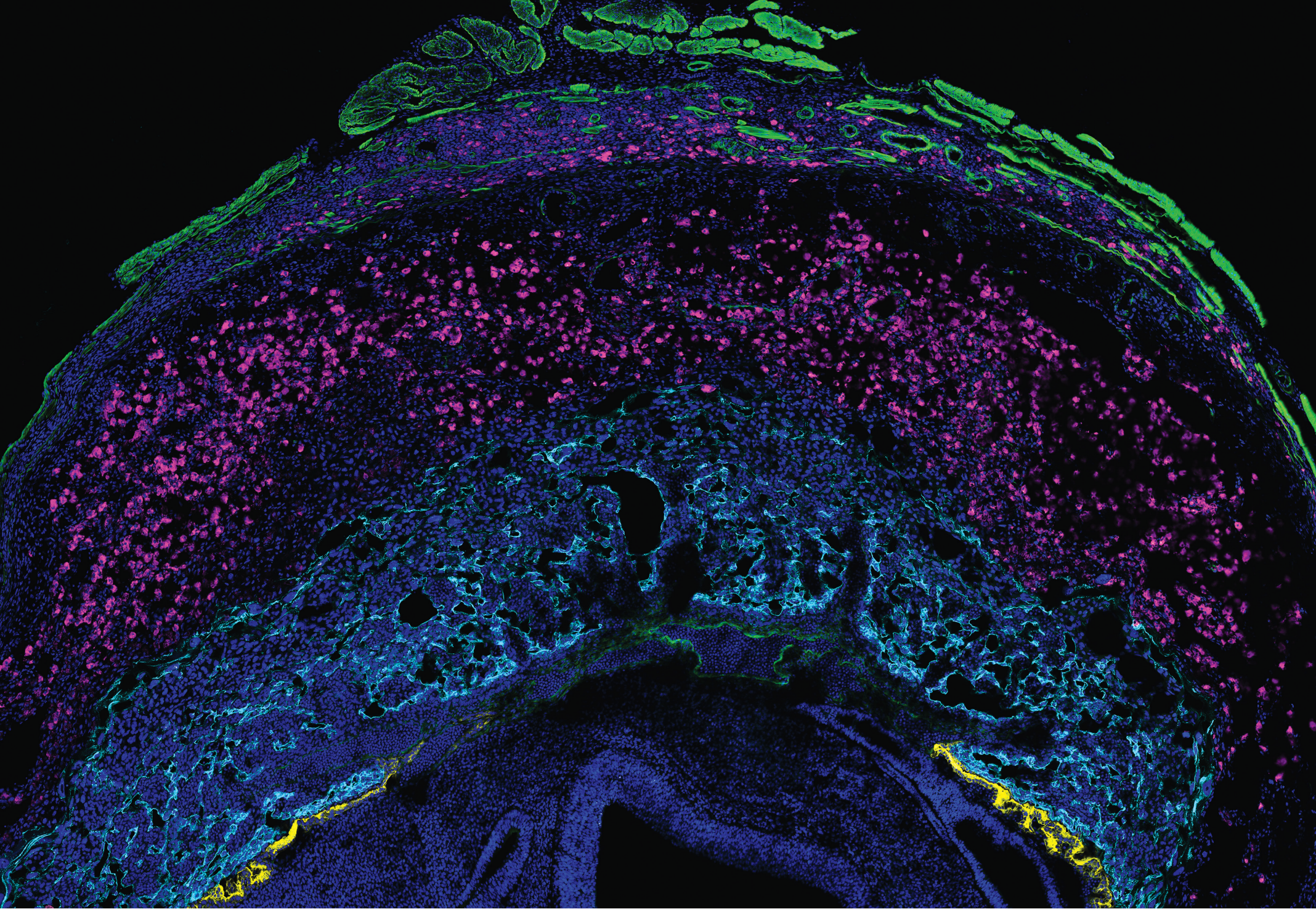
Linyi Chen, PhD

National Tsing Hua University

Embryonic rat brain tissue was sectioned and stained with anti-TUJ1 (green), anti-Tau (red), anti-GFAP (cyan), and DAPI (blue; Cat. No. 422801).

MAY 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																																	
		APRIL <table border="1"> <thead> <tr> <th>S</th> <th>M</th> <th>T</th> <th>W</th> <th>T</th> <th>F</th> <th>S</th> </tr> </thead> <tbody> <tr> <td></td> <td>1</td> <td>2</td> <td>3</td> <td>4</td> <td>5</td> <td>6</td> </tr> <tr> <td>7</td> <td>8</td> <td>9</td> <td>10</td> <td>11</td> <td>12</td> <td>13</td> </tr> <tr> <td>14</td> <td>15</td> <td>16</td> <td>17</td> <td>18</td> <td>19</td> <td>20</td> </tr> <tr> <td>21</td> <td>22</td> <td>23</td> <td>24</td> <td>25</td> <td>26</td> <td>27</td> </tr> <tr> <td>28</td> <td>29</td> <td>30</td> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S		1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30					1	2	3	4							
S	M	T	W	T	F	S																																																	
	1	2	3	4	5	6																																																	
7	8	9	10	11	12	13																																																	
14	15	16	17	18	19	20																																																	
21	22	23	24	25	26	27																																																	
28	29	30																																																					
5	6	7	8	9	10 AAI IMMUNOLOGY2024™ Chicago, IL (until 5/7)	11 CYTO Edinburgh, Scotland (until 5/8)																																																	
12 Mother's Day	13	14	15 CIMT Mainz, Germany (until 5/17)	16	17	18																																																	
19	20	21	22	23	24	25																																																	
26	27 Memorial Day	28	29 ISCT Vancouver, Canada (until 6/1)	30	31	JUNE <table border="1"> <thead> <tr> <th>S</th> <th>M</th> <th>T</th> <th>W</th> <th>T</th> <th>F</th> <th>S</th> </tr> </thead> <tbody> <tr> <td></td> <td></td> <td></td> <td></td> <td></td> <td></td> <td>1</td> </tr> <tr> <td>2</td> <td>3</td> <td>4</td> <td>5</td> <td>6</td> <td>7</td> <td>8</td> </tr> <tr> <td>9</td> <td>10</td> <td>11</td> <td>12</td> <td>13</td> <td>14</td> <td>15</td> </tr> <tr> <td>16</td> <td>17</td> <td>18</td> <td>19</td> <td>20</td> <td>21</td> <td>22</td> </tr> <tr> <td>23</td> <td>24</td> <td>25</td> <td>26</td> <td>27</td> <td>28</td> <td>29</td> </tr> <tr> <td>30</td> <td></td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S							1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30						
S	M	T	W	T	F	S																																																	
						1																																																	
2	3	4	5	6	7	8																																																	
9	10	11	12	13	14	15																																																	
16	17	18	19	20	21	22																																																	
23	24	25	26	27	28	29																																																	
30																																																							

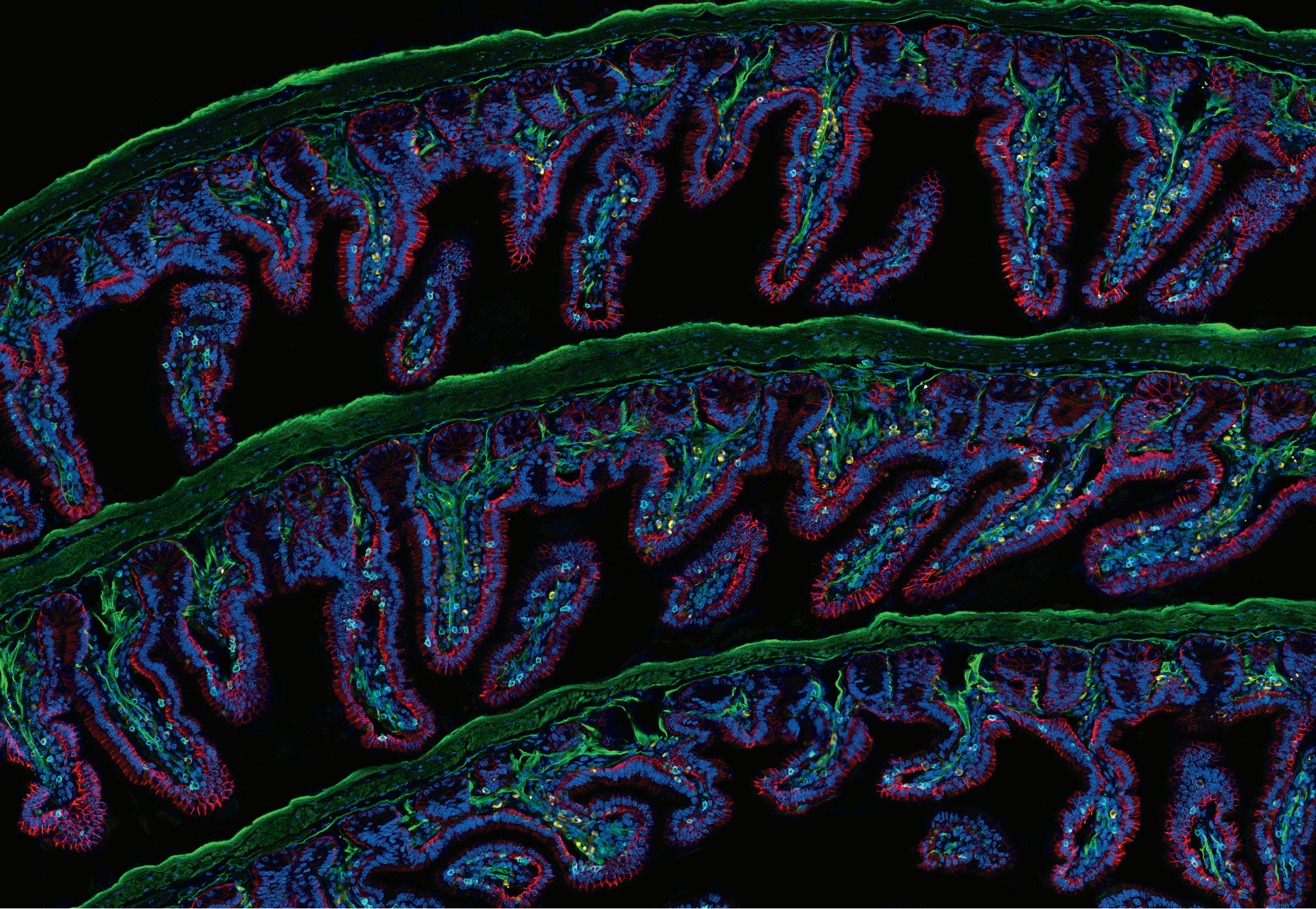


Vitoria Murakami Olyntho
McMaster University

Mouse maternal-fetal interface tissue was stained with Alexa Fluor® 594 anti-mouse CD326 (Ep-CAM)(yellow; Cat. No. 118222), Alexa Fluor® 647 anti-mouse CD138 (Syndecan-1) (cyan; Cat. No. 142526), dolichos biflorus agglutinin (magenta), anti-smooth muscle actin (green), and DAPI (blue).

JUNE 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																										
					<div style="text-align: center; color: #FFC000; font-weight: bold;">MAY</div> <table border="1" style="width: 100%; text-align: center; font-size: small;"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr> <td></td><td></td><td>1</td><td>2</td><td>3</td><td>4</td><td></td> </tr> <tr> <td>5</td><td>6</td><td>7</td><td>8</td><td>9</td><td>10</td><td>11</td> </tr> <tr> <td>12</td><td>13</td><td>14</td><td>15</td><td>16</td><td>17</td><td>18</td> </tr> <tr> <td>19</td><td>20</td><td>21</td><td>22</td><td>23</td><td>24</td><td>25</td> </tr> <tr> <td>26</td><td>27</td><td>28</td><td>29</td><td>30</td><td>31</td><td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S			1	2	3	4		5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31		1
S	M	T	W	T	F	S																																										
		1	2	3	4																																											
5	6	7	8	9	10	11																																										
12	13	14	15	16	17	18																																										
19	20	21	22	23	24	25																																										
26	27	28	29	30	31																																											
2	3	4	5	6	7	8																																										
9	10	11	12	13	14	15																																										
	ISSCR Hamburg, Germany (until 6/13)																																															
16	17	18	19	20	21	22																																										
Father's Day		FOCIS San Francisco, USA (until 6/21)		First Day of Summer																																												
23	24	25	26	27	28	29																																										
30	<div style="text-align: center; color: #FFC000; font-weight: bold;">JULY</div> <table border="1" style="width: 100%; text-align: center; font-size: small;"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr> <td></td><td>1</td><td>2</td><td>3</td><td>4</td><td>5</td><td>6</td> </tr> <tr> <td>7</td><td>8</td><td>9</td><td>10</td><td>11</td><td>12</td><td>13</td> </tr> <tr> <td>14</td><td>15</td><td>16</td><td>17</td><td>18</td><td>19</td><td>20</td> </tr> <tr> <td>21</td><td>22</td><td>23</td><td>24</td><td>25</td><td>26</td><td>27</td> </tr> <tr> <td>28</td><td>29</td><td>30</td><td>31</td><td></td><td></td><td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S		1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31								
S	M	T	W	T	F	S																																										
	1	2	3	4	5	6																																										
7	8	9	10	11	12	13																																										
14	15	16	17	18	19	20																																										
21	22	23	24	25	26	27																																										
28	29	30	31																																													

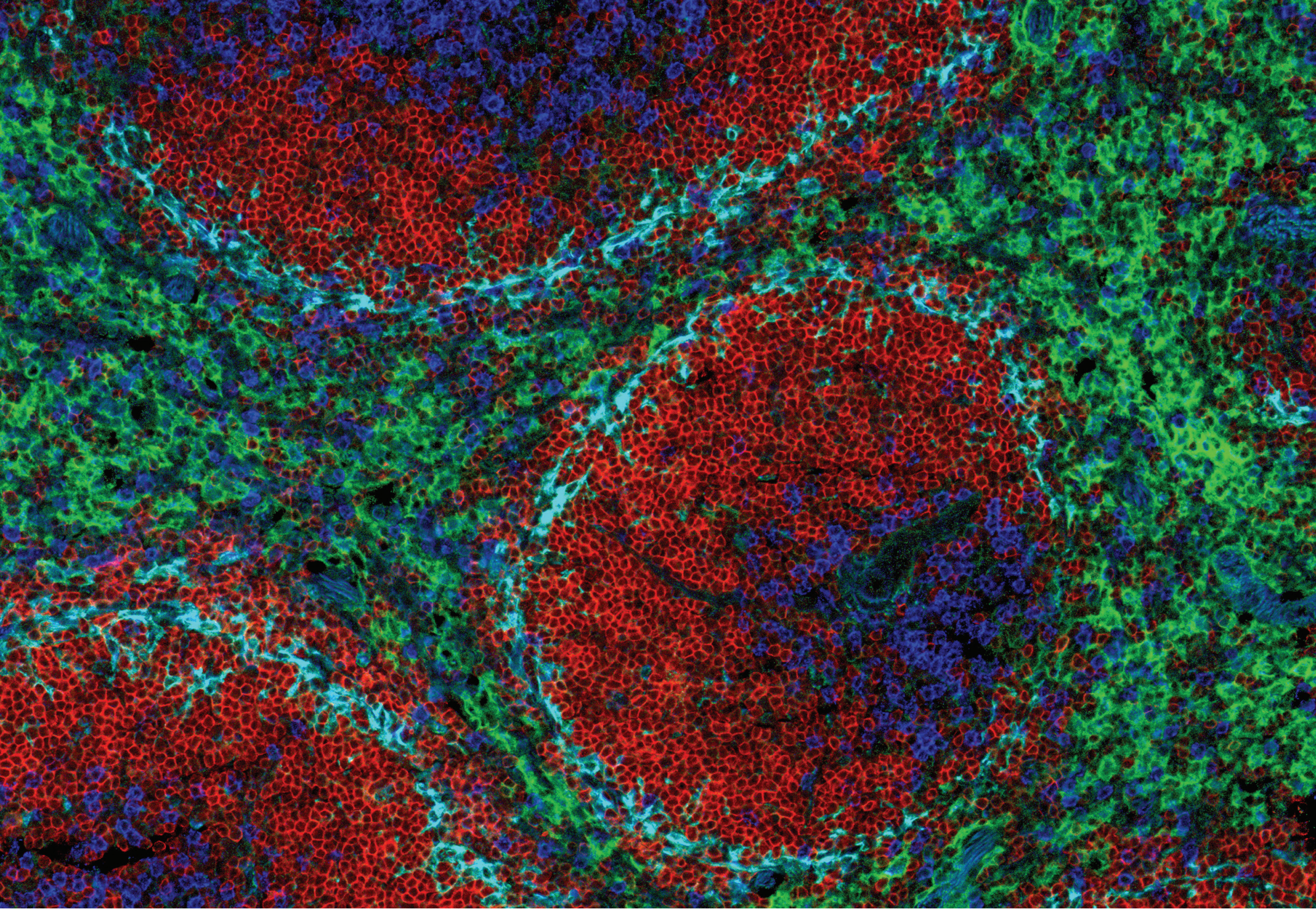


Vitoria Murakami Olyntho
McMaster University

Mouse intestinal tissue in a food-allergic mouse was stained with Alexa Fluor® 488 anti-mouse CD3 (cyan; Cat. No. 100210), Alexa Fluor® 594 anti-mouse CD326 (Ep-CAM)(red; Cat. No. 118222), anti-ST2 (yellow), anti-smooth muscle actin (green), and DAPI (blue).

JULY 2024

							Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday						
JUNE								1	2	3	4	5	6						
S	M	T	W	T	F	S													
						1													
2	3	4	5	6	7	8													
9	10	11	12	13	14	15													
16	17	18	19	20	21	22													
23	24	25	26	27	28	29													
30										Independence Day									
						7	8	9	10	11	12	13							
						14	15	16	17	18	19	20							
						21	22	23	24	25	26	27							
						28	29	30	31	AUGUST									
										S	M	T	W	T	F	S			
													1	2	3				
													4	5	6	7	8	9	10
													11	12	13	14	15	16	17
													18	19	20	21	22	23	24
													25	26	27	28	29	30	31



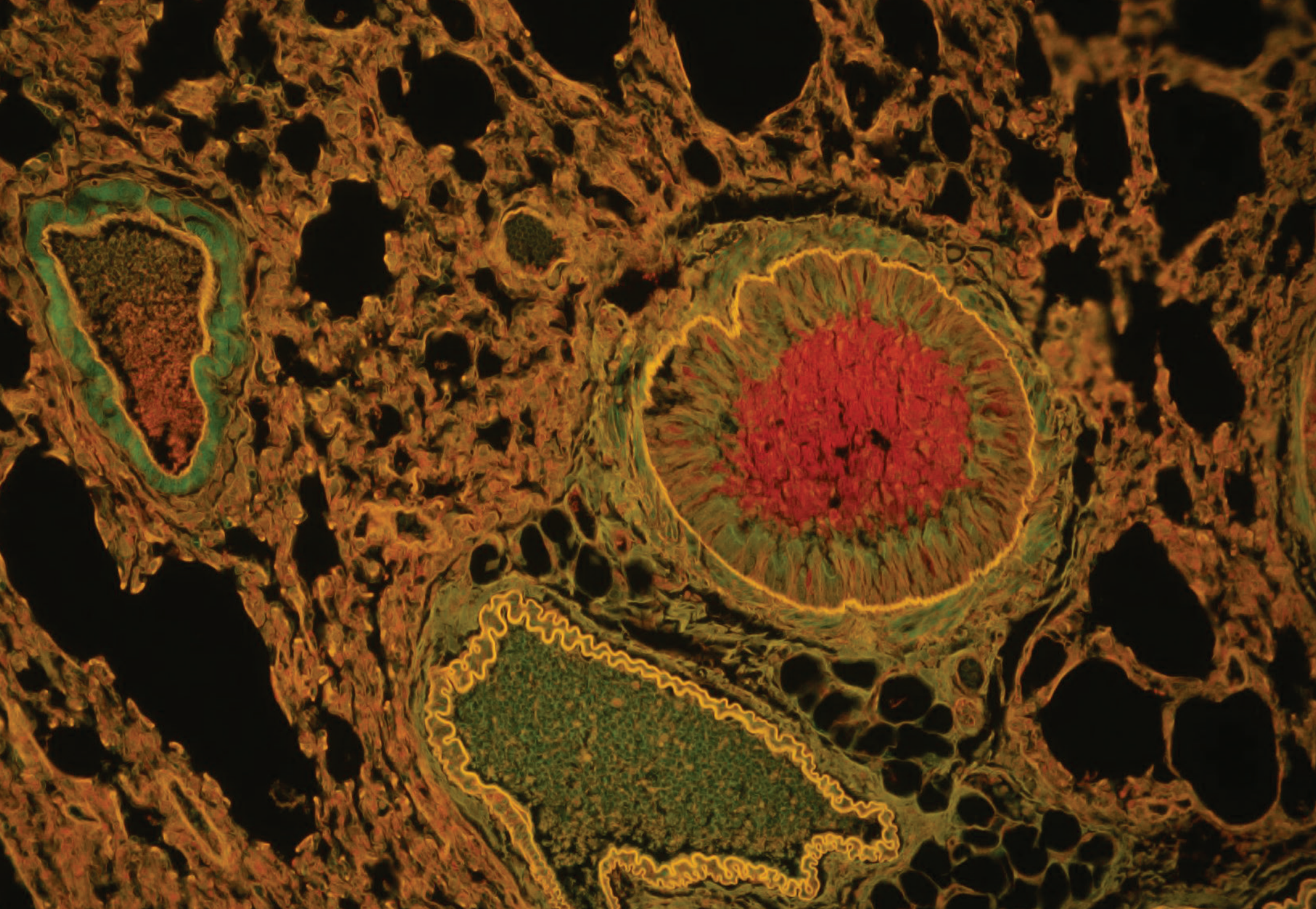
Kianoosh Mempel

University of California, San Diego

Naive mouse spleen tissue was stained with Alexa Fluor® 647 anti-mouse CD169 (Siglec-1)(cyan; Cat. No. 142408), Alexa Fluor® 594 anti-mouse/human CD45R/B220 (red; Cat. No. 103254), FITC anti-mouse F4/80 (green; Cat. No. 123107), and anti-CD8a (blue).

AUGUST 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday
		JULY S M T W T F S 1 2 3 4 5 6 7 8 9 10 11 12 13 14 15 16 17 18 19 20 21 22 23 24 25 26 27 28 29 30 31	SEPTEMBER S M T W T F S 1 2 3 4 5 6 7 8 9 10 11 12 13 14 15 16 17 18 19 20 21 22 23 24 25 26 27 28 29 30	1	2	3
4	5	6	7	8	9	10
11	12	13	14	15	16	17
18	19	20	21	22	23	24
25	26	27	28	29	30	31



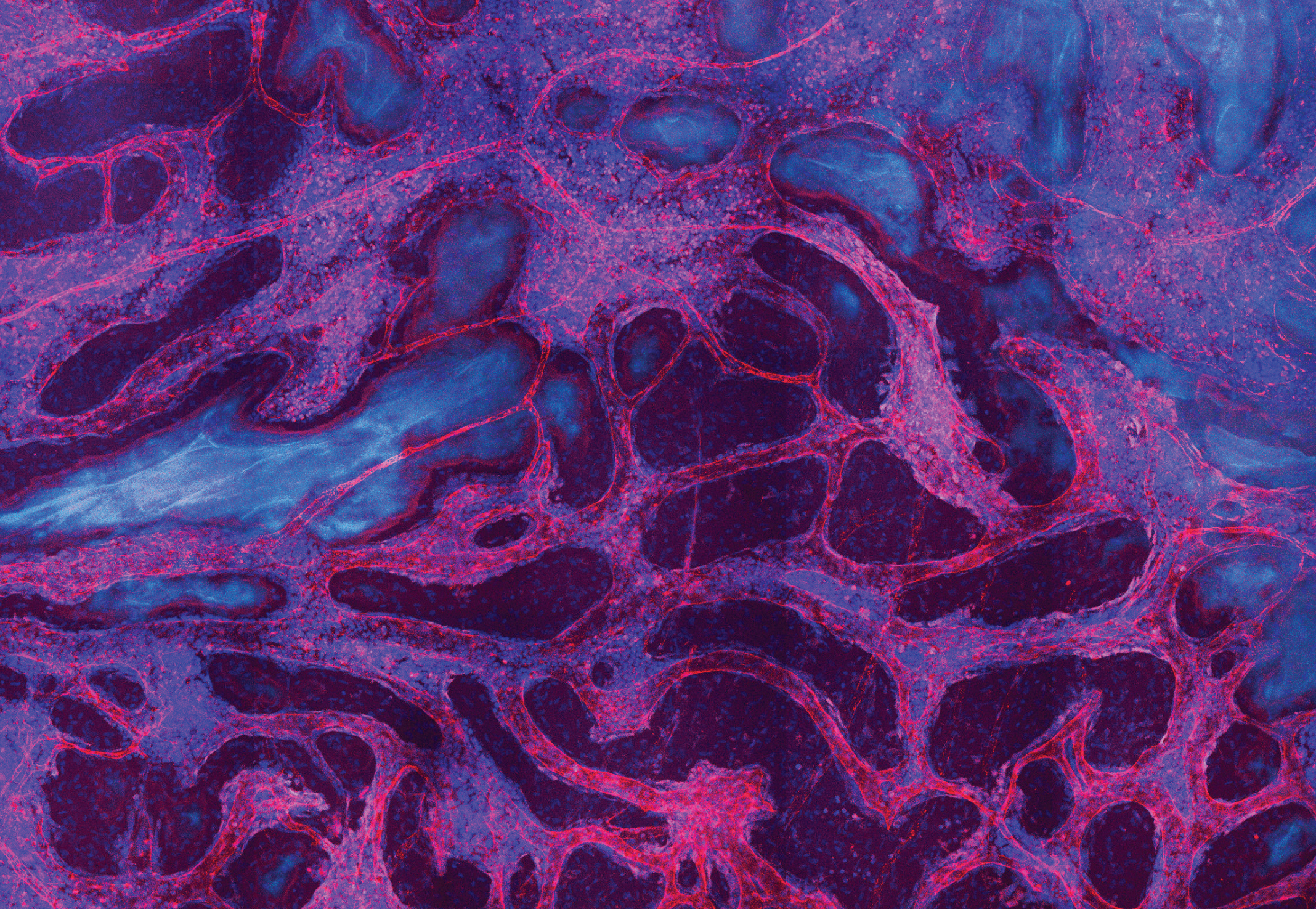
Valerie Oren

The University of Texas MD Anderson Cancer Center

Mucus metaplasia was stimulated in mouse lung tissue using recombinant mouse IL-1 β (Cat. No. 575108). Staining was done with periodic acid fluorescent Schiff reagent (red). Tissue autofluorescence is present in green.

SEPTEMBER 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																																																																																		
1 ECI Dublin, Ireland (until 9/4)	2 Labor Day	3	4	5	6	7																																																																																																		
8	9	10	11 Patriot Day	12	13	14																																																																																																		
15	16	17	18	19	20	21																																																																																																		
22 First Day of Autumn	23	24	25	26	27	28																																																																																																		
29	30	<table border="1"> <thead> <tr> <th colspan="7">AUGUST</th> <th colspan="7">OCTOBER</th> </tr> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr> <td></td><td></td><td></td><td></td><td>1</td><td>2</td><td>3</td> <td></td><td></td><td>1</td><td>2</td><td>3</td><td>4</td><td>5</td> </tr> <tr> <td>4</td><td>5</td><td>6</td><td>7</td><td>8</td><td>9</td><td>10</td> <td>6</td><td>7</td><td>8</td><td>9</td><td>10</td><td>11</td><td>12</td> </tr> <tr> <td>11</td><td>12</td><td>13</td><td>14</td><td>15</td><td>16</td><td>17</td> <td>13</td><td>14</td><td>15</td><td>16</td><td>17</td><td>18</td><td>19</td> </tr> <tr> <td>18</td><td>19</td><td>20</td><td>21</td><td>22</td><td>23</td><td>24</td> <td>20</td><td>21</td><td>22</td><td>23</td><td>24</td><td>25</td><td>26</td> </tr> <tr> <td>25</td><td>26</td><td>27</td><td>28</td><td>29</td><td>30</td><td>31</td> <td>27</td><td>28</td><td>29</td><td>30</td><td>31</td><td></td><td></td> </tr> </tbody> </table>		AUGUST							OCTOBER							S	M	T	W	T	F	S	S	M	T	W	T	F	S					1	2	3			1	2	3	4	5	4	5	6	7	8	9	10	6	7	8	9	10	11	12	11	12	13	14	15	16	17	13	14	15	16	17	18	19	18	19	20	21	22	23	24	20	21	22	23	24	25	26	25	26	27	28	29	30	31	27	28	29	30	31					
AUGUST							OCTOBER																																																																																																	
S	M	T	W	T	F	S	S	M	T	W	T	F	S																																																																																											
				1	2	3			1	2	3	4	5																																																																																											
4	5	6	7	8	9	10	6	7	8	9	10	11	12																																																																																											
11	12	13	14	15	16	17	13	14	15	16	17	18	19																																																																																											
18	19	20	21	22	23	24	20	21	22	23	24	25	26																																																																																											
25	26	27	28	29	30	31	27	28	29	30	31																																																																																													

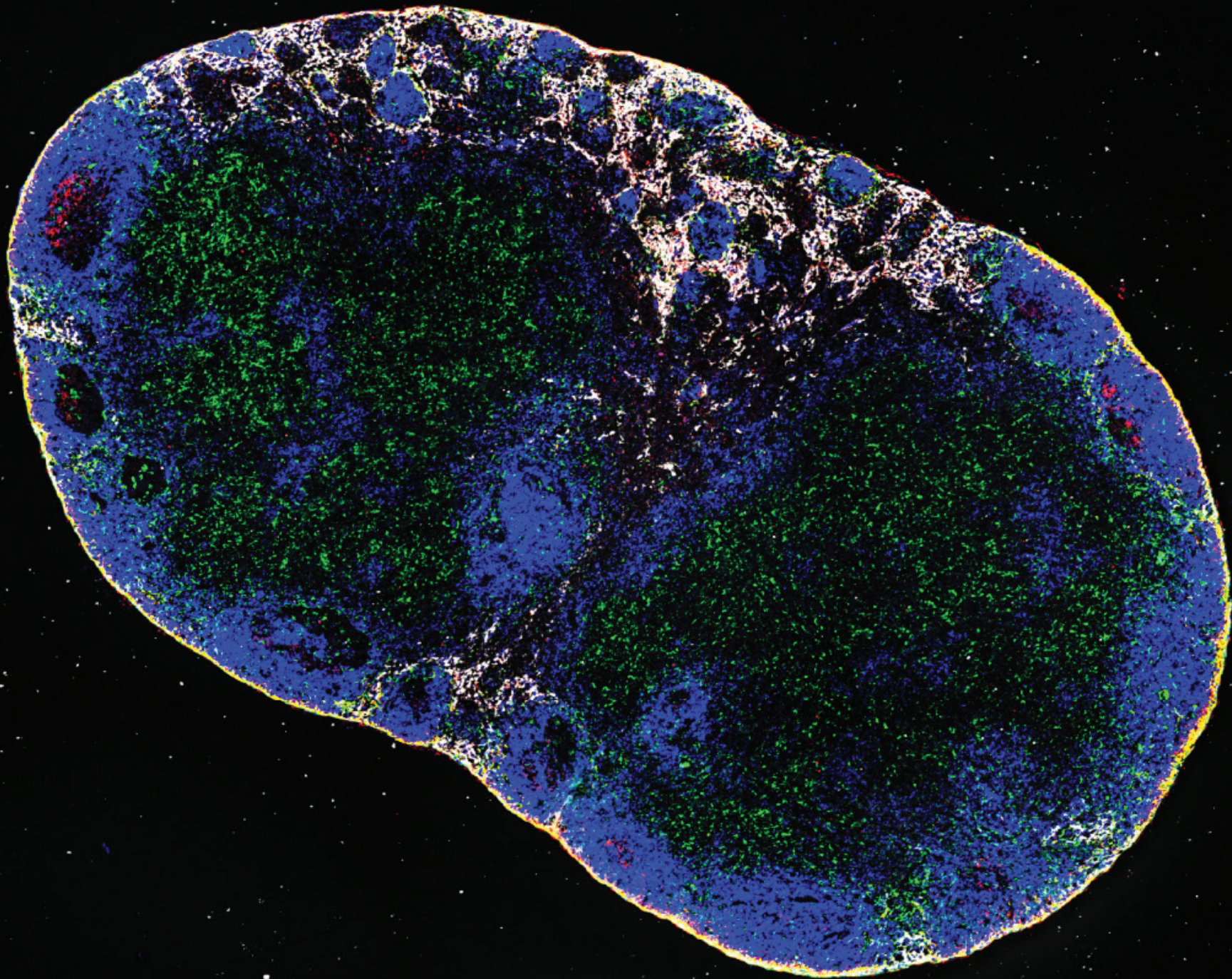


Désirée Zerbst
University of Glasgow

Mouse bone marrow stained with Alexa Fluor® 594 anti-mouse CD31 (magenta; Cat. No. 102432) and DAPI (blue). Bone tissue autofluorescence is shown in cyan.

OCTOBER 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday																																										
	<p>SEPTEMBER</p> <table border="1"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr> <td>1</td><td>2</td><td>3</td><td>4</td><td>5</td><td>6</td><td>7</td> </tr> <tr> <td>8</td><td>9</td><td>10</td><td>11</td><td>12</td><td>13</td><td>14</td> </tr> <tr> <td>15</td><td>16</td><td>17</td><td>18</td><td>19</td><td>20</td><td>21</td> </tr> <tr> <td>22</td><td>23</td><td>24</td><td>25</td><td>26</td><td>27</td><td>28</td> </tr> <tr> <td>29</td><td>30</td><td></td><td></td><td></td><td></td><td></td> </tr> </tbody> </table>	S	M	T	W	T	F	S	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30						1	2	3	4	5
S	M	T	W	T	F	S																																										
1	2	3	4	5	6	7																																										
8	9	10	11	12	13	14																																										
15	16	17	18	19	20	21																																										
22	23	24	25	26	27	28																																										
29	30																																															
6	7	8	9	10	11	12 Neuroscience 2024 Chicago, USA (until 10/9)																																										
13	14	15	16	17	18	19																																										
20	21	22	23	24	25	26																																										
ICIS/Cytokines Seoul, Korea (until 10/23)																																																
27	28	29	30	31	<p>NOVEMBER</p> <table border="1"> <thead> <tr> <th>S</th><th>M</th><th>T</th><th>W</th><th>T</th><th>F</th><th>S</th> </tr> </thead> <tbody> <tr> <td></td><td></td><td></td><td></td><td></td><td>1</td><td>2</td> </tr> <tr> <td>3</td><td>4</td><td>5</td><td>6</td><td>7</td><td>8</td><td>9</td> </tr> <tr> <td>10</td><td>11</td><td>12</td><td>13</td><td>14</td><td>15</td><td>16</td> </tr> <tr> <td>17</td><td>18</td><td>19</td><td>20</td><td>21</td><td>22</td><td>23</td> </tr> <tr> <td>24</td><td>25</td><td>26</td><td>27</td><td>28</td><td>29</td><td>30</td> </tr> </tbody> </table>	S	M	T	W	T	F	S						1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	
S	M	T	W	T	F	S																																										
					1	2																																										
3	4	5	6	7	8	9																																										
10	11	12	13	14	15	16																																										
17	18	19	20	21	22	23																																										
24	25	26	27	28	29	30																																										
				Halloween																																												



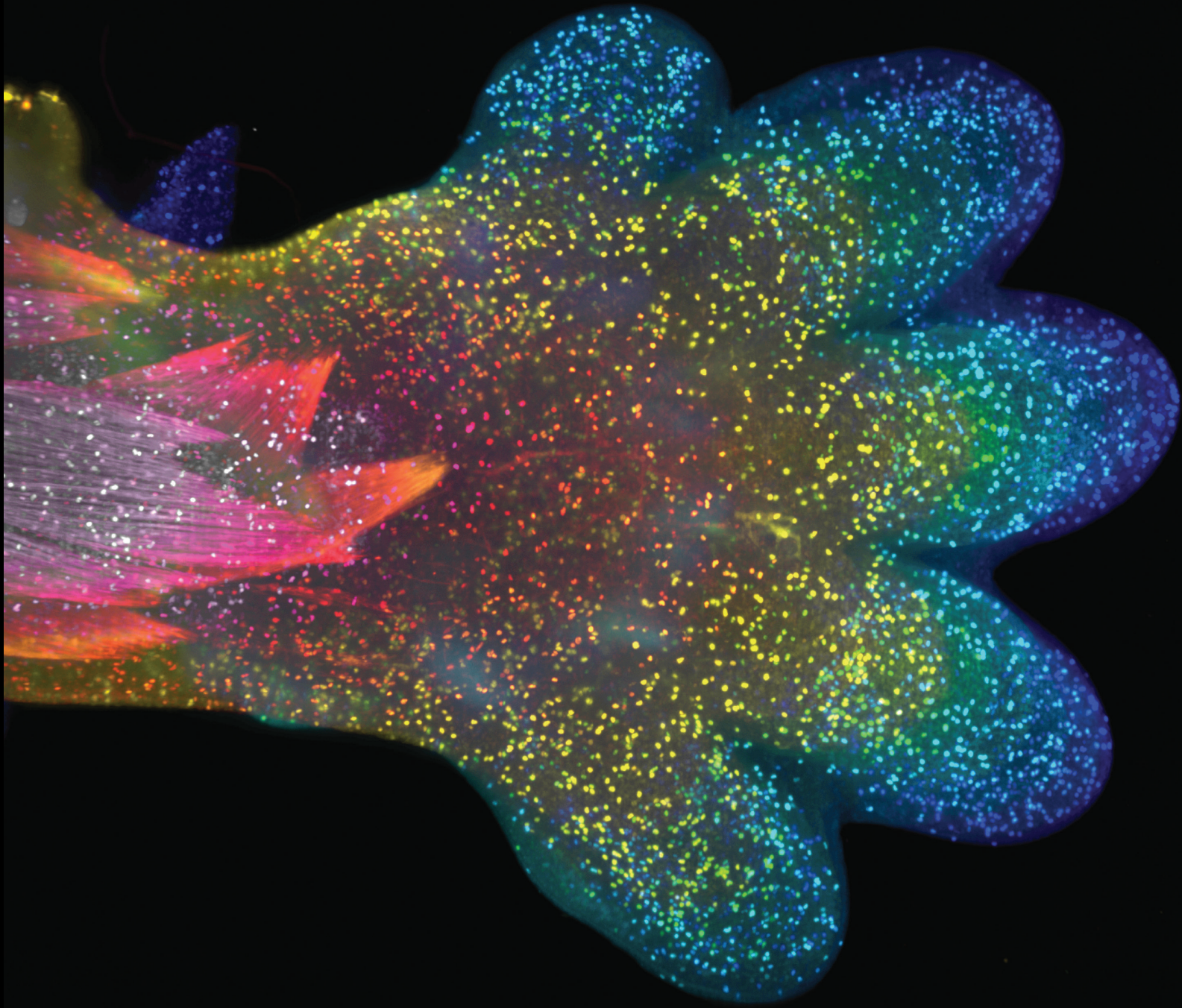
Yuan Wenchen

Chongqing International Institute for Immunology

Mouse mesenteric lymph node tissue stained with Alexa Fluor® 647 anti-mouse CD169 (Siglec-1)(white; Cat. No. 142408), anti-CD209b (red), anti-CD68 (green), and anti-IgD (blue).

NOVEMBER 2024

Sunday	Monday	Tuesday	Wednesday	Thursday	Friday	Saturday
			OCTOBER S M T W T F S 1 2 3 4 5 6 7 8 9 10 11 12 13 14 15 16 17 18 19 20 21 22 23 24 25 26 27 28 29 30 31	DECEMBER S M T W T F S 1 2 3 4 5 6 7 8 9 10 11 12 13 14 15 16 17 18 19 20 21 22 23 24 25 26 27 28 29 30 31	1	2
3	4	5	6	7	8	9
Daylight Saving Time Ends		ASHG Denver, USA (until 11/9)	SITC Houston, USA (until 11/10)			
10	11	12	13	14	15	16
	Veterans Day					
17	18	19	20	21	22	23
24	25	26	27	28	29	30
				Thanksgiving Day		



Oswaldo Contreras, PhD

Victor Chang Cardiac Research Institute

Embryonic mouse paw stained with Alexa Fluor® 594 anti-Histone H3 Phospho (Ser10)(red; Cat. No. 650810) and anti- α smooth muscle actin (cyan). The resulting 3D image was color coded by depth.

