

PerkinElmer NEN[®] Radiochemicals 2019 Fresh Lot Schedules

	JANUARY					FEBRUARY					MARCH					APRIL					MAY					JUNE																
	S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S	S	M	T	W	T	F	S
³²P Radionuclides					3							7							7							4							2							6		
<i>Every Thursday</i>					10							14							14							11							9							13		
(all part numbers with exceptions below)					17							21							21							18							16							20		
					24							28							28							25							23							27		
					31																												30									
³²P Radionuclides					3							7							7							4							2							6		
<i>Every First Thursday</i>																																										
BLU023X, NEG023X																																										
³²P Radionuclides					10							14							14							11							9							13		
<i>Second Thursday of Every Month</i>																																										
BLU004Z, NEG004Z, BLU504Z, NEG504Z, BLU019A, NEG019A																																										
³⁵S Radionuclides					2							5							5							2							7							4		
<i>Every First Tuesday</i>																																										
³⁵S Amino Acids and Glutathione					21	2						11							4	25						15							6	28						17		
<i>Every Third Monday</i>																																										
NEG009A, NEG009C, NEG009H, NEG009L, NEG009T, NEG709A, NEG022T, NEG072, NEG772, NEG029																																										
³²P Radionuclides					2	9						6	13						6	13						3	10						1	8						5	12	
<i>Every Wednesday</i>					16	23						20	27						20	27						17	24						15	22						19	26	
NEX053, NEX054, NEX011, NEX053H, NEX053S, NEX055, NEX060, NEX063					30																												29									
¹²⁵I Radionuclides					9	23						6	20						6	20						3	17						1	15	29					12	26	
<i>Every Other Wednesday</i>																																										
NEZ033, NEZ033A, NEZ033H, NEZ033L																																										
¹³¹I Radionuclides					7	21						4	18						4	18						1	15	29					13	28						10	24	
<i>Every Other Monday</i>																																										
NEZ035A, NEZ035H																																										
⁵¹Cr Radionuclides					4	18						1	15						1	15	29					12	26						10	24						7	21	
<i>Every Other Friday</i>																																										
NEZ030, NEZ030S																																										
⁴⁵Ca Radionuclides					28	2						25							25							22							20							17		
<i>Every Fourth Monday</i>																																										
NEZ013																																										
³⁵S Radionuclides					28	2						25							25							22							20							17		
<i>Every Fourth Monday</i>																																										
NEX041, NEX041H, NEX042																																										

