

PerkinElmer NEN® Radiochemicals 2020 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 Radionucleotides						2							6							5							2							7							4	
<i>Every Thursday</i>						9							13							12							9							14							11	
(all part numbers with exceptions below)						16							20							19							16							21							18	
						23							27							26							23							28							25	
						30																					30															
³²P Radionucleotides						2							6							5							2							7							4	
<i>Every First Thursday</i>																																										
BLU023X, NEG023X																																										
³²P Radionucleotides						9							13							12							9							14							11	
<i>Every Second Thursday</i>																																										
BLU004Z, NEG004Z, BLU504Z, NEG504Z, BLU019A, NEG019A																																										
³⁵S Radionucleotides						7							4							3							7							5							2	
<i>Every First Tuesday</i>																																										
³⁵S Amino Acids						13							3							16							6							18							8	
<i>Every Third Monday</i>													24														27														29	
NEG009A, NEG009C, NEG009H, NEG009L, NEG009T, NEG709A, NEG022T, NEG072, NEG772, and NEG029																																										
³²P Radionuclides						2							5							4							1							6							3	
<i>Every Wednesday</i>						8							12							11							8							13							10	
NEX053, NEX054, NEX011, NEX053H, NEX053S, NEX055, NEX060, NEX063						15							19							18							15							20							17	
						22							26							25							22							27							24	
						29																					29															
¹²⁵I Radionuclides						8							5							4							1							13							10	
<i>Every Other Wednesday</i>						22							19							18							15							27							24	
NEZ033, NEZ033A, NEZ033H, NEZ033L																											29															
¹³¹I Radionuclides						6							3							2							13							11							8	
<i>Every Other Monday</i>						20							18							16							27							26							22	
NEZ035A, NEZ035H																				30																						
⁵¹Cr Radionuclides						3							14							13							10							8							5	
<i>Every Other Friday</i>						17							28							27							24							22							19	
NEZ030, NEZ030S						31																																				
⁴⁵Ca Radionuclides						27							24							23							20							18							15	
<i>Every Fourth Monday</i>																																										
NEZ013																																										
³⁵S Radionuclides						27							24							23							20							18							15	
<i>Every Fourth Monday</i>																																										
NEX041, NEX041H, NEX042																																										

PerkinElmer NEN® Radiochemicals 2020 Fresh Lot Schedules

	JULY							AUGUST							SEPTEMBER							OCTOBER							NOVEMBER							DECEMBER													
	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	S	M	T	W	T	F	S
³²P Radionucleotides					2							6							3							1							5									3							
<i>Every Thursday</i>					9							13							10							8							12									10							
(all part numbers with exceptions below)					16							20							17							15							19									17							
					23							27							24							22							25									-							
					20																		29									31																	
³²P Radionucleotides					2							6							3							1							5									3							
<i>Every First Thursday</i>																																																	
BLU023X, NEG023X																																																	
³²P Radionucleotides					9							13							10							8							12									10							
<i>Every Second Thursday</i>																																																	
BLU004Z, NEG004Z, BLU504Z, NEG504Z, BLU019A, NEG019A																																																	
³⁵S Radionucleotides					7							4							1							6							3							1									
<i>Every First Tuesday</i>																																																	
³⁵S Amino Acids					20							10							21							12							2							14									
<i>Every Third Monday</i>												31																	23							4-Jan													
NEG009A, NEG009C, NEG009H, NEG009L, NEG009T, NEG709A, NEG022T, NEG072, NEG772, and NEG02																																																	
³²P Radionuclides					1							5							2							7							4							2									
<i>Every Wednesday</i>					8							12							9							14							11							9									
NEX053, NEX054, NEX011, NEX053H, NEX053S, NEX055, NEX060, NEX063					15							19							16							21							18							16									
					22							26							23							28							25							23									
					29													30																						30									
¹²⁵I Radionuclides					8							5							2							14							11							9									
<i>Every Other Wednesday</i>					22							19							16							28							25							23									
NEZ033, NEZ033A, NEZ033H, NEZ033L																		30																						6-Jan									
¹³¹I Radionuclides					6							3							14							12							9							7									
<i>Every Other Monday</i>					20							17							28							26							23							21									
NEZ035A, NEZ035H												31																											4-Jan										
⁵¹Cr Radionuclides					6													11							9							6									4								
<i>Every Other Friday</i>																		25							23							20									18								
NEZ030, NEZ030S																																						4-Jan											
⁴⁵Ca Radionuclides					13							10							8							5							2							28									
<i>Every Fourth Monday</i>																																	30							25-Jan									
NEZ013																																																	
³⁵S Radionuclides					13							10							8							5							2							28									
<i>Every Fourth Monday</i>																																	30							25-Jan									
NEX041, NEX041H, NEX042																																																	