



# Revvity NEN<sup>®</sup> Radiochemicals 2025 Fresh Lot Schedule

	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	S	M	T	W	T	F	S							
<b><sup>32</sup>P Radionucleotides</b> <i>Every Thursday</i>					2							6							6							3							10							1							8							5		
					9							13							13							17							22							29							12									
					16							20							20							24							15							19																
					23							27							27														22							26																
					30																																																			
<b><sup>35</sup>S Radionucleotides</b> <i>Every first Tuesday</i>						7						4							4							1							6							3																
<b><sup>35</sup>S Amino Acids</b> <i>Monday, every three weeks</i> NEG009A, NEG009C, NEG009H, NEG009L, NEG009T, NEG709A, NEG022T, NEG072, and NEG772						13						3							17							7							19							9																
<b><sup>32</sup>P Radionuclides</b> <i>Every Wednesday</i> NEX053, NEX053H, NEX060							2					5							5							2							7							4																
							8					12							12							9							14							11																
							15					19							19							16							21							18																
							22					26							26							23							28							25																
							29																	30																																
<b><sup>125</sup>I Radionuclides</b> <i>Every other Wednesday</i> NEZ033, NEZ033A, NEZ033H, NEZ033L							2					12							12							9							7							4																
							15					26							26							23							21							18																
							29																																																	
<b><sup>51</sup>Cr Radionuclides</b> <i>Every other Friday</i> NEZ030, NEZ030S							10					7							7							4							2							13																
							24					21							21							18							16							27																
<b><sup>35</sup>S Radionuclides</b> <i>Every fourth Monday</i> NEX041, NEX041H, NEX042							21					18							17							14							12							9																

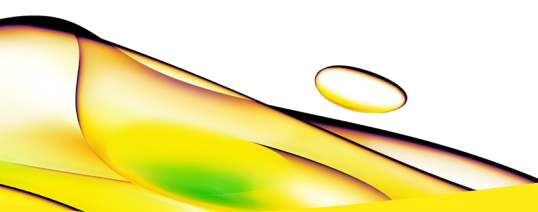
Skipped lot - December 24

■ = Holiday exception

	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
<b><sup>32</sup>P Radionucleotides</b> <i>Every Thursday</i>					3							7							4							2							6							4									
					10							14							11							9							13							11									
					17							20							18							16							20							18									
					24							28							25							23							26							*									
					31																		30												31							08-Jan							
<b><sup>35</sup>S Radionucleotides</b> <i>Every first Tuesday</i>						1							5							2							7							4							2								
																											7								4						06-Jan								
<b><sup>35</sup>S Amino Acids</b> <i>Monday, every three weeks</i> NEG009A, NEG009C, NEG009H, NEG009L, NEG009T, NEG709A, NEG022T, NEG072, and NEG772						21							11							2							13							3							15								
																																	24							05-Jan									
<b><sup>32</sup>P Radionuclides</b> <i>Every Wednesday</i> NEX053, NEX053H, NEX060					2							6							3							1							5							3									
					9							13							10							8							12							10									
					16							20							17							15							19							17									
					23							27							24							22							26							24									
					30																		29																	31									
																																												07-Jan					
<b><sup>125</sup>I Radionuclides</b> <i>Every other Wednesday</i> NEZ033, NEZ033A, NEZ033H, NEZ033L					2							13							10							8							5								3								
					16							27							24							22							19							17									
					30																																	31											
<b><sup>51</sup>Cr Radionuclides</b> <i>Every other Friday</i> NEZ030, NEZ030S						11							8						5							3							14							1		12							
						25							22						19							17												29											
<b><sup>35</sup>S Radionuclides</b> <i>Every fourth Monday</i> NEX041, NEX041H, NEX042						7							4							2							27							24							22								
																																												20-Jan					

Skipped lot - December 24

■ = Holiday exception



revvity