



Revvity NEN[®] Radiochemicals 2024 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																																																																								
³²P Radionucleotides <i>Every Thursday</i>					4								1												7												4													2														9															16																23																	27									
³⁵S Radionucleotides <i>Every first Tuesday</i>							2						6												5													2														7																		4																																																			
³⁵S Amino Acids <i>Monday, every three weeks</i> NEG009A, NEG009C, NEG009H, NEG009L, NEG009T, NEG709A, NEG022T, NEG072, and NEG772													12												4												15														6																		17																																																				
³²P Radionuclides <i>Every Wednesday</i> NEX053, NEX053H, NEX060						3							7												6												3														1															8																15																	22																		29				
¹²⁵I Radionuclides <i>Every other Wednesday</i> NEZ033, NEZ033A, NEZ033H, NEZ033L						3							14												13												10														8															22																		5																																					
⁵¹Cr Radionuclides <i>Every other Friday</i> NEZ030, NEZ030S							12						9												8												5														3															17																		14																																					
³⁵S Radionuclides <i>Every fourth Monday</i> NEX041, NEX041H, NEX042							22							20											18												15														13																		10																																																				

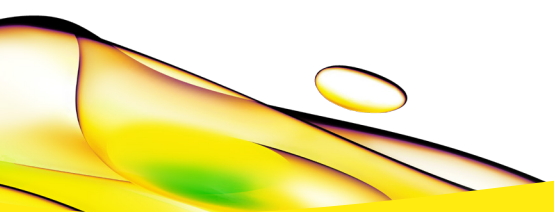
* Skipped lot

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

* Skipped lot

■ = Holiday exception



revvity