

Release Schedule for Select ³H and ¹⁴C Products

Part Number	Description	Availability
¹⁴C Labeled Compounds		
NEC164A050UC	Inulin, Carboxyl-[Carboxyl- ¹⁴ C]-	Third Monday in July
NEC164A250UC		
NEC555010UC	Palmitoyl Coenzyme A, [Palmitoyl-1- ¹⁴ C]-	Third Monday in July
NEC555050UC		
NEC501050UC	Linoleic Acid, [1- ¹⁴ C]-	First Monday in November
NEC501250UC		

Tritium Labeled Compounds		
NET048250UC	Norepinephrine Hydrochloride, DL-[7- ³ H(N)]-	Third Monday in July
NET063001MC	Adenine, [2, 8- ³ H]-	Second Monday in March and September
NET063005MC		
NET1019025UC	Leukotriene D ₄ , [14, 15, 19, 20- ³ H(N)]-	Third Monday in March and September
NET1054025UC	Cytisine HCl, [3, 5- ³ H(N)]-	Second Monday in April
NET1054250UC		
NET1100050UC	Lysophosphatidic Acid, 1-Oleoyl-[Oleoyl-9,10- ³ H]-	Second Monday in November
NET1122001MC	Phenylalanine, L-[2,3,4,5,6- ³ H]-	First Monday in January, May and September
NET1140001MC	Astemizole, [O-Methyl- ³ H]-	Second Monday in April and October
NET1140025UC		
NET1140250UC		
NET1164250UC	WAY100635, [Methoxy- ³ H]-	Second Monday in February and August
NET1165050UC	Quisqualic Acid, [³ H]-	Third Monday in September
NET1167250UC	5-Hydroxy Tryptamine, [³ H]-	First Monday in January, May and September
NET1168001MC	Myo-Inositol, [³ H]-	Second Monday in March and September
NET1177001MC		
NET1169250UC	R(+)-7-HYDROXY DPAT, [³ H]-	Second Monday in April and October
NET1170050UC	α-Bungarotoxin, N-[³ H]-	Second Monday in August
NET1172U100UC	GR125743, [N-Methyl- ³ H]-	Third Monday in February and August
NET1182001MC	Gabapentin, [³ H]-	Third Monday in February and September
NET1198U250UC	SB-269970, [³ H]-	First Monday in June
NET1201001MC	Guanosine 5'-Triphosphate, [8- ³ H]-	Third Monday in April
NET1215001MC	Biotin, D-[³ H(G)]-	First Monday in May and November
NET1215250UC		
NET1227250UC	MK-912, [³ H(G)]-	Second Monday in October
NET127001MC	Tyrosine, L-[RING-3,5- ³ H]-	Second Monday in May
NET127250UC		
NET131001MC	Dihydroxyphenylethylamine, 3,4-[7- ³ H]- (Dopamine)	Second Monday in April and October
NET131250UC		

Release Schedule for Select ³H and ¹⁴C Products

Part Number	Description	Availability
NET309001MC NET309250UC	Cytidine 5'-Triphosphate, [5- ³ H]-, Tetrasodium Salt	Second Monday in November
NET368250UC	Uracil, [5, 6- ³ H]-	Second Monday in February and August
NET376001MC NET376250UC	Lysine, L-[4,5- ³ H(N)]-	Second Monday in May
NET377001MC NET377005MC NET377250UC	Norepinephrine,Levo-[7- ³ H(N)]-	Third Monday in July
NET419250UC	Aldosterone, D-[1,2,6,7- ³ H(N)]-	First Monday in June
NET433025UC NET433050UC NET433250UC	Prostaglandin F _{2α} , [5,6,8,9,11,12,14,15- ³ H(N)]-	Third Monday in April and October
NET443H250UC	Nicotinamide-Adenine Dinucleotide (NAD;DPN), [Adenine-2,8- ³ H]-	First Monday in November
NET517250UC	Estradiol, [2,4,6,7,16,17- ³ H(N)]-	Second Monday in June
NET574250UC	Muscimol, 3-Hydroxy-5-Aminomethylisoxazole, [Methylene- ³ H(N)]-	Second Monday in February and August
NET626005UC	Dihydroxyvitamin D ₃ , 1Alpha, 25-[26,27- ³]	Second Monday in April and October
NET638250UC	Lysergic Acid Diethylamide, [N-Methyl- ³ H]-	Second Monday in April, August and December
NET674250UC	Hydroxycholesterol, 25-[26,27- ³ H]-	First Monday in June
NET869025UC NET869250UC	Paroxetine, [Phenyl-6'- ³ H]-	First Monday in May and November
NET884250UC	Hemicholinium-3 Diacetate Salt, [Methyl- ³ H]-	Second Monday in February and August
NET920100UC NET920250UC	Retinoic Acid, [11, 12- ³ H(N)]-	Third Monday in January and July
NET922050UC NET922250UC	Enkephalin, [Tyrosyl-2,6- ³ H(N)]- (2-D-Penicillamine, 5-D-Penicillamine)- (DPDPE)	Second Monday in February and August
NET927001MC NET927250UC	Retinol, [11, 12- ³ H (N)]-	Third Monday in January
NET966250UC	Guanosine 5'-Diphosphate, Trisodium Salt, [8,5'- ³ H]-	First Monday in April and October
NET975001MC NET975025UC NET975250UC	Raclopride, [Methoxy- ³ H]-	Second Monday in March and September

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2018-2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

014128A_01

PKI