

Release Schedule for Select ³H Products

Tritium Labeled Compounds		
Part Number	Description	Availability
NET048250UC	Norepinephrine Hydrochloride, DL-[7- ³ H(N)]-	Third Monday in January and July
NET063001MC	Adenine, [2, 8- ³ H]-	Second Monday in March and September
NET063005MC		
NET1100050UC	Lysophosphatidic Acid, 1-Oleoyl-[Oleoyl-9,10- ³ H]-	Second Monday in May and November
NET1165050UC	Quisqualic Acid, [³ H]-	Third Monday in March and September
NET1169250UC	R(+)-7-HYDROXY DPAT, [³ H]-	Second Monday in April and October
NET1172U100UC	GR125743, [N-Methyl- ³ H]-	Third Monday in February and August
NET1177001MC	Myo-Inositol, [³ H]-	Second Monday in March and September
NET1182001MC	Gabapentin, [³ H]-	Third Monday in February and September
NET1198U250UC	SB-269970, [³ H]-	First Monday in June
NET1227250UC	MK-912, [3H(G)]-	Second Monday in April and October
NET131001MC	Dihydroxyphenylethylamine, 3,4-[7- ³ H]- (Dopamine)	Second Monday in April and October
NET131250UC		
NET377001MC	Norepinephrine, Levo-[7- ³ H(N)]-	Third Monday in January and July
NET377005MC		
NET377250UC		
NET517250UC	Estradiol, [2,4,6,7,16,17- ³ H(N)]-	Second Monday in June and December
NET574250UC	Muscimol, 3-Hydroxy-5-Aminomethylisoxazole, [Methylene- ³ H(N)]-	Second Monday in February and August
NET638250UC	Lysergic Acid Diethylamide, [N-Methyl- ³ H]-	Second Monday in April, August and December

www.perkinelmer.com

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 © 2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.