

## Release Schedule for Select <sup>3</sup>H Products

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

[www.perkinelmer.com](http://www.perkinelmer.com)

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



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://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.

161377 (94132) PKI