

HTRF® Research Reagents

HTRF Histamine Dyn kit - 96 pts

Product Number : 62HTMDPET Lot Number : 01M

Expiration date : 2024-09-30 Document version : 1

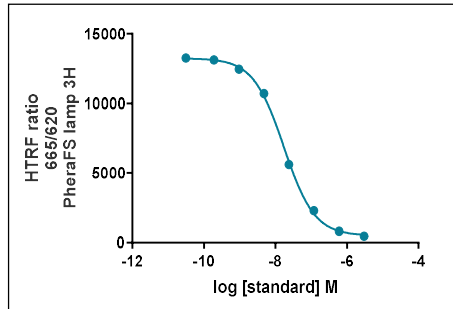
Reagents :	Lot n°
HTRF Diluent 2 - 20 mL	10RA
HTRF Detect. Buf. 2 - 1 mL	13RA
HTRF Histamine Dyn. kit - Acyl. reag.	02A
HTRF Histamine Dyn. kit - Std	02A
HTRF Histamine Dyn. kit - Acyl reagent Dil.	02A
HTRF Histamine Dyn. kit - Acylation Buf.	02A
HTRF Histamine Dyn. kit - DMSO	02A
HTRF Histamine Dyn. kit - 96 pts d2 reagent	01A
HTRF Histamine Dyn. kit - 96 pts Eu Cryptate antibody	01A
HTRF Histamine Dyn. kit - modop	3
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model :

Assay characteristics :

	Conc (nM)	ratio	ratio Accepted Range
Std 0	0	13251	12835-24637
Std 1	0.192	13115	12846-22986
Std 2	0.96	12461	12046-22986
Std 3	4.8	10712	9321 - 17750
Std 4	24	5604	4797 - 8793
Std 5	120	2297	1738 - 3190
Std 6	600	822	740 - 1107
Std 7	3000	470	429 - 644


Assay characteristics :

	Found value	Specifications Rev 01
ratio neg K	351	349-475
DF% max	3672%	3108 - 6530%
Ic 50	18.4 nM	5-28 nM
S/B	28	17-32

HTRF Histamine Dyn kit - 96 pts - Lot n° : 01M

 Accepted

 Rejected

Quality control date : 2023-09-29 Supervisor's initials : FMV

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.