

HTRF® Research Reagents

HTRF Histamine Dyn. kit - 96 pts

Product Number : 62HTMDPET Lot Number : 04B

Expiration date : 2025-12-18 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2303
HTRF Diluent 2 - 20 mL	10RA
HTRF Detect. Buf. 2 - 1 mL	13RA
HTRF Histamine Dyn. kit - Acyl. reag.	03A
HTRF Histamine Dyn. kit - Std	04B
HTRF Histamine Dyn. kit - Acyl reag Dil.	03A
HTRF Histamine Dyn. kit - Acylation Buf.	03A
HTRF Histamine Dyn. kit - DMSO	02RA
HTRF Histamine Dyn. kit - 96 pts d2 reagent	04A
HTRF Histamine Dyn. kit - 96 pts Eu Cryptate antibody	04A

Results :

Incubation :

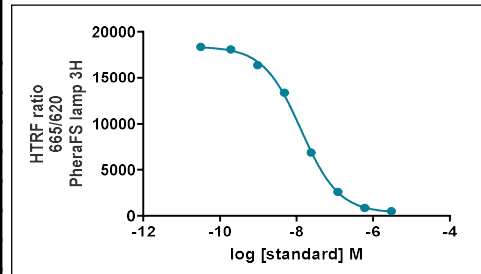
Plate :

Reader : Pherastar Plus

Fitting model :

Assay characteristics :

Histamine	Conc (nM)	ratio	ratio Accepted Range
Std 0	0	18183	12835-24637
Std 1	0.192	17949	12846-22986
Std 2	0.96	17312	12046-22986
Std 3	4.8	13999	9321 - 17750
Std 4	24	7403	4797 - 8793
Std 5	120	2588	1738 - 3190
Std 6	600	977	740 - 1107
Std 7	3000	565	429 - 644



	Found value	Specifications Rev 01
IC50	13.5	5-28 nM
S/B	32	17-32

HTRF Histamine Dyn. kit - 96 pts - Lot n° : 04B

Accepted

Rejected

Quality control date : 2024-07-15	Supervisor's initials : AC
-----------------------------------	----------------------------

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.