

HTRF® Research Reagents

HTRF Histamine Dyn. kit - 96 pts

Product Number : 62HTMDPET Lot Number : 04C

Expiration date : 2026-09-30 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2413
HTRF Diluent 2 - 20 mL	12A
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 reagent Dil.	04A
HTRF Histamine Dyn. kit - Acylation Buf.	04A
HTRF Histamine Dyn. kit - DMSO	03A
HTRF Histamine Dyn. kit - 96 pts d2 reagent	04A
HTRF Histamine Dyn. kit - 96 pts Eu Cryptate antibody	04A

Results :

Incubation : 3h at RT

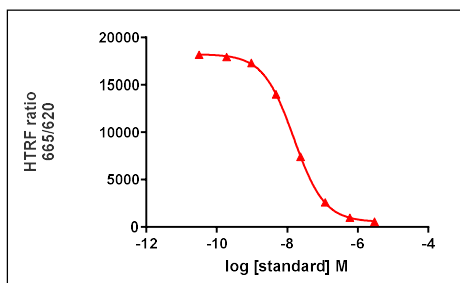
Plate : 384 wells small volume white plate 20 µL

Reader : Envision Lamp

Fitting model : log(inhibitor) vs. response -- Variable slope (four parameters)

Assay characteristics :

Histamine	Conc (nM)	ratio
Std 0	0	18183
Std 1	0,192	17949
Std 2	0,96	17312
Std 3	4,8	13999
Std 4	24	7403
Std 5	120	2588
Std 6	600	977
Std 7	3000	565



	Found value	Specifications Rev 03
Blanc K (ratio)	445	353 - 530
DF% max	3985%	3055% - 4695 %
IC50	13,5 nM	5 - 28 nM
S/B	32,2	26-39

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

☒ Accepted☐ Rejected

Quality control date : 2024-10-02 Supervisor's initials : MP

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.