

HTRF® Research Reagents

HTRF Histamine Dyn. kit - 96 pts

Product Number : 62HTMDPET Lot Number : 05B

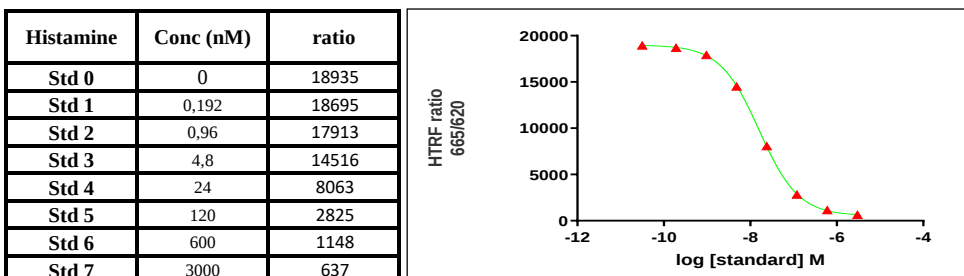
Expiration date : 2026-09-30 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	23
HTRF Diluent 2 - 20 mL	12A
HTRF Detect. Buf. 2 - 1 mL	13RA
HTRF Histamine Dyn. kit - Acyl. reag.	04A
HTRF Histamine Dyn. kit - Std	04B
HTRF Histamine Dyn. kit - Acyl reag Dil.	04A
HTRF Histamine Dyn. kit - Acylation Buf.	03A
HTRF Histamine Dyn. kit - DMSO	03A
HTRF Histamine Dyn. kit - 96 pts d2 reagent	05B
HTRF Histamine Dyn. kit - 96 pts Eu Cryptate antibody	05B

Results :

Incubation : 3h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : EnVision
 Fitting model : log(inhibitor) vs. response -- Variable slope (four parameters)

Assay characteristics :



	Found value	Specifications Rev 03
Blanc K (ratio)	521	353 - 530
DF% max	3534%	3055% - 4695 %
IC50	16nM	5 - 28 nM
S/B	29,7	26-39

HTRF Histamine Dyn. kit - 96 pts - Lot n° : 05A

☒ Accepted☐ Rejected

Quality control date : 2025-04-11 Supervisor's initials : PL

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.