

HTRF® Research Reagents

### HTRF Histamine Dyn. kit - 96 pts

Product Number : 62HTMDPET Lot Number : 03B

Expiration date : 2025-09-11 Document version : 1

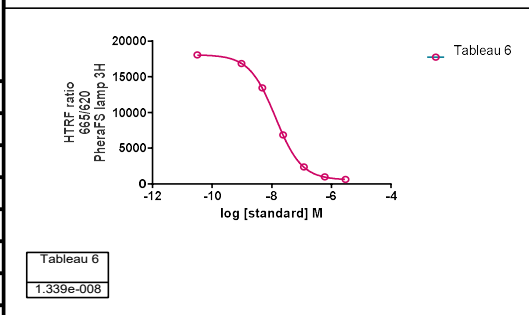
Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF Diluent 2 - 20 mL	10RA
HTRF Detect. Buf. 2 - 1 mL	13RA
HTRF Histamine Dyn. kit - Acyl. reag.	03A
HTRF Histamine Dyn. kit - Std	03A
HTRF Histamine Dyn. kit - Acyl reagent Dil.	02RA
HTRF Histamine Dyn. kit - Acylation Buf.	03A
HTRF Histamine Dyn. kit - DMSO	02RA
HTRF Histamine Dyn. kit - 96 pts d2 reagent	01RA
HTRF Histamine Dyn. kit - 96 pts Eu Cryptate antibody	03A

**Results :**

Incubation : 3h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : /  
 Fitting model : /

**Assay characteristics :**

Histamine	Conc (nM)	ratio	ratio Accepted Range
Std 0	0	18050	12835-24637
Std 1	0.192	17902	12846-22986
Std 2	0.96	16878	12046-22986
Std 3	4.8	13439	9321 - 17750
Std 4	24	6851	4797 - 8793
Std 5	120	2367	1738 - 3190
Std 6	600	976	740 - 1107
Std 7	3000	623	429 - 644



	Found value	Specifications Rev 01
IC50	13.39 nM	5-28 nM
S/B	29	17-32

HTRF Histamine Dyn. kit - 96 pts - Lot n° : 03B

Accepted

Rejected

Quality control date : 2024-05-23 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.