

HTRF® Research Reagents

# HTRF Histamine Dyn kit - 500 pts

Product Number : 62HTMDPEG Lot Number : 02M

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

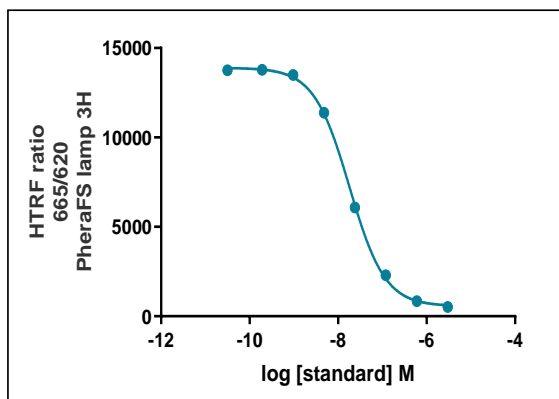
Reagents :		Lot n°
HTRF Diluent 2 - 20 mL		10RA
HTRF Detect. Buf. 2 - 1 mL		16RA
HTRF Histamine Dyn. kit - Acyl. reag.		02A
HTRF Histamine Dyn. kit - Std		02A
HTRF Histamine Dyn. kit - Acyl reag Dil.		02A
HTRF Histamine Dyn. kit - Acylation Buf.		02A
HTRF Histamine Dyn. kit - DMSO		02A
HTRF Histamine Dyn. kit - 500 pts d2 reagent		02A
HTRF Histamine Dyn. kit - 500 pts Eu Cryptate antibody		02A
HTRF Histamine Dyn. kit - modop		3
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND		<b>2203</b>

**Results :**

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

**Assay characteristics :**

Histamine	Conc (nM)	ratio	ratio Accepted Range
Std 0	0	13758	12835-24637
Std 1	0,192	13777	12846 -22986
Std 2	0,96	13488	12046-22986
Std 3	4,8	11381	9321 - 17750
Std 4	24	6081	4797 - 8793
Std 5	120	2291	1738 - 3190
Std 6	600	847	740 - 1107
Std 7	3000	517	429 - 644



Assay characteristics :

	Found value	Specifications Rev 01
ratio neg K	357	349-475
Ic 50	18,4 nM	5-28 nM
S/B	26,7	17-32