

HTRF® Research Reagents

HTRF Human Fc kit - 500 pts

Product Number : 62HFCPEG Lot Number : 13S

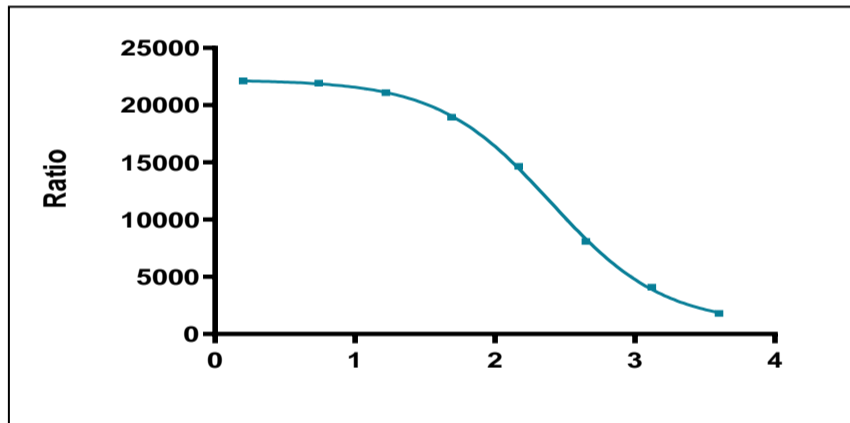
Expiration date : 2025-06-16 Document version : 1

Reagents :		Lot n°
HTRF Diluent 1 - 20 mL		20RA
HTRF Detect. Buf. 2 - 7 mL		16RA
HTRF Human IgG kit - Std		19AR
HTRF Human Fc kit - 500 pts d2 antibody		13RA
HTRF Human Fc kit - 500 pts Eu Cryptate antibody		12RA

Results :

Incubation : **2h30 at RT**
 Plate : **384 wells small volume white plate 20 µL**
 Reader : **Pherastar Plus**
 Fitting model : **nonlinear regression, sigmoïdal dose response (variable slope)**

IgG	Conc. (ng/ml)	Moyenne Ratio
Std 0	0	22105
Std 1	5,5	21928
Std 2	16,5	21095
Std 3	49	18940
Std 4	148	14655
Std 5	444	8085
Std 6	1333	4089
Std 7	4000	1801



Assay characteristics :

	Found value	Specifications Rev 09
Detection limite	3,9	<=9 ng/ml
Df Max	6480	3956 - 6625 %
EC50	249	240-280 ng/ml
S/B	12,3	11-16

HTRF Human Fc kit - 500 pts - Lot n° : 13S

Accepted

Rejected

Quality control date : 2024-02-02	Supervisor's initials : FMV
-----------------------------------	-----------------------------

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.