

HTRF® Research Reagents

HTRF Human Fc kit - 500 pts

Product Number : 62HFCPEG Lot Number : 16A

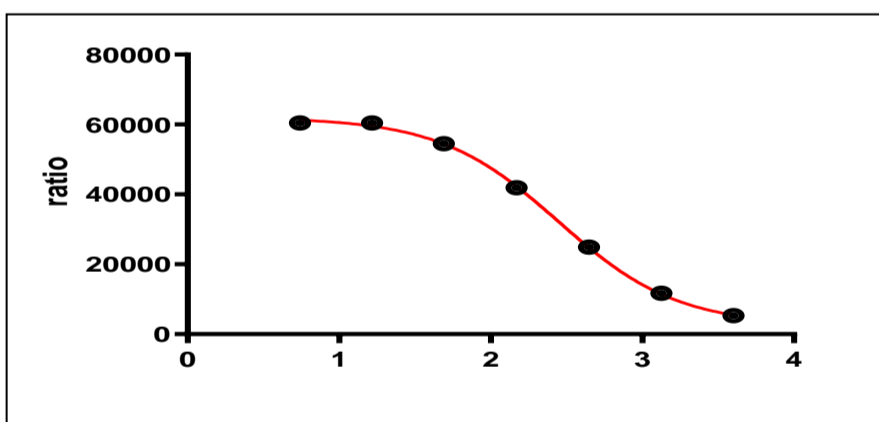
Expiration date : 2026-09-04 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Detect. Buf. 2 - 7 mL	17AR
HTRF Human IgG kit - Std	21A
HTRF Human Fc kit - 500 pts d2 antibody	13RA
HTRF Human Fc kit - 500 pts Eu Cryptate antibody	16A

Results :

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

IgG	Conc. (ng/ml)	Moyenne Ratio
Std 0	0	63444
Std 1	5,5	60401
Std 2	16,5	60462
Std 3	49	54510
Std 4	148	41877
Std 5	444	24880
Std 6	1333	11679
Std 7	4000	5271



	Found value	Specifications Rev 09
Detection limite	0,6	<=9 ng/ml
Df Max	6397	3956 - 6625 %
EC50	279	256-355 ng/ml
S/B	12	11-16

HTRF Human Fc kit - 500 pts - Lot n° : 16A

Accepted

Rejected

Quality control date : 2024-12-06 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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.