

HTRF® Research Reagents

HTRF Human Fc kit - 500 pts

Product Number : 62HFCPEG Lot Number : 13X

Expiration date : 2025-06-16 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	12RA

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	17991
Std 1	5.5	17560
Std 2	16.5	17038
Std 3	49	15014
Std 4	148	11619
Std 5	444	6587
Std 6	1333	3462
Std 7	4000	1508

	Found value	Specifications Rev 09
Detection limite	7	<=9 ng/ml
Df Max	4815	3956 - 6625 %
EC50	246	256-355 ng/ml
S/B	11.9	11-16
	Found value	Specifications Rev 01

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

 Accepted

 Rejected

Quality control date : 2024-04-02 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.