

HTRF® Research Reagents

HTRF Human Fc kit - 10K pts

Product Number : 62HFCPEH Lot Number : 14M

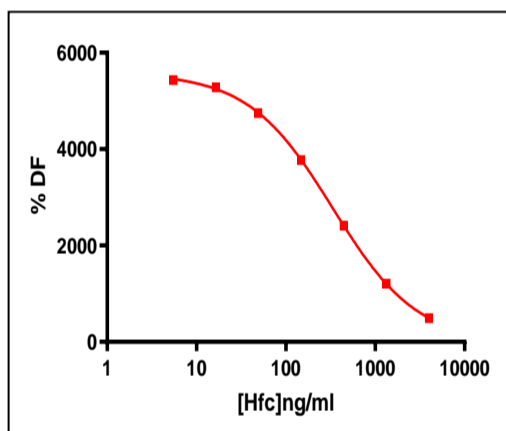
Expiration date : 2025-03-26 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20 A
HTRF Detect. Buf. 3 - 105 mL	29A
HTRF Human IgG kit - Std	17A
HTRF Human Fc kit - 10K pts modop	5
HTRF Human Fc kit - 10K pts d2 antibody	14 A
HTRF Human Fc kit - 10K pts Eu Cryptate antibody	14 A

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)	DF %
Std 0	0	5659%
Std 1	5,5	5430%
Std 2	16,5	5280%
Std 3	49	4749%
Std 4	148	3770%
Std 5	444	2414%
Std 6	1333	1208%
Std 7	4000	491%



Assay characteristics :

	Found value	Specifications Rev 06
Detection limit	7,2	<=9 ng/ml
DF max %	5659	3956-6625 %
fenêtre d'essai sur gamme	11,5	9-16
EC 50	334	246-355 ng/ml

HTRF Human Fc kit - 10K pts - Lot n° : 14M

Accepted

Rejected

Quality control date : 2022-10-18	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.