

HTRF® Research Reagents

HTRF Human IgG kit - 10K pts

Product Number : 6FHIGPEH Lot Number : 17A

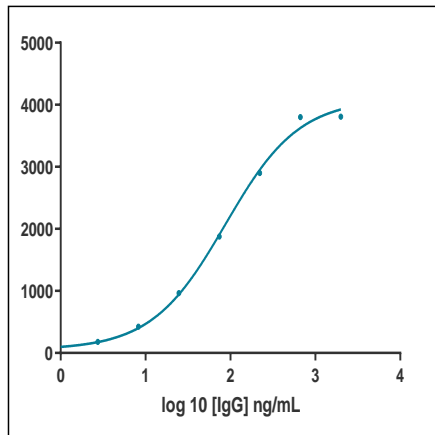
Expiration date : 2025-04-27 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20 A
HTRF Detect. Buf. 3 - 105 mL	28 A
HTRF Human IgG kit - 10K pts modop	3
HTRF Human IgG kit - 10K pts d2 antibody	17A
HTRF Human IgG kit - 10K pts Eu Cryptate antibody	17A
HTRF Human IgG kit - Std	17RA

Results :

Incubation : 2h00 at TA
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL)

human IgG	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	0,91	82%	27-90
Std 2	2,74	148%	81-197
Std 3	8,23	408%	222-429
Std 4	24,7	848%	555-1014
Std 5	74	1780%	1268-2060
Std 6	222	3093%	2314-3355
Std 7	667	4179%	3213-4390
Std 8	2000	4256%	3578-5257



Assay characteristics :

	Found value	Specifications Rev 01
EC 50	128	111-178

Rejected

**Batch validated with EC50 a little bit lower than spécification

HTRF Human IgG kit - 10K pts - Lot n° : 17A

Accepted

Rejected

Quality control date : 2023-10-12	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