

HTRF® Research Reagents

HTRF Human IgG kit - 500 pts

Product Number : 6FHIGPEG Lot Number : 16E

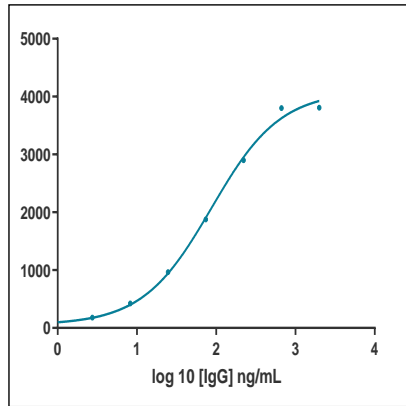
Expiration date : 2025-06-06 Document version : 1

Reagents :		Lot n°
HTRF Diluent 1 - 20 mL		20RA
HTRF Detect. Buf. 3 - 7 mL		30AR
HTRF Human IgG kit - 500 pts d2 antibody		16RA
HTRF Human IgG kit - 500 pts Eu Cryptate antibody		16RA
HTRF Human IgG kit - Std		19AR

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	71%	27-90
Std 2	2,74	179%	81-197
Std 3	8,23	424%	222-429
Std 4	24,7	963%	555-1014
Std 5	74	1875%	1268-2060
Std 6	222	2901%	2314-3355
Std 7	667	3802%	3213-4390
Std 8	2000	3809%	3578-5257



Assay characteristics :

	Found value	Specifications Rev 01
EC 50	87	111-178

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

Rejected

HTRF Human IgG kit - 500 pts - Lot n° : 16E

Accepted

Rejected

Quality control date :	2023-11-17	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.