

HTRF® Research Reagents

HTRF Human Fc kit - 10K pts

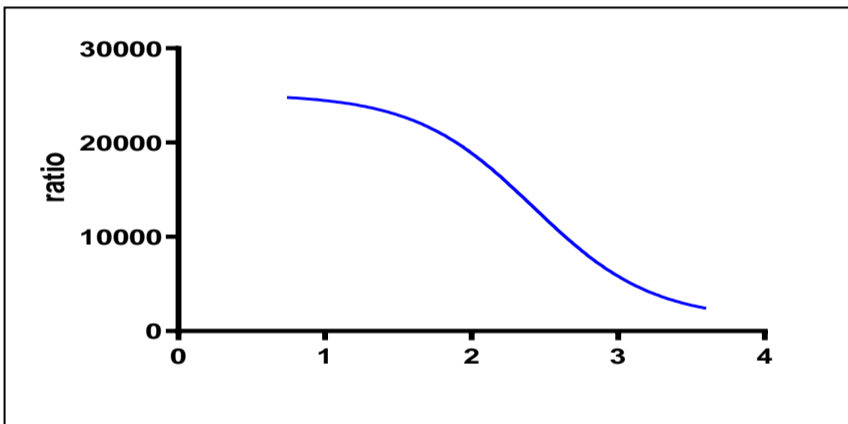
Product Number : 62HFCPEH Lot Number : 16A

Expiration date : 2025-06-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Detect. Buf. 3 - 105 mL	29RA
HTRF Human IgG kit - Std	21A
HTRF Human Fc kit - 10K pts d2 antibody	16AR
HTRF Human Fc kit - 10K pts Eu Cryptate antibody	16AR

Results :
 Incubation :
 Plate :
 Reader : Pherastar Plus
 Fitting model :

hFc	Conc. (ng/ml)	DF %
Std 0	0	0%
Std 1	5,5	24600%
Std 2	16,5	24193%
Std 3	49	21752%
Std 4	148	16795%
Std 5	444	9696%
Std 6	1333	5104%
Std 7	4000	905%



Assay characteristics :

	Found value	Specifications Rev 09
Detection limit	7,2	<=9 ng/ml
DF max %	6476	3958-6625 %
EC 50	263,6	275-355 ng/ml

HTRF Human Fc kit - 10K pts - Lot n° : 16A

Accepted

Rejected

Quality control date : 2024-04-24	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 USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.