

HTRF® Research Reagents

HTRF Human C1q Binding Kit - 100 pts

Product Number : 64C1QPET Lot Number : 05D

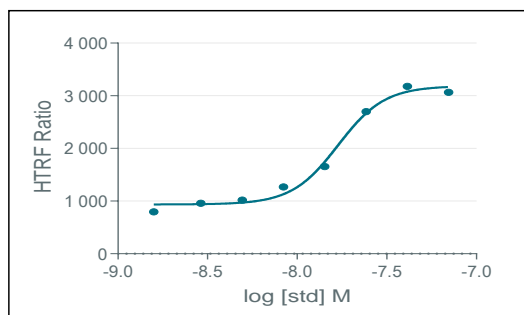
Expiration date : 2027-10-04 Document version : 1

Reagents :	Lot n°
HTRF Diluent 12 - 20 mL	02C
HTRF Human C1q Binding Kit - Std	05A
HTRF Human C1q Binding Kit - 100 pts Streptavidin d2	05B
HTRF Human C1q Binding Kit - 100 pts Anti-Human IgG Fab-biotin	10A
HTRF Human C1q Binding Kit - 100 pts Eu Cryptate antibody	05B
HTRF Human C1q Binding Kit - 100 pts Human C1q	04A
/	/

Results :

Incubation : 3h at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Pherastar Plus
Fitting model : log(inhibitor)vs.response-variable slope (four parameters)

	Conc. (nM)	RATIO
Std 0	0	795
Std 1	2,9	957
Std 2	4,9	1019
Std 3	8,4	1269
Std 4	14,3	1655
Std 5	24	2699
Std 6	41	3178
Std 7	70	3065



Assay characteristics :

	Found value	Specifications Rev 01
S/B	4	4-7
IC50	17	15-23

HTRF Human C1q Binding Kit - 100 pts - Lot n° : 05D

☒ Accepted☐ Rejected

Quality control date : 2025-04-15 Supervisor's initials : MP

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.