

## HTRF® Research Reagents

## HTRF Human C1q Binding Kit - 100 pts

Product Number : 64C1QPET Lot Number : 05B

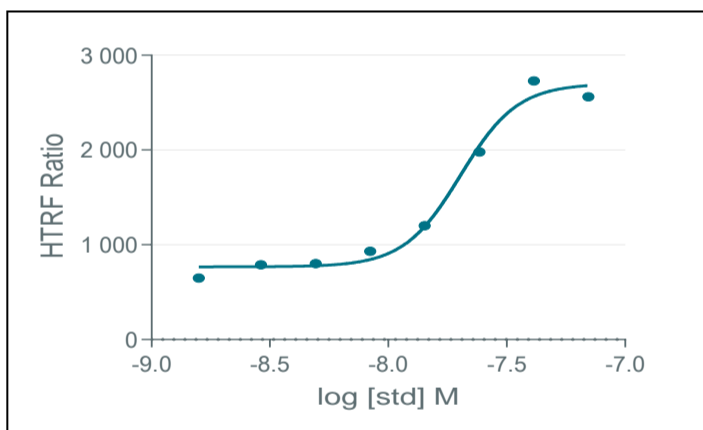
Expiration date : 2027-10-03 Document version : 1

Reagents :		Lot n°
HTRF Diluent 12 - 20 mL		02B
HTRF Human C1q Binding Kit - Std		04A
HTRF Human C1q Binding Kit - 100 pts Streptavidin d2		05A
HTRF Human C1q Binding Kit - 100 pts Anti-Human IgG Fab-biotin		09A
HTRF Human C1q Binding Kit - 100 pts Eu Cryptate antibody		05A
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	648
Std 1	2,9	789
Std 2	4,9	801
Std 3	8,4	931
Std 4	14,3	1201
Std 5	24	1979
Std 6	41	2729
Std 7	70	2562



## Assay characteristics :

	Found value	Specifications Rev 01
S/B	4	4-7
IC50	20	15-23

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

☒ Accepted☐ Rejected

Quality control date : 2024-12-05 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.