

HTRF® Research Reagents

HTRF CD47/SIRP a Bind. kit - 500 pts

Product Number : 64SIRPEG Lot Number : 04M

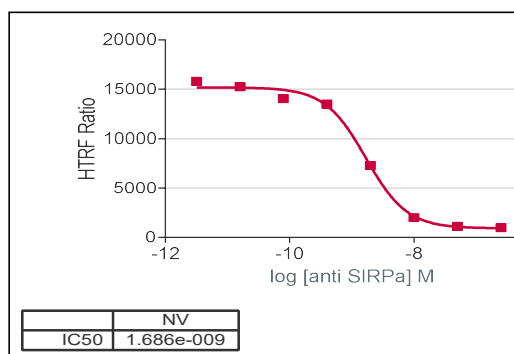
Expiration date : 2026-06-27 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	07RA
HTRF CD47/SIRP a Bind. kit - 500 pts Tag2-SIRPa	04RA
HTRF CD47/SIRP a Bind. kit - Std	04A
HTRF CD47/SIRP a Bind. kit - 500 pts Eu Cryptate reagent	04RA
HTRF CD47/SIRP a Bind. kit - 500 pts Tag1-CD47	04RA
HTRF CD47/SIRP a Bind. kit - 500 pts XL665 antibody	05A

Results :

Incubation : 1h at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Pherastar Plus
Fitting model : log(inhibitor)vs.response-variable slope (four parameters)

anti hSIRPa	Conc. (nM)	RATIO
Std 0	0	15816
Std 1	0,16	15268
Std 2	0,8	14619
Std 3	4	13506
Std 4	20	7288
Std 5	100	2031
Std 6	500	1134
Std 7	2500	1007



Assay characteristics :

	Found value	Specifications Rev 01
S/B	16	15 - 25
IC50	1,7	1.5 - 3.5 nM

HTRF CD47/SIRP a Bind. kit - 500 pts - Lot n° : 04M

☒ Accepted☐ Rejected

Quality control date :	2024-04-24	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.