

HTRF® Research Reagents

HTRF CTLA4/CD80 Bind. kit - 500 pts

Product Number : 64CTLA80PEG Lot Number : 02D

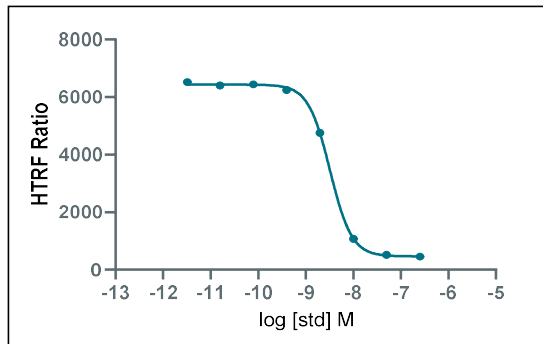
Expiration date : 2026-11-09 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	08A
HTRF CTLA4/CD80 Bind. kit - 500 pts Tag2-CTLA4	02RA
HTRF CTLA4/CD80 Bind. kit - Std	02RA
HTRF CTLA4/CD80 Bind. kit - 500 pts d2 reagent	02RA
HTRF CTLA4/CD80 Bind. kit - 500 pts Eu Cryptate antibody	02RA
HTRF CTLA4/CD80 Bind. kit - 500 pts Tag1-CD80	02RA

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	6524
Std 1	0,016	6401
Std 2	0,08	6445
Std 3	0,4	6251
Std 4	5	4754
Std 5	10	1076
Std 6	50	516
Std 7	250	457



Assay characteristics :

	Found value	Specifications Rev 01
S/B	14	11 17
IC50	3,2	1.5 - 3.5 nM

HTRF CTLA4/CD80 Bind. kit - 500 pts - Lot n° : 02D

Accepted

Rejected

Quality control date : 2024-10-14	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.