

HTRF® Research Reagents

HTRF LAG3/MHCII Bind. kit - 500 pts

Product Number : 64LAG3PEG Lot Number : 02A

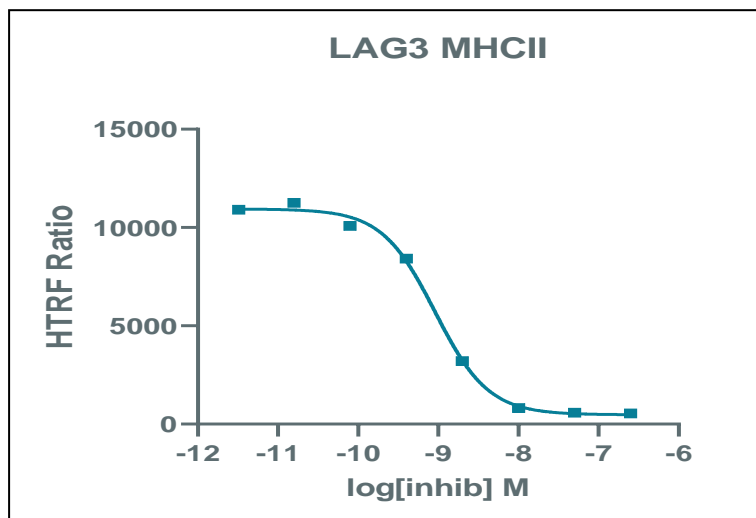
Expiration date : 2025-06-07 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	06A
HTRF LAG3/MHCII Bind. kit - 500 pts Tag2-MHCII	02A
HTRF LAG3/MHCII Bind. kit - Std	02A
HTRF LAG3/MHCII Bind. kit - 500 pts d2 antibody	02A
HTRF LAG3/MHCII Bind. kit - 500 pts Eu Cryptate reagent	02A
HTRF LAG3/MHCII Bind. kit - 500 pts Tag1-LAG3	02A
HTRF LAG3/MHCII Bind. kit - modop	2

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)

std	Conc. (nM)	RATIO
Std 0	0	10911
Std 1	0,16	11258
Std 2	0,8	10086
Std 3	4	8423
Std 4	20	3210
Std 5	100	817
Std 6	500	576
Std 7	2500	538



Assay characteristics :

	Found value	Specifications Rev 01
negatif	514	
S/B	21	15-30
IC50	0,91	0.5 - 2.7 nM

HTRF LAG3/MHCII Bind. kit - 500 pts - Lot n° : 02A

Accepted

Rejected

Quality control date : 2022-08-29	Supervisor's initials : FL
-----------------------------------	----------------------------

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.