

HTRF® Research Reagents

HTRF LAG3/MHCII Bind. kit - 500 pts

Product Number : 64LAG3PEG Lot Number : 03A

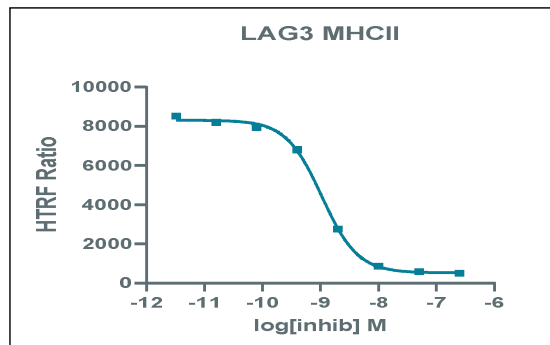
Expiration date : 2025-05-06 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	05RA
HTRF LAG3/MHCII Bind. kit - 500 pts Tag2-MHCII	03BR
HTRF LAG3/MHCII Bind. kit - Std	02RA
HTRF LAG3/MHCII Bind. kit - 500 pts d2 antibody	03A
HTRF LAG3/MHCII Bind. kit - 500 pts Eu Cryptate reagent	03A
HTRF LAG3/MHCII Bind. kit - 500 pts Tag1-LAG3	03BR
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	8528
Std 1	0.16	8202
Std 2	0.8	7965
Std 3	4	6802
Std 4	20	2762
Std 5	100	860
Std 6	500	590
Std 7	2500	504



Assay characteristics :

	Found value	Specifications Rev 01
negatif	482	
S/B	18	15-30
IC50	1.06	0.5 - 2.7 nM

HTRF LAG3/MHCII Bind. kit - 500 pts - Lot n° : 03A

Accepted

Rejected

Quality control date : 2023-11-10	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.