

HTRF® Research Reagents

HTRF CD32a (FcγRIIa,H131) kit - 100 pts

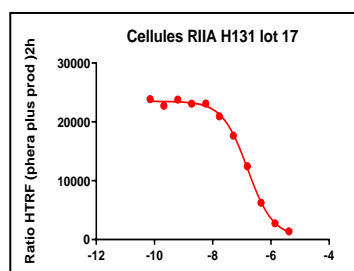
Product Number : 6FRAHPAG Lot Number : 17C

Expiration date : 2024-12-08 Document version : 1

Reagents :	Lot n°
CD32A (H131 prelabelled cells)	17
HTRF Tag-lite Buffer (5X) - 1.5 mL	15RA
HTRF IgG d2 conjugate CD32A H131 - 100 pts	18RB

Results :

Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model :



Assay characteristics :

	Found value
IC50 (nM)	158.8
S/B	17

HTRF CD32a (FcγRIIa,H131) kit - 100 pts - Lot n° : 17C

Accepted

Rejected

Quality control date : 2021-09-14	Supervisor's initials : JB
-----------------------------------	----------------------------

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.