

HTRF® Research Reagents

HTRF CD32b (FcγRIIb) kit - 100 pts

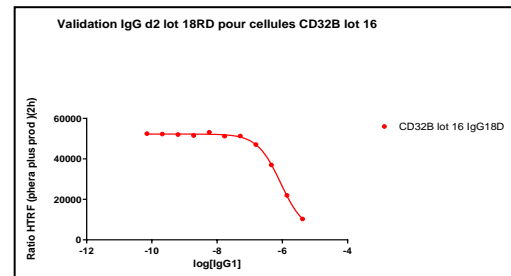
Product Number : 6FR2BPAG Lot Number : 16C

Expiration date : 2026-02-07 Document version : 1

Reagents :	Lot n°
CD32B (prelabelled cells)	16
HTRF Tag-lite Buffer (5X) - 1.5 mL	15RA
HTRF IgG d2 conjugate CD32B - 100 pts	18RD

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)	886
HTRF CD32b (FcγRIIb) kit - 100 pts - Lot n° : 16C	13

HTRF CD32b (FcγRIIb) kit - 100 pts - Lot n° : 16C

Accepted

Rejected

Quality control date :	2024-02-07	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.