

HTRF® Research Reagents

HTRF CD64 (FcγRI) - 5 x 6FC64PAG

Product Number : 6FC64PAE Lot Number : 21A

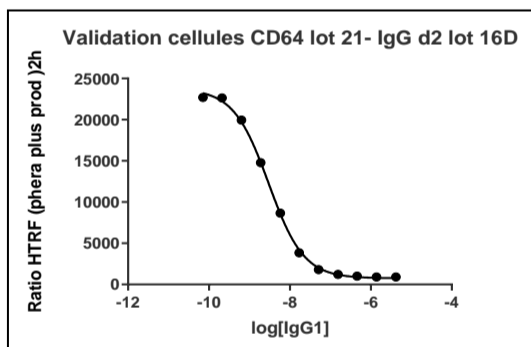
Expiration date : 2026-03-19 Document version : 1

Reagents :	Lot n°
HTRF IgG d2 conjugate CD64 - 100 pts	16D
CD64 (prelabelled cells)	21
HTRF CD64 cellular kit 100 pts - modop	2
HTRF Tag-lite Buffer (5X) - 1.5 mL	15RA

Results :

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

Assay characteristics :



	Found value
Window	31
IC50 (nM)	2.988

HTRF CD64 (FcγRI) - 5 x 6FC64PAG - Lot n° : 21A

Accepted

Rejected

Quality control date : 2023-09-29	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.