

HTRF® Research Reagents

**HTRF CD64 (FcγRI) - 5 x 6FC64PAG**

Product Number : 6FC64PAE Lot Number : 21C

Expiration date : 2026-06-19 Document version : 1

Reagents :		Lot n°
HTRF Tag-lite Buffer (5X) - 1.5 mL		20AR
CD64 (prelabelled cells)		21
HTRF IgG d2 conjugate CD64 - 100 pts		16RD

Results :

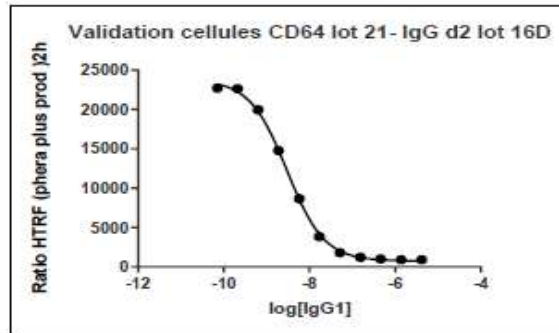
Incubation :

Plate :

Reader : Pherastar Plus

Fitting model :

Assay characteristics :



	Found value
Window	31
IC50 (nM)	2.988

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

Accepted

Rejected

Quality control date : 2023-09-29	Supervisor's initials : AC
-----------------------------------	----------------------------

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