

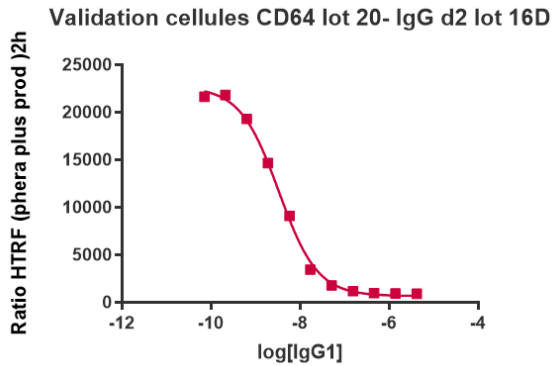
Quality Control report
 CD64 (FcγRI) - cellular kit - 100 tests
 (6FC64PAG)
 Lot n°20D
 Expiration date : 2026-03-19

Reagents :	Lot n°
Tag-lite buffer 1.5 ml (5X)	15RA
CD64 (prelabelled cells)	20
CD64 cellular kit 100 tests modop	2
IgG d2 conjugate CD64 - 100 tests	16D

Results :

Incubation : 2H
 Plate :
 Reader : Pherastar Plus
 Fitting model :

Assay characteristics :



	Value
Window	34
IC50 (nM)	3.322

CD64 (FcγRI) cellular kit - 100 tests Lot n°20D

Accepted

Rejected

Quality control date : 19/06/2023	Supervisor's initials : JB
-----------------------------------	----------------------------

