

HTRF® Research Reagents

Tag-lite Glucagon GIPR labeled cells -200 tests

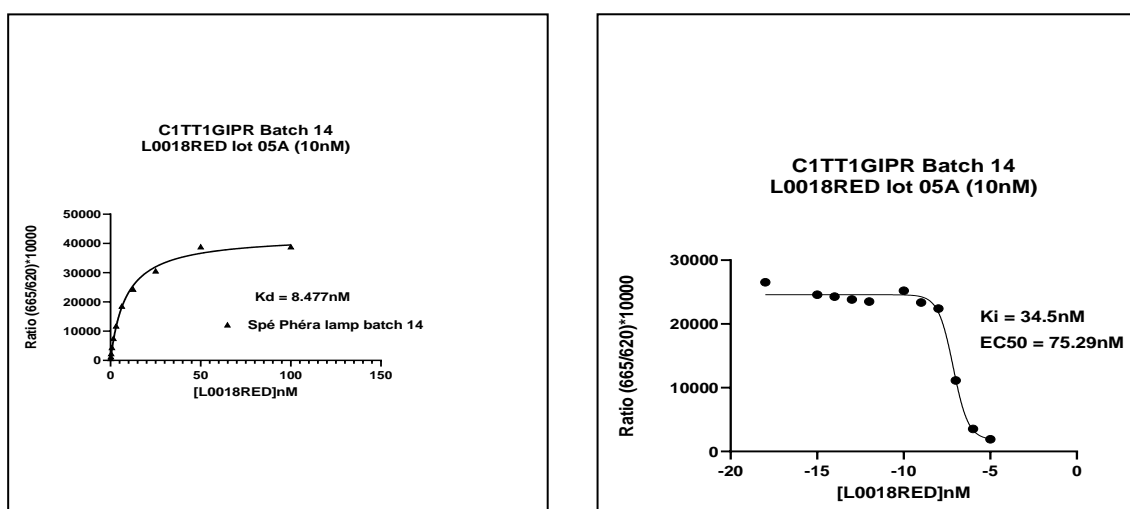
Product Number : C1TT1GIPR Lot Number : 14

Expiration date : 2025-12-31 Document version : 1

Reagents :		Lot n°
Tag-lite Glucagon GIPR labaled cells-200 tests		14
LABMED	SNAP/CLIP lab medium 5X - 100 ml	17A
L0018RED	GIPR receptor red agonist	05A

Results :

Incubation : 1H  
 Plate : 384 wells small volume white plate 20µL  
 Reader : Pherastar FS (flash lamp)  
 Fitting model :



Assay characteristics :

	Found value	Specifications Rev 01
Kd (nM)	8.477	5 - 22 nM
Ki (nM)	34.5	9 - 55 nM
Window (Max/Min)	13.5	6 -18
620 signal (cps)	26 692	> 10 000

Tag-lite Glucagon GIPR labeled cells - 200 tests - Lot n°14

☒ Accepted

☐ Rejected

Quality control date : 2023-12-07 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.