

Quality Control report

Tag-lite Dopamine D2 labeled cells-200 tests

(C1TT1D2)

Lot n°14

Expiration date : 2026-04-23

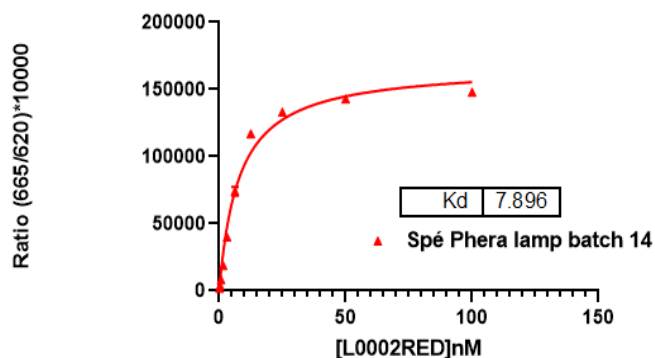
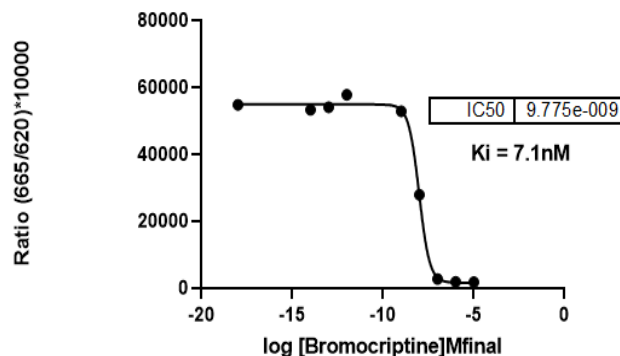
Reagents used for the Quality Control :	Lot n°
Tag-lite Dopamine D2 labeled cells-200 tests	14
LABMED SNAP/CLIP lab medium 5X - 100 mL	16A
L0002RED Dopamine D2 receptor red antagonist	07B

Results :

Incubation : 2h at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Pherastar FS (flash lamp)

C1TT1D2 Batch 14
L0002RED lot 07B (3nM)C1TT1D2 Batch 14
L0002RED lot 07B (3nM)

Assay characteristics :

	Found value	Specifications Rev 01
Kd (nM)	7.896	1.9 - 7.6 nM
Ki (nM)	7.1	3.4 - 21 nM
Window (Max/Min)	33	17 - 49
620 signal (cps)	22 728	> 10 000

Tag-lite Dopamine D2 labeled cells-200 tests Lot n°14

☒ Accepted

x

☐ Rejected

Quality control date :	2023-10-23	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.