

HTRF® Research Reagents

HTRF HIF-1 alpha kit - 500 pts

Product Number : 64HIFPEG Lot Number : 02Q

Expiration date : 2026-04-03 Document version : 1

Reagents :	Lot n°
HTRF HIF-1 alpha kit - 500 pts d2 antibody	02RA
HTRF HIF-1 alpha kit - 500 pts Eu Cryptate antibody	02RA
HTRF HIF-1 alpha kit - Ctrl lysate	02RA
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28B
HTRF P-T prot. - Detect. Buf. 2 mL	32B

Results :

Incubation :

Plate :

Reader : Pherastar Plus

Fitting model :

Assay characteristics :

	Found value	Specifications Rev 01
Negative	402	365-548
Control Lysate	16460	13415-20123

HTRF HIF-1 alpha kit - 500 pts - Lot n° : 02Q

Accepted

Rejected

Quality control date : 2023-12-21	Supervisor's initials : AC
-----------------------------------	----------------------------

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