

HTRF® Research Reagents

HTRF EIF2 alpha p-S52 kit - 10K pts

Product Number : 64EF2PEH Lot Number : 08D

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

Reagents :	Lot n°
HTRF EIF2 alpha p-S52 kit - 10K pts d2 antibody	08A
HTRF EIF2 alpha p-S52 kit - 10K pts Eu Cryptate antibody	08A
HTRF EIF2 alpha p-S52 kit - Ctrl lysate	08A
HTRF P-T prot. - Lysis Buf.4 (4X) 130 mL	14RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	29A
HTRF P-T prot. - Detect. Buf. 50 mL	33A

Results :

Incubation : Overnight at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Pherastar Plus
Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	395	/
Control Lysate	4249	/
S/B	10,7	>5

HTRF EIF2 alpha p-S52 kit - 10K pts - Lot n° : 08D

☒ Accepted☐ Rejected

Quality control date : 2025-03-07	Supervisor's initials : LM
-----------------------------------	----------------------------

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.