

## HTRF® Research Reagents

**HTRF Human c-Myc cell. kit - 10K pts**

Product Number : 63ADK053PEH Lot Number : 06E

Expiration date : 2026/01/15 Document version : 1

Reagents :	Lot n°
HTRF Lysis Buffer 4 - 130 mL	14AR
HTRF Detect. Buf. 3 - 105 mL	29RA
HTRF Human c-Myc cell. kit - Ctrl lysate	11AR
HTRF Human c-Myc cell. kit - 10K pts d2 antibody	06RA
HTRF Human c-Myc cell. kit - 10K pts Eu Cryptate antibody	06RA
HTRF Human c-Myc cell. kit - modop	4

**Results :**

 Incubation : **3h at RT**  
 Plate : **384 wells small volume white plate 20 µL**  
 Reader : **Pherastar Plus**  
 Fitting model : /

**Assay characteristics :**

	Found value	Specifications Rev 01
<b>Control Lysate</b>	3451	> 3000
<b>Ratio control lysate / negative</b>	5.9	≥ 5
	/	/
	/	/

HTRF Human c-Myc cell. kit - 10K pts - Lot n° : 06E

 Accepted

 Rejected

Quality control date : 2023-09-01	Supervisor's initials : PL
-----------------------------------	----------------------------

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.