

HTRF® Research Reagents

HTRF Human c-Myc cell. kit - 500 pts

Product Number : 63ADK053PEG Lot Number : 07G

Expiration date : 2026-06-15 Document version : 1

Reagents :	Lot n°
HTRF Lysis Buffer 4 - 16 mL	14RA
HTRF Detect. Buf. 3 - 7 mL	30B
HTRF Human c-Myc cell. kit - Ctrl lysate	11AR
HTRF Human c-Myc cell. kit - 500 pts d2 antibody	07A
HTRF Human c-Myc cell. kit - 500 pts Eu Cryptate antibody	07A

Results :

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

Assay characteristics :

	Found value	Specifications Rev 01
Negative	638	462-694
Control Lysate	5536	>3000
S/B	8,7	>7

HTRF Human c-Myc cell. kit - 500 pts - Lot n° : 07G

 Accepted

 Rejected

Quality control date : 2025-01-27	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.