

## AlphaLISA<sup>®</sup> SureFire<sup>®</sup> Ultra<sup>™</sup>

# Human and Mouse p-RNA POL II (Ser2) Detection Kit

Product number:

ALSU-PRNAP-A500, ALSU-PRNAP-A10K,

ALSU-PRNAP-A50K, ALSU-PRNAP-A-HV



## Kit specificity:

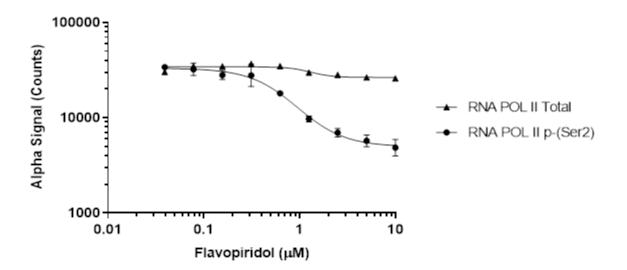
This assay kit contains antibodies which recognize the phospho-Ser2 epitope and a distal epitope on RNA POL II. The protein detected by this kit corresponds to UniProt ID P24928. RNA POL II is also known as DNA-directed RNA polymerase II subunit RPB1 and RNA polymerase II subunit B1. These antibodies recognize RNA POL II of human and mouse origin. Other species should be tested on a case-by-case basis.

#### **Control lysate information:**

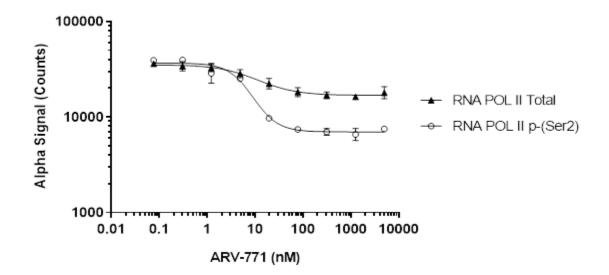
Positive Control Lysate: Prepared from A549 cells, cultured to confluence in T175 flasks in 10% FBS containing medium, lysed with 4 mL of Lysis Buffer.

#### **Representative data**

Data obtained with a 2-plate, 2-incubation protocol. A431 cells were seeded at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with Flavopiridol at the indicated concentrations for 2 hours, lysed with Lysis Buffer and assayed separately for Phospho (Ser2) and Total RNA POL II using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



Data obtained with a 2-plate, 2-incubation protocol. LNCaP cells were seeded at 20K cells/well in a 96 well plate and incubated for 48 hours. Cells were treated with ARV-771 at the indicated concentrations for 20 hours. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser2) and Total RNA POL II using respective *SureFire Ultra* kits. Equivalent to approximately 4,000 cells/datapoint.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. 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"

Revvity, Inc.

940 Winter Street

Waltham, MA 02451 USA

(800) 762-4000 www.revvity.com For a complete listing of our global offices, visit <u>www.revvity.com</u> Copyright ©2023, Revvity, Inc. All rights reserved.



VersionTGRSU531.1