

AlphaLISA™ SureFire® Ultra™

Human and Mouse PAX8 Total Detection Kit

Product number: ALSU-TPAX8-A500, ALSU-TPAX8-A10K,
 ALSU-TPAX8-A50K, ALSU-TPAX8-A-HV



Kit specificity:

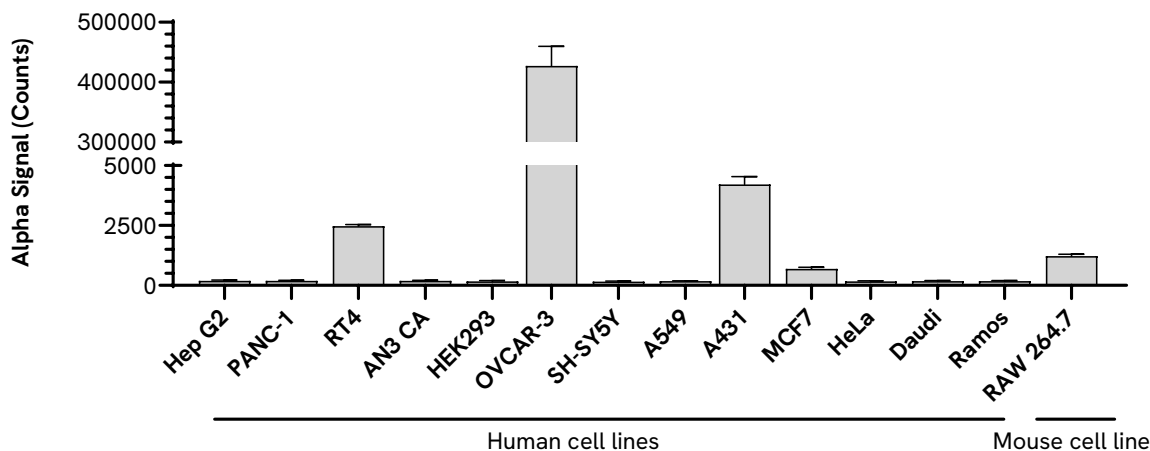
This assay kit contains antibodies which recognize distinct epitopes on PAX8. The protein detected by this kit corresponds to UniProt ID Q02962. PAX8 is also known as Paired box protein Pax-8. These antibodies recognize PAX8 of human and mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from OVCAR-3 cells, cultured to confluence in T175 flasks in 0.1% insulin and 20% FBS containing medium, and lysed with 4 mL of Lysis Buffer.

Representative data:

Data obtained from measurement of PAX8 Total in various cell types. Adherent cell lines were seeded at 40K cells/well in a 96-well plate and incubated overnight. Cells were lysed with 100 µL of Lysis Buffer. Suspension cell lines (Daudi, Ramos) were seeded at 100K cells/well in a 96-well plate in HBSS + 0.1% BSA, cells were spun down and lysed with 100 µL of Lysis Buffer. Suspension cell lysates were further diluted 1:2.5 with Lysis Buffer. Approximately 4,000 cells/datapoint for the various cell lines.



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"