

Research use only. Not for use in diagnostic procedures.

AlphaLISA® SureFire® Ultra™

Human JAK1 Total Detection Kit

Product number: ALSU-TJAK1-A500, ALSU TJAK1-A10K,

ALSU-TJAK1-A50K, ALSU-TJAK1-A-HV



Kit specificity:

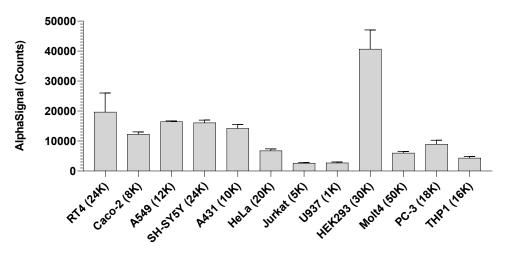
This assay kit contains antibodies which recognize distinct epitopes on JAK1. The protein detected by this kit corresponds to UniProt ID P23458. JAK1 is also known as Janus kinase 1. These antibodies recognize JAK1 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from HEK293 cells grown to confluence in a T175mL flask and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained from measurement of JAK1 Total in various cell types lysed with Lysis Buffer. Approximate number of cells/datapoint is indicated for the various cell lines.



Cell Line (Approximate Number of 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"



