

Research use only. Not for use in diagnostic procedures.

AlphaLISA™ SureFire® Ultra™

Human IRF5 Aggregate Detection Kit

Product number: ALSU-AIRF5-A500, ALSU-AIRF5-A10K,

ALSU-AIRF5-A50K, ALSU-AIRF5-A-HV



Kit specificity:

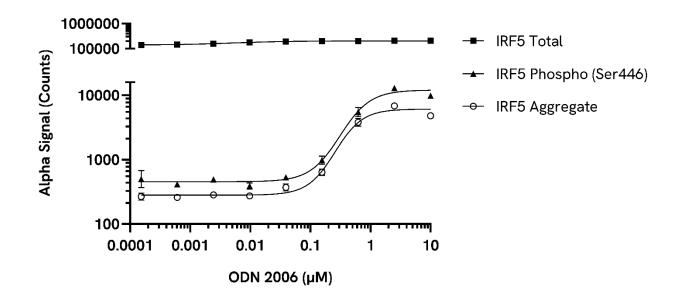
This assay kit contains antibodies which recognize aggregated IRF5. The protein detected by this kit corresponds to UniProt ID Q13568. IRF5 is also known as Interferon regulatory factor 5. These antibodies recognize IRF5 of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

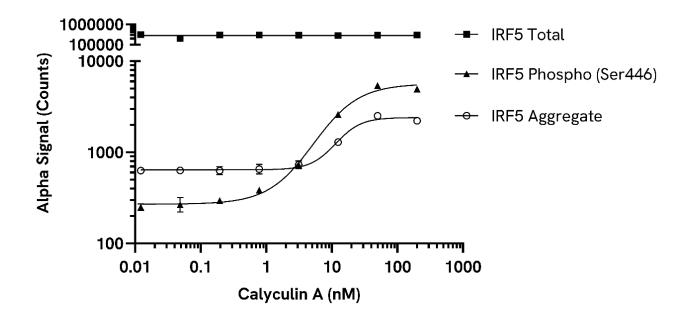
Positive Control Lysate: Prepared from THP-1 cells seeded at 500K cells/mL and incubated overnight in 10% FBS containing medium. Cells were harvested, adjusted to 5×10^6 cells/mL and treated with 100 nM Calyculin A for 3 hours. Following treatment, cells were washed with HBSS + 0.1% BSA and lysed with Lysis Buffer at a final concentration of 10×10^6 cells/mL.

Representative data:

Data obtained using a 2-plate, 2-incubation protocol. RPMI 8226 cells were seeded at 500K cells/well in a 96-well plate. Cells were treated with CpG-B ODN 2006 at the indicated concentrations for 3 hours. Cells were washed with HBSS + 0.1% BSA, then lysed with Lysis Buffer and assayed separately for Aggregate, Phospho (Ser446) and Total IRF5 using respective *SureFire Ultra* assay kits. Equivalent to approximately 50,000 cells/datapoint.



Data obtained using a 2-plate, 2-incubation protocol. THP-1 cells were seeded at 200K cells/well in a 96-well plate. Cells were treated with Calyculin A at the indicated concentrations for 3 hours. Cells were washed with HBSS + 0.1% BSA, then lysed with Lysis Buffer and assayed separately for Aggregate, Phospho (Ser446) and Total IRF5 using respective *SureFire Ultra* assay kits. Equivalent to approximately 50,000 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"



