

AlphaLISA® SureFire® Ultra™

Human p-Estrogen Receptor α (Ser167) Detection Kit

Product number: ALSU-PERA-B500, ALSU-PERA-B10K,
 ALSU-PERA-B50K, ALSU-PERA-B-HV



Kit specificity:

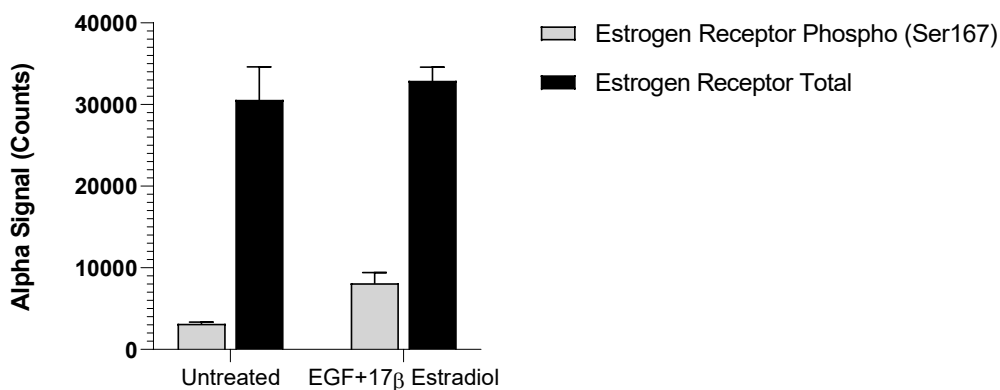
This assay kit contains antibodies which recognize the phospho-Ser167 epitope and a distal epitope on Estrogen Receptor α . The protein detected by this kit corresponds to UniProt ID P03372. Estrogen Receptor is also known as ER-alpha and Estradiol Receptor. These antibodies recognize Estrogen Receptor of human origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from MCF7 cells, cultured to confluence in T175 flasks in 10% FBS containing media, treated with 2 $\mu\text{g}/\text{mL}$ EGF for 10 minutes and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. MCF7 cells were seeded at 10K cells/well in a 96 well plate and incubated for 72 hours. Cells were treated with EGF and 17 β -estradiol at 50 $\mu\text{g}/\text{mL}$ and 0.2nM, respectively for 10 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Ser167) or Total Estrogen Receptor α 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"