

Research use only. Not for use in diagnostic procedures.

AlphaLISA[®] SureFire[®] Ultra[™]

Human Insulin Receptor β Total Detection Kit

Product number:

ALSU-TINR-A500, ALSU-TINR-A10K,

ALSU-TINR-A50K, ALSU-TINR-A-HV



Kit specificity:

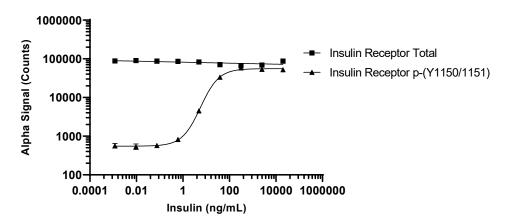
This assay kit contains antibodies which recognize distinct epitopes, on Insulin Receptor β . The protein detected by this kit corresponds to UniProt ID P06213. These antibodies recognize Insulin Receptor β of human origin and are predicted to recognize Insulin Receptor of mouse origin. Other species should be tested on a case-by-case basis.

Control lysate information:

Positive Control Lysate: Prepared from CHO cells transfected with the insulin receptor (CHO-IR cells), cultured to confluence in T175 flasks in 10% FBS containing media, then treated with 2 μ g/mL Insulin and 0.1% BSA for 15 minutes and lysed with 4 mL of Lysis Buffer.

Representative data

Data obtained with a 2-plate, 2-incubation protocol. CHO-IR cells were plated at 40K cells/well in a 96 well plate and incubated overnight. Cells were treated with insulin in serum-free medium at the indicated concentrations for 15 minutes. Cells were lysed with Lysis Buffer and assayed separately for Phospho (Tyr1150/1151) and Total Insulin Receptor β 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.