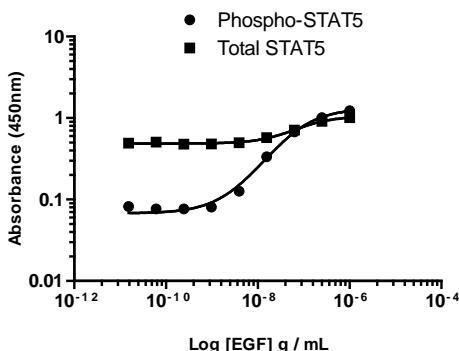


STAT5A/B (Total/Phospho) InstantOne ELISA™ Kit

Catalog Number: 85-86113

Also known as: Signal transducer and activator of transcription 5

RUO: For Research Use Only. Not for use in diagnostic procedures.



Induction of STAT5 A/B (pY694/699) phosphorylation in A431 cells in response to EGF treatment. A431 cells were cultured in 96-well tissue culture plates and treated (15 min) with a dose-range of EGF before cell lysis. Lysates were analysed by InstantOne ELISA™. Data from quadruplicate measurements of STAT5 A/B (pY694/699) are plotted and compared against Total STAT5 A/B levels.

Product Information

Contents: STAT5A/B (Total/Phospho)
InstantOne ELISA™



Catalog Number: 85-86113



Temperature Limitation: Store at 2-8°C.

Batch Code: Refer to vial

Use By: Refer to vial

Description

The InstantOne ELISA™ is specifically engineered for accurate measurement of total and phosphorylated human STAT5A/B (Tyr694/Tyr699) in cell lysates. Detection of mouse and rat STAT5A/B is expected. The InstantOne ELISA™ kit allows for fast analysis of samples in approximately one hour. All reagents used in a traditional sandwich ELISA are added in solution to a plate followed by a wash step and detection with the TMB colorimetric substrate.

STAT5, which exists as STAT5A and STAT5B, is a member of the Signal Transducers and Activators of Transcription family of transcription factors that are activated via phosphorylation by JAK tyrosine kinases in response to various stimuli, including cytokines and growth factors. STAT5A is activated by phosphorylation at Tyr694, and STAT5B is phosphorylated at Tyr699. STAT5A and STAT5B show differential, cell-specific regulation. STAT5A is expressed predominantly in mammary tissue, while STAT5B is more abundant in muscle and liver tissues. The STAT5 transcription factor plays an integral role in immune cell development and regulation, and is an important mediator of IL-2 and IL-15 signaling in regulatory T cells.

Components

This (Total/Phospho) kit includes sufficient reagents to perform 48 wells of the total and 48 wells of the phospho-specific detection for a total of 96 wells.

User Manual

STAT5A/B (Total) Capture Antibody Reagent (Part No. IOSTAT5CAR1M) – 1 x 1.5 mL
STAT5A/B (Total) Detection Antibody Reagent (Part No. IOSTAT5DAR1M) – 1 x 1.5 mL
phospho-STAT5A/B (Tyr694/Tyr699) Capture Antibody Reagent (Part No. IOPSTAT5CAR1M) – 1 x 1.5 mL
phospho-STAT5A/B (Tyr694/Tyr699) Detection Antibody Reagent (Part No. IOPSTAT5DAR1M) – 1 x 1.5 mL
InstantOne ELISA™ Assay Plate (Part No. IOAP96) - 96 well (sealed)
Cell Lysis Buffer (5X) (Part No. IOCLB1) - 12 mL
Enhancer Solution (Part No. IOES1) - 1 mL
Wash Buffer (10X) (Part No. IOWB1) - 15 mL
Detection Reagent (TMB Substrate) (Part No. IODR1) - 12 mL
Stop Solution (Part No. IOSS1) - 12 mL
Positive Control Cell Lysate (lyophilized) (Part No. IOSTAT5PC1) – 0.25 mL
Plate seal - 1 ea

Not for further distribution without written consent. Copyright © 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Tel: 1 800 9556288 • thermofisher.com/askaquestion • thermofisher.com/support

STAT5A/B (Total/Phospho) InstantOne ELISA™ Kit

Catalog Number: 85-86113

Also known as: Signal transducer and activator of transcription 5

RUO: For Research Use Only. Not for use in diagnostic procedures.

Related Products

85-86095 STAT Family Activation InstantOne ELISA™

85-86111 STAT5 A/B (Total) InstantOne ELISA™

85-86112 Phospho-STAT5A/B (Tyr694/Tyr699) InstantOne ELISA™

IOCLB InstantOne ELISA Cell Lysis Buffer (5X) – 100 mL

IOES InstantOne ELISA Enhancer Solution – 100 mL

IOWB InstantOne ELISA Wash Buffer (10X) – 500 mL