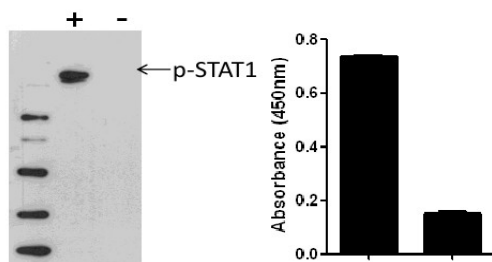


phospho-STAT1 (Tyr701) InstantOne ELISA™ Kit

Catalog Number: 85-86092

Also known as: Signal transducer and activator of transcription 1-alpha/beta

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



A431 cells were either unstimulated (-) or treated with EGF (1 ug/mL for 10 min) (+). Lysates were prepared as described in the InstantOne™ User Manual. Phosphorylated STAT1 (Tyr701) was analyzed in cell lysates by immunoblotting (left) and InstantOne ELISA™ (right) 5 ug/analysis. The intensity of the band as detected by immunoblot correlates with InstantOne ELISA™ absorption values.

Product Information

Contents: phospho-STAT1 (Tyr701)
InstantOne ELISA™



Catalog Number: 85-86092

Clone: ELISA



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 phosphorylated human STAT1 (Tyr701) in cell lysates. The InstantOne ELISA™ kit allows for fast analysis of samples in approximately two hours. 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.

STAT1 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. More specifically, numerous studies have demonstrated the involvement of STAT1 in controlling gene expression in response to interferon (IFN) stimulation. Depending on the stimulus, tyrosine phosphorylation leads to STAT1 homodimerization (e.g., IFN γ) or heterodimerization with STAT2 (e.g., IFN α and IFN β) via its SH2 domains. After dimerization, STAT1 translocates into the nucleus to regulate transcription. Research shows that STAT1 binds interferon-stimulated response elements within the promoters of genes involved in inducing the cellular antiviral response.

Components

User Manual

phospho-STAT1 (Tyr701) Capture Antibody Reagent (Part No. IOPSTAT1CAR1) - 3 mL

phospho-STAT1 (Tyr701) Detection Antibody Reagent (Part No. IOPSTAT1DAR1) - 3 mL

InstantOne™ Assay Plate (Part No. IOAP96) - 96 well (sealed)

Cell Lysis Buffer (5X) (Part No. IOCLB1) - 10 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. IOSTAT1PC1) - 0.25 mL

Plate seal - 1 ea

Related Products

85-86095 STAT Family Activation 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

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