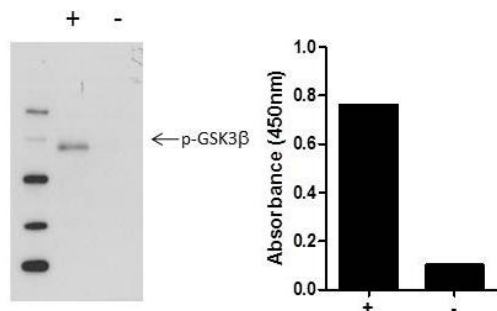


phospho-GSK3 beta (Ser9) ELISA Kit

Catalog Number: 85-86172

Also known as: Glycogen synthase kinase-3 beta

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



MCF7 cells were treated with either the PI3 Kinase inhibitor Wortmannin (10uM) (-) or insulin (10ng/mL for 10 min.) (+). Cell lysates were prepared as described in the InstantOne ELISA User Manual. The cell lysates were analyzed by immunoblotting (left) and phospho-GSK3 beta (Ser9) InstantOne™ ELISA (right). The intensity of the band at 50 kDa as detected by immunoblot correlates with InstantOne™ ELISA absorption values.

Product Information

Contents: phospho-GSK3 beta (Ser9) ELISA Kit



Catalog Number: 85-86172

Clone: ELISA



Batch Code: Refer to vial



Use By: Refer to vial

Description

The InstantOne™ ELISA is specifically engineered for accurate measurement of phosphorylated human and mouse GSK3b (Ser9) in cell lysates. Detection of phosphorylated rat GSK3b (Ser9) is also 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.

Glycogen Synthase Kinase 3 (GSK3) is a Ser/Thr kinase that exists as two isoforms, GSK3a and GSK3b. Unlike other kinases, phosphorylation at Ser9 and Ser21 for GSK3a and GSK3b, respectively, leads to inactivation of this enzyme. GSK3 has been implicated in several diseases, including cancer and diabetes, through its phosphorylation of glycogen synthase. GSK3 has also been associated with neurodegenerative diseases such as Alzheimer's disease due to its phosphorylation of the neuronal Tau protein.

Components

The kit includes sufficient reagents to perform 96 wells of the kit specific target.

User Manual

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

Cell Lysis Buffer (5X) (Part No. IOCLB1) - 8 mL

Enhancer Solution (Part No. IOES1) - 0.5 mL

Wash Buffer (10X) (Part No. IOWB1) - 8 mL

phospho-GSK3b (Ser9) Capture Antibody Reagent (Part No. IOPGSKCAR1) - 3 mL

phospho-GSK3b (Ser9) Detection Antibody Reagent (Part No. IOPGSKDAR1) - 3 mL

Positive Control Cell Lysate (lyophilized) (Part No. IOGSKPC1) - 1 ea

Detection Reagent (TMB Substrate) (Part No. IODR1) - 10 mL

Stop Solution (Part No. IOSS1) - 10 mL

Plate Seal - 1 ea

Applications Reported

This ELISA antibody has been reported for use in ELISA.

Applications Tested

This assay has been validated for the detection of total human GSK3b in stimulated MCF7 cell lysates. This assay was evaluated for specificity by immunoblot analysis using the same lysates and antibody (data shown above).

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience • info@ebioscience.com

phospho-GSK3 beta (Ser9) ELISA Kit

Catalog Number: 85-86172

Also known as: Glycogen synthase kinase-3 beta

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

Related Products

85-86171 GSK3 beta (Total) ELISA Kit (ELISA)

85-86173 GSK3 beta (Total/Phospho) ELISA Kit (ELISA)

IOCLB InstantOne ELISA Cell Lysis Buffer (5X)

IOES InstantOne ELISA Enhancer Solution

IOWB InstantOne ELISA Wash Buffer(10X)

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience •
info@ebioscience.com