



ClariCELL™ Mouse ITK Kinase Assay Service

4-Step Assay Validation

Description



The **ClariCELL™ Mouse ITK Kinase Assay** quantifies auto-phosphorylation of mouse full-length ITK in human cells. The assay is useful to determine potencies of small-molecule inhibitors against the specified kinase in the context of a cellular environment. Compound testing services are available utilizing the assay.

Overview

Human Embryonic Kidney (HEK 293) cells transiently expressing sequence verified mouse full-length ITK are exposed to test compound or control, then lysed to release cellular proteins. ITK is captured onto an assay plate, and the extent of autophosphorylation is quantified by ELISA using an antibody specific for the phosphorylation event. Cells expressing kinase deficient mouse ITK [K390M] are also utilized as controls to calculate the % inhibition of test compounds.

サービスの詳細、最新情報はHPをご覧ください

カルナバイオサイエンス株式会社

〒650-0047 神戸市中央区港島南町1-5-5 BMA3F

TEL: 078-302-7091 (営業部直通) / FAX: 078-302-7086

E-mail: info@carnabio.com

Mouse ITK Expression in Cells

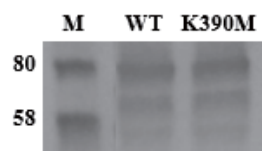


Figure 1: Wild type (wt) or kinase dead (K390M) mouse ITK was expressed transiently in 293 cells. Following cell lysis, an IP Western was performed with appropriate antibodies to capture and detect total ITK protein.

Mouse ITK Autophosphorylation in Cells

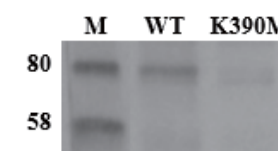


Figure 2: Wild type (wt) or kinase dead (K390M) mouse ITK was expressed transiently in 293 cells. Following cell lysis, an IP Western was performed with appropriate antibodies to capture and detect phospho-ITK protein.

Quantification of Phosphorylation

ELISA Activity

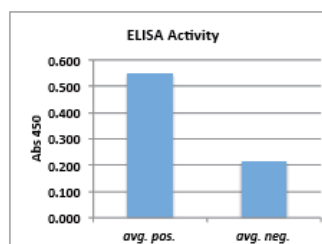


Figure 3: Wild type (wt) or kinase dead (K390M) mouse ITK was expressed transiently in 293 cells. Following cell lysis, an ELISA was performed to quantify the extent of autophosphorylation of ITK.

Reference Inhibitor Data

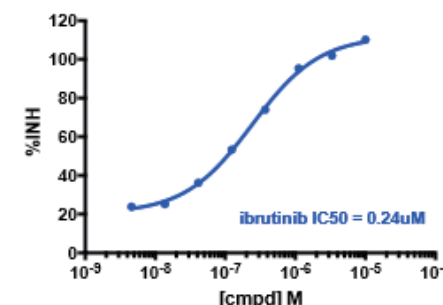


Figure 4: An autophosphorylation assay was performed in the presence of ibrutinib, an ITK inhibitor. Percent inhibition data were plotted to determine the IC₅₀s.