

Anti-phospho-MET (Tyr1349 Tyr1356)

Rabbit Polyclonal Antibody

Catalog # **M52-65R**

Lot # B3216-35

Cited Applications

ELISA, WB

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the human c-MET phosphorylated at Tyrosine 1349 and 1356

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to residues surrounding Tyr1349 and Tyr1356 of human c-MET protein.

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

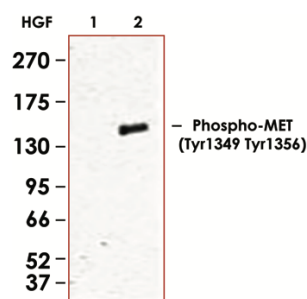
Scientific Background

c-MET is a receptor tyrosine kinase that binds hepatocyte growth factor (also known as scatter factor, HGF/SF). Stimulation by HGF causes c-MET to autophosphorylate at multiple tyrosines, including Y1234 and Y1235. Activated c-MET binds to various substrates, such as GAB1, GRB2, phosphatidylinositol 3-kinase (PI3K) and others. GAB1 is scaffold protein, which enables c-MET to elicit downstream signaling cascades, which include the well-characterized ERK/MAPK, PI3K-Akt/PKB, Crk-Rap and Rac-Pak pathways; leading to cell proliferation and cell survival.

References:

1. Birchmeier, C., et al.: Met, metastasis, motility and more. Nat. Rev. Mol. Cell Bio. 2003; 4; 915-925.
2. Fan, S., et al.: The multisubstrate adapter Gab1 regulates hepatocyte growth factor (scatter factor)-c-Met signaling for cell survival and DNA repair. Mol. Cell Biol. 2001; 21; 4968-84.
3. Longati, P., et al. Receptor tyrosine kinases as therapeutic targets: the model of the MET oncogene. Curr. Drug Targets. 2001; 2; 41-55.

Sample Data



Western Blot using Anti-phospho-MET (Tyr1349 Tyr1356) antibody shows detection of phosphorylated c-Met in Human mammary B5/589 epithelial cells serum-deprived and treated with or without HGF.

Anti-Phospho-MET (Tyr1349 Tyr1356)

Rabbit Polyclonal Antibody

Catalog #	M52-65R
Lot #	B3216-35
Purification	Immunoaffinity chromatography
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.