

Anti-Phospho-Rel A (ser276)

Rabbit Polyclonal Antibody

Catalog # R29-365R

Lot # B3216-59

Cited Applications

ELISA, IHC, WB

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

Specificity

Recognizes the human p65 protein phosphorylated at Serine 276

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, Antiserum

Immunogen

The antibody was produced against synthesized peptide corresponding to a region near phosphorylated Ser276 of the protein conjugated to KLH

Formulation

0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

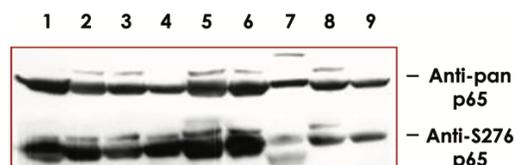
Scientific Background

Transcription factors of the nuclear factor κB (NF-κB)/Rel family play a pivotal role in inflammatory and immune responses. In unstimulated cells, NF-κB is sequestered in the cytoplasm by its inhibitory proteins, the IκBs (1). NF-κB-activating agents can induce the phosphorylation of IκB's, targeting them for rapid degradation through an ubiquitin-proteasome pathway, releasing NF-κB to enter the nucleus, where it regulates gene expression.

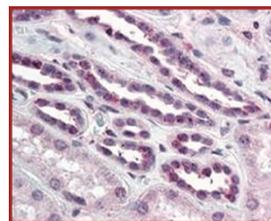
References:

1. Baeuerle, P. et al: Function and activation of NF-kappa B in the immune system. Annu Rev Immunol. 1994;12:141-79.

Sample Data



Western Blot of total protein lysates from head and neck tumors using anti-Phospho-Rel A (Ser276) antibody. Lanes 1-6: squamous carcinoma cells. Lane 7: primary culture of human keratinocytes. Lane 8: ATCC SCC9 cells (also a head and neck squamous carcinoma). Lane 9: EGF-induced human derived A431 cells



Anti-pS529 shows phospho p65 staining in carcinoma cells.

Anti-phospho-Rel A (Ser276)

Rabbit Polyclonal Antibody

Catalog #	R29-365R
Lot #	B3216-59
Purification	Delipidation and defibrination
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.