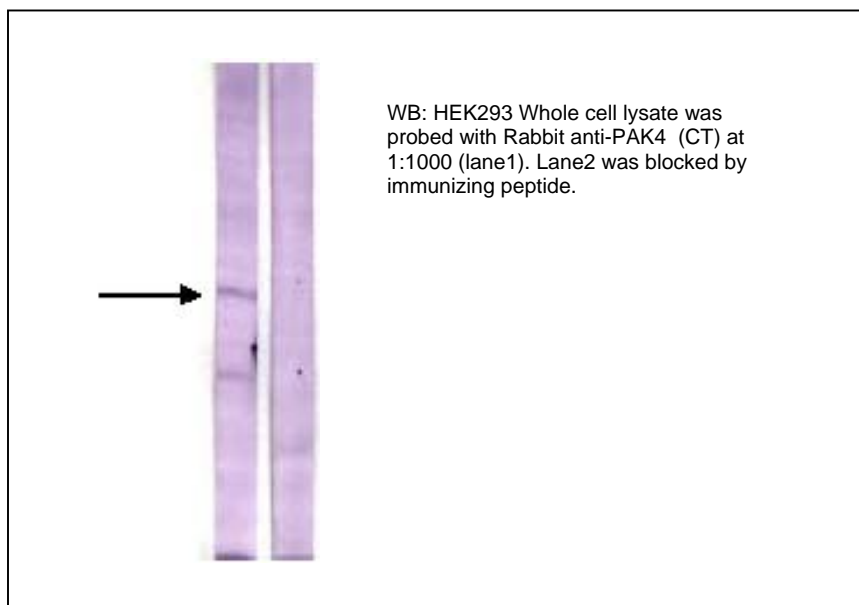




Product Data Sheet

Product Name:	Anti - PAK4 (CT)
Catalog Number:	29663
Lot Number:	See label on vial
Product Description:	The antibody is supplied as an epitope affinity purified rabbit IgG, 50 µg in 250 µl (0.2 mg/ml) of 1X PBS (pH 7.4) containing 0.05% Sodium Azide.
Immunogen:	Rabbit anti-PAK4 (CT) polyclonal antibody was raised against a synthetic peptide corresponding to the amino acids 470-485 of human PAK4.
Species Reactivity:	Species reactivity includes human, while others remain unknown.
Application Notes:	The following are recommended starting points for different applications: WB: 0.5 to 2 µg/ml IP: 3.0-5.0 µg/extract from 10 ⁷ cells



Background:

P21-activated kinases (PAKs) belong to the family of serine/threonine kinases involved in the control of various cellular processes, including the cell cycle, dynamics of the cytoskeleton, apoptosis, oncogenic transformation, and transcription. All PAK family members are characterized by the presence of p21-binding domain. p21-activated kinases are regulated by the small GTP-binding proteins Rac and Cdc42, and lipids, which stimulate autophosphorylation and phosphorylation of exogenous substrates. Serine (Ser-474) is the likely autophosphorylation site in the kinase domain of PAK4 *in vivo*. Phosphospecific antibodies directed against serine 474 detect activated PAK4 on the Golgi membrane when PAK4 is co-expressed with activated Cdc42. Current data strongly implicates PAK-4 in oncogenesis. PAK4 is frequently overexpressed in human tumor cell lines of various tissue origins.

Storage:

Store at 2-8 °C for up to one year. Avoid repeated freezing and thawing.

References:

1. Qu, J. et al. *Mol Cell Biol* **23**, 7122 (2003).
2. Callow, M. et al. *J Biol Chem* **277**, 550 (2002).
3. Zenke, F. et al *J Biol Chem* **274**, 32565 (1999).
4. Arie, A. et al. *The EMBO Journal* **17**, 6527 (1998).
5. Sells, M. et al. *Trends in Cell Biol* **7**, 162 (1997).

This product is for *in vitro* research use only.