

IEX-1 antibody (pAb)

Rabbit Anti-Human/Mouse IEX-1

Instruction Manual

Catalog Number	PK-AB718-4247
Synonyms	IEX-1 Antibody: Immediate early response 3, IER3, DIF2, PRG1, GLY96
Description	IEX-1 is a stress inducible gene that is induced by ionizing radiation, ultraviolet radiation, and a variety of growth factors, i.e., FAS and TNF- α . IEX-1 is widely expressed in epithelial and endocrine tissues, as well as in vascular endothelium. It plays an important role in the regulation of cellular growth, cell death and oncogenesis. IEX-1 is precisely regulated by multiple transcription factors such as p53, NF-kappaB/rel, Sp1 and c-Myc, to ensure rapid and transient expression of IEX-1 in cells under a variety of stress conditions. IEX-1 is expressed as both a longer form (IEX1L) and a splice variant, designated IEX1S. It is localized to the nucleus and perinuclear region. Overexpression of IEX-1 facilitates apoptosis and cell cycle progression, whereas disruption of IEX-1 expression is associated with decreases in both apoptosis and cell cycle progression.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	IEX1 antibody was raised in rabbits against a 24 amino acid peptide near the center of human IEX1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 μ l or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. IEX1 antibody can be used for detection of IEX1 by Western blot at 2 - 4 μ g/mL. Antibody can also be used for immunohistochemistry starting at 2.5 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	Kondratyev AD, Chung K-N and Jung MO. Identification and characterization of a radiation-inducible glycosylated human early-response gene. <i>Cancer Res.</i> 1996; 56:1498-502. Feldmann KA, Piddelkow MR, Roche PC, et al. Expression of an immediate early gene, IEX-1, in human tissues. <i>Histochem. Cell Biol.</i> 2001; 115:489-97. Wu MX, Zhaohui A, Prasad KVS, et al. IEX-1L, an apoptosis inhibitor involved in NFkappaB mediated cell survival. <i>Science</i> 1998; 281:998-1001. Arlt A, Grobe O, Sieke A, et al. Expression of the NF-kappa B target gene IEX-1 (p22/PRG1) does not prevent cell death but instead triggers apoptosis in Hela cells. <i>Oncogene</i> 2001; 20:69-76.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4247P; IEX-1 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.