

IFN-beta (IN) antibody (pAb)

Rabbit Anti-Human/Mouse IFN-beta (IN)

Instruction Manual

Catalog Number	PK-AB718-4243
Synonyms	IFN-beta Antibody; Interferon beta, beta-interferon
Description	Type I Interferons (IFN-a/b) are produced primarily in response to viral infection by "Natural IFN-producing cells" (NPCs) as part of the host immune response and can also inhibit the development of tumors. IFN-b binding by its receptor results in the activation of the tyrosine kinases Jak1 and Tyk2 and phosphorylation of members of the STAT family of transcription factors, leading to the transcription and expression of the immune response genes. More recently, several members of the toll-like receptor (TLR) family were found to stimulate the production IFN-b. IFN-b is currently used clinically for treatment of tumors, infections and multiple sclerosis.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	IFN-beta antibody was raised in rabbits against an 17 amino acid peptide from near the center of human IFN-b.
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 µl 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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. IFN-b antibody can be used for detection of IFN-b by Western blot at 5 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Gresser I. Wherefore interferon? J. Leuk. Biol. 1997; 61:567-74. Colonna M. TLR pathways and IFN-regulatory factors: to each its own. Eur. J. Immunol. 2007; 37:306-9. Kock M, Mostert J, De Keyser J, et al. Interferon-beta treatment and the natural history of relapsing-remitting multiple sclerosis. Ann. Neurol. 2007; epub.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4243P; IFN-beta Peptide Cat.No. PK-AB718-1288; A20 Cell Lysate

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