IFN-β polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

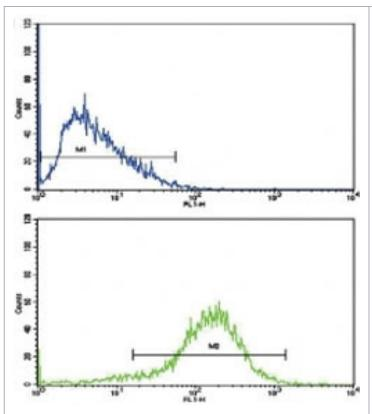
Order Online »

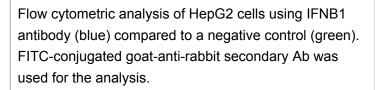
ENZ-ABS901-0400

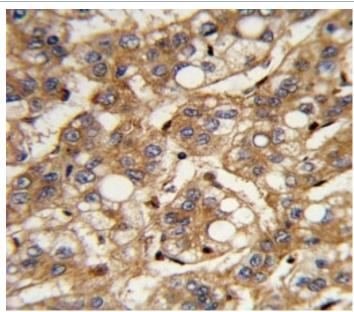
0.4ml

Manuals, SDS & CofA

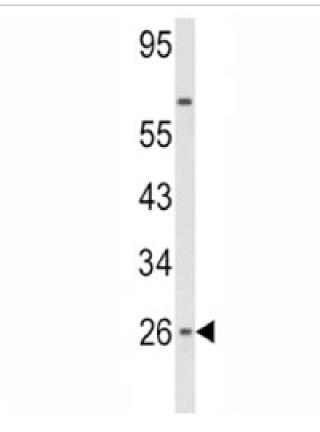
View Online »







IHC analysis of FFPE human hepatocarcinoma with IFNB1 antibody



Western blot analysis of IFNB1 antibody and HepG2 lysate.

Handling & Storage

Handling Aliquot the Interferon beta antibody and store frozen at -20°C or colder. Avoid

freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interferon-β

Application ELISA, Flow Cytometry, IHC, WB

Formulation Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.

Host Rabbit

Immunogen Human IFN-β (aa 39-66).

Isotype Ig

Purity Detail Antigen affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:10-1:50)Immunohistochemistry (paraffin

sections) (1:10-1:50)Western blot (1:1,000)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P01574

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com