IFN-γ recombinant monoclonal antibody (IFNG/3996R)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

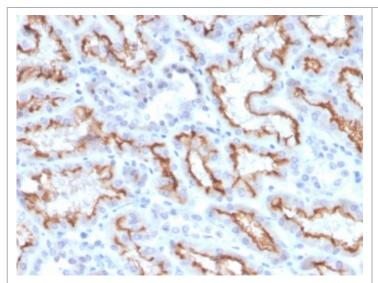
Order Online »

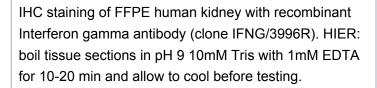
ENZ-ABS902-0100

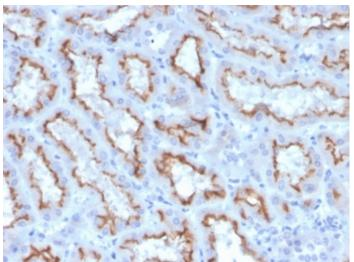
100µg

Manuals, SDS & CofA

View Online »







IHC staining of FFPE human kidney with recombinant Interferon gamma antibody (clone IFNG/3996R). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 10-20 min and allow to cool before testing.

Handling & Storage

Handling Store the recombinant Interferon gamma antibody at 2-8°C (with azide) or aliquot and

store at -20°C or colder (without azide).

-20°C **Long Term Storage**

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interferon-y

Application IHC (PS)

Clone IFNG/3996R

Formulation Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA

(US sourced) and 0.05% sodium azide.

Host Mouse

Purified recombinant human IFN-y. **Immunogen**

Isotype lgGκ

Purity Detail Protein A affinity chromatography purified.

Recommendation Dilutions/Conditions Immunohistochemistry (FFPE) (1-2µg/ml)Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P01579

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

Last modified: July 12, 2024



info-

eu@enzolifesciences.com