## IL-5 monoclonal antibody (IL5/4161)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

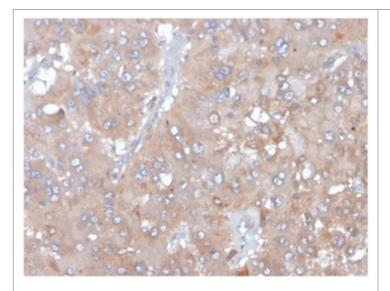
Order Online »

ENZ-ABS903-0100

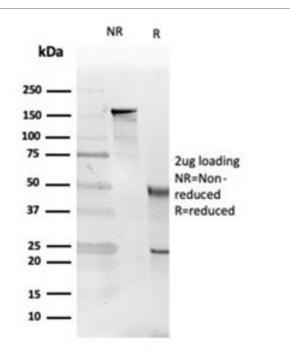
100µg

Manuals, SDS & CofA

**View Online** »



IHC staining of FFPE human adrenal gland tissue with Interleukin 5 antibody (clone IL5/4161). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



SDS-PAGE analysis of purified, BSA-free Interleukin 5 antibody (clone IL5/4161) as confirmation of integrity and purity.

## **Handling & Storage**

Handling Aliquot the Interleukin 5 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 Interleukin-5

Application IHC (PS)

Clone IL5/4161

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

(US sourced), 0.05% sodium azide.

**Host** Mouse

Immunogen Recombinant fragment of human IL-5 protein

corresponding to the mature protein sequence.

Isotype IgG2bk

Purity Detail Protein A/G affinity 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 P05113

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

-

Last modified: July 12, 2024



info-

eu@enzolifesciences.com