

EPX Polyclonal Antibody

(Catalog # A69483)

Background

Mediates tyrosine nitration of secondary granule proteins in mature resting eosinophils. Shows significant inhibitory activity towards *Mycobacterium tuberculosis* H37Rv by inducing bacterial fragmentation and lysis.

Description

EPX Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P11678

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Eosinophil peroxidase protein (140-300AA)

Storage

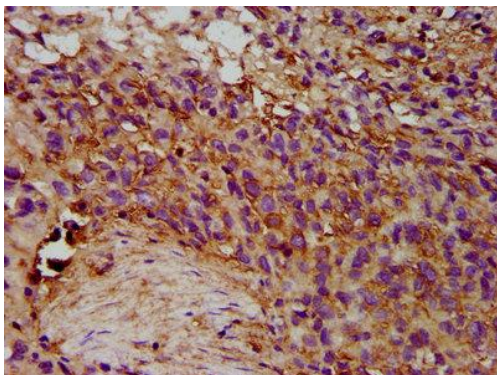
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

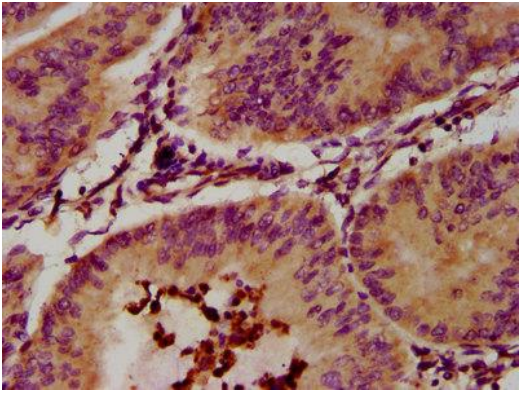
Eosinophil peroxidase, EPO, Eosinophil peroxidase light chain, Eosinophil peroxidase heavy chain, EPX, EPER, EPO, EPP

Application

ELISA, IHC; Recommended dilution: IHC: 1:200-1:500



IHC image of EPX Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human ovarian cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of EPX Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human endometrial cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.