

Cxcl11 Polyclonal Antibody

(Catalog # A54097)

Description

Cxcl11 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

Q9JHH5

Purification

Protein G purified

Immunogen

Recombinant Mouse C-X-C motif chemokine 11 protein (22-98AA)

Storage

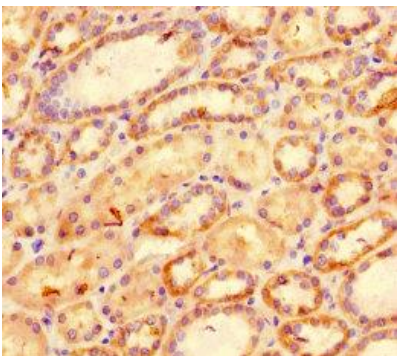
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Scyb11, Cxcl11, Interferon-inducible T-cell alpha chemoattractant, I-TAC, Small-inducible cytokine B11

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human kidney tissue at dilution of 1:100