

[pTyr³⁹⁷]Focal adhesion kinase polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	FAK
Application	ICC, IHC, WB
Formulation	Liquid. In PBS, pH 7.4. with 50% glycerol, 1 mg/mL BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of FAK (focal adhesion kinase) phosphorylated at Tyr ³⁹⁷ . The sequence is conserved in mouse, rat, chicken and frog.
Isotype	IgG
Purity Detail	Purified by antigen affinity chromatography.
Recommendation Dilutions/Conditions	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	Chicken, Human, Mouse, Xenopus
Specificity	Recognizes FAK phosphorylated at Tyr ³⁹⁷ .
UniProt ID	Q05397
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee