[pTyr³⁹⁷]Focal adhesion kinase polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

BML-SA433-0020

20µl

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 Tyr397. 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 Tyr³⁹⁷.

UniProt ID Q05397

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

.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com