

[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 Tyr ³⁹⁷ .
UniProt ID	Q05397
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES,
INC.
Phone: +1 949 946 9120
info@enzolifesciences.com
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

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

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

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