

FLAG polyclonal antibody

FLAG-tag stands for an octapeptide with a sequence of DYKDDDDK. It is a polypeptide protein tag that can be added to a protein using recombinant DNA technology. It is widely used for affinity chromatography to purify recombinant proteins from variety of systems. It can also be used in the isolation of protein complexes with multiple subunits. A FLAG-tag can be used in many different assays that require recognition by an antibody. If there is no antibody against the studied protein, adding a FLAG-tag to this protein allows one to follow the protein with an antibody against the FLAG sequence, such as in ELISA, Western Blot, etc.

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

Ordering Information

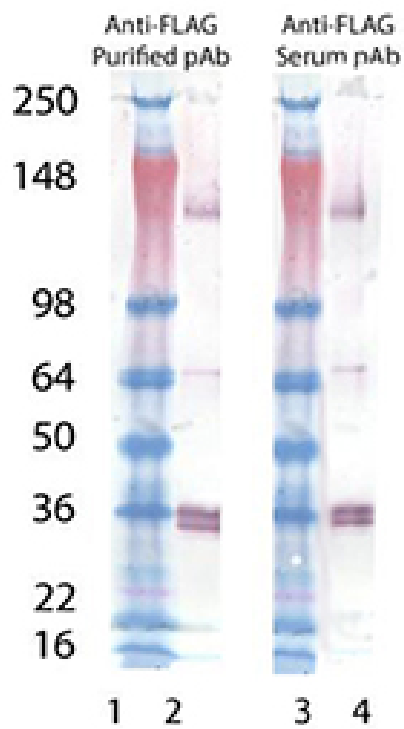
[Order Online »](#)

ADI-SAB-410-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

ADI-SAB-410_WB



Western blot analysis of anti-FLAG, pAb (Prod. No. ADI-SAB-410): Lane 1: MW marker, Lane 2: SuperFasLigand™ (soluble) (human), (recombinant) (Prod. No. ALX-522-020), Lane 3: MW marker, Lane 4: SuperFasLigand™ (soluble) (human), (recombinant) (Prod. No. ALX-522-020).

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, WB
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide (CDYKDDDDK) coupled to KLH.
Isotype	IgG
Purity Detail	Immunoaffinity purified.
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.
Source	Purified from rabbit serum.
Species Reactivity	Species independent
Specificity	Recognizes FLAG-tagged proteins.
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co. LLC
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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