

α -Fetoprotein Recombinant monoclonal antibody (AF5) (Rabbit IgG κ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS608-0200	200 μ g
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AFP
Clone	AF5
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	AF5 was prepared by immunizing female BALB/c mice with human amniotic fluid containing AFP, from a pregnancy affected by a neural tube defect, diluted and emulsified in Freund's complete adjuvant.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Technical Info / Product Notes	This chimeric rabbit antibody was made using the variable domain sequences of the original murine IgG1 format, for improved compatibility with existing reagents, assays and techniques.
UniProt ID	P02771



Worry-Free Guarantee

ENZO LIFE SCIENCES

INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee

European Sales Office

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

Belgium, The Netherlands & Luxembourg

Phone: +32 3 466 0420
info-be@enzolifesciences.com

France

Phone: +33 1 72 440 1335
info-fr@enzolifesciences.com

Germany

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

UK & Ireland

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