

Anti-Pig Interferon Alpha, Clone K9 (MAb)

Binds to and neutralizes pig IFN alpha

Ordering Information

[Order Online »](#)

PBL-27100-1	500µg
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	After opening, prepare aliquots and store at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	IFN- α , Interferon- α
Clone	K9
Formulation	Liquid. 25 mM Tris; 75 mM Sodium Chloride; 50 % glycerol, pH 8.0.
Host	Mouse
Immunogen	Pig IFN- α (interferon- α).
Isotype	IgG1
Species Reactivity	Porcine



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

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: +33 472 440 655
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