

# Glucagon like peptide 1 polyclonal antibody

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

Citations: 1

[View Online »](#)

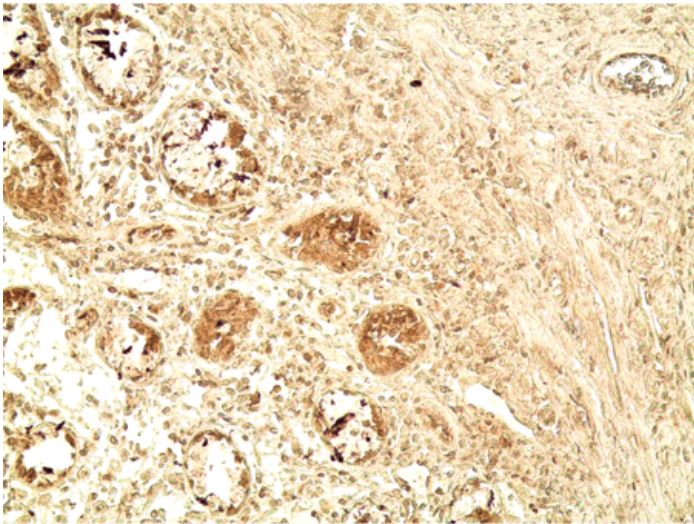
## Ordering Information

[Order Online »](#)

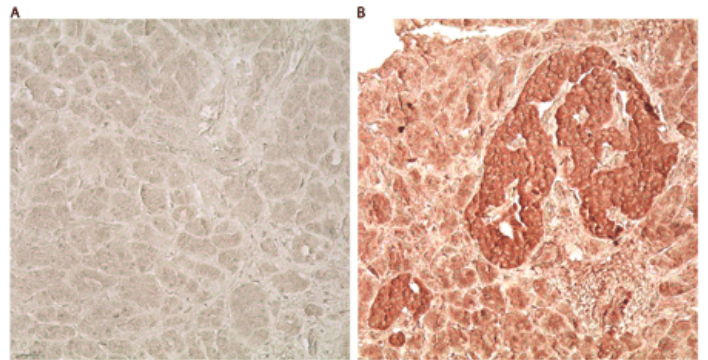
BML-GA1176-0100	100µl
-----------------	-------

Manuals, SDS & CofA

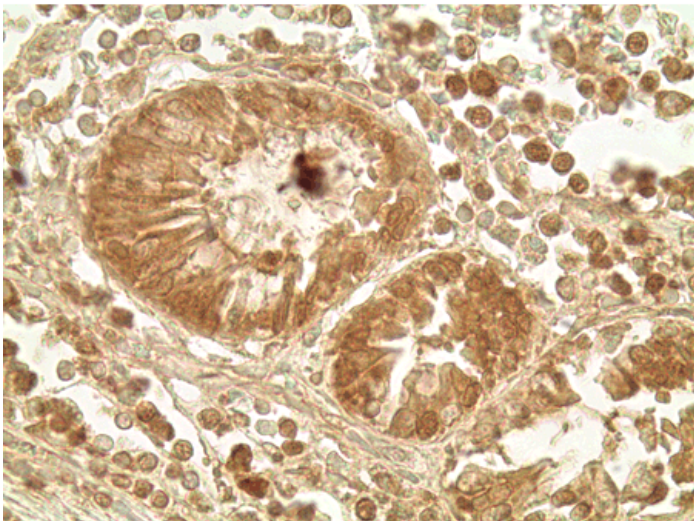
[View Online »](#)



Immunohistochemistry analysis of human colon tumor tissue using Glucagon like peptide 1 pAb, (Prod. No. BML-GA1176) at a dilution of 1:500.



Immunohistochemistry analysis of human pancreas tissue using Glucagon like peptide 1 pAb, (Prod. No. BML-GA1176) at a dilution of 1:1000. A: Negative control: normal rabbit serum, B: GLP 1, rabbit pAb (Prod. No. BML-GA1176).



Immunohistochemistry analysis of human colon tumor tissue using Glucagon like peptide 1 pAb, (Prod. No. BML-GA1176) at a dilution of 1:500.

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

Alternative Name	GLP 1
Application	IHC
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 1-19 of human GLP-1.
Source	Purified from rabbit serum.
Species Reactivity	Mammalian
UniProt ID	P01275
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-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)