

Glucagon like peptide 1 polyclonal antibody

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

Citations: 1

[View Online »](#)

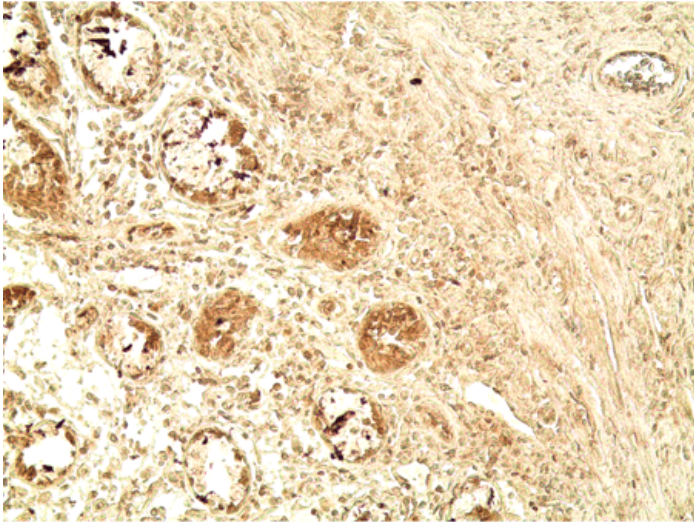
Ordering Information

[Order Online »](#)

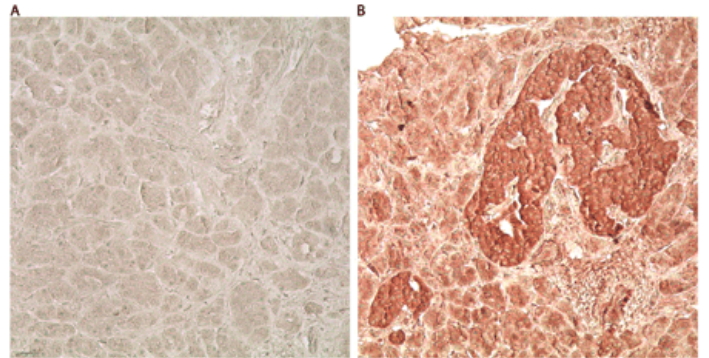
BML-GA1176-0100	100µl
-----------------	-------

Manuals, SDS & CofA

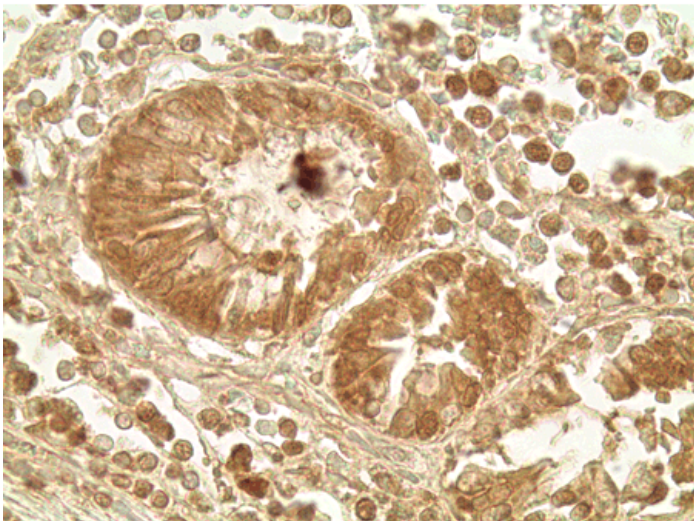
[View Online »](#)



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 [Worry-Free Guarantee](#).



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)